

**CKIP SERVICES**

**Version 10MY1**

# **CUSTOM SERVICES**

## **KINASE-INHIBITOR PROFILING CUSTOMER INFORMATION PACKAGE**

**Toll free: 1-866-KINEXUS or 604-323-2547**

**Facsimile: 604-323-2548**

**E-mail: [info@kinexus.ca](mailto:info@kinexus.ca)**

**[www.kinexus.ca](http://www.kinexus.ca)**



## TABLE OF CONTENTS

<b>Sample Preparation</b>	<b>PDF Page No.</b>
1. Introduction .....	3
2. Quantity of compound required .....	6
3. Kinase assay conditions .....	6
<b>Shipping &amp; Pricing</b>	
4. Shipping information .....	6
5. Pricing information .....	7
<b>Description of Follow Up Services</b>	
6. Follow up services .....	7
7. Forms to be completed .....	9
<b>Kinex™ Antibody Microarray Screening Services</b>	
8. Appendix A – List of available active protein kinases for compound profiling .....	13
<b>Forms to Complete and Return with your Samples</b>	
9. Service Order Form (KICP-SOF-01) .....	40
10. Service Identification Forms (KICP-SIF-01A and KICP-SIF-01B) .....	41
11. Confidential Sample Description Form (KICP-CSDF-01) .....	43
12. Commercial Invoice – No ice (required for customers outside of Canada) .....	44
13. Commercial Invoice – Dry ice (required for customers outside of Canada) .....	45
14. Kinexus Service Agreement (first time KICP customers only) .....	46

# KINASE INHIBITOR COMPOUND PROFILING SERVICE

## 1. INTRODUCTION

The Kinase Inhibitor Compound Profiling (KICP) Service is a convenient and cost-effective solution to assist scientists in ascertaining of the specificity of lead compounds and their mechanisms of action for drug discovery. This information is critical for the selection of better drug candidates for clinical testing. Approximately a third of all pharmaceutical R&D is now focused on protein kinases as drug targets. Kinexus currently has over 265 human protein kinases available for screening with our KICP Service. This number will continue to increase in the near future. At least 515 human protein kinases target the phosphorylation of apparently more than 500,000 phosphorylation sites in the proteome. In view of this, it is critical to establish the specificity of any kinase drug candidate for clinical studies. The more specific the kinase inhibitors, the lower the chances of off targets that could compromise on the utility of the drug from toxicity and other undesired side-effects. This service presently relies on the use of gamma phosphate-radiolabeled ATP to phosphorylate peptide and recombinants protein substrates with purified and active preparations of human protein kinases. Kinexus performs the KICP Service under strict confidentiality, and all materials, information and results are used as directed by the client. We recommend the KICP Service as a useful follow on from our Kinex™ kinase microarray services, which is an alternative method to counter screen for kinase drug specificity by assaying the ability of test compounds to prevent the binding of an ATP analogue probe to over 150 human protein kinases. These *in vitro* services from Kinexus compliments the *in vivo* compound profiling analyses that are offered with our Kinex™ antibody microarray and Kinetworks™ multi-immunoblotting services. With these proteomics services, it is feasible to determine the effects of compounds and other treatments of animals and isolated cells on the expression levels of over 193 known protein kinases and 174 other proteins as well as the status of at least 270 different phospho-sites to monitor indirectly the activities of their upstream kinases.

The KICP Service uses the most reliable direct assay of protein kinase phosphotransferase activity that is known. The methodology is based on the direct quantification of radio-labeled phosphate from ATP (gamma-<sup>32</sup>P or gamma-<sup>33</sup>P) on to a peptide or protein substrate of a target protein kinase. This provides for a high signal to noise detection of phosphorylation, high reproducibility, and reduces the opportunity for artifacts inherent in other methods, such as the measurement of the production of ADP or disappearance of ATP. Furthermore, the assay provides a direct measure of the effect of a compound on the enzymatic phosphotransferase activity of a target protein kinase, rather than a measure of the ability of a compound to bind near the active site of the kinase, as is performed with some other approaches to compound screening.

The preparations of recombinant protein kinases that we use in our KICP Service possess high specific activities, and generally represent full-length human clones. In some instances, we use kinases that feature activating mutations that may occur *in vivo*. But generally, the kinases are activated by endogenous phosphorylation in the baculovirus-infected insect cells or by the addition of the purified and activate upstream protein kinase. For each kinase used in the KICP Service, the assay conditions have been carefully optimized to ensure the highest levels of phosphotransferase activity. An abbreviated list of the protein kinases that are presently available with our KICP Service is provided on the next two pages in Table 1, and more information is provided in Appendix A. You can also download an MS-Excel spreadsheet with very detailed information on each protein kinase along with active hyperlinks to other websites from our website at [http://www.kinexus.ca/ourServices/inhibitor\\_profiling/index.html](http://www.kinexus.ca/ourServices/inhibitor_profiling/index.html).

Table 1. Active Human Protein Kinases Available in the KICP Service.

Protein Kinase Name	ID Code	Protein Kinase Name	ID Code	Protein Kinase Name	ID Code
Abl (Abl1)	ABL1	CDK6/CyclinD1	C6D1	GRK7	GRK7
ACK (TNK2)	ACK1	CDK6/CyclinD3	C6D3	GSK3 $\alpha$	GS3A
Akt1/PKB $\alpha$	AKT1	CDK7/Cyclin H1/MAT1	C7H1	GSK3 $\beta$	GS3B
Akt2/PKB $\beta$	AKT2	CDK9/Cyclin K	C9CK	Hck	HCK1
Akt3/PKB $\gamma$	AKT3	CHK1 (CHEK1)	CHK1	HER2 (ErbB2, Neu)	ERB2
ALK1	ALK1	CHK2 (CHEK2)	CHK2	HER4 (ErbB4)	ERB4
ALK2	ALK2	CK2 $\alpha$ 1 (CSNK2A1)	CK21	HIPK1 (Myak )	HIP1
ALK4 (ACVR1B)	ALK4	CK2 $\alpha$ 2 (CSNK2A2)	CK22	HIPK3	HIP3
AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 (PRKAA1/B1/G1)	AM11	CLK1	CLK1	HIPK4	HIP4
AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 2 (PRKAA1/B1/G2)	AM12	CLK2	CLK2	IGF1R	IGF1
AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 3 (PRKAA1/B1/G3)	AM13	CLK3	CLK3	InsR	INSR
AMPK $\alpha$ 1/ $\beta$ 2/ $\gamma$ 1 (PRKAA1/B2/G1)	AM21	COT (MAP3K8)	COT1	INSRR (IRR )	ISRR
AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1 (PRKAA2/B1/G1)	AP11	Csk	CSK1	IRAK2	IRA2
AMPK $\alpha$ 2/ $\beta$ 2/ $\gamma$ 2 (PRKAA2/B2/G2)	AP22	DAPK1	DAP1	IRAK4	IRA4
Arg (Abl2)	ABL2	DAPK2	DAP2	Itk	ITK1
Ask1 (MAP3K5)	ASK1	DAPK3 (ZIPK )	DAP3	JAK3	JAK3
Aurora A (AurA, STK6)	AUR1	DCAMKL2	DCA2	JNK1 (MAPK8) - mouse	JNK1
Aurora B (AurB, AURKB)	AUR2	DDR2	DDR2	JNK2 (MAPK9)	JNK2
Aurora C (AurC, AURKC)	AUR3	DMPK	DMPK	JNK3 (MAPK10)	JNK3
Axl	AXL1	DYRK1A	DY1A	KDR (VEGFR2)	KDR1
BARK1 (ADRBK1, GRK2)	GRK2	DYRK3	DYR3	KHS1 (MAP4K5)	KHS1
Blk	BLK1	eEF2K	EF2K	Kit	KIT1
Bmx	BMX1	EPHA1	EPA1	Kit [V654A]	KITM
B-Raf	RAF2	EPHA2	EPA2	Lck	LCK1
B-Raf [V599E]	RAF4	EPHA3	EPA3	LIMK1	LIM1
Brk (PTK6)	BRK1	EPHA4	EPA4	LKB1	LKB1
Btk	BTK1	EPHA6	EPA6	LOK (STK10)	LOK1
CaMK1 $\alpha$ (CAMK1)	CM1A	EphB1	EPB1	LRRK2	LRR2
CaMK1 $\beta$ (PNCK)	CM1B	EphB2	EPB2	Lyn A	LYNA
CaMK1 $\delta$ (CAMK1D)	CM1D	EphB3	EPB3	Lyn B	LYNB
CAMK1 $\gamma$	CM1G	EphB4	EPB4	MAPKAPK2	MKA2
CAMK2 $\alpha$ (CAMK2A)	CM2A	Erk1 (MAPK3)	ERK1	MAPKAPK3	MKA3
CaMK2 $\beta$ (CAMK2B)	CM2B	Erk2 (MAPK1)	ERK2	MAPKAPK5	MKA5
CAMK2 $\delta$	CM2D	FAK (PTK2)	FAK1	MARK1	MAR1
CAMK2 $\gamma$	CM2G	Fer	FER1	MARK2	MAR2
CAMK4	CMK4	Fes (Fps)	FES1	MARK3	MAR3
CAMKK1 (CAMKKA )	CKK1	FGFR1 (Flt2)	FGR1	MARK4	MAR4
CAMKK2	CKK2	FGFR1 (Flt2) [V561M]	FGRM	MEK1 [S218E, S222E]	MEKA
CDK1 (CDC2)/CyclinA1	C1A1	FGFR2	FGR2	MEK1 (MKK1, MAP2K1)	MEK1
CDK1 (CDC2)/CyclinA2	C1A2	FGFR3	FGR3	MEK2 (MAP2K2)	MEK2
CDK1 (CDC2)/CyclinB1	C1B1	FGFR4	FGR4	MEKK2 (MAP3K2 )	MKK2
CDK2/CyclinA1	C2A1	Fgr	FGR1	MEKK3 (MAP3K3)	MKK3
CDK2/CyclinA2	C2A2	Flt1	FLT1	Mer (MERTK )	MER1
CDK2/CyclinE1	C2E1	Flt3	FLT3	Met	MET1
CDK3/CyclinE1	C3E1	Fms (CSF1R)	FMS1	MLK3 (MAP3K11)	MLK3
CDK4/Cyclin D1	C4D1	FRK	FRK1	MNK1 (MKNK1)	MNK1
CDK4/CyclinD3	C4D3	FynA	FYN1	MNK2 (MKNK2)	MNK2
CDK5/p25	C5P2	GCK	GCK1	MRCK $\alpha$ (CDC42BPA)	MRCA
CDK5/p29	C5P9	GPRK5 (GRK5)	GRK5	MRCK $\beta$ (CDC42BPB)	MRCB
CDK5/p35	C5P3	GRK6	GRK6	MSK1 (PRS6KA5)	MSK1

Table 1. Active Human Protein Kinases Available in the KICP Service. (Cont'd)

Protein Kinase Name	ID Code	Protein Kinase Name	ID Code	Protein Kinase Name	ID Code
MSSK1 (STK23)	MSS1	PKC $\epsilon$ (PRKCE)	PKCE	TGF $\beta$ R1	TGF1
MST1 (STK4)	MST1	PKC $\gamma$ (PRKCG)	PKCG	TGF $\beta$ R2	TGF2
MST2 (STK3)	MST2	PKC $\eta$ (PRKCH)	PKCH	Tie2 (Tek)	TIE2
MST3 (STK24)	MST3	PKC $\iota$ (PRKCI)	PKCI	TLK2	TLK2
MST4	MST4	PKC $\theta$ (PRKCQ)	PKCQ	TrkA (NTRK1)	TRKA
MUSK	MUSK	PKC $\zeta$ (PRKCZ)	PKCZ	TrkB (NTRK2)	TRKB
MYO3 $\beta$	MYO3	PKD1 (PRKD1, PKC $\mu$ )	PKD1	TrkC	TRKC
MYT1	MYT1	PKD2 (PRKD2)	PKD2	TSSK1 (STK22D)	TSS1
NDR1	NDR1	PKD3 (PRKD3)	PKD3	TSSK2 (STK22B)	TSS2
NEK11	NK11	PKG1 $\alpha$ (PRKG1A)	PKG1	TTK	TTK1
NEK2	NK02	PKN1	PKN1	TXK	TXK1
NEK3	NK03	PKN2	PKN2	TYK2	TYK2
NEK6	NK06	PKR (EIF2AK2)	PKR1	TYRO3 (RSE )	TYR3
NEK7	NK07	PLK1	PLK1	ULK1	ULK1
NEK9	NK09	PLK2	PLK2	ULK2	ULK2
NUAK2	NUA2	PLK4	PLK4	WEE1	WEE1
p38 $\alpha$ (MAPK14)	P38A	PYK2 (PTK2B)	PYK2	Yes (Yes1)	YES1
p38 $\beta$ (MAPK11)	P38B	QIK (SNF1LK2)	QIK1	YSK1 (STK25, SOK1)	YSK1
p38 $\delta$ (MAPK13)	P38D	Raf1	RAF1	ZAK	ZAK1
p38 $\gamma$ (MAPK12)	P38G	Raf1 [Y340E, Y341E]	RAFM	ZAP70	ZA70
p70S6K (RPS6KB1)	P70A	Ret	RET1	ZC3 (MINK1)	ZC03
p70S6Kb (RPS6KB2)	P70B	RIPK2	RIP2		
PAK1/CDC42	PAK1	RIPK5 (SGK496)	RIP5		
PAK2	PAK2	ROCK1 (ROK $\beta$ )	ROKB		
PAK3	PAK3	ROCK2 (ROK $\alpha$ )	ROKA		
PAK4	PAK4	RON (MST1R)	RON1		
PAK5 (PAK7)	PAK5	ROS	ROS1		
PAK6	PAK6	RSK1 (RPS6KA2)	RSK1		
PASK	PASK	RSK2 (PRS6KA3)	RSK2		
PBK (TOPK)	PBK1	RSK3 (RPS6KA1)	RSK3		
PDGFR $\alpha$	PDGA	RSK4 (RPS6KA6)	RSK4		
PDGFR $\beta$	PDGB	SGK1	SGK1		
PDK1 (PDPK1)	PDK1	SGK2	SGK2		
PEK (EIF2AK3)	PEK1	SGK3	SGK3		
PHKG2	PHK2	SIK	SIK1		
PI3K (p110 $\alpha$ /p65 $\alpha$ )	PIKT	skMLCK (MYLK2)	SKMK		
PI3K (p110 $\alpha$ /p85 $\alpha$ )	PIK1	SLK	SLK1		
PI3K (p110 $\alpha$ (E545K)/p85 $\alpha$ )	PIKM	smMLCK (MLCK2 )	SMMK		
PI3K (p110 $\delta$ /p85 $\alpha$ )	PIK4	Src	SRC1		
PI3K (p110 $\beta$ /p85 $\alpha$ )	PIK2	SRPK1	SRP1		
PI3K (p120 $\gamma$ )	PIK3	SRPK2	SRP2		
Pim1	PIM1	STK33	ST33		
Pim2	PIM2	SYK	SYK1		
PKA $\alpha$ (PRKACA)	PKAA	TAK1-TAB1	TAK1		
PKA $\beta$ (PRKACB)	PKAB	TAO1 (TAOK1)	TAO1		
PKA $\gamma$ (PRKACG)	PKAG	TAO2 (TAOK2)	TAO2		
PKC $\alpha$ (PRKCA)	PKCA	TAO3 (TAOK3, JIK )	TAO3		
PKC $\beta$ 1 (PRKCB1)	PKC1	TBK1	TBK1		
PKC $\beta$ 2 (PRKCB2)	PKC2	TEC	TEC1		
PKC $\delta$ (PRKCD)	PKCD	TESK1	TES1		

We provide a wide range of options to our clients with the KICP Service. Individual compounds may be profiled against a panel of protein kinase targets to establish the specificity of the compound. Alternatively, a panel of compounds may be tested against a single kinase target to identify a lead compound with the highest potency. Compounds may be tested either using a single dose or at multiple concentrations in order to allow in-depth IC determinations. In addition, the protein kinase assays can be performed under varying ATP concentrations to evaluate competition with respect to ATP. Compounds can be supplied by the client as DMSO stocks of known concentration, as solid material in vials, or in 96-well plates.

A large body of information and instruction is provided with this Kinase Inhibitor Compound Profiling Service Customer Information Package. Your careful review of this package will ensure that we can offer the highest level of quality in providing this and our other unique proteomics services to you. Should you have any questions or concerns, we would be pleased to hear from you. Thank you in advance for letting Kinexus become one of your trusted service providers.

## 2. QUANTITY OF COMPOUND REQUIRED

The amount of each compound required for the KICP Service depends on how many kinase activity measures are to be performed and the concentrations at which each compound will be tested. The final volume of the KICP assays are 25  $\mu$ l, and the stock concentration of the compound to be tested should be at least 10-times the final concentration of the highest dose desired for KICP analysis. For example, if 100  $\mu$ M is the single concentration of a compound to be tested against one kinase in triplicate, then (3 x 2.5  $\mu$ l =) 7.5  $\mu$ l of a 1 mM compound solution would be the minimum amount required. **However, we recommend that a minimum volume of 50  $\mu$ l of 10X concentrated compound stock solution in water or 2% DMSO is provided in a 1.5 ml Eppendorf vial; please Paraffin wax wrap the closed lid for further protection. If the compound is supplied in powder form, please provide sufficient material so that the compound can be prepared at as 10X concentrated solution with a volume of at least 500  $\mu$ l.**

## 3. KINASE ASSAY CONDITIONS

Due to the distinct protein/peptide substrate and other assay conditions for the different protein kinases, the components of the various assays are optimized for each enzyme and are not described here. If left unspecified by the client, most assays are performed for 15 minutes duration, at 30°C, with 50  $\mu$ M [ $\gamma$ -<sup>33</sup>P]ATP in a final volume of 25  $\mu$ l. The assays are typically terminated by spotting 20  $\mu$ l of the reaction mixture onto a phosphocellulose P81 plate. The phosphocellulose P81 plate is washed 3 times for approximately 15 minutes each in a 1% phosphoric acid solution to remove unreacted [ $\gamma$ -<sup>33</sup>P]ATP. The radioactivity in the captured <sup>33</sup>P-labeled peptide/protein substrate on the P81 plate is quantified in a scintillation counter.

## 4. SHIPPING INFORMATION

The aforementioned procedure has been designed to reduce the use of shipping materials and courier costs, and to ensure that your precious compound samples arrive in a safe and stable form at our laboratory facilities. Note that clients are responsible for payment of courier costs. The sample vials should be sent to the address listed below by any express courier that accepts dry ice shipments if necessary. We recommend Federal Express for shipments



originating in North America, and World Express is the preferred courier choice outside of North America. Ship the samples to the following address:

Kinase Inhibitor Compound Profiling Services  
Kinexus Bioinformatics Corporation  
Suite 1, 8755 Ash Street  
Vancouver, B.C. Canada V6P 6T3  
Telephone: (604) 323-2547  
Facsimile: (604) 323-2548  
E-mail [info@kinexus.ca](mailto:info@kinexus.ca)

Please ensure 3 copies of a signed commercial invoice accompany your shipment which specifies your samples are non hazardous. Since the samples are not for resale, the value of your shipment should be priced at approximately \$1.00 per sample in order to avoid paying additional duties and taxes on entry into Canada. **It is highly recommended that customers e-mail their courier airway bill number and the date of departure to [info@kinexus.ca](mailto:info@kinexus.ca) so we can track your shipment in transit and ensure it arrives in a timely manner.** We will send a confirmation e-mail once your shipment arrives at our facility.

## 5. PRICING INFORMATION

In view of the customized nature and the opportunity for volume discounts for the KICP Service, it is necessary to obtain a quotation from Kinexus for the desired work. To obtain a quotation, clients should complete the Kinase Inhibitor Compound Profiling - Service Order Form (KICP-SOF-01) in the sections under “Customer Information” and “Requested Work and Pricing Information” up to the point indicated on this form. In addition, clients should complete the appropriate Kinase Inhibitor Compound Profiling - Service Information Forms (KICP-SIF-01A and KICO-SIF-01B). These forms should be transmitted by facsimile to 604-323-2548 or by e-mail to [info@kinexus.ca](mailto:info@kinexus.ca) attention to “KICP Service.” Completion and submission of the KICP-SOF01 and KICP-SIF-01A/B forms will permit us to determine your specific needs and provide an accurate estimate of the costs.

Issued quotations are valid for 30 days from issuance. In the event that a special promotion on the KICS Service is offered at a lower price within 30 days of the receipt of Samples for KICP Service from a client, the lower promotional price will be applied at the time of billing the client.

## 6. FOLLOW UP SERVICES

Once you receive your Kinase Inhibitor Compound Profiling results back, you may find that your compound(s) targets additional protein kinases as perhaps even greater potency and selectivity than the intended target(s). This could offer new opportunities for development of intellectual property and clinical exploitation of your compounds. Only about 75 protein kinases are presently being targeted with drug leads in clinical trials. But, over 400 different diseases have been correlated with defects in cell signalling pathways that involve protein kinases. However, for the majority of the human protein kinases, it is unclear how they participate within cell signalling pathways, and no physiological substrates have been identified. With our unique suite of integrated systems proteomics and

bioinformatics services, Kinexus can further assist you in your disease biomarker and therapeutic drug discovery programs in so many ways in a cost-effective and timely fashion.

For example, with our Custom Kinase Substrate Profiling Service that uses our Kinex™ 800 Antibody Microarray (KAM-1.2 chip), we can identify a panel of physiological phosphoprotein substrates for over 360 different human protein kinases. This service can further identify the actual site of phosphorylation by the target kinase as well as commercial phospho-site antibodies for its detection in follow up studies. This is extremely useful for development of even more specific *in vitro* kinase activity assays. Moreover, this aids in the identification of a panel of biomarkers for the activity of the target kinase in the tissues of animals and isolated cells treated with compounds *in vivo*. Using our free access on-line PhosphoNET services ([www.phosphonet.ca](http://www.phosphonet.ca)), you can identify the protein kinases that we have determined with our proprietary algorithms are most likely to target more than 90,000 known human phospho-sites.

Our Kinex™ 800 Antibody Microarray services with the KAM-1.2 chip provide both qualitative and semi-quantitative analyses of the expression and phosphorylation states of cell signalling proteins in cell and tissue lysates. This can be useful to ascertain whether a lead compound actually targets the intended kinase *in vivo*, has potentially toxic effects in living systems, and how wide spread its actions may be. The qualitative analyses from the Kinex™ services include TIFF and JPEG files of the scanned Kinex™ antibody microarray that features the detected target signalling proteins in control and experimental (e.g. drug-treated) samples artificially labeled in two distinct colours by Adobe Photoshop and presented side-by-side in a coloured overlay. The quantitative analysis of the strength of the fluorescence signals for each target protein is provided in duplicate in a Microsoft Excel spreadsheet and includes the (average) percent change from the control sample, the percent range in error, and fold-changes ratios. Because the Kinex™ services use lysates that contain non-denatured proteins, the interpretation of results with the KAM-1.2 microarray chip is complicated by the possibilities of changes in protein-protein interactions and possible cross-reactivity with off target proteins.

Kinexus offers two types of Western blotting follow-up services to cost-effectively validate the results from the KAM-1.2 microarray chip. Clients can choose the Kinetworks™ Custom KCPS 1.0 (Multi-Antibody) Protein Screen, where any 18 antibodies can be selected and we will optimize it to their model system. Alternatively, with our Kinetworks™ Custom KCSS 1.0 (Multi-Sample) Protein Screen, customers send up to 8 different samples and choose up to 3 different antibodies (provided the molecular weights are significantly separated by SDS-PAGE).

Once the results are confirmed by Western blotting, clients can correlate their own data with hundreds of other data points from hundreds of different model systems using our on-line KiNET database. For more information about this and our other bioinformatics services, please contact one of our customer service representatives at [info@kinexus.ca](mailto:info@kinexus.ca).



## 7. FORMS TO BE COMPLETED

**All customers are required to complete the following forms for each order placed:**

- A. Kinexus Kinase Inhibitor Compound Profiling Services Agreement - Customers are required to complete and sign our standard Kinexus KICP Services Agreement before their first order can be processed. Unless otherwise specified, this Agreement is valid for all future orders with a standard term of 15 years.
- B. Service Order Form (KICP-SOF-01). The Service Order Form (SOF) allows us to track all of the requested work to be performed with your order. And your contact and billing information.
- C. Service Identification Form (KICP-SIF-01A and KICP-01B). The Service Identification Forms (SIF) permits us to determine which kinases and compounds are to be assayed. Two possible forms are provided depending on whether multiple kinases or multiple compounds are to be analyzed.
- D. Confidential Sample Description Form (KICP-CSDF-01). The Confidential Sample Description Form (CSDF) provides information on the specific compounds to be tested for safety and efficiency purposes. It is not necessary to disclose the name or structure of the compound to be analyzed.
- E. Commercial Invoice (required for all customers located outside of Canada). Two possible forms are provided depending on the stability of the compounds to be tested for shipping without freezing.

All orders with compounds should have as a minimum: 1 SOF, 1 SIF, and 1 CSDF forms completed, along with a courier airway bill and commercial invoice. A new KICP Service Agreement is not necessary if the client has previously placed an order with Kinexus and submitted a signed KICP Service Agreement at that time.

### FOR ALL CUSTOMERS

#### **A. Kinexus Kinase Inhibitor Compound Profiling Services Agreement**

A Kinexus KICP Services Agreement is required to be signed before the first order can be processed.

- This Agreement is required to be signed and dated by an authorized representative, typically a Senior Officer, Senior Scientist, or Principal Investigator, before the first order can be processed, but does not have to be signed again for repeat orders. The Kinexus Service Agreement is typically valid for 15 years. If you require changes or modifications to be made to our standard Kinexus KICP Services Agreement, please email us at [sales@kinexus.ca](mailto:sales@kinexus.ca) to request a Microsoft Word version of the document so your requested changes can be made directly into the agreement and emailed to us for our final approval.

#### **B. Service Order Form (KICP-SOF-01)**

*Please ensure:*

- Shipping address and contact name and numbers are specified
- Billing information is completed
- Any promotional vouchers or quotations are listed in the “Requested Works and Pricing Information” section

- Include a Purchase Order, Visa or MasterCard number for payment
- The form is signed and dated

### C. Service Identification Form (KICP-SIF-01A and KICP-SIF-01B)

*Please ensure:*

- You have selected the most appropriate form for your needs. With multiple compounds and multiple kinases, it will be necessary to use multiple copies of these forms.
- In Section A, you must assign a unique KICP-SIF ID name for each separate KICP-SIF form submitted and these are all recorded on the KICP-SOF form in the “Requested Works and Pricing Information” section
- In Section B, you must calculate the total number of individual assays requested for each separate KICP-SIF form submitted and these are all recorded on the KICP-SOF form in the “Requested Works and Pricing Information” section
- For Section B of the **KICP-SIF-01A** form, you must list the “Compound ID” name on each line in the first blank column that you provided in Section A of the KICP-CSDF-01 forms that have been completed
- For Section B of the **KICP-SIF-01B** form, you must list the “Kinase Code and Name from Table 1 or Appendix A together on each line in the first blank column that you provided in Section A of the KICP-CSDF-01 forms that have been completed
- When completing Section B of either the KICP-SIF-01A or KICP-SIF-01B forms for different concentrations of compound, make sure that you include a zero concentration control. Also ensure that you specify whether the concentrations shown in the table in Section B are in  $\mu\text{M}$  or  $\mu\text{g/ml}$ .
- Complete Section D if you require an assay time different from 15 minutes
- Complete Section F if you have any special instructions, such as a different temperature from  $30^{\circ}\text{C}$  for the assay incubation temperature or a concentration of ATP different from  $50\ \mu\text{M}$  for use in the assays
- The form is certified correct and signed and dated

For an example, in the completion of the Section B, consider the case of where 3 kinases (i.e. CDK1/CyclinB1, ERK2 and p38 $\alpha$ ) are to be tested against four concentrations of a compound in duplicate as well as in the absence of the compound. Section B of the KICP-SIF-01B form would be completed as exemplified on the next page:

### C. KINASE SELECTION:

Use Appendix A of the Kinase Inhibitor Compound Profiling Customer Information Package to obtain the code (Column A) and name (Column B) of the protein kinase that will be tested for inhibition by the compounds listed in Section B. A more detailed MS-Excel spreadsheet with information about each kinase can be downloaded from the Kinexus website at [www.kinexus.ca/services/kinase\\_inhibitor\\_profiling.php](http://www.kinexus.ca/services/kinase_inhibitor_profiling.php). For each concentration of the compound to be tested (up to 6) indicate the amount and whether this represents "µM" or "µg/ml" by ticking the appropriate box. Also indicate whether each assay is to be performed in single ("Once"), duplicate ("Dupl.") or triplicate ("Tripl.") measurements. Under "Subtotal" for each compound, indicate the number of individual assays that you intend to be performed.

Kinase Code + Name	Compound Final Conc. (µM) <input checked="" type="checkbox"/> or (µg/ml) <input type="checkbox"/>						Replicates			Subtotal
	Dil. 1	Dil. 2	Dil. 3	Dil. 4	Dil. 5	Dil. 6	Once	Dupl.	Tripl.	
1. C1B1-CDK1 /CyclinB1	0	0.01	0.1	1.0	10		<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	10
2. ERK2 - ERK2	0	0.01	0.1	1.0	10		<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	10
3. P38A - p38α	0	0.01	0.1	1.0	10		<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	10
4. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
5. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
6. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
7. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
8. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
9. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
10. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
11. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
12. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
13. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
14. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
15. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	

Total Number of Assays 30

### D. Confidential Sample Description Form (KICP-CSDF-01)

For each compound submitted, please ensure the following:

- Each sample tube is labeled and properly identified with the "Client ID Name" on the form in Section A, including final concentration and volume if it is a solution or weight if it is a powder.
- Up to two different compounds can be listed on the same page. Use additional copies of this form for more compounds to be analyzed
- You have indicated whether there are any safety concerns with the compounds on this form
- The form is certified correct and signed and dated

### E. Airway bill for Federal Express or any courier that you choose to ship your compounds to Kinexus

Complete the airway bill and specify:

- Priority overnight delivery
- Bill transportation charges to your institute

- Place sufficient dry ice if necessary to last several days into a Styrofoam shipping container
- Seal the edges of the Styrofoam container with tape to preserve dry ice longer
- Dry ice is a “*hazardous*” item, so ensure proper labels are attached to the outside of the box
- Do not specify Saturday delivery or hold at courier location
- For Federal Express shipments telephone 1-800-GO-FEDEX or visit them on-line at [www.fedex.com](http://www.fedex.com) or [www.fedex.ca](http://www.fedex.ca) to schedule a pick up or complete your forms
- For shipments coming from within Canada or the United States, please ship any day from Monday to Wednesday. **Do not ship on a Thursday or Friday.**
- For international shipments coming from outside of North America, the best day to ship is on a Monday to ensure arrival in Canada for delivery later the same week
- It is recommended that customers e-mail the date of your shipment and the courier airway bill number with the number of samples to Kinexus at [info@kinexus.ca](mailto:info@kinexus.ca) to ensure we can track your package should it get held up in Canadian Customs
- For any customer located outside of Canada, 3 copies of a commercial invoice is required to accompany your shipment (see below)

## FOR U.S AND INTERNATIONAL CUSTOMER ONLY

### F. Commercial Invoice (not required by Canadian customers)

*Please complete the attached commercial invoice with the following information:*

- Date of exportation
- Shipper/Exporter name, address, phone number
- Country of export/Country of origin
- Name of courier and the airway bill number
- Number, type and total weight of package(s)
- Total declared value of shipment (number of samples x \$1.00 per sample) and please specify currency
- Date, name, signature, and title of authorized person
- Include three (3) copies of the commercial invoice with the airway bill

**NOTE: Do not change the value of your shipment to more than \$1.00 per sample as this will prompt the custom brokers to charge Kinexus with a duty and GST fee on your package. Since the samples are processed internally and not returned to the customer or resold, there is no real commercial value.**

The international air waybill is required for all international shipments between Canada and the rest of the world. It is also your customs declaration, which can possibly be used to clear your shipment through customs at the destination. The customs clearance process begins with the description of the air waybill. If the description is too vague or missing, customs authorities may select the shipment for further inspection. All customs paperwork, such as the commercial invoice, must have detailed commodity descriptions. A detailed description on the air waybill and other customs documentation will help speed up the clearance time and reduce your delivery time. In the event that Kinexus must go to a Canada Customs facility to claim the package of samples for client order due to inadequate completion of the commercial invoice, additional charges will apply.

## Appendix A - List of Available Active Protein Kinases for Compound Profiling - 2010 May 1

This list may change depending on availability.

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
ABL1	Abl (Abl1)	Proto-oncogene tyrosine-protein kinase Abl1	ABL; JTK7; p150; c-ABL; v-abl; bcr/abl	P00519	NP_005148	Recombinant human ABL1 (27-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~135
ACK1	ACK (TNK2)	Activated CDC42 kinase 1	ACK, ACK1, FLJ44758, FLJ45547, p21cdc42Hs	Q07912	NP_005772	Recombinant human ACK (110-476) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~66
AKT1	Akt1/PKB $\alpha$	RAC-alpha serine/threonine-protein kinase	PKB; RAC; PRKBA; MGC99656; RAC-ALPH	P31749	NP_005154	Recombinant full-length human AKT1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~85
AKT2	Akt2/PKB $\beta$	RAC-beta serine/threonine-protein kinase	PRKBB; PKBBETA; RAC-BETA	P31751	NP_001617	Recombinant full-length human AKT2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~85
AKT3	Akt3/PKB $\gamma$	Rac-gamma serine/threonine-protein kinase	PKBG; PRKBG; STK-2; RAC-gamma; RAC-PK-gamma; DKFZP434N0250	Q9Y243	NP_005456	Recombinant full-length human AKT3 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~85
ALK1	ALK1	Serine/threonine-protein kinase receptor R3	ACVRL1, ACVRLK1, ALK1, HHT, HHT2, ORW2, SKR3, ALK1, TSR-I	P37023	NP_000011	Recombinant human ALK1 (144-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~64
ALK2	ALK2	Activin receptor type I	ACVR1, ACTRI, ACVR1A, ACVRLK2, FOP, SKR1, TSRI	Q04771	NP_001096	Recombinant human ALK2 (147-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~67
ALK4	ALK4 (ACVR1B)	Serine/threonine-protein kinase receptor R2	ACVR1B, SKR2, ACTRIB, ACVRLK4	P36896	NP_004293	Recombinant human ALK4 (150-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~64

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
AM11	AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 (PRKAA1/B1/G1)	5'-AMP-activated protein kinase, catalytic alpha-1 chain	Subunit A1: PRKAA1, MGC33776, MGC57364 Subunit B1: PRKAB1, AMPK, HAMPKb, MGC17785 Subunit G1: PRKAG1, AMPKG, MGC8666	Q13131	NP_996790	Recombinant full-length human AMPK (combination of A1/B1/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~68k (A1), ~38 (B1), ~4 (G1)
AM12	AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 2 (PRKAA1/B1/G2)	5'-AMP-activated protein kinase, catalytic alpha-1 chain	Subunit A1: PRKAA1, MGC33776, MGC57364 Subunit B1: PRKAB1, AMPK, HAMPKb, MGC17785 Subunit G2: PRKAG2, AAKG, CMH6, WPWS, AAKG2, H91620p	Q13131	NP_996790	Recombinant full-length human AMPK (combination of A1/B1/G2 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~68k Da (A1), ~38 (B1), ~65 (G2)
AM13	AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 3 (PRKAA1/B1/G3)	5'-AMP-activated protein kinase, catalytic alpha-1 chain	Subunit A1: PRKAA1, MGC33776, MGC57364 Subunit B1: PRKAB1, AMPK, HAMPKb, MGC17785 Subunit G3: PRKAG3	Q13131	NP_996790	Recombinant full-length human AMPK (combination of A1/B1/G3 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~68 (A1), ~38 (B1), ~51 (G3)
AM21	AMPK $\alpha$ 1/ $\beta$ 2/ $\gamma$ 1 (PRKAA1/B2/G1)	5'-AMP-activated protein kinase, catalytic alpha-1 chain	Subunit A1: PRKAA1, MGC33776, MGC57364 Subunit B2: PRKAB2, MGC61468 Subunit G1: PRKAG1, AMPKG, MGC8666	Q13131	NP_996790	Recombinant full-length human AMPK (combination of A1/B2/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~68 (A1), ~36 (B2), ~40 (G1)
AP11	AMPK $\alpha$ 2/ $\beta$ 1/ $\gamma$ 1 (PRKAA2/B1/G1)	5'-AMP-activated protein kinase, catalytic alpha-2 chain	Subunits A2: PRKAA2, AMPK, AMPK2, PRKAA Subunit B1: PRKAB1, AMPK, HAMPKb, MGC17785 Subunit G1: PRKAG1, AMPKG, MGC8666	P54646	NP_006243	Recombinant full-length human AMPK (combination of A2/B1/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~69 (A2), ~38 (B1), and ~40 (G1)



Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
AP22	AMPK $\alpha$ 2/ $\beta$ 2/ $\gamma$ 2 (PRKAA2/B2/G2)	5'-AMP-activated protein kinase, catalytic alpha-2 chain	Subunits A2: PRKAA2, AMPK, AMPK2, PRKAA Subunit B2: PRKAB2, AMPK, HAMPKb, MGC17785 Subunit G2: PRKAG2, AMPKG, MGC8666	P54646	NP_006243	Recombinant full-length human AMPK (combination of A2/B1/G1 subunits) was expressed by baculovirus in Sf9 insect cells using a C-terminal His tags.	~69 (A2), ~38 (B1), and ~40 (G1)
ABL2	Arg (Abl2)	Tyrosine-protein kinase Abl2	ARG; ABL	P42684	NP_005149	Recombinant human ABL2 (38-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~132
ASK1	Ask1 (MAP3K5)	Mitogen-activated protein kinase kinase 5	MAP3K5; MEKK5; MAPKKK5	Q99683	NP_005914	Recombinant human ASK1(649-946) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~60
AUR1	Aurora A (AurA, STK6)	Serine/threonine-protein kinase 6	AURKA, AIK, ARK1, AURA, BTAK, STK6, STK7, STK15, AURORA2, MGC34538	O14965	NP_940835	Recombinant full-length human AURORA A was expressed by baculovirus in Sf9 cells using an N-terminal GST tag.	~72
AUR2	Aurora B (AurB, AURKB, INCENP)	Serine/threonine-protein kinase 12	AURKB, AIK2; AIM1; ARK2; AurB; IPL1; AIM-1; STK12	Q96GD4	NP_004208	Recombinant full-length human AURORA B was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~68
AUR3	Aurora C (AurC, AURKC)	Serine/threonine-protein kinase 13	AURKC, AIE2, AIK3, AurC, STK13, aurora-C	Q9UQB9	NP_001015878	Recombinant full-length human AURORA C was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~61
AXL1	Axl	Tyrosine-protein kinase receptor UFO	UFO, JTK11	P30530	NP_068713	Recombinant human AXL (473-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~55
GRK2	BARK1 (ADRBK1, GRK2)	Adrenergic, beta, receptor kinase 1	BETA-ARK1, FLJ16718, GRK2	P25098	NP_001610	Recombinant full length human GRK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~102
BLK1	Blk	B lymphoid tyrosine kinase	MGC10442	P51451	NP_001706	Recombinant full-length human BLK was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~84

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
BMX1	Bmx	Cytoplasmic tyrosine-protein kinase BMX	ETK; PSCTK2; PSCTK3	P51813	NP_001712	Recombinant full-length human BMX was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~110
RAF2	B-Raf	B-Raf proto-oncogene serine/threonine-protein kinase	BRAF1, RAFB1, B-raf 1, MGC126806, MGC138284	P15056	NP_004324	Recombinant human BRAF (416-766) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~63
RAF4	B-Raf [V599E]	B-Raf proto-oncogene serine/threonine-protein kinase	BRAF1, RAFB1, B-raf 1, MGC126806, MGC138284	P15056	NP_004324	Recombinant human BRAF (416-end; V599E) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
BRK1	Brk (PTK6)	Tyrosine-protein kinase 6	PTK6; FLJ42088	Q13882	NP_005966	Recombinant full-length human BRK was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~80
BTK1	Btk	Tyrosine-protein kinase BTK	AT; ATK; BPK; XLA; IMD1; AGMX1; PSCTK	Q06187	NP_000052	Recombinant full-length human BTK was expressed by baculovirus in Sf9 insect cells using a N-terminal His tag.	~78
CM1A	CaMK1 $\alpha$ (CAMK1)	Calcium/calmodulin-dependent protein kinase type 1	AI505105; D6ErtD263e	Q14012	NP_003647	Recombinant full-length mouse CAMK1 $\alpha$ was expressed in E-coli cells using an N-terminal GST tag.	~70
CM1B	CaMK1 $\beta$ (PNCK)	Calcium/calmodulin-dependent protein kinase 1beta	Punc; Bstk3; CaMKIb2; caMKIb1; Pnck	Q6P2M8	AAH64422	Recombinant full-length mouse CAMK1 $\beta$ was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~64
CM1D	CaMK1 $\delta$ (CAMK1D)	Calcium/calmodulin-dependent protein kinase I delta	RP11-462F15.1; CKLiK; CaM-K1; CaMKID	Q8IU85	NP_705718	Recombinant full-length human CAMK1 $\delta$ was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~68
CM1G	CAMK1 $\gamma$	Calcium/calmodulin-dependent protein kinase 1 gamma	VWS1; CLICKIII; dJ272L16.1	Q96NX5	NP_065172	Recombinant human CAMK1 $\gamma$ (C-terminal truncation) protein was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag. Recombinant , GST-tagged human CAMK1 $\gamma$ (1-330) is also available.	~60
CM2A	CAMK2 $\alpha$ (CAMK2A)	Calcium/calmodulin-dependent protein kinase type II alpha	CAMKA; KIAA0968	Q9UQM7	NP_057065	Recombinant full-length human CAMK2 $\alpha$ was expressed in Sf9 cells using an N-terminal GST tag.	~74
CM2B	CaMK2 $\beta$ (CAMK2B)	Calcium/calmodulin-dependent protein kinase type II beta	CAMKB, CAM2, CAMK2, MGC29528	Q13554	NP_001211	Recombinant full-length human CAMK2 $\beta$ was expressed by baculovirus in Sf9 cells using an N-terminal His tag.	~58

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
CM2D	CAMK2 $\delta$	Calcium/calmodulin-dependent protein kinase type delta	CAMKD, MGC44911	Q8IU85	NP_742125	Recombinant full length human CAMK2delta was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~77
CM2G	CAMK2 $\gamma$	Calcium/calmodulin-dependent protein kinase type II gamma	CAMKG, CAMK, CAMK-II, MGC26678	Q13555	NP_751911	Recombinant human CAMK2 $\gamma$ (C-terminal truncation) was expressed by baculovirus in Sf9 cells using an N-terminal GST tag.	~60
CMK4	CAMK4	Calcium/calmodulin-dependent protein kinase type IV	CaMK-GR; MGC36771	Q16566	NP_001735	Recombinant full-length human CAMK4 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~79
CKK1	CAMKK1 (CAMKKA )	Calcium/calmodulin-dependent protein kinase kinase 1	CAMKKA, MGC34095, DKFZp761M0423	Q8N5S9	NP_115670	Recombinant full-length human CAMKK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~94
CKK2	CAMKK2	Calcium/calmodulin-dependent protein kinase kinase 2	CAMKK; CAMKKB; KIAA0787; MGC15254	Q96RR4	NP_006540	Recombinant full-length human CAMKK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~88
C1A1	CDK1 (CDC2)/CyclinA1	Cell division control protein 2 homolog	CDK1: CDC2 CDC28A; MGC111195; DKFZp686L20222	P06493	NP_001777	Recombinant full-length human CDK1 and CyclinA1 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and CyclinA1 ~78
C1A2	CDK1 (CDC2)/CyclinA2	Cell division control protein 2 homolog	CDK1: CDC2 CDC28A; MGC111195; DKFZp686L20222 CyclinA2: CCN1; CCNA	P06493	NP_001777	Recombinant full-length human CDK1 and CyclinA2 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and CyclinA2 ~78
C1B1	CDK1 (CDC2)/CyclinB1	Cell division control protein 2 homolog	CDK1: CDC2 CDC28A; MGC111195; DKFZp686L20222	P06493	NP_001777	Recombinant full-length human CDK1 and CyclinB1 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and CyclinB1 ~62
C2A1	CDK2/CyclinA1	Cell division protein kinase 2	CDK2: p33(CDK2)	P24941	NP_001789	Recombinant full-length human CDK2 and CyclinA1were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~58

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
C2A2	CDK2/CyclinA2	Cell division protein kinase 2	CDK2: p33(CDK2) CyclinA2: CCN1; CCNA	P24941	NP_001789	Recombinant full-length human CDK2 and CyclinA2 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~58 and CyclinA2 ~78
C2E1	CDK2/CyclinE1	Cell division protein kinase 2	CDK2: p33(CDK2) CyclinE1: CCNE1; CCNE	P24941	NP_001789	Recombinant full-length human CDK2 and CyclinA2 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~58 and CyclinE1 ~78
C3E1	CDK3/CyclinE1	Cell division protein kinase 3	CDK3: None; CyclinE1: CCNE1, CCNE	Q00526	NP_001249	Recombinant full-length human CDK3 and Cyclin E1 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~60 and CyclinE1 ~73
C4D1	CDK4/Cyclin D1	Cell division protein kinase 4	CDK4: CMM3; PSK-J3; MGC14458 CyclinD1: BCL1, PRAD1, U21B31, D11S287E	P11802	NP_000066	Recombinant full-length human CDK4 and CyclinD1 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~57 and CyclinD1 ~61
C4D3	CDK4/CyclinD3	Cell division protein kinase 4	CDK4: CMM3; PSK-J3; MGC14458 Cyclin D3: CCND3	P11802	NP_000066	Recombinant full-length human CDK4 and CyclinD3 were co-expressed by baculovirus in Sf9 cells using an N-terminal GST tag on both proteins.	~58 and CyclinD3 ~58
C5P2	CDK5/p25	Cell division protein kinase 5	CDK5: PSSALRE; p25: CDK5R1; CDK5P35, CDK5R, NCK5A, p23, p35, p35nck5a	Q00535	NP_004926	Recombinant full-length human CDK5 and p25 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and p25 ~49
C5P9	CDK5/p29	Cell division protein kinase 5	CDK5: PSSALRE	Q00535	NP_004926	Recombinant full-length human CDK5 and p29 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and p25 ~53
C5P3	CDK5/p35	Cell division protein kinase 5	CDK5: PSSALRE	Q00535	NP_004926	Recombinant full-length human CDK5 and p29 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~59 and p35 ~63
C6D1	CDK6/CyclinD1	Cell division protein kinase 6	PLSTIRE, MGC59692 /CCND1	Q00534	NP_001250	Recombinant full-length human CDK6 and CyclinD3 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~40 and CyclinD1 ~35

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
C6D3	CDK6/CyclinD3	Cell division protein kinase 6	PLSTIRE, MGC59692 /CCND3	Q00534	NP_001250	Recombinant full-length human CDK6 and CyclinD3 were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~40 and CyclinD3 ~35
C7H1	CDK7/Cyclin H1/MAT1	Cell division protein kinase 7	CDK7: CAK1, STK1, CDKN7, p39MO15; Cyclin H1: CCNH, CAK, p34, p37; MNAT1: MAT1, RNF66	P50613	NP_001790	Recombinant full-length human CDK7, Cyclin H1 and MNAT1 were co-expressed by baculovirus in Sf9 insect cells using N-terminal His tags.	~40, Cyclin H1 ~39, and MNAT1 ~37
C9CK	CDK9/Cyclin K	Cell division protein kinase 9	CDC2L4, C-2k, TAK, PITALRE; CCNK, CPR4, MGC9113	P50750	NP_001252	Recombinant full-length human CDK9 and CyclinK were co-expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag on both proteins.	~68 and CyclinK~67
CHK1	CHK1 (CHEK1)	Serine/threonine-protein kinase Chk1	CHEK1	O14757	NP_001265	Recombinant full-length human CHK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag. Also available with N-terminal His tag.	~82
CHK2	CHK2 (CHEK2)	Serine/threonine-protein kinase Chk2	RP11-436C9.1, CDS1, CHEK2, HuCds1, LFS2, PP1425, RAD53	O96017	NP_009125	Recombinant full-length human CHK2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~88
CK21	CK2 $\alpha$ 1 (CSNK2A1)	Casein kinase II, alpha chain	CKII; CK2A1; CKII $\alpha$	P68400	NP_808227	Recombinant full-length human CK2 $\alpha$ 1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~70
CK22	CK2 $\alpha$ 2 (CSNK2A2)	Casein kinase 2, alpha prime	CKII; CK2A2; CKII $\alpha$ , CSNK2A1, FLJ43934	P19784	NP_001887	Recombinant full-length human CK2 $\alpha$ 2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~70
CLK1	CLK1	Dual specificity protein kinase CLK1	CLK, STY, CLK/STY	P49759	NP_004062	Recombinant human CLK1 (129-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~66
CLK2	CLK2	Dual specificity protein kinase CLK2	hCLK2, MGC61500	P49760	NP_003984	Recombinant human CLK2 (137-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~68
CLK3	CLK3	Dual specificity protein kinase CLK3	CDC-like kinase 3	P49761	NP_003983	Recombinant full-length human CLK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~86

KinexUS ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
COT1	COT (MAP3K8)	Mitogen-activated protein kinase kinase kinase 8	MAP3K8, EST, ESTF, TPL2, Tpl-2, c-COT, FLJ10486	P41279	NP_005195	Recombinant human COT (30-397) was expressed by baculovirus in Sf9 insect cells.	~70
CSK1	Csk	Tyrosine-protein kinase CSK	MGC117393	P41240	NP_004374	Recombinant full-length human CSK was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~78
DAP1	DAPK1	Death-associated protein kinase 1	DAPK, DKFZp781I035	P53355	NP_004929	Recombinant human DAPK1 (1-363) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
DAP2	DAPK2	Death-associated protein kinase 2	DRP-1; MGC119312	Q9UIK4	NP_055141	Recombinant human full-length DAPK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~67
DAP3	DAPK3 (ZIPK )	Death-associated protein kinase 3	ZIP, ZIPK, FLJ36473	O43293	NP_001339	Full-length recombinant human DAPK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~79
DCA2	DCAMKL2	Serine-threonine-protein kinase DCAMKL2	DCLK2; DCK2; DCDC3; DCDC3B; DKFZp761I032; MGC45428	Q8N568	NP_689832	Recombinant human DCAMKL2 (1-690) was expressed by baculovirus in Sf9 cells using an N-terminal GST tag.	~105
DDR2	DDR2	Discoidin domain receptor 2	TKT; MIG20a; NTRKR3; TYRO10	Q16832	NP_001014796	Recombinant human DDR2 (467-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~70
DMPK	DMPK	Myotonic Dystrophy Protein Kinase	DM; DM; DMK; MDPK; DM1PK; MT-PK	Q09013	NP_004400	Recombinant full-length human DMPK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~105
DY1A	DYRK1A	Dual-specificity tyrosine-(Y)-phosphorylation regulated kinase 1A	DYRK, DYRK1, HP86, MNB, MNBH	Q13627	NP_569120		
DYR3	DYRK3	Dual-specificity tyrosine-(Y)-phosphorylation regulated kinase 3	DYRK5, RED, REDK, hYAK3-2	O43781	NP_003573		
EF2K	eEF2K	Eukaryotic elongation factor-2 kinase	HSU93850, MGC45041, eEF-2K	O00418	NP_037434	Recombinant full-length human EEF2K was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~125
EPA1	EPHA1	Ephrin type-A receptor 1	EPH, EPHT, EPHT1	P21709	NP_005223	Recombinant mouse EPHA1 (569-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~71



Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
EPA2	EPHA2	Ephrin type-A receptor 2	ECK	P29317	NP_004422	Recombinant human EPHA2 (561-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~72
EPA3	EPHA3	Ephrin type-A receptor 3	ETK, HEK, ETK1, HEK4, TYRO4	P29320	NP_005224	Recombinant human EPHA3 (571-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~72
EPA4	EPHA4	Ephrin type-A receptor 4	SEK, HEK8, TYRO1	P54764	NP_004429	Recombinant human EPHA4 (610-887) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~58
EPA6	EPHA6	Ephrin type-A receptor 6	EPA6, FLJ35246, PRO57066, DKFZp434C1418	Q9UF33	NP_775926	Recombinant human EPHA6 (561-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
EPB1	EphB1	Ephrin type-B receptor 1	Elk, Net, Cek6, Elkh, Hek6, EPHT2, AW488255, 9330129L11	P54762	NP_004432	Recombinant mouse EPHB1 (591-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~62
EPB2	EphB2	Ephrin type-B receptor 2	DRT, ERK, CAPB, Hek5, PCBC, EPHT3, Tyro5, MGC87492	P29323	NP_059145	Recombinant human EPHB2 (570-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~66
EPB3	EphB3	Ephrin type-B receptor 3	ETK2, HEK2, Cek10, Mdk5, TYRO6	P54753	NP_004434	Recombinant human EPHB3 (585-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~68
EPB4	EphB4	Ephrin type-B receptor 4	HTK, MYK1, TYRO11, Mdk2	P54760	NP_004435	Recombinant human EPHB4 (561-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~50
ERK1	Erk1 (MAPK3)	Mitogen-activated protein kinase 3	MAPK3; PRKM3; P44ERK1; P44MAPK; HS44KDAP; HUMKER1A; MGC20180	P27361	AAA36142.1	Recombinant full-length, tag-free human ERK1 was expressed in Sf9 cells and activated by active MEK1 in vitro.	~44
ERK2	Erk2 (MAPK1)	Mitogen-activated protein kinase 1	MAPK1, P42MAPK, MAPK2, PRKM2, PRKM1, p41mapk, ERT1, p40, p41, ERK	P28482	NP_002736	Recombinant full-length human ERK2 was expressed by E. coli cells using an N-terminal GST tag and activated by MEK1 in vitro.	~68
FAK1	FAK (PTK2)	Focal adhesion kinase 1	PTK2, FADK, FAK1, pp125FAK	Q05397	NP_722560	Recombinant human FAK (393-698) was expressed in Sf9 insect cells using an N-terminal His tag.	~35

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
FER1	Fer	Proto-oncogene tyrosine-protein kinase Fer	Fert, Fert2, AV082135, C330004K01Rik	P16591	NP_005237	Recombinant mouse FER (542-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~59
FES1	Fes (Fps)	Proto-oncogene tyrosine-protein kinase Fes/Fps	Fps	P07332	NP_001996	Recombinant full-length human FES was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~125
FGR1	FGFR1 (Flt2)	Basic fibroblast growth factor receptor 1	CEK; FLG; FLT2; KAL2; BFGFR; CD331; FGFR; HBGFR; N-SAM; C-FGR	P11362	NP_000595	Recombinant human FGFR1 (399-822) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
FGRM	FGFR1 (Flt2) [V561M]	Basic fibroblast growth factor receptor 1	FLT2, CEK, FLG, KAL2, BFGFR, C-FGR, CD331, N-SAM	P11362	NP_000595	Recombinant human FGFR1 (399-822; V561M) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~75
FGR2	FGFR2	Basic fibroblast growth factor receptor 2	K-SAM, BFR-1, CEK3, ECT1, TK14, TK25, CD332, JWS, TK14	P21802	NP_000132	Recombinant human FGFR2 (285-end) was expressed by baculovirus in Sf9 cells using an N-terminal GST tag.	~72
FGR3	FGFR3	Fibroblast growth factor receptor 3	ACH, CEK2, JTK4, CD333, HSFGFR3EX	P22607	NP_000133	Recombinant human FGFR3 (397-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
FGR4	FGFR4	Fibroblast growth factor receptor 4		P22455	NP_002002	Recombinant human FGFR4 (460-end) was expressed by baculovirus in Sf9 cells using an N-terminal GST tag.	~65
FGR1	Fgr	Proto-oncogene tyrosine-protein kinase Fgr	SRC2, c-fgr, c-src, FLJ43153, MGC75096, p55c-fgr, p58c-fgr	P09769	NP_005239	Recombinant full-length human FGR was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~86
FLT1	Flt1	Vascular endothelial growth factor receptor 1 [Precursor]	FLT, VEGFR1	P17948	NP_002010	Recombinant human FLT1 (784-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~94
FLT3	Flt3	FL cytokine receptor	FLK2, STK1, CD135	P36888	NP_004110	Recombinant human FLT3 (571-993) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~73
FMS1	Fms (CSF1R)	Macrophage colony stimulating factor 1 receptor	CSF1R, CSFR, FIM2, C-FMS, CD115	P07333	NP_005202	Recombinant human FMS (539-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~76

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
FRK1	FRK	Fyn-related kinase	GTK; RAK; PTK5	P42685	NP_002022	Recombinant human FRK (208-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~60
FYN1	FynA	Fyn oncogene related to Src, Fgr, Yes	RP1-66H14.1; MGC45350; SLK; SYN	P06241	NP_002028	Recombinant full length human FYN A was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~85
GCK1	GCK	Mitogen-activated protein kinase kinase kinase 2	MAP4K2, BL44, RAB8IP	Q12851	NP_004570	Recombinant full-length human GCK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~125
GRK5	GPRK5 (GRK5)	G protein-coupled receptor kinase 5	GPRK5	P34947	NP_005299	Recombinant full-length human GRK5 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~95
GRK6	GRK6	G protein-coupled receptor kinase 6	GPRK6	P43250	NP_002073	Recombinant full length human GRK6 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~93
GRK7	GRK7	G protein-coupled receptor kinase 7	GPRK7	Q8WTQ7	NP_631948	Recombinant full length human GRK7 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~89
GS3A	GSK3 $\alpha$	Glycogen synthase kinase-3 alpha	GSK3 alpha	P49840	NP_063937	Recombinant full-length human GSK3? was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~81
GS3B	GSK3 $\beta$	Glycogen synthase kinase-3 beta	GSK3 beta	P49841	NP_002084	Recombinant full-length human GSK3? was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~73
HCK1	Hck	Hemopoietic cell kinase	JTK9	P08631	NP_002101	Recombinant human HCK (230-497) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~57
ERB2	HER2 (ErbB2, Neu)	v-erb-b2 erythroblastic leukemia viral oncogene homologue 2	CD340, HER-2, HER-2/neu, NEU, NGL, TKR1	P04626	NP_004439	Recombinant human HER2 (676-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~116
ERB4	HER4 (ErbB4)	Receptor tyrosine-protein kinase erbB4	HER4, MGC138404, p180erbB4	Q15303	NP_005226	Recombinant human ErbB4 (682-993) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~57

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
HIP1	HIPK1 (Myak )	Homeodomain-interacting protein kinase 1	Myak, Nbak2, KIAA0630, MGC26642, MGC33446, MGC33548	Q86Z02	NP_938009	Recombinant human HIPK1 (156-555) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
HIP3	HIPK3	Homeodomain-interacting protein kinase 3	PKY, YAK1, DYRK6, FIST3	Q9H422	NP_005725	Recombinant human HIPK3 (163-562) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~49
HIP4	HIPK4	Homeodomain interacting protein kinase 4	FLJ32818	Q8NE63	NP_653286	Recombinant full-length human HIPK4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
IGF1	IGF1R	Insulin-like growth factor I receptor	CD221, IGFIR, JTK13, MGC142170, MGC142172	P08069	NP_000866	Recombinant human IGF1R (960-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~53
INSR	InsR	Insulin receptor [Precursor]	HHF5, CD220	P06213	NP_000199	Recombinant human InsR (1011-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~70
ISRR	INSRR (IRR )	Insulin receptor-related protein	IRR	P14616	NP_055030	Recombinant human IRR (945-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~65
IRA2	IRAK2	Interleukin-1 receptor-associated kinase-like 2	IRAK-2, MGC150550	O43187	NP_001561	Recombinant full-length human IRAK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~103
IRA4	IRAK4	Interleukin-1 receptor-associated kinase 4	IPD1, REN64, NY-REN-64	Q9NWZ3	NP_057207	Recombinant full-length human IRAK4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~81
ITK1	Itk	Tyrosine-protein kinase Itk/Tsk	EMT, LYK, PSCTK2, MGC126257, MGC126258	Q08881	NP_005537	Recombinant human ITK (352-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~53
JAK3	JAK3	Tyrosine-protein kinase JAK3	JAKL, LJAK, JAK-3, L-JAK, JAK3_HUMAN	P52333	NP_000206	Recombinant human JAK3 (781-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~64
JNK1	JNK1 (MAPK8) - mouse	Mitogen-activated protein kinase 8	JNK, PRKM8, SAPK1, AI849689, JNK1A2, JNK21B1/2	P45983	NP_002741	Recombinant full-length mouse JNK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~71

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
JNK2	JNK2 (MAPK9)	Mitogen-activated protein kinase 9	JNK-55, JNK2, JNK2A, JNK2ALPHA, JNK2B, JNK2BETA, PRKM9, SAPK, p54a, p54aSAPK	P45984	NP_002743	Recombinant full-length human JNK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~70
JNK3	JNK3 (MAPK10)	Mitogen-activated protein kinase 10	FLJ12099, FLJ33785, JNK3, JNK3A, MGC50974, PRKM10, p493F12, p54bSAPK	P53779	NP_002744	Recombinant full-length human JNK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
KDR1	KDR (VEGFR2)	Kinase insert domain receptor	FLK1, VEGFR, VEGFR2	P35968	NP_002244	Recombinant human KDR (789-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
KHS1	KHS1 (MAP4K5)	Mitogen-activated protein kinase kinase kinase 5	MAP4K5, KHS, GCKR, MAPKKKK5	Q9Y4K4	NP_006566	Recombinant full-length human KHS1 was expressed by baculovirus in Sf9 insect cells using an N-terminal tag.	~135
KIT1	Kit	Mast/stem cell growth factor receptor	PBT, SCFR, CD117	P10721	NP_000213	Recombinant human c-KIT (544-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~73
KITM	Kit [V654A]	Mast/stem cell growth factor receptor	PBT, SCFR, CD117	P10721	NP_000213	Recombinant human c-KIT (544-end, V654A) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
LCK1	Lck	Proto-oncogene tyrosine-protein kinase Lck	YT16, p56lck, pp58lck	P06239	NP_005347	Recombinant full-length human LCK was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~84
LIM1	LIMK1	LIM domain kinase 1	LIMK	P53667	NP_002305		~73
LKB1	LKB1	Serine-threonine-protein kinase 11	STK11; MO25a; STRADa	Q15831	NP_000446		~49
LOK1	LOK (STK10)	Serine/threonine-protein kinase 10	STK10, PRO2729	O94804	NP_005981	Recombinant human LOK (1-348) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~65
LRR2	LRRK2	Leucine-rich repeat kinase 2	AURA17, DARDARIN, PARK8, RIPK7, ROCO2	Q5S007	NP_940980.3		~290

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
LYNA	Lyn A	Tyrosine-protein kinase Lyn	JTK8, FLJ26625	P07948	NP_002341	Recombinant full-length human LYN A was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~81
LYNB	Lyn B	v-yes-1 Yamaguchi sarcoma viral related oncogene homolog	JTK8, FLJ26625, p65	P07948	NP_002341	Recombinant full-length human LYN B was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~85
MKA2	MAPKAPK2	Mitogen-activated protein kinase-activated protein kinase 2		P49137	NP_004750	Recombinant human MAPKAPK2 (46-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~68
MKA3	MAPKAPK3	Mitogen activated protein kinase activated protein kinase 3	3PK, MAPKAP3, 3pK	Q16644	NP_004626	Recombinant full-length human MAPKAPK3 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~69
MKA5	MAPKAPK5	Mitogen-activated protein kinase-activated protein kinase 5	PRAK	Q8IW41	NP_620777	Recombinant full-length human MAPKAPK5 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~79
MAR1	MARK1	MAP/microtubule affinity-regulating kinase 1	MARK, KIAA1477, MGC126512, MGC126513	Q9P0L2	NP_061120	Recombinant full-length human MARK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~125
MAR2	MARK2	MAP/microtubule affinity-regulating kinase 2	EMK1, MGC99619, PAR-1, Par1b, ELKL	Q7KZI7	NP_059672		~88
MAR3	MARK3	Microtubule affinity-regulating kinase 3	KP78, CTAK1, PAR1A	P27448	NP_002367	Recombinant full-length human MARK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
MAR4	MARK4	MAP-microtubule affinity-regulating kinase 4	MAP/microtubule affinity-regulating kinase-like 1	Q96L34	NP_113605		~83
MEKA	MEK1 (MAP2K1) [S218E, S222E]	Dual specificity mitogen-activated protein kinase kinase 1	MAP2K1, MKK1, MAPKK1, PRKMK1	Q02750	NP_002746	Recombinant full-length mouse MEK1 (S218E, S222E) was expressed in E. coli cells using a N-terminal GST tag.	~69
MEK1	MEK1 (MKK1, MAP2K1)	Dual specificity mitogen-activated protein kinase kinase 1	MAP2K1, MKK1, MAPKK1, PRKMK1	Q02750	NP_002746	Recombinant full-length human MEK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~72



Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
MEK2	MEK2 (MAP2K2)	Dual specificity mitogen-activated protein kinase kinase 2	MAP2K2, MKK2, PRKMK2, MAPKK2	P36507	NP_109587	Recombinant full-length human MEK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
MKK2	MEKK2 (MAP3K2)	Mitogen-activated protein kinase (MAPK) kinase kinase 2	MAP3K2, MEKK2B	Q9Y2U5	NP_006600	Recombinant full-length human MEKK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~115
MKK3	MEKK3 (MAP3K3)	Mitogen-activated protein kinase kinase kinase 3	MAP3K3, MAPKKK3	Q99759	NP_002392	Recombinant full-length human MEKK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~105
MER1	Mer (MERTK)	Proto-oncogene tyrosine-protein kinase Mer	MER, MERTK, MGC133349	Q12866	NP_006334	Recombinant human MERTK (578-872) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~58
MET1	Met	Hepatocyte growth factor receptor	HGFR	P08581	NP_000236	Recombinant rat MET (958-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~74
MLK3	MLK3 (MAP3K11)	Mitogen-activated protein kinase kinase kinase 11	MGC17114; Mixed lineage kinase 3; MLK-3; MLK3; PTK1; SPRK	Q16584	NP_002410		~93
MNK1	MNK1 (MKNK1)	MAP kinase interacting serine/threonine kinase 1	RP11-49P4.3	Q9BUB5	NP_003675		~51
MNK2	MNK2 (MKNK2)	MAP kinase-interacting kinase 2	MKNK2, GPRK7	Q9HBH9	NP_060042	Recombinant full-length human MNK2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~74
MRCA	MRCK $\alpha$ (CDC42BPA)	Myotonic dystrophy kinase-related Cdc42-binding kinase alpha	CDC42BPA, MRCK, PK428, FLJ23347, KIAA0451, DKFZp686L1738, DKFZp686P1738	Q5VT25	NP_003598	Recombinant human MRCK $\alpha$ (1-473) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag	~73
MRCB	MRCK $\beta$ (CDC42BPB)	Myotonic dystrophy kinase-related Cdc42-binding kinase beta	CDC42BPB, KIAA1124	Q9Y5S2	NP_006026	Recombinant human MRCK $\beta$ (1-473) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
MSK1	MSK1(PRS6KA5)	Mitogen- and stress-activated protein kinase 1	RPS6KA5, RLPK, MSPK1, MGC1911	O75582	NP_004746	Recombinant full-length human MSK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~120

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
MSS1	MSSK1 (STK23)	Serine/threonine-protein kinase 23	STK23, SRPK3, SFRS protein kinase 3, MGC102944	Q9UPE1	NP_055185	Recombinant full-length human MSSK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~98
MST1	MST1 (STK4)	Mammalian Sterile 20-like kinase 1	KRS2, YSK3, DKFZp686A2068, STK4	Q13043	NP_006273	Recombinant full-length human MST1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~83
MST2	MST2 (STK3)	Serine/threonine kinase 3	KRS1; MST2; FLJ90748	Q13188	NP_006272	Recombinant full-length human STK3 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~87
MST3	MST3 (STK24)	Mammalian Sterile 20-like kinase 3	STK24, MST-3, STK3, MST3B, RP11-111L24.5	Q9Y6E0	NP_003567	Recombinant human MST3 (1-311) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
MST4	MST4	Mammalian Sterile 20-like kinase 4	MASK, RP6-213H19.1	Q9P289	NP_057626	Recombinant full-length human MST4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~72
MUSK	MUSK	Muscle, skeletal receptor tyrosine protein kinase	MGC126323, MGC126324	O15146	NP_005583	Recombinant human MUSK (519-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
MYO3	MYO3 $\beta$	Myosin IIIB	Myosin IIIB	Q8WXR4	NP_620482	Recombinant Human MYO3 $\beta$ (1-326) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
MYT1	MYT1	Membrane-associated tyrosine-and threonine-specific cdc2-inhibitory kinase	MYT1 kinase; PKMYT1	O14731	NP_004194		~55
NDR1	NDR1	Serine-threonine-protein kinase 38	STK38	Q15208	NP_009202	Recombinant full length human NDR was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~54
NK11	NEK11	Serine-threonine-protein kinase Nek11	BC009414; FLJ23495	Q8TBY1	NP_079076	Recombinant full-length human NEK11 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
NK02	NEK2	Serine/threonine-protein kinase Nek2	NLK1, HsPK21, NEK2A	P51955	NP_002488	Recombinant full-length human NEK2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~76
NK03	NEK3	Serine-threonine-protein kinase Nek3	NimA-related protein kinase 3; HSPK 36; MGC29949	P51956	NP_002489	Recombinant full-length human NEK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~86

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
NK06	NEK6	NimA-related protein kinase 6	SID6-1512	Q9HC98	NP_055212	Recombinant full-length human NEK6 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~63
NK07	NEK7	NimA-related protein kinase 7		Q8TDX7	NP_598001	Recombinant full-length human NEK7 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~63
NK09	NEK9	Serine-threonine-protein kinase Nek9	NEK8, NERCC, NERCC1, MGC16714, MGC138306, DKFZp434D0935	Q8TD19	NP_149107	Recombinant human NEK9 (1-979; deleted fragment of 346-731) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~115
NUA2	NUAK2	NUAK family, SNF1-like kinase, 2	DKFZp434J037, DKFZp686F01113, FLJ90349, SNARK	Q9H093	NP_112214		
P38A	p38 $\alpha$ (MAPK14)	Mitogen-activated protein kinase 14	CSBP1, CSBP2, CSPB1, PRKM14, PRKM15, SAPK2A, MAPK14, RK, p38, EXIP, Mxi2	Q16539	NP_001306	Recombinant full-length human p38alpha was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~67
P38B	p38 $\beta$ (MAPK11)	Mitogen-activated protein kinase 11	MAPK11, SAPK2, p38-2, PRKM11, SAPK2B, p38b, P38b2	Q15759	NP_002742.3	Recombinant full-length human p38beta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~71
P38D	p38 $\delta$ (MAPK13)	Mitogen-activated protein kinase 13	SAPK4, PRKM13, MAPK13, MGC99536	O15264	NP_002745	Recombinant full-length human p38delta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~71
P38G	p38 $\gamma$ (MAPK12)	Mitogen-activated protein kinase 12	ERK3, ERK6, SAPK3, PRKM12, MAPK12	P53778	NP_002960	Recombinant full-length human p38gamma was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~71
P70A	p70S6K (RPS6KB1)	Ribosomal protein S6 kinase 1	S6K, PS6K, S6K1, STK14A, RPS6KB1, p70-alpha, p70(S6K)-alpha	P23443	NP_003152	Recombinant full-length human p70S6K was expressed by baculovirus in Sf9 insect cells using a N-terminal His tag.	~76

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
P70B	p70S6Kb (RPS6KB2)	Ribosomal protein S6 kinase, 70kDa, polypeptide 2	RPS6KB2, STK14B, KLS, P70-beta, P70-beta-1, P70-beta-2, S6K2, p70(S6K)-beta, S6K-beta2, SRK	Q9UBS0	NP_003943	Recombinant full-length human p70S6Kb was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~85
PAK1	PAK1/CDC42	Serine/threonine-protein kinase PAK 1	PAKalpha, MGC130000, MGC130001	Q13153	NP_002567	Recombinant full-length human PAK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. Combined with Cdc42 in assays.	~96
PAK2	PAK2	p21 protein (Cdc42/Rac)-activated kinase 2	PAK65, PAKgamma, S6/H4 kinase	Q13177	NP_002568	Recombinant full-length mouse PAK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~58
PAK3	PAK3	Serine/threonine-protein kinase PAK3	Stk4, mPAK-3, Pak65beta, Pak65alpha	O75914	NP_002569	Recombinant full-length mouse PAK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~89
PAK4	PAK4	Serine/threonine-protein kinase PAK4	p21-activated kinase 4; PAK-4	O96013	NP_005875	Recombinant full-length human PAK4 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~90
PAK5	PAK5 (PAK7)	Serine/threonine-protein kinase PAK 7	PAK5, KIAA1264, MGC26232	Q9P286	NP_065074	Recombinant full-length human PAK7 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~130
PAK6	PAK6	Serine-threonine-protein kinase PAK 6	p21-activated kinase 6; PAK-6; PAK-5	Q9NQU5	NP_064553	Recombinant full-length human PAK6 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
PASK	PASK	PAS domain containing serine-threonine-protein kinase	STK37; PASKIN; KIAA0135; DKFZp434O051; DKFZp686P2031	Q96RG2	NP_055963	Recombinant human PASK (981-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~66
PBK1	PBK (TOPK)	T-lymphokine-activated killer cell-originated protein kinase	SPK, Nori-3, FLJ14385	Q96KB5	NP_060962	Recombinant full-length human TOPK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~68
PDGA	PDGFR $\alpha$	Platelet-derived growth factor receptor alpha	CD140A, PDGFR2, MGC74795, Rhe-PDGFR $\alpha$	P16234	NP_006197	Recombinant human PDGFR $\alpha$ (550-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~100

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
PDGB	PDGFR $\beta$	Platelet-derived growth factor receptor beta	JTK12, PDGFR, CD140B, PDGFR1	P09619	NP_002600	Recombinant human PDGFRbeta (557-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~104
PDK1	PDK1 (PDPK1)	3-Phosphoinositide-dependent protein kinase 1	PRO0461, PDPK1, MGC20087, MGC35290, PkB-like, PkB-like 1	O15530	NP_002604	Recombinant full-length human PDK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal His tag.	~67
PEK1	PEK (EIF2AK3)	Eukaryotic translation initiation factor 2-alpha kinase 3	DKFZp781H1925, HRI, PEK, PERK, WRS	Q9NZJ5	NP_004827		
PHK2	PHKG2	Phosphorylase b kinase gamma catalytic chain, testis-liver isoform	GSD9C; PHK-gamma-T; Phosphorylase kinase gamma subunit 2; PSK-C3	P15735	NP_000285	Recombinant full-length human PHKG2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~70
PIKT	PI3K (p110 $\alpha$ /p65 $\alpha$ )	Phosphatidylinositol 3-kinase-alpha	p110 $\alpha$ ; PIK3CA; MGC142161; MGC142163; p110-alpha			Recombinant full-length human p110alpha and truncated human p85alpha were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	
PIK1	PI3K (p110 $\alpha$ /p85 $\alpha$ )	Phosphatidylinositol 3-kinase-alpha	p110 $\alpha$ ; PIK3CA; MGC142161; MGC142163; p110-alpha			Recombinant full-length mouse p110alpha and human p85alpha were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~111
PIKM	PI3K (p110 $\alpha$ (E545K)/p85 $\alpha$ )	Phosphatidylinositol 3-kinase-alpha	p110 $\alpha$ ; PIK3CA; MGC142161; MGC142163; p110-alpha			Recombinant full-length human p110 $\alpha$ (E545K) mutant subunit and human p85-alpha wild-type subunit were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~111
PIK4	PI3K (p110 $\delta$ /p85 $\alpha$ )	Phosphatidylinositol 3-kinase-delta	p110 $\delta$			Recombinant full-length human p110 $\delta$ and human p85 $\alpha$ were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~111

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
PIK2	PI3K (p110 $\beta$ /p85 $\alpha$ )	Phosphatidylinositol 3-kinase-beta	p110 $\alpha$ ; PIK3CB; PI3K; PIK3C1; PI3Kbeta; MGC133043; p110-BETA; DKFZp779K1237;			Recombinant full-length human p110beta and human p85alpha were co-expressed by baculovirus in Sf9 insect cells using an N-terminal His tag on both proteins.	~111
PIK3	PI3K (p120 $\gamma$ )	Phosphatidylinositol 3-kinase-gamma	p110 $\gamma$ ; PI3K; PIK3; PI3CG; PI3Kgamma			Recombinant full-length human PI3K (p120 $\gamma$ ) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~121
PIM1	Pim1	Proto-oncogene serine/threonine-protein kinase Pim-1	PIM	P11309	NP_002639	Recombinant full-length human PIM1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~62
PIM2	Pim2	Serine/threonine-protein kinase Pim2		Q9P1W9	NP_006866	Recombinant full-length human PIM2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~61
PKAA	PKA $\alpha$ (PRKACA)	cAMP-dependent protein kinase, alpha-catalytic subunit	PKAa, cAPKa, MGC48865, MGC102831	P17612	NP_002721	Recombinant full-length human PKAalpha was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~69
PKAB	PKA $\beta$ (PRKACB)	cAMP-dependent protein kinase, beta-catalytic subunit	PKAb, cAPKb, MGC9320, MGC41879, DKFZp781I2452	P22694	NP_891993	Recombinant full-length human PKAcbeta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~65
PKAG	PKA $\gamma$ (PRKACG)	cAMP-dependent protein kinase, gamma-catalytic subunit	KAPG, PKAr, cAPKr	P22612	NP_002723	Recombinant full-length human PKAcgamma was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~65
PKCA	PKC $\alpha$ (PRKCA)	Protein kinase C, alpha type	AAG6, PKCA, PRKCA, MGC129900, MGC129901	P17252	NP_002728	Recombinant full-length human PKCalpha was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~103
PKC1	PKC $\beta$ 1 (PRKCB1)	Protein kinase C, beta	PKCB; PRKCB; PRKCB2; MGC41878; PKC-beta	P05771	NP_002729	Recombinant full-length human PKCbeta I was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~102



Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
PKC2	PKC $\beta$ 2 (PRKCB2)	Protein kinase C, beta type	PRKCB1, PKCB2, PRKCB1, PRKCB (X07109)	P05771	NP_002729	Recombinant full-length human PKCbeta II was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~105
PKCD	PKC $\delta$ (PRKCD)	Protein kinase C, delta	PRKCD, MAY1, MGC49908, nPKC-delta	Q05655	NP_006245	Recombinant full-length human PKCdelta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~104
PKCE	PKC $\epsilon$ (PRKCE)	Protein kinase C, epsilon type	PRKCE, MGC125656, MGC125657, nPKC-epsilon	Q02156	NP_005391	Recombinant full-length human PKCepsilon was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~110
PKCG	PKC $\gamma$ (PRKCG)	Protein kinase C, gamma type	PKCC, PKCG, SCA14, MGC57564, PKC-gamma, PRKCG	P05129	NP_002730	Recombinant full-length human PKCgamma was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~105
PKCH	PKC $\eta$ (PRKCH)	Protein kinase C, eta	PKCL, PKC-L, PRKCL, MGC5363, MGC26269, nPKC-eta, PRKCH	P24723	NP_006246	Recombinant full-length human PKCeta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~103
PKCI	PKC $\iota$ (PRKCI)	Protein kinase C, iota type	PRKCI, DXS1179E, MGC26534, nPKC-iota	P41743	NP_002731	Recombinant full-length human PKCiota was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~98
PKCQ	PKC $\theta$ (PRKCQ)	Protein kinase C, theta	PRKCQ, PRKCT, MGC126514, MGC141919, nPKC-theta	Q04759	NP_006248	Recombinant full-length human PKC-theta was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~110
PKCZ	PKC $\zeta$ (PRKCZ)	Protein kinase C, zeta type	PRKCZ, PRKCZ, PKC2	Q05513	NP_002735	Recombinant full-length human PKCzeta was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~93
PKD1	PKD1 (PRKD1, PKC $\mu$ )	Protein kinase C, mu	PKD, PKCM, PRKCM	Q15139	NP_002733	Recombinant full-length human PKCmu was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~131
PKD2	PKD2 (PRKD2)	Protein kinase C, D2	HSPC187, DKFZp586E0820, PRKD2	Q9BZL6	NP_057541	Recombinant full-length human PKD2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~130

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
PKD3	PKD3 (PRKD3, PKCv)	Protein kinase C, nu	PKCv, EPK2, PRKCN, PKD3, nPKC-NU	O94806	NP_005804	Recombinant full-length human PKCnu was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~142
PKG1	PKG1 $\alpha$ (PRKG1A)	cGMP-dependent protein kinase 1, alpha	PGK, CGKI, PRKG1B, PRKGR1B, FLJ36117, MGC71944, cGKI-BETA, cGKI-alpha, DKFZp686K042	Q13976	NP_006249	Recombinant full-length human PRKG1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~100
PKN1	PKN1	Protein kinase N1	DBK, MGC46204, PAK1, PKN, PKN-ALPHA, PRK1, PRKCL1	Q16512	NP_998725		
PKN2	PKN2	Protein kinase N2	PRK2; PRKCL2, PKN2, PRK2, PAK2, Pak-2, PRO2042, MGC71074, MGC150606; Protein kinase C-like 2; Protein-kinase C-related kinase 2	Q16513	NP_006247	Recombinant full-length human PKN2/PRK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~145
PKR1	PKR (EIF2AK2)	Interferon-induced, double-stranded RNA-activated protein kinase	PKR, PRKR, EIF2AK1	P19525	NP_002750	Recombinant human EIF2AK2 (252-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~64
PLK1	PLK1	Serine/threonine-protein kinase Plk1	STPK13	P53350	NP_005021	Recombinant full-length human PLK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal His tag.	~70
PLK2	PLK2	Polo-like kinase 2	SNK	Q9NYY3	NP_006613		
PLK4	PLK4			O00444	NP_055079		
PYK2	PYK2 (PTK2B)	Protein tyrosine kinase 2 beta	FAK2, PTK2B, CADTK, FADK2, CAKB, PKB, PTK, RAFTK, PYK2	Q14289	NP_775266	Recombinant human PYK2 (360-690) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~39

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
QIK1	QIK (SNF1LK2)	SNF1-like kinase 2	DKFZp434K1115, KIAA0781, LOH11CR1I, SIK2	Q9H0K1	NP_056006		
RAF1	Raf1	Raf proto-oncogene serine/threonine-protein kinase		P04049	NP_002871	Recombinant human Raf1(306-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~63
RAFM	Raf1 [Y340E, Y341E]	Raf proto-oncogene serine/threonine-protein kinase		P04049	NP_002871	Recombinant human RAF1(Y340E Y341E, 306-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~63
RET1	Ret	Ret proto-oncogene-encoded kinase	CDHF12, HSCR1, MEN2A, MEN2B, MTC1, PTC, RET-ELE1	P07949	NP_066124		
RIP2	RIPK2	Receptor-interacting serine/threonine-protein kinase 2	RICK, RIP2, CARD3, CARDIAK, CCK, GIG30	O43353	NP_003812	Recombinant human RIPK2 (1-299) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~59
RIP5	RIPK5 (SGK496)	Receptor interacting protein kinase 5, isoform 1	DSTYK; DustyPK; HDCMD38P; KIAA0472; RIP5; SqK496	Q6XUX3	NP_056190	Recombinant full-length human RIPK5 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~140
ROKB	ROCK1 (ROK $\beta$ )	Rho-associated protein kinase 1	P160ROCK, ROCK-I, ROKbeta	Q13464	NP_005397	Recombinant human ROCK1 (17-535) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~85
ROKA	ROCK2 (ROK $\alpha$ )	Rho-associated protein kinase 2	KIAA0619, ROCK-II, ROKalpha	O75116	NP_004841	Recombinant human ROCK2 (5-554) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~88
RON1	RON (MST1R)	Macrophage-stimulating protein receptor	MST1R, PTK8, CDw136	Q04912	NP_002438	Recombinant human RON (983-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
ROS1	ROS	Proto-oncogene tyrosine-protein kinase ROS	c-ros-1; MCF3	P08922	NP_002935	Recombinant human ROS1 (1883-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~82
RSK1	RSK1 (RPS6KA2)	Ribosomal protein S6 kinase alpha 2	RPS6KA1, HU-1, MAPKAPK1A, S6K-alpha 1	Q15418	NP_066958	Recombinant full-length human RSK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~108

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
RSK2	RSK2 (PRS6KA3)	Ribosomal S6 kinase 2 (alpha 3)	RPS6KA3, HU-3, MAPKAPK1B, CLS, MRX19, ISPK-1, p90-RSK2, pp90RSK2, S6K-alpha3	P51812	NP_004577	Recombinant full-length human RSK2 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~112
RSK3	RSK3 (RPS6KA1)	Ribosomal protein S6 kinase alpha 1	RSK, HU-2, RPS6KA2, p90-RSK3, pp90RSK3, MAPKAPK1C, S6K-alpha, S6K-alpha2	Q15349	NP_002944	Recombinant full-length human RSK3 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~112
RSK4	RSK4 (RPS6KA6)	Ribosomal protein S6 kinase alpha 6	RPS6KA6	Q9UK32	NP_055311	Recombinant full-length human RSK4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~112
SGK1	SGK1	Serine-threonine-protein kinase Sgk1	Serum/glucocorticoid-regulated kinase 1	O00141	NP_005618	Recombinant human SGK1 (60-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~73
SGK2	SGK2	Serine-threonine-protein kinase Sgk2	Serum/glucocorticoid-regulated kinase 2; H-SGK2; dJ138B7.2	Q9HBY8	NP_057360	Recombinant full-length human SGK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~71
SGK3	SGK3	Serine-threonine-protein kinase Sgk3	Serum/glucocorticoid-regulated kinase 3; CISK; SGK1	Q96BR1	NP_037389	Recombinant full-length human SGK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~82
SIK1	SIK	Serine-threonine-protein kinase SNF1-like kinase 1	BG719047; SNF1LK; MSK; SIK1	P57059	NP_775490	Recombinant human SIK (1-303) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~36
SKMK	skMLCK (MYLK2)	Myosin light chain kinase 2, skeletal/cardiac muscle	skMLCK, KMLC, MLCK, MLCK2	Q9H1R3	NP_149109	Recombinant full-length human MYLK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~74
SLK1	SLK	CTCL tumor antigen se20-9	Ste20-related serine/threonine kinase; LOSK; STK2; se20-9; KIAA0204; MGC133067; bA16H23.1	Q9H2G2	NP_055535	Recombinant full-length human SLK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~180
SMMK	smMLCK (MLCK2)	Myosin light chain kinase, smooth muscle and non-muscle isozymes	MYLK, MLCK2, KRP, MLCK108, MLCK210, MSTP083, FLJ12216, DKFZp686I10125	Q15746	NP_444253	Recombinant human MLCK (1425-1776) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~70

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
SRC1	Src	Proto-oncogene tyrosine-protein kinase Src	ASV, SRC1, c-SRC, p60-Src	P12931	NP_005408	Recombinant full-length human SRC was expressed in E.coli cells using an N-terminal GST tag.	~83
SRP1	SRPK1	SFRS protein kinase 1	SFRSK1	Q96SB4	NP_003128	Recombinant full-length human SRPK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~130
SRP2	SRPK2	SFRS protein kinase 2	FLJ36101, SFRSK2	P78362	NP_872633	Recombinant full-length human SRPK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~135
ST33	STK33	Serine/threonine-protein kinase 33		Q8NEF5	NP_112168	Recombinant full-length human STK33 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~94
SYK1	SYK	Tyrosine-protein kinase SYK	Spleen tyrosine kinase	P43405	NP_003168	Recombinant full-length human SYK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~100
TAK1	TAK1-TAB1	Mitogen-activated protein kinase kinase kinase 7	Transforming growth factor-beta-activated kinase 1; TGF-beta-activated kinase 1; MAP3K7	O43318	NP_003179	Recombinant human proteins TAK1 (1-303) and TAB1 (437-end), linked by a small peptide (DFGGGGG), were expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~74
TAO1	TAO1 (TAOK1)	Serine/threonine-protein kinase TAO1	PSK2, MARKK, MAP3K16, FLJ14314, KIAA1361	Q7L7X3	NP_065842	Recombinant human TAOK1 (1-314) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
TAO2	TAO2 (TAOK2)	Serine/threonine-protein kinase TAO2	PSK, PSK1, TAO1, TAO2, MAP3K17, KIAA0881	Q9UL54	NP_004774	Recombinant human TAOK2 (1-314) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~63
TAO3	TAO3 (TAOK3, JIK )	TAO kinase 3	DKFZp666H245, DPK, FLJ31808, JIK, MAP3K18, TAOK3	Q9H2K8	NP_057365		
TBK1	TBK1	TANK-binding kinase 1	NFkB-activating kinase, NAK, T2K, FLJ11330	Q9UHD2	NP_037386	Recombinant full-length human TBK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~105
TEC1	TEC	Tec protein-tyrosine kinase	MGC126760, MGC126762, PSCTK4	P42680	NP_003206		
TES1	TESK1	Dual-specificity testis-specific protein kinase 1	Testicular protein kinase 1	Q15569	NP_006276		

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
TGF1	TGFβR1	Transforming growth factor, beta receptor 1	AAT5, ACVRLK4, ALK-5, ALK5, LDS1A, LDS2A, SKR4; TGFR1	P36897	NP_004603	Recombinant human TGFβR1 (80-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~65
TGF2	TGFβR2	TGF-beta receptor type II	AAT3, FAA3, MFS2, RIIC, HNPCC6, TGFR-2, TGFbeta-RII, TAAD2	P37173	NP_003233	Recombinant human TGFβR2 (190-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~68
TIE2	Tie2 (Tek)	Angiopoietin 1 receptor	TEK, VMCM, VMCM1, CD202B	Q02763	NP_000450	Recombinant human TIE 2 (771-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~61
TLK2	TLK2	Serine-threonine-protein kinase tousled-like 2	MGC44450; PKU-ALPHA; Tousled-like kinase 2	Q86UE8	NP_006843	Recombinant human TLK2 (388-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~67
TRKA	TrkA (NTRK1)	High affinity nerve growth factor receptor	NTRK1, MTC, TRK, TRK1, p140-TrkA, DKFZp781I14186	P04629	NP_002520	Recombinant human TRKA (440-end) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~66
TRKB	TrkB (NTRK2)	BDNF/NT-3 growth factors receptor	Neurotrophic tyrosine kinase receptor type 2; NTRK2, GP145-TrkB	Q16620	NP_001018074	Recombinant human TRKB (455-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~67
TRKC	TrkC	NT-3 growth factor receptor	Neurotrophic tyrosine kinase receptor type 3; NTRK3, gp145(trkC)	Q16288	NP_002521	Recombinant human TRKC (507-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag.	~38
TSS1	TSSK1 (STK22D)	Testis-specific serine kinase 1B	FKSG81, SPOGA4, STK22D, TSSK1	Q9BXA7	NP_114417		
TSS2	TSSK2 (STK22B)	Testis-specific serine/threonine protein kinase 2	DGS-G, SPOGA2, STK22B, FLJ38613	Q96PF2	NP_443732	Recombinant full-length human TSSK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~66
TTK1	TTK	Dual specificity protein kinase TTK	Phosphotyrosine picked threonine kinase; ESK; PYT; MPS1; MPS1L1; FLJ38280	P33981	NP_003309	Recombinant full-length human TTK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~130
TXK1	TXK	Tyrosine-protein kinase TXK	RLK, TKL, BTKL, PTK4, PSCTK5, MGC22473	P42681	NP_003319	Recombinant human TXK (239-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~53

Kinexus ID Code	Protein Kinase Name	Full Name	Alias	UniProt	Refseq	Description	Recombinant Mol. Mass (KDa)
TYK2	TYK2	Tyrosine kinase 2	JTK1	P29597	NP_003322		~125
TYR3	TYRO3 (RSE )	Tyrosine-protein kinase receptor TYRO3	BYK, RSE, Dtk, Sky, Tif, Brt	Q06418	NP_006284	Recombinant human TYRO3 (455-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~77
ULK1	ULK1	Serine-threonine-protein kinase ULK1	ATG1; FLJ38455; Unc-51-like kinase 1; UNC51; Unc51.1	O75385	NP_003556	Recombinant human ULK1 (1-649) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~125
ULK2	ULK2	Serine-threonine-protein kinase ULK2	KIAA0623; Unc-51-like kinase 2; Unc51.2	Q8IYT8	NP_055498	Recombinant human ULK2 (1-631) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~98
WEE1	WEE1	WEE1 homologue	DKFZp686I18166, FLJ16446, WEE1A, WEE1hu	P30291	NP_003381		
YES1	Yes (Yes1)	v-Yes-1 Yamaguchi sarcoma viral oncogene homologue 1	HsT441, P61-YES, Yes1, c-yes	P07947	NP_005424	Recombinant full length human YES1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~91
YSK1	YSK1 (STK25, SOK1)	Serine/threonine kinase 25	DKFZp686J1430, SOK1, STK25	O00506	NP_006365	Recombinant full length human YSK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~76
ZAK1	ZAK	Sterile alpha motif and leucine zipper containing kinase AZK	AZK, MLK7, MLT, MLTK, MRK, mlklak	Q9NYL2	NP_057737	Recombinant full-length human ZAK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag.	~82
ZA70	ZAP70	Tyrosine-protein kinase ZAP-70	SRK, STD, ZAP-70	P43403	NP_001070	Recombinant full-length human ZAP70 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~96
ZC03	ZC3 (MINK1)	Mitogen-activated protein kinase kinase kinase 6	B55, ZC3, MINK, YSK2, hMIN, MAP4K6, MGC21111, hMINK $\beta$	Q8N4C8	NP_056531	Recombinant human MINK1 (1-320) was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag.	~61

Form: **KICP-SOF-01****KINEXUS ORDER NUMBER****KINASE INHIBITOR  
COMPOUND PROFILING SERVICE ORDER FORM****CUSTOMER INFORMATION** ☐ REPEAT CUSTOMER **OR** ☐ NEW CUSTOMER☐ Dr. ☐ Mr. ☐ Ms.

Name of Authorized Representative or Principal Investigator

Title/Position

Company Name or Institute

Department

Street Address

City

State or Province

Country

Zip or Postal Code

Email Address

(Area Code)

Telephone Number

(Area Code)

Facsimile Number

Contact Person (if different from Authorized Representative)

Email Address

(Area Code)

Telephone Number

**KINASE INHIBITOR COMPOUND PROFILING REPORTS**RESULTS SENT BY EMAIL TO: ☐ AUTHORIZED REPRESENTATIVE/INVESTIGATOR AND/OR ☐ CONTACT PERSON**REQUESTED WORK AND PRICING INFORMATION**

Refer to Service Identification Forms (KICP-SIF-01A and KICP-SIF-01B)

All prices in U.S. Funds

Sheet No.	KICP-SIF ID Name (from Section A of KICP-01 Forms)	Total Number of Assays (from Section B of KICP-01 Forms)
1.	_____	_____
2.	_____	_____
3.	_____	_____
4.	_____	_____
5.	_____	_____
6.	_____	_____
7.	_____	_____
8.	_____	_____

Total Number Assays (Sum 1-8) \_\_\_\_\_

(Complete up to this point for quotation purposes only)

Quotation or Reference Number: \_\_\_\_\_

TOTAL COST AS PER QUOTATION = \$ \_\_\_\_\_

FOR CANADIAN CUSTOMERS ONLY:

Add an additional 5% to the above total for GST (No. 893907329 RT0001): + \$ \_\_\_\_\_ = \$ \_\_\_\_\_

TOTAL AMOUNT PAYABLE IN U.S FUNDS

**PAYMENT METHOD**☐ PURCHASE ORDER ACCEPTED FROM COMPANIES AND INSTITUTES WITH APPROVED CREDIT. P.O. NUMBER: \_\_\_\_\_☐ VISA **OR** ☐ MASTERCARD

Print Cardholder Name

Visa Number

Expires (M/Y)

Cardholder Signature

**BILLING INFORMATION**☐ SEND INVOICE TO CUSTOMER AT ABOVE ADDRESS **OR** ☐ SEND INVOICE TO ACCOUNTS PAYABLE CONTACT:☐ Dr. ☐ Mr. ☐ Ms.

Accounts Payable Contact Name

Company Name or Institute

Street Address

City

State or Province

Country

Zip or Postal Code

(Area Code)

Telephone Number

**AUTHORIZATION**

CUSTOMER HAS READ THE KINEXUS SERVICE AGREEMENT AND AGREES TO BE BOUND BY THE TERMS AND CONDITIONS:

Print Name of Authorized Representative or Principal Investigator

Authorized Signature

Date (m/d/y)

How did you originally hear about the KICP Services?

☐ Direct Mail☐ Email☐ Web Site☐ Advertisement☐ Referral☐ Conference or Trade Show☐ Other





Form: KICP-SIF-01A

**KINASE INHIBITOR  
COMPOUND PROFILING****SERVICE INFORMATION FORM**

Subject to terms of the Kinexus Service Agreement

**KINEXUS ORDER NUMBER**NAME: \_\_\_\_\_ COMPANY/INSTITUTE: \_\_\_\_\_  
(Authorized Representative or Principal Investigator)**Confidential Service Requested: One Kinase; Multiple Numbers and/or Dilutions of Compounds**

Use the KICP-SIF-01B Form if you wish to have many kinases tested against only one compound. Please refer to the Kinase Inhibitor Compound Profiling Customer Information Package for further details about this in vitro custom compound screening service. Initially use this form to obtain a quotation from Kinexus for pricing of your custom order. Please also provide a copy of this completed form at the time of submission of your compounds for analysis. Use additional copies of this form if you have more compounds or a larger number of kinases to be tested. If you need assistance completing this form, contact a technical service representative by calling toll free in North America 1-866-KINEXUS (866-546-3987) or by email at [info@kinexus.ca](mailto:info@kinexus.ca).

<b>CUSTOM SERVICE REQUESTED: KICP-1.0</b> Custom multi-compound (1-15) and single target kinase (1) screen  <i>Depending on the scale of your order, make sure that you are supplying at least five-times as much as minimally necessary of the compound to be tested..</i>		<b>KINEXUS ID NUMBER</b> (Bar Code Identification Number)  For Kinexus Internal Use Only.		<b>A. KICP-SIF IDENTIFICATION NAME:</b>  Client ID: _____  <i>Use this ID name of your choice for your internal reference and completion of the KICP-SOF form. This is useful when multiple copies of this KICP-SIF forms are to be used in your order.</i>						
<b>B. COMPOUND SELECTION:</b>  <i>For each separate compound, under "Compound ID" please provide the assigned client name that has been entered into Box A from the completed and attached "Confidential Sample Description" form (KICP-CSDP-01). For each concentration of the compound to be tested (up to 6) indicate the amount and whether this represents "µM" or "µg/ml" by ticking the appropriate box. Also indicate whether each assay is to be performed in single ("Once"), duplicate ("Dupl.") or triplicate ("Tripl.") measurements. Under "Subtotal" for each compound, indicate the number of individual assays that you intend to be performed.</i>				<b>C. KINASE SELECTION:</b>  <i>Use Appendix A of the Kinase Inhibitor Compound Profiling Customer Information Package to obtain the code (Column A) and name (Column B) of the protein kinase that will be tested for inhibition by the compounds listed in Section B. A more detailed MS-Excel spreadsheet with information about each kinase can be downloaded from the Kinexus website at <a href="http://www.kinexus.ca/services/kinase_inhibitor_profiling.php">www.kinexus.ca/services/kinase_inhibitor_profiling.php</a>.</i>						
Compound ID.	Final Conc. (µM) <input type="checkbox"/> or (µg/ml) <input type="checkbox"/>						Replicates			Subtotal
	Dil. 1	Dil. 2	Dil. 3	Dil. 4	Dil. 5	Dil. 6	Once	Dupl.	Tripl.	
1. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
2. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
3. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
4. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
5. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
6. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
7. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
8. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
9. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
10. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
11. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
12. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
13. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
14. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
15. _____							<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Total Number of Assays _____										
<b>D. ASSAY TIME:</b>  <i>If you have a recommended or required incubation time, please provide this here.</i> Assay time (minutes): _____										
<b>E. PRICING:</b>  <i>Pricing varies with the number of kinases and compounds to be tested in each custom order. It is also feasible for Kinexus to test compounds against protein kinases that are provided by clients. Therefore, it is necessary to complete this form and transmit it by facsimile to 1-604-323-2548 with your e-mail address, facsimile and telephone number to obtain a quotation for each custom analysis. We will endeavor to issue a quotation that is valid for four weeks within 24 hours of receipt.</i>										
<b>F. SPECIAL INSTRUCTIONS:</b>  _____ _____ _____ _____ _____										

Name of person completing this form

Email Address/Facsimile Number/ Phone Number

Date (m/d/y)



Form: KICP-SIF-01B

**KINASE INHIBITOR  
COMPOUND PROFILING****SERVICE INFORMATION FORM**

Subject to terms of the Kinexus Service Agreement

**KINEXUS ORDER NUMBER**NAME: \_\_\_\_\_ COMPANY/INSTITUTE: \_\_\_\_\_  
(Authorized Representative or Principal Investigator)**Confidential Service Requested: One Compound at One or More Dilutions; Multiple Kinases**

Use the KICP-SIF-01A Form instead if you wish to have many compounds tested against only one kinase. Please refer to the Kinase Inhibitor Compound Profiling Customer Information Package for further details about this in vitro custom compound screening service. Initially, use this form to obtain a quotation from Kinexus for pricing of your custom order. Please also provide a copy of this completed form at the time of submission of your compounds for analysis. Use additional copies of this form if you have more compounds or a larger number of kinases to be tested. If you need assistance completing this form, contact a technical service representative by calling toll free in North America 1-866-KINEXUS (866-546-3987) or by email at [info@kinexus.ca](mailto:info@kinexus.ca).

<b>CUSTOM SERVICE REQUESTED: KICP-1.0</b> Custom multi-kinase (1-15) and single compound (1) screen  <i>Depending on the scale of your order, make sure that you are supplying at least three-times as much as minimally necessary of the compound to be tested.</i>	<b>KINEXUS ID NUMBER</b> (Bar Code Identification Number)  For Kinexus Internal Use Only.	<b>A. KICP-SIF IDENTIFICATION NAME:</b>  Client ID: _____  <i>Use this ID name of your choice for your internal reference and completion of the KICP-SOF form. This is useful when multiple copies of this KICP-SIF forms are to be used in your order.</i>																																																																																																																																																															
<b>C. KINASE SELECTION:</b>  <i>Use Appendix A of the Kinase Inhibitor Compound Profiling Customer Information Package to obtain the code (Column A) and name (Column B) of the protein kinase that will be tested for inhibition by the compounds listed in Section B. A more detailed MS-Excel spreadsheet with information about each kinase can be downloaded from the Kinexus website at <a href="http://www.kinexus.ca/services/kinase_inhibitor_profiling.php">www.kinexus.ca/services/kinase_inhibitor_profiling.php</a>. For each concentration of the compound to be tested (up to 6) indicate the amount and whether this represents "µM" or "µg/ml" by ticking the appropriate box. Also indicate whether each assay is to be performed in single ("Once"), duplicate ("Dupl.") or triplicate ("Tripl.") measurements. Under "Subtotal" for each compound, indicate the number of individual assays that you intend to be performed.</i>		<b>B. COMPOUND SELECTION:</b>  <i>Please provide the assigned client name that has been entered into Box A from the completed and attached "Confidential Sample Description" form (KICP-CSDF-01).</i>  Compound ID: _____																																																																																																																																																															
<table border="1"><thead><tr><th rowspan="2">Kinase Code + Name</th><th colspan="6">Compound Final Conc. (µM) <input type="checkbox"/> or (µg/ml) <input type="checkbox"/></th><th rowspan="2">Replicates Once Dupl. Tripl.</th><th rowspan="2">Subtotal</th></tr><tr><th>Dil. 1</th><th>Dil. 2</th><th>Dil. 3</th><th>Dil. 4</th><th>Dil. 5</th><th>Dil. 6</th></tr></thead><tbody><tr><td>1. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>2. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>3. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>4. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>5. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>6. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>7. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>8. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>9. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>10. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>11. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>12. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>13. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>14. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td>15. _____</td><td></td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/></td><td></td></tr><tr><td colspan="8">Total Number of Assays _____</td><td></td></tr></tbody></table>		Kinase Code + Name	Compound Final Conc. (µM) <input type="checkbox"/> or (µg/ml) <input type="checkbox"/>						Replicates Once Dupl. Tripl.	Subtotal	Dil. 1	Dil. 2	Dil. 3	Dil. 4	Dil. 5	Dil. 6	1. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		2. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		3. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		4. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		5. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		6. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		7. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		8. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		9. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		10. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		11. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		12. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		13. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		14. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		15. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>		Total Number of Assays _____									<b>D. ASSAY TIME:</b>  <i>If you have a recommended or required incubation time, please provide this here.</i>  Assay time (minutes): _____
Kinase Code + Name	Compound Final Conc. (µM) <input type="checkbox"/> or (µg/ml) <input type="checkbox"/>						Replicates Once Dupl. Tripl.	Subtotal																																																																																																																																																									
	Dil. 1	Dil. 2	Dil. 3	Dil. 4	Dil. 5	Dil. 6																																																																																																																																																											
1. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
2. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
3. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
4. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
5. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
6. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
7. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
8. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
9. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
10. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
11. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
12. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
13. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
14. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
15. _____							<input type="checkbox"/> <input type="checkbox"/> <input type="checkbox"/>																																																																																																																																																										
Total Number of Assays _____																																																																																																																																																																	
		<b>E. PRICING:</b>  <i>Pricing varies with the number of kinases and compounds to be tested in each custom order. It is also feasible for Kinexus to test compounds against protein kinases that are provided by clients. Therefore, it is necessary to complete this form and transmit it by facsimile to 1-604-323-2548 with your e-mail address, facsimile and telephone number to obtain a quotation for each custom analysis. We will endeavor to issue a quotation that is valid for four weeks within 24 hours of receipt.</i>																																																																																																																																																															
		<b>F. SPECIAL INSTRUCTIONS:</b>  _____ _____ _____ _____ _____ _____ _____																																																																																																																																																															



Form: KICP-CSDf-01

**KINASE INHIBITOR  
COMPOUND PROFILING****CLIENT SUPPLIED  
CONFIDENTIAL SAMPLE DESCRIPTION FORM***Subject to terms of the Kinexus Service Agreement***KINEXUS ORDER NUMBER**NAME: \_\_\_\_\_ COMPANY/INSTITUTE: \_\_\_\_\_  
(Authorized Representative or Principal Investigator)**Compound Details:**

Please refer to the Kinase Inhibitor Compound Profiling Customer Information Package for details on how to prepare and ship your compounds to Kinexus for testing. Clients are required to complete Sections A and B for a confidential analysis with this in vitro screening service. Use additional copies of this form if you have more than two compounds for testing. For each compound to be tested, please send enough material to complete the analysis. If you need assistance completing this form, contact a technical service representative by calling toll free in North America 1-866-KINEXUS (866-546-3987) or by email at [info@kinexus.ca](mailto:info@kinexus.ca).

<b>A. CLIENT IDENTIFICATION NAME</b>  CLIENT ID: _____  <i>Use the Client ID Name that you entered here in Box A on the Custom Kinase Inhibitor Compound Profiling Service Identification Form" (KICP-SIF-01A or KICP-SIF-01B). Provide a separate Client ID number for each different compound to be tested.</i>	<b>KINEXUS ID NUMBER</b> (FOR INTERNAL USE ONLY)  <i>(Bar Code Identification Number)</i>
<b>B. COMPOUND DETAILS:</b> Compound name (not required): _____ Supplied form of compound <input type="checkbox"/> Solid <input type="checkbox"/> Liquid  <b>FOR SOLIDS PROVIDE:</b> Mass (mg): _____ Formula Weight: _____ Is the compound or solution toxic? <input type="checkbox"/> Yes <input type="checkbox"/> No  <b>FOR LIQUIDS PROVIDE:</b> Molarity: _____ or Concentration: _____ Volume: _____ MSDS or safety sheets provided <input type="checkbox"/> Yes <input type="checkbox"/> No  For solids, what is the recommended solvent? _____ Is the compound soluble in water or 2% DMSO? _____  Provide safety instructions and storage details for handling, or any other special information: _____ _____ _____ _____	

<b>A. CLIENT IDENTIFICATION NAME</b>  CLIENT ID: _____  <i>Use the Client ID Name that you entered here in Box A on the Custom Kinase Inhibitor Compound Profiling Service Identification Form" (KICP-SIF-01). Provide a separate Client ID number for each different compound to be tested.</i>	<b>KINEXUS ID NUMBER</b> (FOR INTERNAL USE ONLY)  <i>(Bar Code Identification Number)</i>
<b>B. COMPOUND DETAILS:</b> Compound name (not required): _____ Supplied form of compound <input type="checkbox"/> Solid <input type="checkbox"/> Liquid  <b>FOR SOLIDS PROVIDE:</b> Mass (mg): _____ Formula Weight: _____ Is the compound or solution toxic? <input type="checkbox"/> Yes <input type="checkbox"/> No  <b>FOR LIQUIDS PROVIDE:</b> Molarity: _____ or Concentration: _____ Volume: _____ MSDS or safety sheets provided <input type="checkbox"/> Yes <input type="checkbox"/> No  For solids, what is the recommended solvent? _____ Is the compound soluble in water or 2% DMSO? _____  Provide safety instructions and storage details for handling, or any other special information: _____ _____ _____ _____	

I hereby certify that all the sample information provided in this order is correct and accurate to the best of my knowledge. I further acknowledge that I may be contacted by a Kinexus representative for additional details if the information provided is unclear.

\_\_\_\_\_  
Name of person completing this form\_\_\_\_\_  
Signature\_\_\_\_\_  
Date (m/d/y)

# COMMERCIAL INVOICE

<b>DATE OF EXPORTATION</b>	<b>EXPORT REFERENCES</b> (not required)
<b>SHIPPER/EXPORTER</b>	<b>CONSIGNEE</b>  Kinexus Bioinformatics Corporation Suite 1, 8755 Ash Street Vancouver, B.C. Canada V6P 6T3  Telephone: (604) 323-2547 Ext. 1 or 10 Facsimile: (604) 323-2548 Email: info@kinexus.ca
<b>COUNTRY OF EXPORT</b>	<b>TERMS OF SALE</b> Not for resale, sample for analysis
<b>COUNTRY OF ORIGIN</b>	<b>PURPOSE</b> Research and development
<b>COUNTRY OF ULTIMATE DESTINATION</b> Canada	<b>EXPORTING CARRIER</b>
<b>INTERNATIONAL AIR WAYBILL NUMBER</b>  Courier Number:	

NO. OF PKGS	TYPE OF PACKAGING	QUANTITY OF SAMPLES	COMPLETE AND ACCURATE COMMODITY DESCRIPTION	UNIT VALUE
	<input type="checkbox"/> FedEx Letter <input type="checkbox"/> FedEx Pak <input type="checkbox"/> Box <input type="checkbox"/> Other	Total number of 1.5 ml Eppendorf tubes:	Non-hazardous, compounds for research and development for therapeutic purposes. Samples are not for resale and there is no commercial value.	\$1.00 per sample
<b>TOTAL NO. OF PACKAGES</b>		<b>TOTAL WEIGHT OF PACKAGES</b>		<b>TOTAL DECLARED VALUE</b>
				\$

*These commodities were exported from the Country indicated above in accordance with the Export Administration Regulations and are licensed for the ultimate designation shown. It is hereby certified that this commercial invoice shows the actual price of the goods described, that no other invoice has been or will be issued for these goods, and that all particulars are true and correct.*

## SIGNATURE AND STATUS OF AUTHORIZED PERSON

_____	_____
<i>Print Name</i>	<i>Title</i>
_____	_____
<i>Authorized Signature</i>	<i>Date (month/day/year)</i>

INCLUDE THREE (3) COPIES OF THIS INVOICE WITH YOUR SHIPMENT

# COMMERCIAL INVOICE

DATE OF EXPORTATION	EXPORT REFERENCES
SHIPPER/EXPORTER	<b>CONSIGNEE</b>  Kinexus Bioinformatics Corporation Suite 1, 8755 Ash Street Vancouver, B.C. Canada V6P 6T3  Telephone: (604) 323-2547 Ext. 1 or 10 Facsimile: (604) 323-2548 Email: info@kinexus.ca
COUNTRY OF EXPORT	<b>TERMS OF SALE</b> Not for resale, sample for analysis
COUNTRY OF ORIGIN	<b>PURPOSE</b> Research and development
COUNTRY OF ULTIMATE DESTINATION Canada	<b>EXPORTING CARRIER</b>
<b>INTERNATIONAL AIR WAYBILL NUMBER</b> Courier Number:	

NO. OF PKGS	TYPE OF PACKAGING	QUANTITY OF SAMPLES	COMPLETE AND ACCURATE COMMODITY DESCRIPTION	UNIT VALUE
	— FedEx Letter — FedEx Pak — Box — Other	<i>Total number of 1.5 ml Eppendorf tubes:</i>	Non-hazardous, compounds for research and development for therapeutic purposes. Samples are not for resale and there is no commercial value.  Samples are packaged on Dry Ice, Class 9, UN 1845, Group 3 (____ X ____ kgs).	<b>\$1.00</b> <i>per sample</i>
<b>TOTAL NO. OF PACKAGES</b>			<b>TOTAL WEIGHT OF PACKAGES</b>	<b>TOTAL DECLARED VALUE</b>
				\$

*These commodities were exported from the Country indicated above in accordance with the Export Administration Regulations and are licensed for the ultimate designation shown. It is hereby certified that this commercial invoice shows the actual price of the goods described, that no other invoice has been or will be issued for these goods, and that all particulars are true and correct.*

## SIGNATURE AND STATUS OF AUTHORIZED PERSON

_____	_____
<i>Print Name</i>	<i>Title</i>
_____	_____
<i>Authorized Signature</i>	<i>Date (month/day/year)</i>

INCLUDE THREE (3) COPIES OF THIS INVOICE WITH YOUR SHIPMENT



## PROTEOMICS SERVICES AGREEMENT

SERVICE AGREEMENT NO.

This Agreement is entered into effective as of the Effective Date by and between Kinexus Bioinformatics Corporation (“**Kinexus**”), a Canadian corporation with a principal place of business at Suite 1, 8755 Ash Street, Vancouver, British Columbia, Canada, V6P 6T3 **AND** the corporation or other entity (“**Customer**”) having the following name and business or institution address: \_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

### RECITALS

**WHEREAS** Kinexus is a bioinformatics company employing proprietary proteomics and bioinformatics services to create and interpret data to map protein signalling networks and compile databases with this knowledge to enable disease biomarker and therapeutics discovery.

**WHEREAS** the Customer desires to have Kinexus perform standard and/or customized proteomics services with materials and/or information provided by the Customer.

**WHEREAS** Kinexus is willing to provide these proteomics services under the terms and conditions set forth herein.

**THEREFORE**, in consideration of the premises and covenants and agreements contained herein, and other good and valuable consideration the receipt and sufficiency of which is hereby acknowledged, Kinexus and the Customer agree as follows:

#### 1. DEFINITIONS

1.1 “Academic Collaborator” means a principal investigator, employed at a university or other not-for-profit academic research institution.

1.2 “Affiliate” means any corporation or other entity that directly or indirectly controls, is controlled by or is under common control with a party to this Agreement. A corporation or other entity shall be regarded as in control of another corporation or entity if it owns or directly or indirectly controls more than fifty percent (50%) of the outstanding voting stock or other ownership interest of the other corporation or entity.

1.3 “Corporate Partner” means any Third Party which enters into an agreement with the Customer or its Affiliates involving the grant to such Third Party of rights for the development or commercialization of a product that was discovered, identified, selected, characterized or determined to have therapeutic or diagnostic use through use of the Proteomics Analyses provided to the Customer pursuant to this Agreement.

1.4 “Confidential Information” means any information or data received by a party (the “Receiving Party”) from the other party (the “Disclosing Party”) in connection with the performance of this Agreement that, if

disclosed in writing, is marked or otherwise identified by the Disclosing Party as confidential or, if disclosed orally is identified in writing by the Disclosing Party as confidential within ten (10) days following the disclosure. Confidential Information shall not include any information or data that the Receiving Party can demonstrate:

- (a) was generally available to the public before its disclosure to the Receiving Party or became generally available to the public after its disclosure to the Receiving Party, provided that such information or data did not become generally available to the public by means of an unauthorized act or omission of the Receiving Party;
- (b) was already in the possession of the Receiving Party before its disclosure under this Agreement, as demonstrated by Receiving Party's written records, provided that such information or data was not obtained directly or indirectly from the Disclosing Party under an obligation of confidentiality;
- (c) was disclosed to the Receiving Party, whether before or after its disclosure under this Agreement, by a Third Party, provided that such information or data was not obtained directly or indirectly from the Disclosing Party under an obligation of confidentiality; or
- (d) was independently developed or discovered by employees or agents of the Receiving Party without any use of Confidential Information of the Disclosing Party as demonstrated by Receiving Party's written records.

All of the Proteomics Services technologies provided by Kinexus will be deemed to have been identified as proprietary and considered the Confidential Information of Kinexus.

1.5 "Contact" means the contact person of the Customer that is designated on the Service Order Forms, who is deemed to have the authority to deliver Samples, Service Order Forms, Service Information Forms, and Sample Description Forms to Kinexus, on behalf of the Customer, under this Agreement.

1.6 "Proteomics Analyses" means one or more of the Custom and Standard Proteomics Services offered by Kinexus that may permit the identification and/or quantification of proteins, their phosphorylation states, their interactions with proteins, peptides, and other compounds, and the regulation of their functional activities by these agents.

1.7 "Proteomics Products" means the products of the Custom Proteomics Services offered by Kinexus to manufacture one or more proteins using recombinant DNA technology, and designer peptides by chemical synthesis.

1.8 "Sample" means a lysate or semi-purified fraction from cells and tissues, a protein, and/or a compound provided to Kinexus by the Customer, which the Customer has prepared and shipped in a manner that it can be properly used by Kinexus for the Proteomics Analyses. Samples for Proteomics Analyses may also be provided by Kinexus at the request of the Customer.

1.9 "Sample Description Form" means the Kinexus form to be completed by the Customer to provide information on the nature of each Sample submitted for the Proteomics Analyses. It is included in the Proteomics Services Customer Information Package with this Agreement, and may be amended from time to time as updated on the Kinexus website.

1.10 "Antibody" means the immunoglobulin reagent that permits detection of a target protein or phosphorylation site.

1.11 "Antibody Description Form" means the Kinexus form to be completed by the Customer to provide information on the nature of each Antibody submitted by the Customer for the Proteomics Analyses. It is included

in the Proteomics Services Customer Information Package with this Agreement, and may be amended from time to time as updated on the Kinexus website.

1.12 "Service Order Form" means the Kinexus form to be completed by the Customer to provide Kinexus with the Customer's contact and billing information for the Proteomics Analyses or Proteomics Products. This form indicates the level of confidentiality requested by the Customer. It is included in the Proteomics Services Customer Information Package with this Agreement, and may be amended from time to time as updated on the Kinexus website.

1.13 "Service Information Form" means the Kinexus form to be completed by the Customer to provide Kinexus with a specific listing of the Samples to be tested for the Proteomics Analysis or a specific description of the Proteomics Products that are requested. It is included in the Proteomics Services Customer Information Package with this Agreement, and may be amended from time to time as updated on the Kinexus website.

1.14 "Report" means the underlying raw data and the report provided to The Customer hereunder consisting of the Proteomic Analyses of Samples, including, but not limited to tables of the experimental results. For Proteomics Products, the Report may include raw data confirming the composition and purity of the Proteomics Products.

1.15 "Field of Use" means use by Kinexus and its Affiliates and Academic Collaborators of data from the Report for research and commercial purposes relating to the creation and interpretation of knowledge about the composition, architecture and operation of cell signalling networks, improving its Proteomics Services, and the compilation of databases that may become accessible to Third Parties on-line over the Internet.

1.16 "Third Party" means any entity other than Kinexus', Kinexus' Affiliates, the Customer and the Customer's Affiliates.

1.17 "Effective Date" means the date of the last signature on this Agreement.

## **2. REQUEST FOR AND DELIVERY OF PROTEOMICS SERVICES**

2.1 Request for Proteomics Services. From time to time, over the Term of this Agreement (as defined in Section 6.1 herein), the Customer can engage Kinexus to provide its Proteomics Analyses or Proteomics Products. After submission of a quotation from Kinexus to the Customer, by delivery to Kinexus of a Service Order Form, a Service Information Form and a Sample Description Form with Samples as appropriate, the Customer hereby requests and authorizes Kinexus to perform Proteomics Services and deliver the results of these services to the Customer, pursuant to the terms and conditions in this Agreement. In the case of Customer requested Proteomics Analyses, this would include the delivery of a Report. In the case of Customer requested Proteomics Products, this would include the delivery of the Proteomics Products and a Report.

2.2 Representation and Warranty. The Customer represents and warrants that: (a) it has all right and authority to provide the Sample to Kinexus for analysis under the terms and conditions of this Agreement, (b) it collected the Sample lawfully and with all necessary consents and approvals, and (c) that the collection, use and disclosure of the Sample by Kinexus pursuant to this Agreement will not violate the rights of any Third Party.

2.3 Delivery Conditions for Customer Sample. The Customer shall be responsible for making shipping arrangements to deliver Samples to Kinexus. The Customer shall also be responsible for complying with all applicable laws and regulations (including but not limited to customs requirements and relevant handling procedures and protocols) and obtaining any and all permits, forms or permissions that may be required by all regulatory authorities to ship and deliver the Sample; to Kinexus and for Kinexus to accept delivery of the Sample.



2.4 Processing and Delivery of Report and Proteomics Products. Subject to the terms of this Agreement, Kinexus shall analyze Samples with the Customer-specified Proteomics Services or produce Customer-specified Proteomics Products, and deliver a Report to the Customer as requested on the Service Order Form and Service Information Form.

2.5 Quality of Samples for Proteomics Analyses. Kinexus shall not deliver a Report on any Sample that Kinexus, in its sole discretion, believes has not been prepared and delivered in a manner that would compromise its ability to provide a reliable result. Under such a circumstance, the Sample will be destroyed by Kinexus after ten (10) days notification by e-mail to the Customer or at the request of the Customer prior to the scheduled destruction of the Sample, it will be returned to the Customer provided that the Customer agrees to reimburse Kinexus for the courier costs for its delivery.

### 3. PAYMENTS

3.1 Payments for Proteomics Services. For each Proteomics Analyses and Proteomics Product requested under this Agreement, the Customer shall pay to Kinexus a fee in accordance with the amount specified on the Service Order Form and the Service Identification Form for the requested service, which may be amended from time to time as updated on Kinexus' website. This amount will be based on a formal quotation issued by Kinexus to the Customer. In the absence of a formal quotation, the pricing will be based on the pricing specified in the latest versions of the Customer Information Packages for Proteomics Services that are downloadable from the Kinexus website ([www.kinexus.ca](http://www.kinexus.ca)). The category of pricing depends on the level of requested confidentiality for analysis:

- (a) Non-Confidential Analyses. If the Samples are provided by the Customer, then all of the Sample information on the Client Supplied **Non-Confidential** Sample Description Form is completed and **is not** designated as Confidential Information on the Service Identification Form. If Antibodies are supplied by the Customer, then all of the Antibody information on the Client Supplied Antibody Description Form (see example in Appendix) must be completed and **is not** designated as Confidential Information on the Service Identification Form.
- (b) Confidential Analyses. If the Samples are provided by the Customer, then all of the Sample information on the Client Supplied **Confidential** Sample Description Form must be completed and **is** designated as Confidential Information on the Service Identification Form.

3.2 The Customer shall issue a purchase order or provide a charge account at the time the Customer sample arrives at Kinexus' offices at Suite 1, 8755 Ash Street, Vancouver, British Columbia, Canada, V6P 6T3. Kinexus will invoice Customer when the Proteomics Analyses or Proteomics Products are complete and delivered to Customer. Payment terms are net 30 days from date of invoice.

3.3 Interest on Late Payments. Any overdue payments by the Customer to Kinexus under this Agreement shall bear interest, to the extent permitted by applicable law at 18% per annum, calculated on the total number of days payment is delinquent; provided, however, that interest shall not accrue pursuant to this Section 3.3 on any amounts payable under this Agreement with respect to which payment is disputed in good faith; provided, further that interest shall accrue pursuant to this Section 3.3 once such dispute has been resolved if payment is not made promptly thereafter.

## 4. INTELLECTUAL PROPERTY RIGHTS

4.1 Ownership of Sample Information. The Customer owns all rights to the Sample information provided to Kinexus. For Non-Confidential Proteomics Analyses, the Customer grants Kinexus a non-exclusive, royalty-free fully paid up worldwide perpetual license to use, copy, publish, compile, display, communicate, modify, translate and otherwise exploit (and authorize Third Parties to do any of the foregoing) to use the information on the Client Supplied **Non-Confidential** Sample Description Form in the Field of Use, provided that the Customer's identity is not linked to, or otherwise disclosed with respect to, such data.

4.2 Ownership of Report. The Customer shall own the data in the Report. For Non-Confidential Proteomics Analyses, the Customer grants Kinexus a non-exclusive, royalty-free fully paid up worldwide perpetual license to use, copy, publish, compile, display, communicate, modify, translate and otherwise exploit (and authorize Third Parties to do any of the foregoing) data from the Report in the Field of Use.

4.3 Confidentiality of Sample Information. Kinexus will have no rights with respect to the Confidential Sample information until the Sample information is published or otherwise enters the public domain. Thereafter, Kinexus can use the results of the Proteomics Analyses of the Customer Samples for its internal research and development programs.

4.4 Ownership of Proteomics Products. The Customer owns the Proteomics Products that have been delivered to the Customer in the amounts specified in the Service Order Form and the Service Information Form. Kinexus owns any excess Proteomics Products and may dispose of these in its best interests.

4.5 Ownership of New Intellectual Property.

- (a) The Customer shall own and have rights to all inventions, discoveries, improvements, know-how, technical information, data or other technology discovered, conceived, made, developed and/or reduced to practice through the use of the data in the Report and Proteomics Products solely by employees of the Customer or jointly with its Affiliates;
- (b) Kinexus shall own and have rights to all inventions, discoveries, improvements, know-how, technical information, data or other technology discovered, conceived, made, developed and/or reduced to practice through the use of the data in the Report and Proteomics Products solely by employees of Kinexus or jointly with its Affiliates.

4.6 Non-Exclusive License to Preserve Kinexus Proteomics Services Freedom of Operation. In the event one or more claims of an issued patent arising from the use of a Report by the Customer, its Affiliates, Academic Collaborators or Corporate Partners would, absent a license from the Customer or its Affiliates, prevent Kinexus from using or permitting others to use the Kinexus Proteomics Services or any data therein, then the Customer and/or its Affiliates (as applicable) shall grant to Kinexus a non-exclusive, royalty-free fully-paid up perpetual license, including the right to grant sublicenses, under any such patent claim to use and permit others to use the Proteomics Services.

## 5. CONFIDENTIALITY

5.1 Confidentiality. Each Receiving Party shall treat the Confidential Information of the Disclosing Party as strictly confidential and (a) take reasonable precautions to protect such Confidential Information (including, without limitation, all precautions such as the Receiving Party employs with respect to its own confidential information), (b) not disclose or make available to any Third Party such Confidential Information without the express prior written consent of the Disclosing Party and (c) use such Confidential Information only for purposes specifically authorized under this Agreement. Each Receiving Party may disclose Confidential

Information to its employees, consultants, Affiliates and agents, and to licensees or prospective licensees of its rights to any invention, on a need-to-know basis and on the condition that such employees, Affiliates, agents, licensees and prospective licensees are obligated to maintain the confidentiality of the Confidential Information under written agreements that contain terms and conditions no less restrictive than the terms and conditions of this Section 5. Each Receiving Party may disclose Confidential Information of the Disclosing Party pursuant to a demand issued by a court or governmental agency or as otherwise required by law, provided, however, that the Receiving Party notifies the Disclosing Party promptly upon receipt thereof, giving the Disclosing Party sufficient advance notice to permit it to seek a protective order or other similar order with respect to such Confidential Information, and provided, further, that the Receiving Party furnishes only that portion of the Confidential Information which it is advised by counsel is legally required whether or not a protective order or other similar order is obtained by the Disclosing Party.

5.2 Publication. The Customer may publish and/or present the Report, abstracts or manuscripts generated utilizing the Report, and any data and/or results generated by the Customer utilizing the Report. The Customer is encouraged to disclose in scientific publications any Proteomics Analyses that were performed by Kinexus and any Proteomics Products were produced by Kinexus that meaningfully contributed to the described work. Please refer to “Kinexus Bioinformatics Corporation (Vancouver, Canada).” For all Samples submitted for analysis and identified as Non-Confidential by the Customer, Kinexus will not use, copy, publish, compile, display, communicate, modify, or translate the Sample Information or the data from the Report for a period of 180 days (6 months) following the return of the Report to the Customer. At any time, the Customer may opt to pay the difference in price between the Non-Confidential pricing level to the Confidential pricing level for each applicable Sample, to ensure the confidentiality status of such sample is changed.

5.3 Confidential Sample Information. All parties agree that the term of confidentiality pertaining to that Sample information will expire when the Sample information is published or otherwise enters public domain through no fault of Kinexus.

5.4 Use of Customer Name. Except as expressly provided in Section 9.5, no right or license is granted hereunder by Customer for Kinexus to use the Customer’s name in relation to data from a Report to a third party.

## **6. TERM AND TERMINATION**

6.1 Term. The term of this Agreement (“**Term**”) shall commence on the Effective Date and shall remain in effect for fifteen (15) years or until the termination of this Agreement pursuant to the terms hereof.

6.2 Early Termination. Each party shall have the right to terminate this Agreement at any time prior to Kinexus' delivery of a Report or Proteomics Product to the Customer hereunder, upon ten (10) days written notice to the other party, if such party reasonably determines that the production, or use of such Sample infringes intellectual property rights of any Third Party, and the Customer elects not to obtain a license under the necessary Third Party intellectual property rights at its sole expense. If this Agreement is terminated by either party pursuant to this Section 6.2, neither party shall have any obligation to the other with respect to payments under this Agreement regarding the Sample or Proteomics Product at issue.

Kinexus shall have the right to terminate any work order for any Proteomics Services upon ten (10) days written notice to the Customer, upon the identification of a technical difficulty related to the Sample or Proteomics Product which would prevent it from delivering the Report or Proteomics Product using reasonable efforts. If Kinexus terminates a work order as a result of a technical difficulty related to a Customer Sample that is the fault of Kinexus, Kinexus shall provide for the reanalysis of the same number of problematic Customer Samples for the Proteomics Analyses at the original agreed upon price without any additional expenses incurred by the Customer, or Kinexus shall repay any prepayment fee paid by the Customer for such a Customer Sample and neither party shall have any further obligation to the other with respect to that Customer Sample.

If Kinexus terminates a work order for Proteomics Analyses as a result of a technical difficulty related to the Customer Sample (including insufficient material or other problems associated with the quality of the Sample) that is the fault of the Customer, then Kinexus shall provide for the reanalysis of the problematic Customer Samples at the original agreed upon price without any additional expenses incurred by the Customer, provided Kinexus completes the full Proteomics Analyses for all Samples. For any subsequent resubmission of Customer Samples for Proteomics Analyses due to technical difficulty that is again the fault of the Customer, Kinexus shall provide for the reanalysis of the problematic Customer Samples at an additional charge per sample at a price mutually agreed by the Customer and Kinexus. If the Customer elects not to resubmit Samples for Proteomics Analyses, then the Customer will pay Kinexus an amount equivalent to 50% of the quoted price for the work performed by Kinexus to this point.

6.3 Events of Default. An event of default (an “Event of Default”) shall be deemed to occur upon a material breach of this Agreement by a party (including, without limitation, any breach of the provisions of Section 5) if the breaching party fails to remedy such breach within thirty (30) days after written notice thereof by the non-breaching party.

6.4 Effect of an Event of Default.

- (a) Remedies Available to Kinexus. If an Event of Default occurs relating to a material breach by the Customer, then Kinexus shall have the right, at its option exercisable in its sole discretion, in addition to any other rights or remedies available to it at law or in equity, to immediately terminate this Agreement upon notice thereof to the Customer, in which case the Customer shall return to Kinexus, or, upon Kinexus' written instruction, destroy any Report, Proteomics Products, and all information, other materials or documentation provided or made available by Kinexus pursuant to this Agreement, and any copies thereof (including electronic copies).
- (b) Remedies Available to the Customer. If an Event of Default occurs relating to a material breach by Kinexus, then the Customer shall have the right, at its option exercisable in its sole discretion, in addition to any other rights or remedies available to it at law or in equity and subject to the limitations set forth in Section 7, to terminate this Agreement upon notice thereof to Kinexus.

6.5 Effect of Expiration or Termination of Agreement. The expiration or termination of this Agreement shall not relieve the parties of any obligation accruing prior to such expiration or termination. Kinexus will not be required to continue Custom Immunohistochemistry Analyses on a Sample after termination, and the Customer will be required to pay for work done prior to termination. The provisions of Sections 4, 5, 6, 7, 8, and 9 hereof shall survive any expiration or termination of this Agreement.

## **7. DISCLAIMER OF WARRANTIES AND LIMITATION OF LIABILITY**

7.1 Disclaimer of Warranties. THE PROTEOMICS SERVICES ARE BEING SUPPLIED TO CUSTOMER WITH NO EXPRESS, IMPLIED, STATUTORY OR OTHER WARRANTIES, REPRESENTATIONS, CONDITIONS OR GUARANTEES, INCLUDING THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, TITLE AND DURABILITY. WITHOUT LIMITING THE FOREGOING, KINEXUS MAKES NO REPRESENTATION OR WARRANTY THAT THE USE OF THE REPORT, ANY PROTEOMICS PRODUCTS OR THE DATA THEREIN OR THE PERFORMANCE OF THIS AGREEMENT WILL NOT INFRINGE ANY INTELLECTUAL PROPERTY OR OTHER RIGHTS OF ANY THIRD PARTY.

7.2 Limitation of Liability. Kinexus shall not be liable for any use by the Customer, its Affiliates, Corporate Partners, or Academic Collaborators of the Report and any Proteomics Products or any loss, claim,

damage or liability, of whatever kind or nature, which may arise from or in connection with the use of the Report or the data therein, and any Proteomics Products. NOTWITHSTANDING ANYTHING ELSE IN THIS AGREEMENT OR OTHERWISE TO THE CONTRARY, NEITHER KINEXUS NOR CUSTOMER WILL BE LIABLE TO EACH OTHER WITH RESPECT TO ANY MATTER ARISING UNDER THIS AGREEMENT UNDER ANY CONTRACT, NEGLIGENCE, STRICT LIABILITY OR OTHER LEGAL OR EQUITABLE THEORY FOR (I) ANY PUNITIVE, EXEMPLARY, INCIDENTAL OR CONSEQUENTIAL DAMAGES OR LOST PROFITS OR (II) COST OF PROCUREMENT OF SUBSTITUTE GOODS, TECHNOLOGY OR SERVICES. WITHOUT IN ANY WAY LIMITING THE FOREGOING, KINEXUS SHALL NOT, IN ANY EVENT, HAVE ANY LIABILITY WHATSOEVER IN CONNECTION WITH THIS AGREEMENT IN EXCESS OF AN AMOUNT EQUAL TO THE FEES PAID TO KINEXUS BY CUSTOMER HEREUNDER IN RESPECT OF THE PROTEOMICS SERVICES AT ISSUE.

## **8. INDEMNIFICATION**

Except to the extent prohibited by law, the Customer shall assume all liability for, and shall defend, indemnify and hold Kinexus, its Affiliates and their respective directors, officers, employees and agents harmless from, all claims, losses, damages or expenses (including reasonable attorneys' fees) arising directly or indirectly as a result of: (a) the use of the Report or the data therein and any Proteomics Products by the Customer or its Affiliates, Corporate Partners or Academic Collaborators, or (b) the breach, untruthfulness or inaccuracy of any of the Customer's representations and warranties in this Agreement.

## **9. MISCELLANEOUS**

9.1 Entire Agreement. The Appendices to this Agreement, together with all terms and conditions contained within this Agreement constitute the entire understanding between the parties with respect to the subject matter hereof and, with respect to any conflicting terms from prior agreements between the parties, supersedes and cancels such conflicting sections from all previous registrations, agreements, commitments and writings in respect thereof. This Agreement may be amended, or any term hereof modified, only by a written instrument duly executed by both parties hereto.

9.2 Assignment and Waiver. This Agreement may not be assigned or otherwise transferred by either party without the written consent of the other party, such consent will not be unreasonably withheld. Notwithstanding the foregoing, Kinexus may, without such consent, assign its rights and obligations under this Agreement (a) to any Affiliate or (b) to a Third Party in connection with a merger, consolidation or sale of such portion of its assets that includes rights under this Agreement provided, however, that Kinexus' rights and obligations under this Agreement shall be assumed by its successor in interest in any such transaction. In the event of such a transaction with Third Party, notwithstanding the other provisions of this Agreement, the intellectual property rights of such Third Party shall not be subject to the licenses granted by Kinexus under this Agreement. Any purported assignment in violation of the provisions of this Section 9.2 shall be void. Any permitted assignee shall assume all obligations of its assignor under this Agreement. The waiver by either party hereto of any right hereunder or the failure to perform or of a breach by the other party shall not be deemed a waiver of any other right hereunder or of any other breach or failure by said other party whether of a similar nature or otherwise.

9.3 Force Majeure. Neither party shall be held liable or responsible to the other party nor be deemed to have defaulted under or breached this Agreement for failure or delay in fulfilling or performing any obligation under this Agreement when such failure or delay is caused by or results from causes beyond the reasonable control of the affected party, including but not limited to fire, floods, embargoes, war, acts of war (whether war is declared or not), insurrections, riots, civil commotions, strikes, lockouts or other labor or supply disturbances, acts of God or acts, omissions or delays in acting by any governmental authority or the other party; provided, however, that the party so affected shall use reasonable commercial efforts to avoid or remove such causes of nonperformance, and

shall continue performance hereunder with reasonable dispatch whenever such causes are removed. Either party shall provide the other party with prompt written notice of any delay or failure to perform that occurs by reason of force majeure. The parties shall mutually seek a resolution of the delay or the failure to perform as noted above.

9.4 Notices. Any consent, notice, or report required or permitted to be given or made under this Agreement by one of the notification parties hereto to the other shall be in writing, delivered personally, by email or by facsimile (and promptly confirmed by telephone, personal delivery or courier) or courier, postage prepaid (where applicable), addressed to such other party at its address indicated below, or to such other address as the addressee shall have last furnished in writing to the addressor and shall be effective upon receipt by the addressee.

*If to Kinexus:*

Kinexus Bioinformatics Corporation  
Suite 1, 8755 Ash Street  
Vancouver, British Columbia, Canada V6P 6T3  
Attention: Dr. Steven Pelech  
President & C.S.O.  
Telephone: (604) 323-2547 extension 10  
Facsimile: (604) 323-2548

*If to the Customer:*

To the Customer at the address designated at the front of this Agreement and to the attention of the duly authorized representative signing this Agreement.

9.5 Publicity. Except as required by law, the terms of this Agreement shall be treated as Confidential Information and shall not be disclosed to anyone (except for the parties' respective directors, officers, employees, consultants, agents and attorneys assisting in the review and negotiation of this Agreement and/or who have a need to know the terms of this Agreement) without the written consent of the other party, such consent which will not be unreasonably withheld. Notwithstanding the foregoing, (a) Kinexus may, without such consent, publicly announce the execution of this Agreement with the Customer and may reference the Customer as a Kinexus client.

9.6 No Partnership. It is expressly agreed that the relationship between Kinexus and the Customer shall not constitute a partnership, joint venture or agency. Neither Kinexus nor the Customer shall have the authority to make any statements, representations or commitments of any kind, or to take any action, which shall be binding on the other, without the prior consent of the other party to do so.

9.7 Applicable Law. This Agreement shall be governed by, construed, interpreted and enforced in accordance with, the laws of the province of British Columbia and the laws of Canada, without reference to conflict of laws principles.

9.8 Dispute Resolution.

- (a) The parties hereby agree that they will attempt in good faith to resolve any controversy or claim arising out of or relating to this Agreement promptly by negotiations. If a controversy or claim should arise hereunder, the matter shall be referred to an individual designated by the Chief Executive Officer or President of Kinexus and an individual designated by the Chief Executive Officer (or the equivalent position) of the Customer (the "Representatives"). If the matter has not been resolved within twenty-one (21) days of the first meeting of the Representatives of the parties (which period may be extended by mutual agreement) concerning such matter, subject to rights to injunctive relief and specific performance, and unless otherwise specifically provided for herein, any controversy or claim arising out of or relating to this Agreement, or the breach thereof, will be settled as set forth in Section 9.8(b).

- (b) All disputes arising in connection with this Agreement that are not resolved pursuant to Section 9.8(a) above shall be finally settled in Vancouver, British Columbia, by a single arbitrator appointed pursuant to the provisions of the *Commercial Arbitration Act* (British Columbia). Notwithstanding the above, either party has the right to bring an action in a court of competent jurisdiction against the other party for (i) any breach of such other party's duties of confidentiality pursuant to Section 5 of this Agreement; (ii) any infringement of its proprietary rights by the other party; and (iii) for interim protection such as, by way of example, an interim injunction. Judgment upon the arbitrator's award may be entered in any court of competent jurisdiction. The award of the arbitrator may include compensatory damages against either party, but under no circumstances will the arbitrator be authorized to, nor shall he/she, award punitive, consequential or incidental damages against either party. The parties agree not to institute any litigation or proceedings against each other in connection with this Agreement except as provided in this Section 9.8.

9.9 Severability. Each party hereby agrees that it does not intend to violate any public policy, statutory or common laws, rules, regulations, treaty or decision of any government agency or executive body thereof of any country or community or association of countries. Should one or more provisions of this Agreement be or become invalid, the parties hereto shall substitute, by mutual consent, valid provisions for such invalid provisions which valid provisions in their economic effect are sufficiently similar to the invalid provisions that it can be reasonably assumed that the parties would have entered into this Agreement with such valid provisions. In case such valid provisions cannot be agreed upon, the invalidity of one or several provisions of this Agreement shall not affect the validity of this Agreement as a whole, unless the invalid provisions are of such essential importance to this Agreement that it is to be reasonably assumed that the parties would not have entered into this Agreement without the invalid provisions.

9.10 Counterparts. This Agreement may be executed in counterparts, each of which when executed and delivered is an original, but both of which together shall constitute one and the same instrument.

9.11 Fax Delivery. This Agreement may be executed by the parties and transmitted by facsimile and if so executed and transmitted this Agreement will be for all purposes as effective as if the parties had delivered an executed original Agreement.

**IN WITNESS WHEREOF**, the parties have caused their duly authorized officer to execute and deliver this Agreement as of the Effective Date.

\_\_\_\_\_  
*Printed Name of Institute or Company*

Per: \_\_\_\_\_  
*Signature of Authorized Representative*

Name: \_\_\_\_\_  
*Printed Name of Authorized Representative*

Title: \_\_\_\_\_  
*Printed Title of Authorized Representative*

Date signed: \_\_\_\_\_

**KINEXUS BIOINFORMATICS CORPORATION**

Per: \_\_\_\_\_  
*Signature of Dr. Steven Pelech*

**Dr. Steven Pelech**

**President and Chief Scientific Officer**

Date signed: \_\_\_\_\_