

CUSTOM KINASE SUBSTRATE PROFILING (CKSP) SERVICE

List of Available Active Protein Kinases

This list may change with respect to availability and pricing.

Protein Kinase Name	Code	U.S. Price	Protein Kinase Name	Code	U.S. Price
Abl1	AB01	\$200.00	BUBR1(BUB1B)	BU01	\$600.00
Abl1 [E255K]	AB02	\$400.00	CaMK1δ (CAMK1D)	CA03	\$200.00
Abl1 [G250E]	AB03	\$400.00	CAMK1γ	CA04	\$200.00
Abl1 [H369P]	AB04	\$600.00	CAMK2α (CAMK2B)	CA05	\$200.00
Abl1 [T315I]	AB05	\$400.00	CaMK2β (CAMK2B)	CA06	\$200.00
Abl1 [Y253F]	AB06	\$400.00	CaMK2δ (CAMK2D)	CA07	\$400.00
Abl1 [M351T]	AB07	\$600.00	CaMK2γ (CAMK2G)	CA08	\$600.00
Abl1 [Q252H]	AB08	\$600.00	CAMK3γ	CA09	\$200.00
Abl2 (Arg)	AB09	\$200.00	CAMK4	CA10	\$200.00
ACK	AC01	\$200.00	CAMK4 (CaMKIV)	CA11	\$400.00
ACVR1 (ALK2)	AC02	\$400.00	CAMKK1 (CAMKKA)	CA12	\$200.00
ACVRL1	AC03	\$400.00	CAMKK2	CA13	\$200.00
ADRBK1 (GRK2)	AD01	\$200.00	CAMKK2 (CaMKK beta)	CA14	\$400.00
ADRBK2 (GRK3)	AD02	\$400.00	CDC42 BPA (MRCKA)	CD01	\$400.00
Akt1/PKBα	AK01	\$200.00	CDC42 BPB (MRCKB)	CD02	\$400.00
Akt1/PKBα [δPH, S473D]	AK02	\$600.00	CDC7/ASK	CD03	\$600.00
Akt1/PKBα [δPH]	AK03	\$600.00	CDK1/cyclin B1	CD04	\$400.00
Akt2/PKBβ	AK04	\$200.00	CDK1/CyclinA2	CD05	\$200.00
Akt2/PKBβ [δPH, S474D]	AK05	\$600.00	CDK2/cyclin A	CD06	\$400.00
Akt3/PKBγ	AK06	\$200.00	CDK2/Cyclin E1	CD07	\$600.00
Akt3/PKBγ [S472D]	AK07	\$600.00	CDK2/CyclinA2	CD08	\$200.00
ALK1	AL01	\$200.00	CDK3/Cyclin E1	CD09	\$600.00
ALK4 (ACVR1B)	AL02	\$200.00	CDK4//Cyclin D3	CD10	\$600.00
AMPKα1/β1/γ1 (PRKAA1/B1/G1)	AM01	\$200.00	CDK4/Cyclin D1	CD11	\$200.00
AMPKα1/β1/γ2 (PRKAA1/B1/G2)	AM02	\$200.00	CDK5	CD12	\$600.00
AMPKα1/β1/γ3 (PRKAA1/B1/G3)	AM03	\$200.00	CDK5/p25	CD13	\$200.00
AMPKα1/β2/γ1 (PRKAA1/B2/G1)	AM04	\$200.00	CDK5/p35	CD14	\$400.00
AMPKα2/β1/γ1 (PRKAA2/B1/G1)	AM05	\$200.00	CDK6//Cyclin D3	CD15	\$600.00
Ark5	AR01	\$600.00	CDK6/cyclin D1	CD16	\$400.00
Ask1 (MAP3K5)	AS01	\$200.00	CDK7/Cyclin H1/MAT1	CD17	\$200.00
Aurora A (AurA, AURKA)	AU01	\$200.00	CDK8/Cyclin C	CD18	\$400.00
Aurora B (AurB, AURKB, INCENP)	AU02	\$200.00	CDK9/Cyclin K	CD19	\$200.00
Aurora C (AurC, AURKC)	AU03	\$200.00	CDK9/Cyclin T1	CD20	\$400.00
Axl	AX01	\$200.00	CGK2 (PRKG2)	CG01	\$600.00
BARK1 (ADRBK1)	BA01	\$600.00	CHK1 (CHEK1)	CH01	\$200.00
BARK2 (ADRBK2)	BA02	\$600.00	CHK2 (CHEK2)	CH02	\$200.00
Blk	BL01	\$200.00	CK1α (CSNK1A1)	CK01	\$400.00
BMPR1A (ALK3)	BM01	\$600.00	CK1δ (CSNK1D)	CK02	\$400.00
BMPR1B (ALK6)	BM02	\$600.00	CK1δ (CSNK1D) [1-294]	CK03	\$600.00
Bmx	BM03	\$200.00	CK1ε (CSNK1E)	CK04	\$400.00
B-Raf	BR01	\$200.00	CK1γ1 (CSNK1G1)	CK05	\$400.00
B-Raf [δ1-415]	BR02	\$600.00	CK1γ2 (CSNK1G2)	CK06	\$400.00
B-Raf [V599E]	BR03	\$200.00	CK1γ3 (CSNK1G3)	CK07	\$400.00
Brk (PTK6)	BR04	\$200.00	CK1γ3 (CSNK1G3)	CK08	\$400.00
BrSK1 (SAD1)	BR05	\$400.00	CK2α1 (CSNK2A1)	CK09	\$200.00
BrSK2	BR06	\$600.00	CK2α2 (CSNK2A2)	CK10	\$400.00
Btk	BT01	\$200.00	CLK1	CL01	\$200.00
Btk [E41K]	BT01	\$600.00	CLK2	CL02	\$200.00
Btk [R28H]	BT02	\$600.00	CLK3	CL03	\$400.00

List of Available Active Protein Kinases (Cont'd) - 2

Protein Kinase Name	Code	U.S. Price	Protein Kinase Name	Code	U.S. Price
CLK4	CL04	\$400.00	FGFR4	FG08	\$200.00
COT (MAP3K8)	CO01	\$200.00	Fgr	FG09	\$600.00
CRIK	CR01	\$600.00	Flt1 (VEGFR1)	FL01	\$400.00
Csk	CS02	\$200.00	Flt3	FL02	\$200.00
C-TAK1	CT01	\$600.00	Flt3 [D835Y]	FL03	\$400.00
CTK	CT02	\$600.00	Flt4 (VEGFR3)	FL04	\$200.00
DAPK1	DA01	\$200.00	Fms (CSF1R)	FM01	\$400.00
DAPK2	DA02	\$400.00	FRAP1 (mTOR)	FR01	\$600.00
DAPK3 (ZIPK)	DA03	\$200.00	FRK (PTK5)	FR02	\$400.00
DCAMKL1	DC01	\$600.00	Fyn	FY01	\$200.00
DCAMKL2 (DCK2)	DC02	\$400.00	GRK4	GR01	\$400.00
DDR1	DD01	\$600.00	GRK5	GR02	\$200.00
DDR2	DD02	\$200.00	GRK6	GR03	\$400.00
DLK (MAP3K12)	DL01	\$600.00	GRK7	GR04	\$600.00
DMPK1 (DMPK)	DM01	\$400.00	GSK3 α	GS01	\$200.00
DRAK1 (STK17A)	DR01	\$600.00	GSK3 β	GS02	\$400.00
DYRK1A	DY01	\$400.00	Haspin (GSG2)	HA01	\$200.00
DYRK1B	DY02	\$400.00	Hck	HC01	\$200.00
DYRK2	DY03	\$600.00	Hgk (MAP4K4)	HG01	\$400.00
DYRK3	DY04	\$400.00	HIPK1 (Myak)	HI01	\$200.00
DYRK4	DY05	\$400.00	HIPK2	HI02	\$200.00
eEF2K	EE01	\$400.00	HIPK3	HI03	\$200.00
EGFR (ErbB1)	EG01	\$400.00	HIPK3 (YAK1)	HI04	\$200.00
EGFR (ErbB1) [L858R]	EG02	\$400.00	HIPK4	HI05	\$600.00
EGFR (ErbB1) [L861Q]	EG03	\$400.00	IGF1R	IG01	\$400.00
EGFR (ErbB1) [T790M, L858R]	EG04	\$400.00	IGF1R [δ 1-958]	IG02	\$600.00
EGFR (ErbB1) [T790M]	EG05	\$400.00	IKK α (CHUK)	IK01	\$200.00
EIF2AK3 (PERK)	EI06	\$400.00	IKK β (IKBKB)	IK02	\$200.00
EphA1	EP01	\$200.00	IKK ϵ (IKBKE)	IK03	\$400.00
EphA2	EP02	\$400.00	INSRR (IRR)	IN01	\$200.00
EphA3	EP03	\$400.00	Insulin Rec. (INSR)	IN02	\$400.00
EphA4	EP04	\$400.00	IRAK1	IR01	\$600.00
EphA5	EP05	\$400.00	IRAK2	IR02	\$200.00
EphA6	EP06	\$600.00	IRAK4	IR03	\$400.00
EphA7	EP07	\$600.00	Itk	IT01	\$200.00
EphA8	EP08	\$600.00	JAK1	JA01	\$600.00
EphA8	EP09	\$400.00	JAK2	JA02	\$200.00
EphB1	EP10	\$600.00	JAK2 [JH1, JH2, V617F]	JA03	\$600.00
EphB2	EP11	\$400.00	JAK2 [JH1, JH2]	JA04	\$400.00
EphB3	EP12	\$200.00	JAK3	JA05	\$400.00
EphB4	EP13	\$200.00	JNK1a1 (MAPK8)	JN01	\$400.00
ErbB2 (HER2, Neu)	ER01	\$200.00	JNK2 α 2 (MAPK9)	JN02	\$600.00
ErbB4 (HER4)	ER02	\$400.00	JNK3 (MAPK10)	JN03	\$600.00
Erk1 (MAPK3)	ER03	\$600.00	KDR (VEGFR2)	KD01	\$400.00
Erk2 (MAPK1)	ER04	\$600.00	KHS1(MAP4K5)	KH01	\$400.00
Erk5 (MAPK7)	ER05	\$200.00	Kit	KI01	\$200.00
FAK (PTK2)	FA01	\$400.00	Kit [D816H]	KI02	\$600.00
Fer	FE01	\$200.00	Kit [D816V]	KI03	\$600.00
Fes (Fps)	FE02	\$200.00	Kit [T670I]	KI04	\$400.00
FGFR1 (FLT2)	FG01	\$200.00	Kit [V560G]	KI05	\$200.00
FGFR1 (FLT2) [V561M]	FG02	\$200.00	Kit [V654A]	KI06	\$200.00
FGFR2	FG03	\$400.00	Lck	LC01	\$200.00
FGFR2 [N549H]	FG04	\$600.00	LIMK1	LI01	\$200.00
FGFR3	FG05	\$200.00	LIMK2	LI02	\$400.00
FGFR3 [K650E]	FG06	\$400.00	LKB1 (MO25 α , STRAD α , STK11)	LK01	\$600.00
FGFR3 [K650M]	FG07	\$600.00	LOK (STK10)	LO01	\$600.00

List of Available Active Protein Kinases (Cont'd) - 3

Protein Kinase Name	Code	U.S. Price	Protein Kinase Name	Code	U.S. Price
LRRK2	LR01	\$600.00	MST4	MS07	\$200.00
LRRK2 [G2019S]	LR02	\$400.00	MUSK	MU01	\$200.00
LTK (TYK1)	LT01	\$400.00	MYLK2 (skMLCK)	MY01	\$400.00
Lyn A	LY01	\$200.00	MYO3 β	MY02	\$400.00
Lyn B	LY02	\$200.00	MYT1 (PKMYT1)	MY03	\$200.00
MKK5 (MEK5, MAP2K5)	MK01	\$600.00	NDR1 (STK38)	ND01	\$200.00
MAP3K14 (NIK)	MA01	\$600.00	NEK1	NE01	\$200.00
MEKK3 (MAP3K3)	ME01	\$200.00	NEK11 (FLJ23495)	NE02	\$200.00
MAP3K4	MA02	\$600.00	NEK2	NE03	\$200.00
MAP3K7-MAP3K7IP1 (TAK1-TAB1)	MA03	\$400.00	NEK3	NE04	\$200.00
GCK (MAP4K2)	GC01	\$400.00	NEK4	NE05	\$400.00
MAP4K4 (HGK)	MA04	\$600.00	NEK6	NE06	\$200.00
MAPKAPK2	MA05	\$400.00	NEK7	NE07	\$600.00
MAPKAPK3	MA06	\$400.00	NEK9	NE08	\$600.00
MAPKAPK5 (PRAK)	MA07	\$400.00	NLK	NL01	\$400.00
MARK1 (MARK)	MA08	\$400.00	NUAK1 (ARK5)	NU01	\$400.00
MARK2	MA09	\$400.00	p38 α (MAPK14)	MA13	\$200.00
MARK3	MA10	\$400.00	p38 α (MAPK14) [T106M]	MA14	\$400.00
MARK4	MA11	\$400.00	p38 β (MAPK11)	MA15	\$400.00
MATK (HYL)	MA12	\$400.00	p38 δ (MAPK13)	MA16	\$200.00
MEK1 (MKK1, MAP2K1)	ME01	\$200.00	p38 γ (MAPK12)	MA17	\$200.00
MEK1 (MKK1, MAP2K1) [S218E, S222E]	ME02	\$200.00	p70S6K (RPS6KB1)	RS01	\$400.00
MEK2 (MKK2, MAP2K2)	ME03	\$200.00	p70S6K (RPS6KB1) [T412E]	RS02	\$400.00
MEKK2 (MAP3K2)	ME04	\$200.00	p70S6K β (RPS6KB2)	RS03	\$400.00
MEKK3 (MAP3K2)	ME05	\$200.00	PAK1/CDC42	PA01	\$200.00
MELK	ME06	\$400.00	PAK2 (PAK65)	PA02	\$600.00
MERTK (Mer)	ME07	\$200.00	PAK3	PA03	\$200.00
Met	ME08	\$400.00	PAK4	PA04	\$200.00
Met [M1250T]	ME09	\$200.00	PAK6	PA05	\$200.00
Met [Y1235D]	ME10	\$400.00	PAK7 (KIAA1264, PAK5)	PA06	\$600.00
MGC42105	MG01	\$200.00	PAR-1Ba/MARK2	PA07	\$200.00
MINK1 (MINK)	MI01	\$200.00	PASK	PA08	\$400.00
MKK3 (MEK3, MAP2K3)	MK01	\$200.00	PBK (TOPK)	PB01	\$200.00
MKK4 (MEK4, MAP2K4)	MK02	\$400.00	PCTAIRE1 (PCTK1)	PC01	\$400.00
MKK6 (MEK6, MAP2K6)	MK03	\$200.00	PDGFR α	PD01	\$200.00
MKK6 (MEK6, MAP2K6) [S599D, T603D]	MK04	\$600.00	PDGFR α [550-end, D842V]	PD02	\$200.00
MKK7 α 1 (MEK7 α 1, MAP2K7A1)	MK05	\$400.00	PDGFR α [550-end, V561D]	PD03	\$400.00
MKK7 β 1 (MKK7 β 1, MAP2K7B1)	MK06	\$600.00	PDGFR α [D842V]	PD04	\$200.00
MLCK (MLCK2, MYLK)	ML01	\$200.00	PDGFR α [T674I]	PD05	\$600.00
MLK1 (MAP3K9)	ML02	\$200.00	PDGFR α [V561D]	PD06	\$400.00
MLK2 (MAP3K10)	ML03	\$200.00	PDGFR β	PD07	\$400.00
MLK3 (MAP3K11)	ML04	\$200.00	PDHK2 (PDK2)	PD08	\$600.00
MNK1 (MKNK1)	MN01	\$600.00	PDHK3 (PDK3)	PD09	\$200.00
MNK2 (MKNK2)	MN02	\$600.00	PDHK4 (PDK4)	PD10	\$600.00
MOS	MO01	\$600.00	PDK1	PD11	\$600.00
MRCK α (CDC42BPA)	MR01	\$600.00	PEK (EIF2AK3)	PE01	\$400.00
MRCK β (CDC42BPB)	MR02	\$200.00	PhK γ 1 (PHKG1)	PH01	\$400.00
MSK1 (RPS6KA5)	MS01	\$600.00	PhK γ 2 (PHKG2)	PH02	\$400.00
MSK2 (RPS6KA4)	MS02	\$600.00	Pim1	PI01	\$200.00
MSSK1 (STK23)	MS03	\$200.00	Pim2	PI02	\$600.00
MST1 (STK4)	MS04	\$600.00	Pim3	PI03	\$600.00
MST2 (STK3)	MS05	\$200.00	PKA α (PRKACA)	PK01	\$600.00
MST3 (STK24)	MS06	\$600.00	PKA β (PRKACB)	PK02	\$200.00

List of Available Active Protein Kinases (Cont'd) - 4

Protein Kinase Name	Code	U.S. Price	Protein Kinase Name	Code	U.S. Price
PKA γ (PRKACG)	PK03	\$600.00	SGK2	SG03	\$400.00
PKC μ (PRKD1)	PK04	\$400.00	SGK3 (SGKL)	SG04	\$600.00
PKC α (PRKCA)	PK05	\$400.00	SGT220	SG05	\$200.00
PKC β 1 (PRKCB1)	PK06	\$200.00	SGT222-25UG	SG06	\$200.00
PKC β 2 (PRKCB2)	PK07	\$200.00	SIK	SI01	\$200.00
PKC δ (PRKCD)	PK08	\$600.00	skMLCK (MYLK2)	SK01	\$400.00
PKC ϵ (PRKCE)	PK09	\$200.00	SLK (STK2)	SL01	\$200.00
PKC γ (PRKCG)	PK10	\$200.00	smMLCK (MYLK)	SM01	\$600.00
PKC η (PRKCH)	PK11	\$200.00	SNF1LK2 (QIK)	SN01	\$200.00
PKC ι (PRKCI)	PK12	\$200.00	SOK	SO01	\$200.00
PKC λ (PRKCL)	PK13	\$200.00	Src	SR01	\$200.00
PKC θ (PRKCQ)	PK14	\$200.00	Src [T341M]	SR02	\$600.00
PKC ζ (PRKCZ)	PK15	\$200.00	SRMS (Srm)	SR03	\$600.00
PKD1 (PRKD1)	PK16	\$200.00	SRPK1	SR04	\$600.00
PKD2 (PRKD2)	PK17	\$200.00	SRPK2	SR05	\$600.00
PKD3 (PRKD3, PRKCN)	PK18	\$200.00	STK16 (PKL12)	ST01	\$400.00
PKG1 α (PRKG1A)	PK19	\$200.00	STK25 (YSK1)	ST02	\$400.00
PKG1 β (PRKG1B)	PK20	\$200.00	STK33	ST03	\$200.00
PKG2 (PRKG2)	PK21	\$600.00	Syk	SY01	\$600.00
PKN1 (PRK1)	PK22	\$600.00	TAO2	TA01	\$400.00
PKN2 (PRK2)	PK23	\$200.00	TAO3	TA02	\$200.00
PKR (EIF2AK2)	PK24	\$200.00	TAOK1	TA03	\$400.00
PLK1	PL01	\$200.00	TAOK2 (TAO1)	TA04	\$400.00
PLK2	PL02	\$600.00	TAOK3 (JIK)	TA05	\$200.00
PLK3	PL03	\$200.00	TBK1	TB01	\$200.00
PLK4	PL04	\$400.00	TEC	TE01	\$400.00
PRKX	PR01	\$600.00	TESK1	TE02	\$200.00
PTK2B (FAK2)	PT01	\$600.00	TGF β R1 (TGFB1, ALK5)	TG01	\$600.00
PTK5	PT02	\$400.00	TGF β R2	TG02	\$600.00
Pyk2	PY01	\$400.00	Tie2 (Tek)	TI01	\$200.00
Raf1 [Y340E, Y341E]	RA01	\$400.00	Tie2 (Tek) [R849W]	TI02	\$200.00
Raf1 (truncated)	RA02	\$600.00	Tie2 (Tek) [Y1108F]	TI03	\$400.00
Raf1 [Y340D, Y341D]	RA03	\$200.00	Tie2 (Tek) [Y897S]	TI04	\$400.00
Ret	RE01	\$600.00	TLK2	TL01	\$400.00
Ret [V804L]	RE02	\$400.00	TNK1	TN01	\$600.00
Ret [V804M]	RE03	\$400.00	TNK2 (ACK)	TN02	\$600.00
Ret [Y791F]	RE04	\$600.00	TrkA (NTRK1)	TR01	\$200.00
RIPK2	RI01	\$200.00	TrkB (NTRK2)	TR02	\$200.00
RIPK5	RI02	\$400.00	TrkC (NTRK3)	TR03	\$600.00
ROCK1 (ROK β)	RO01	\$600.00	TSSK1 (STK22D)	TS01	\$200.00
ROCK2 (ROK α)	RO02	\$200.00	TSSK2 (STK22B)	TS02	\$200.00
RON (MST1R)	RO03	\$200.00	TTBK1	TT01	\$600.00
ROR1	RO04	\$600.00	TTK	TT02	\$600.00
ROR2	RO05	\$400.00	TXK	TX01	\$400.00
Ros	RO06	\$400.00	TYK2	TY01	\$200.00
Rse	RO07	\$200.00	TYRO3 (RSE)	TY02	\$200.00
RSK1 (RPS6KA1)	RS04	\$400.00	VRK1	VR01	\$600.00
RSK2 (RPS6KA3)	RS05	\$200.00	Wee1	WE01	\$600.00
RSK3 (RPS6KA2)	RS06	\$200.00	WNK1	WN01	\$200.00
RSK4 (RPS6KA6)	RS07	\$200.00	WNK2	WN02	\$600.00
SGK1	SG01	\$600.00	WNK3	WN03	\$600.00
SGK1 [δ 1-59, S422D]	SG02	\$200.00	WNK4	WN04	\$200.00

List of Available Active Protein Kinases (Cont'd) - 5

Protein Kinase Name	Code	U.S. Price
YES1	YE01	\$600.00
ZAK	ZA01	\$200.00
ZAP70	ZA02	\$600.00
ZIPK	ZI01	\$600.00



Suite 1, 8755 Ash Street
Vancouver, B.C
Canada V6P 6T3

Phone: 1-866-KINEXUS
Phone: 1-604-323-2547
Facsimile: 1-604-323-2548
E-Mail: info@kinexus.ca
Internet: www.kinexus.ca