

Appendix 1 - List of 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	CaMK1 δ (CAMK1D)	CA03	\$200.00
Abl1 [G250E]	AB03	\$400.00	CAMK1 γ	CA04	\$200.00
Abl1 [H369P]	AB04	\$600.00	CAMK2 α (CAMK2B)	CA05	\$200.00
Abl1 [T315I]	AB05	\$400.00	CaMK2 β (CAMK2B)	CA06	\$200.00
Abl1 [Y253F]	AB06	\$400.00	CaMK2 δ (CAMK2D)	CA07	\$400.00
Abl1 [M351T]	AB07	\$600.00	CaMK2 γ (CAMK2G)	CA08	\$600.00
Abl1 [Q252H]	AB08	\$600.00	CAMK3 γ	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/PKB α	AK01	\$200.00	CDC42 BPB (MRCKB)	CD02	\$400.00
Akt1/PKB α [δ PH, S473D]	AK02	\$600.00	CDC7/ASK	CD03	\$600.00
Akt1/PKB α [δ PH]	AK03	\$600.00	CDK1/cyclin B1	CD04	\$400.00
Akt2/PKB β	AK04	\$200.00	CDK1/CyclinA2	CD05	\$200.00
Akt2/PKB β [δ PH, S474D]	AK05	\$600.00	CDK2/cyclin A	CD06	\$400.00
Akt3/PKB γ	AK06	\$200.00	CDK2/Cyclin E1	CD07	\$600.00
Akt3/PKB γ [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
AMPK α 1/ β 1/ γ 1 (PRKAA1/B1/G1)	AM01	\$200.00	CDK4/Cyclin D1	CD11	\$200.00
AMPK α 1/ β 1/ γ 2 (PRKAA1/B1/G2)	AM02	\$200.00	CDK5	CD12	\$600.00
AMPK α 1/ β 1/ γ 3 (PRKAA1/B1/G3)	AM03	\$200.00	CDK5/p25	CD13	\$200.00
AMPK α 1/ β 2/ γ 1 (PRKAA1/B2/G1)	AM04	\$200.00	CDK5/p35	CD14	\$400.00
AMPK α 2/ β 1/ γ 1 (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	CK1 α (CSNK1A1)	CK01	\$400.00
BMPR1A (ALK3)	BM01	\$600.00	CK1 δ (CSNK1D)	CK02	\$400.00
BMPR1B (ALK6)	BM02	\$600.00	CK1 δ (CSNK1D) [1-294]	CK03	\$600.00
Bmx	BM03	\$200.00	CK1 ϵ (CSNK1E)	CK04	\$400.00
B-Raf	BR01	\$200.00	CK1 γ 1 (CSNK1G1)	CK05	\$400.00
B-Raf [δ 1-415]	BR02	\$600.00	CK1 γ 2 (CSNK1G2)	CK06	\$400.00
B-Raf [V599E]	BR03	\$200.00	CK1 γ 3 (CSNK1G3)	CK07	\$400.00
Brk (PTK6)	BR04	\$200.00	CK1 γ 3 (CSNK1G3)	CK08	\$400.00
BrSK1 (SAD1)	BR05	\$400.00	CK2 α 1 (CSNK2A1)	CK09	\$200.00
BrSK2	BR06	\$600.00	CK2 α 2 (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

Appendix 1 - List of 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

Appendix 1 - List of 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

Appendix 1 - List of 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

Appendix 1 - List of 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