

Search below to match your KINASE with our SENSOR PEPTIDE substrate!

Bulk Sensor Peptide: Designed for higher throughput applications. 1mg is sufficient for ~1,200 assays in a low-volume 384-well plate (10µM sensor peptide in 25µL reaction volume). Available from 1mg to +200mg aliquots. Optimized ancillary reagents sold separately.

Assay Kits: Designed for streamlined technology evaluation. Each kit enables 96 reactions in a low-volume 384-well plate (10µM sensor peptide in 25µL reaction volume). Includes all assay components (excluding your kinase).

See your target of interest? Contact us here. Or email us at hello@assayquant.com!

Don't see your target of interest? Contact us. Our PhosphoSens Technology enables rapid assay development.

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
ABL1	ABL proto-oncogene 1, non-receptor tyrosine kinase	NM_005157	ABL1	CSKS-AQT0031B CSKS-AQT0031K	Y	Carna (08-001)	VS
ABL1 [E255K]	ABL proto-oncogene 1, non-receptor tyrosine kinase	NM_005157	ABL1	CSKS-AQT0031B CSKS-AQT0031K	Y	Carna (08-094)	VS
ABL1 [T315I]	ABL proto-oncogene 1, non-receptor tyrosine kinase	NM_005157	ABL1	CSKS-AQT0031B CSKS-AQT0031K	Y	Carna (08-093)	VS
ABL2 (ARG)	ABL proto-oncogene 2, non-receptor tyrosine kinase	NM_005157	ARG	CSKS-AQT0031B CSKS-AQT0031K	Y	Carna (08-102)	VS
AKT1 (PKB)	AKT Serine/Threonine Kinase 1	NM_005163	AKT1	CSKS-AQT0536B CSKS-AQT0536K	S/T	Carna (01-101)	VS
AKT2 (PKB β)	AKT Serine/Threonine Kinase 2	NM_001626	AKT2	CSKS-AQT0536B CSKS-AQT0536K	S/T	Carna (01-102)	VS
AKT3 (PKB γ)	AKT Serine/Threonine Kinase 3	NM_005465	AKT3	CSKS-AQT0536B CSKS-AQT0536K	S/T	Carna (01-103)	VS
ALK	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-518)	VS
ALK [C1156Y]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-530)	VS
ALK [F1174L]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-519)	VS
ALK [G1269A]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-537)	VS
ALK [G1202R]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-544)	VS
ALK [L1196M/G1202R]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	SignalChem (A19-12NG)	VS
ALK [L1196M]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-529)	VS
ALK [R1275Q]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-520)	VS
ALK [T1151-L1152 INS T]	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-539)	VS
ALK:NPM1 (NPM1-ALK)	ALK receptor tyrosine kinase	NM_004304	ALK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-517)	S
AURKA (AURA)	Aurora kinase A	NM_003600	AurA	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (05-101)	VS
AURKB (AURB)	aurora kinase B	NM_004217	AurB	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (05-102)	VS
AURKC (AURC)	aurora kinase C	NM_004217	AurB	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (05-103)	VS
AXL	AXL receptor tyrosine kinase	NM_001699	AXL	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-107)	VS
BLK	BLK proto-oncogene, Src family tyrosine kinase	NM_001715	BLK	CSKS-AQT0032B CSKS-AQT0032K	Y	Carna (08-164)	VS
BMX	BMX non-receptor tyrosine kinase	NM_001721	BMX	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-179)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
BRSK1	BR serine/threonine kinase 1	NM_032430	BRSK1	CSKS-AQT0123B CSKS-AQT0123K	S/T	Carna (02-115)	S
BRSK2	BR serine/threonine kinase 2	NM_003957	BRSK2	CSKS-AQT0123B CSKS-AQT0123K	S/T	Carna (02-116)	S
BTK	Bruton tyrosine kinase	NM_000061	BTK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-180)	VS
BTK [C481S]	Bruton tyrosine kinase	NM_000061	BTK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-547)	VS
BUB1	BUB1 mitotic checkpoint serine/threonine kinase	NM_004336	BUB1	CSKS-AQT0848B CSKS-AQT0848K	S/T	Carna (05-187)	M
CAMK1A (CAMK1 α)	calcium/calmodulin dependent protein kinase I	NM_003656	CaMK1a	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-104)	VS ¹
CAMK1D (CAMK1 δ)	calcium/calmodulin dependent protein kinase ID	NM_020397	CaMK1d	CSKS-AQT0420B CSKS-AQT0420K	S/T	SignalChem (C09-10G)	VS ¹
CAMK1G (CAMK1 γ)	calcium/calmodulin dependent protein kinase IG	NM_020439	CaMK1g	CSKS-AQT0420B CSKS-AQT0420K	S/T	SignalChem (C10-11G)	VS ¹
CAMK2A (CAMK2 α)	calcium/calmodulin dependent protein kinase II alpha	NM_003656	CaMK2a	CSKS-AQT0425B CSKS-AQT0425K	S/T	Carna (02-109)	VS ¹
CAMK2B (CAMK2 β)	calcium/calmodulin dependent protein kinase II beta	NM_001220	CaMK2b	CSKS-AQT0425B CSKS-AQT0425K	S/T	Carna (02-110)	VS
CAMK2D (CAMK2 δ)	calcium/calmodulin dependent protein kinase II delta	NM_001221	CaMK2d	CSKS-AQT0425B CSKS-AQT0425K	S/T	Carna (02-111)	VS ¹
CAMK2G (CAMK2 γ)	calcium/calmodulin dependent protein kinase II gamma	NM_001222	CaMK2g	CSKS-AQT0425B CSKS-AQT0425K	S/T	Carna (02-112)	VS ¹
CAMK4	calcium/calmodulin dependent protein kinase IV	NM_001744	CaMK4	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-108)	VS ¹
CAMKK1	alcium/calmodulin dependent protein kinase kinase 2	NM_006549	CaMKK2	CSKS-AQT0699B CSKS-AQT0699K	S/T	Carna (05-157)	S ¹
CDC42BPB (MRCKB)	CDC42 binding protein kinase beta	NM_006035	MRCKb	CSKS-AQT0713B CSKS-AQT0713K	S/T	Carna (01-108)	M
CDC42BPG (DMPK2)	CDC42 binding protein kinase gamma	NM_017525	DMPK2	CSKS-AQT0324B CSKS-AQT0324K	S/T	SignalChem (D17-11G)	S
CDK1/CycB1(CDC2)	cyclin dependent kinase 1	NM_001786	CDC2	CSKS-AQT0297B CSKS-AQT0297K	S/T	Carna(04-102)	VS
CDK1/CycE1	cyclin dependent kinase 1	NM_001786	CDC2	CSKS-AQT0297B CSKS-AQT0297K	S/T	Proqinase (0134-0055-1)	VS
CDK2/CycA2	cyclin dependent kinase 2	NM_001798	CDK2	CSKS-AQT0297B CSKS-AQT0297K	S/T	Carna (04-103)	VS
CDK2/CycE1	cyclin dependent kinase 2	NM_001798	CDK2	CSKS-AQT0297B CSKS-AQT0297K	S/T	Carna (04-165)	VS
CDK3/CycE1	cyclin dependent kinase 3	NM_001258	CDK3	CSKS-AQT0297B CSKS-AQT0297K	S/T	Carna (04-104)	VS
CDK4 /CycD3	cyclin dependent kinase 4	NM_000075	CDK4	CSKS-AQT0258B CSKS-AQT0258K	S/T	Carna (04-105)	S
CDK5/p25	cyclin dependent kinase 5	NM_004935	CDK5	CSKS-AQT0255B CSKS-AQT0255K	S/T	Carna (04-106)	VS
CDK5/p35	cyclin dependent kinase 5	NM_004935	CDK5	CSKS-AQT0255B CSKS-AQT0255K	S/T	SignalChem (C33-10BG)	VS
CDK6/CycD1	cyclin dependent kinase 6	NM_001259	CDK6	CSKS-AQT0258B CSKS-AQT0258K	S/T	Carna (04-114)	VS
CDK6/CycD3	cyclin dependent kinase 6	NM_001259	CDK6	CSKS-AQT0258B CSKS-AQT0258K	S/T	Carna (04-107)	S
CDK7/CycH/MAT1	cyclin dependent kinase 7	NM_001799	CDK7	CSKS-AQT0733B CSKS-AQT0733K	S/T	Carna (04-108)	S
CDK9 (TAK) w/CycK	cyclin dependent kinase 9	NM_001261.2	CDK9	CSKS-AQT0449B CSKS-AQT0449K	S/T	Carna (04-111)	M
CDK9 (TAK) w/CycT1	cyclin dependent kinase 9	NM_001261.2	CDK9	CSKS-AQT0449B CSKS-AQT0449K	S/T	Carna (04-110)	VS
CDK14 (PFTAIRE1)/CycY	cyclin dependent kinase 14	NM_012395	PFTAIRE1	CSKS-AQT0730B CSKS-AQT0730K	S/T	SignalChem (P15-10G)	M
CDK16/CycY	cyclin dependent kinase 16	NM_006201	PCTAIRE1	CSKS-AQT0258B CSKS-AQT0258K	S/T	Carna (04-116)	S

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
CDK17 (PCTAIRE2)/CycY	cyclin dependent kinase 17	NM_002595	PCTAIRE2	CSKS-AQT0730B CSKS-AQT0730K	S/T	SignalChem (P10-10G)	S
CDK18/CycY (PCTAIRE3)	cyclin dependent kinase 18	NM_002596	PCTAIRE3	CSKS-AQT0255B CSKS-AQT0255K	S/T	Thermo (A30492)	S
CDPK1 (toxoplasma gondii)				CSKS-AQT0124B CSKS-AQT0124K	S/T	Academic Lab	VS
CDPK3 (toxoplasma gondii)				CSKS-AQT0124B CSKS-AQT0124K	S/T	Academic Lab	VS
CHEK1 (CHK1)	Check point kinase 1	NM_001274	CHK1	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-117)	VS
CHEK2 (CHK2)	Check point kinase 2	NM_007194	CHK2	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-162)	VS
CHUK (IKK α , IKK α)	conserved helix-loop-helix ubiquitous kinase	NM_001278	IKK α	CSKS-AQT0603B CSKS-AQT0603K	S/T	Carna (05-112)	M
CIT (CRIK)	citron rho-interacting serine/threonine kinase	NM_007174	CRIK	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (01-104)	M
CLK1	CDC like kinase 1	NM_004071	CLK1	CSKS-AQT0360B CSKS-AQT0360K	S/T	SignalChem (C57-11G)	S
CLK2	CDC like kinase 2	NM_001291	CLK2	CSKS-AQT0360B CSKS-AQT0360K	S/T	Carna (04-127)	VS
CLK3	CDC like kinase 3	NM_001292	CLK3	CSKS-AQT0360B CSKS-AQT0360K	S/T	Carna (04-128)	VS
CSF1R (FMS)	Colony stimulating factor 1 receptor	NM_005211	FMS	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-155)	VS
CSK	C-terminal Src kinase	NM_004383	CSK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-111)	M
CSNK1A1 (CK1A1, CK1 α)	casein kinase 1 alpha 1	NM_001892	CK1a	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-101)	VS
CSNK1A1L (CK1A1L, CK1 α 2)	casein kinase 1 alpha 1 like	NM_145203	CK1a2	CSKS-AQT0709B CSKS-AQT0709K	S/T	SignalChem (C75-10G)	VS
CSNK1D (CK1D, CK1 δ)	casein kinase 1 delta	NM_001893	CK1d	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-103)	VS
CSNK1E (CK1E, CK1 ϵ)	casein kinase 1 epsilon	NM_001894	CK1e	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-104)	VS
CSNK1G1 (CK1G1, CK1 γ 1)	casein kinase 1 gamma 1	NM_022048	CK1g1	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-105)	VS
CSNK1G2 (CK1G2, CK1 γ 2)	casein kinase 1 gamma 2	NM_001319	CK1g2	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-106)	VS
CSNK1G3 (CK1G3, CK1 γ 3)	casein kinase 1 gamma 3	NM_004384	CK1g3	CSKS-AQT0709B CSKS-AQT0709K	S/T	Carna (03-107)	VS
CSNK2A1 (CK2A1, CK2 α 1)	casein kinase 2 alpha 1	NM_001895	CK2a1	CSKS-AQT0696B CSKS-AQT0696K	S/T	Thermo (PV3248)	VS
CSNK2A2 (CK2A2, CK2 α 2)	casein kinase 2 alpha 2	NM_001896	CK2a2	CSKS-AQT0696B CSKS-AQT0696K	S/T	Thermo (PV3624)	S
DAPK1	Death-associated protein kinase 1	NM_004938	DAPK1	CSKS-AQT0173B CSKS-AQT0173K	S/T	SignalChem (D01-11G)	VS ¹
DAPK2 (DRP-1)	Death-associated protein kinase 2	NM_014326	DAPK2	CSKS-AQT0236B CSKS-AQT0236K	S/T	SignalChem (D02-10G)	VS ¹
DAPK3 (ZIPK)	Death-associated protein kinase 3	NM_001348	DAPK3	CSKS-AQT0237B CSKS-AQT0237K	S/T	Carna (02-136)	VS ¹
DCLK1 (DCAMKL1)	doublecortin like kinase 1	NM_004734	DCAMKL1	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-139)	S ¹
DCLK2 (DCAMKL2)	doublecortin like kinase 2	NM_152619	DCAMKL2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-140)	VS ¹
DDR1 (SRC treated)	discoidin domain receptor tyrosine kinase 1	NM_001954	DDR1	CSKS-AQT0235B CSKS-AQT0235K	Y	Carna (08-132)	M
DDR2	discoidin domain receptor tyrosine kinase 2	NM_006182	DDR2	CSKS-AQT0235B CSKS-AQT0235K	Y	Carna (08-114)	VS ²
DYRK1A	dual specificity tyrosine phosphorylation regulated kinase 1A	NM_001396	DYRK1A	CSKS-AQT0260B CSKS-AQT0260K	S/T	Carna (04-130)	VS
DYRK1B	dual specificity tyrosine phosphorylation regulated kinase 1B	NM_004714	DYRK1B	CSKS-AQT0260B CSKS-AQT0260K	S/T	Carna (04-131)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
DYRK2	dual specificity tyrosine phosphorylation regulated kinase 2	NM_006482	DYRK2	CSKS-AQT0260B CSKS-AQT0260K	S/T	Carna (04-132)	VS
DYRK3	dual specificity tyrosine phosphorylation regulated kinase 3	NM_003582	DYRK3	CSKS-AQT0260B CSKS-AQT0260K	S/T	Carna (04-133)	VS
DYRK4	dual specificity tyrosine phosphorylation regulated kinase 4	NM_003845	DYRK4	CSKS-AQT0260B CSKS-AQT0260K	S/T	Carna (04-434)	VS
EEF2K (eEF2K)	eukaryotic elongation factor 2 kinase	NM_013302	eEF2K	CSKS-AQT0349B CSKS-AQT0349K	S/T	Carna (10-113)	VS ¹
EGFR (ERBB1)	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	BPS (40187)	VS
EGFR (D770_N771insNPG)	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-553)	VS
EGFR [C797S/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122ZG)	VS
EGFR [d746-750/C797S/T790M]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122UG)	VS
EGFR [d746-750/C797S]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122TG)	VS
EGFR [d746-750/T790M]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122KG)	VS
EGFR [d746-750]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-527)	VS
EGFR [L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122BG)	VS
EGFR [L861Q]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-513)	VS
EGFR [T790M/C797S/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	SignalChem (E10-122VG)	VS
EGFR [T790M/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-510)	VS
EGFR [T790M]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-194)	VS
EIF2AK1 (HRI)	eukaryotic translation initiation factor 2 alpha kinase 1	NM_014413	HRI	CSKS-AQT0748B CSKS-AQT0748K	S/T	Carna (05-154)	M
EIF2AK2 (PKR), Full length	eukaryotic translation initiation factor 2 alpha kinase 2	NM_002759	PKR	CSKS-AQT0355B CSKS-AQT0355K	S/T	Academic Lab	VS ³
EIF2AK3 (PERK/PEK)	eukaryotic translation initiation factor 2 alpha kinase 3	NM_004836	PEK	CSKS-AQT0356B CSKS-AQT0356K	S/T	Carna (05-155)	VS
EIF2AK4 (GCN2)	eukaryotic translation initiation factor 2 alpha kinase 4	NM_001013703	GCN2	CSKS-AQT0349B CSKS-AQT0349K	S/T	Carna (05-153)	VS ⁴
EPHA1	EPH receptor A1	NM_005232	Epha1	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-119)	VS
EPHA2	EPH receptor A2	NM_004431	Epha2	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-121)	VS
EPHA3	EPH receptor A3	NM_005233	Epha3	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-122)	VS
EPHA4	EPH receptor A4	NM_004438	Epha4	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-123)	VS
EPHA5	EPH receptor A5	NM_004439	Epha5	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-124)	VS
EPHA6	EPH receptor A6	XM_114973	Epha6	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-125)	VS
EPHA7	EPH receptor A7	NM_004440	Epha7	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-126)	VS
EPHA8	EPH receptor A8	NM_020526	Epha8	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-127)	VS
EPHB1	EPH receptor B1	NM_004441	Ephb1	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-128)	VS
EPHB2	EPH receptor B2	NM_004442	Ephb2	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-129)	VS
EPHB3	EPH receptor B3	NM_004443	Ephb3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-130)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
EPHB4	EPH receptor B4	NM_004444	Ephb4	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-131)	VS
ERBB2 (HER2)	Erb-B2 Receptor Tyrosine Kinase 2	NM_004448	HER2/ErbB2	CSKS-AQT0794B CSKS-AQT0794K	Y	BPS (40230)	VS
ERBB2 (HER2) [A775/G776insYVMA]	Erb-B2 Receptor Tyrosine Kinase 2	NM_004448	HER2/ErbB2	CSKS-AQT0794B CSKS-AQT0794K	Y	SignalChem (E27-13BG)	VS
ERBB4 (HER4)	Erb-B2 Receptor Tyrosine Kinase 4	NM_005235	Erb-B2 Receptor Tyrosine Kinase 4	CSKS-AQT0794B CSKS-AQT0794K	Y	Carna (08-118)	VS
ERN1 (IRE1)	endoplasmic reticulum to nucleus signaling 1	NM_001433	IRE1	CSKS-AQT0715B CSKS-AQT0715K	S/T	Thermo (A30971)	S
ERN2 (IRE2)	endoplasmic reticulum to nucleus signaling 2	NM_033266	IRE2	CSKS-AQT0715B CSKS-AQT0715K	S/T	SignalChem (E32-11G)	M
FAM20C	FAM20C golgi associated secretory pathway kinase	NM_020223		CSKS-AQT0698B CSKS-AQT0698K	S/T	SignalChem (F18-11G)	S
FER	Fer tyrosine kinase	NM_005246	FER	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-139)	VS
FES	FES proto-oncogene, tyrosine kinase	NM_002005	FES	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-140)	VS
FGFR1	Fibroblast growth factor receptor 1	NM_000604	FGFR1	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-133)	VS
FGFR2	Fibroblast growth factor receptor 2	NM_022970.3	FGFR2	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-134)	VS
FGFR3	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-135)	VS
FGFR3 [G697C]	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Proqinase (1071-0000-1)	S
FGFR3 [K650E]	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-501)	VS
FGFR3 [K650M]	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-199)	VS
FGFR3 [V555L]	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-548)	VS
FGFR3 [V555M]	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-543)	VS
FGFR4	Fibroblast growth factor receptor 4	NM_002011	FGFR4	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-136)	VS
FGR	FGR proto-oncogene, Src family tyrosine kinase	NM_005248	FGR	CSKS-AQT0032B CSKS-AQT0032K	Y	Carna (08-166)	VS
FLT1	FMS-related tyrosine kinase 1	NM_002019	FLT1	CSKS-AQT0001B CSKS-AQT0001K	Y	Carna (08-189)	VS
FLT3	FMS-related tyrosine kinase 3	NM_004119	FLT3	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-154)	VS
FLT4	FMS-related tyrosine kinase 4	NM_002020	FLT4	CSKS-AQT0001B CSKS-AQT0001K	Y	Carna (08-190)	VS
FRK	Fyn related Src family tyrosine kinase	NM_002031	FRK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-167)	VS
FYNa	FYN proto-oncogene, Src family tyrosine kinase	NM_002037	FYN	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-168)	VS
FYNb	FYN proto-oncogene, Src family tyrosine kinase	NM_002037	FYN	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-531)	VS
GRK1 (RHOK)	G protein-coupled receptor kinase 1	NM_002929	RHOK	CSKS-AQT0603B CSKS-AQT0603K	S/T	SignalChem (R03-10G)	S
GRK2 (ADRBK1, BARK1)	G protein-coupled receptor kinase 2	NM_001619	BARK1	CSKS-AQT0641B CSKS-AQT0641K	S/T	Thermo (PV3361)	VS
GRK3 (ADRBK2, BARK2)	G protein-coupled receptor kinase 3	NM_005160	BARK2	CSKS-AQT0641B CSKS-AQT0641K	S/T	Carna (01-112)	S
GRK4 (GPRK4)	G protein-coupled receptor kinase 4	NM_005307	GPRK4	CSKS-AQT0641B CSKS-AQT0641K	S/T	Carna (01-313)	M
GRK5 (GPRK5)	G protein-coupled receptor kinase 5	NM_005308	GPRK5	CSKS-AQT0641B CSKS-AQT0641K	S/T	SignalChem (G03-10G)	M

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
GRK7 (GPRK7)	G protein-coupled receptor kinase 7	NM_139209	GPRK7	CSKS-AQT0641B CSKS-AQT0641K	S/T	Carna (01-316)	S
GSK3A (GSK3 α)	glycogen synthase kinase 3 alpha	NM_019884	GSK3A	CSKS-AQT0157B CSKS-AQT0157K	S/T	Carna (04-140)	VS
GSK3B (GSK3 β)	glycogen synthase kinase 3 beta	NM_001146156	GSK3B	CSKS-AQT0157B CSKS-AQT0157K	S/T	Carna (04-141)	VS
HASPIN	histone H3 associated protein kinase	NM_031965	Haspin	CSKS-AQT0798B CSKS-AQT0798K	S/T	Carna (05-111)	VS
HCK	HCK proto-oncogene, Src family tyrosine kinase	NM_002110	HCK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-169)	VS
HIPK1	homeodomain interacting protein kinase 1	NM_152696	HIPK1	CSKS-AQT0849B CSKS-AQT0849K	S/T	Carna (04-135)	M
HIPK2	Homeodomain interacting protein kinase 2	NM_022740	HIPK2	CSKS-AQT0134B CSKS-AQT0134K	S/T	Carna (04-136)	VS
HIPK3	homeodomain interacting protein kinase 3	NM_005734	HIPK3	CSKS-AQT0849B CSKS-AQT0849K	S/T	Carna (04-137)	M
HIPK4	homeodomain interacting protein kinase 4	NM_144685	HIPK4	CSKS-AQT0849B CSKS-AQT0849K	S/T	Carna (04-138)	M
ICK	intestinal cell kinase	NM_014920	ICK	CSKS-AQT0257B CSKS-AQT0257K	S/T	Carna (04-156)	S
IGF1R	Insulin-like growth factor 1 receptor	NM_000875	IGF1R	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-141)	VS
IKKB (IKK β)	Inhibitor of nuclear factor kappa-B kinase subunit beta	NM_001556	IKKb	CSKS-AQT0215B CSKS-AQT0215K	S/T	Carna (05-084)	VS
IKKE (IKK ϵ)	Inhibitor of nuclear factor kappa-B kinase subunit epsilon	NM_014002	IKKe	CSKS-AQT0215B CSKS-AQT0215K	S/T	Thermo (PV4875)	VS
INSR (IR)	Insulin receptor	NM_000208	INSR	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-142)	VS
INSRR (IRR)	Insulin receptor-related receptor	NM_014215	IRR	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-143)	VS
IRAK1	interleukin 1 receptor associated kinase 1	NM_001569	IRAK1	CSKS-AQT0326B CSKS-AQT0326K	S/T	Carna (09-101)	VS
IRAK4	interleukin 1 receptor associated kinase 4	NM_016123	IRAK4	CSKS-AQT0326B CSKS-AQT0326K	S/T	Carna (09-145)	VS
ITK	IL2-inducible T-cell kinase	NM_005546	ITK	CSKS-AQT0101B CSKS-AQT0101K	Y	Thermo (PV3875)	VS
JAK1 [JH1]	Janus kinase 1	NM_002227	JAK1	CSKS-AQT0663B CSKS-AQT0663K	Y	Thermo (PV4774)	S
JAK1 [JH1-JH2]	Janus kinase 1	NM_002227	JAK1	CSKS-AQT0663B CSKS-AQT0663K	Y	SignalChem (J01-11G)	S
JAK2 [JH1]	Janus kinase 2	NM_004972	JAK2	CSKS-AQT0661B CSKS-AQT0661K	Y	Carna (08-045)	VS
JAK2 [JH1-JH2]	Janus kinase 2	NM_004972	JAK2	CSKS-AQT0661B CSKS-AQT0661K	Y	Carna (08-514)	VS
JAK3 [JH1]	Janus kinase 3	NM_000215	JAK3	CSKS-AQT0658B CSKS-AQT0658K	Y	Carna (08-046)	VS
KDR (VEGFR2)	Kinase insert domain receptor	NM_002253	KDR	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-191)	VS
KIT	KIT proto-oncogene receptor tyrosine kinase	NM_000222	KIT	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-156)	VS
KSR1	kinase suppressor of ras 1	NM_014238	KSR1	CSKS-AQT0694B CSKS-AQT0694K	S/T	SignalChem (J01-11G)	M
KSR2	kinase suppressor of ras 2	NM_173598	KSR2	CSKS-AQT0694B CSKS-AQT0694K	S/T	SignalChem (J01-11G)	S
LATS1	large tumor suppressor kinase 1	NM_004690	LATS1	CSKS-AQT0476B CSKS-AQT0476K	S/T	Carna (01-123)	VS
LATS2	large tumor suppressor kinase 2	NM_014572	LATS2	CSKS-AQT0476B CSKS-AQT0476K	S/T	Carna (01-124)	VS
LCK	LCK proto-oncogene, Src family tyrosine kinase	NM_005356	LCK	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-170)	VS
LRRK2	leucine rich repeat kinase 2	NM_198578	LRRK2	CSKS-AQT0615B CSKS-AQT0615K	S/T	Thermo (PV48783)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
LRRK2 [G2019S]	leucine rich repeat kinase 2	NM_198578	LRRK2	CSKS-AQT0615B CSKS-AQT0615K	S/T	Thermo (PV4881)	VS
LTK	Leukocyte receptor tyrosine kinase	NM_002344	LTK	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-106)	VS
LYNa	LYN proto-oncogene, Src family tyrosine kinase	NM_002350	LYN	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-171)	VS
LYNb	LYN proto-oncogene, Src family tyrosine kinase	NM_002350	LYN	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-172)	VS
MAK	male germ cell associated kinase	NM_005906	MAK	CSKS-AQT0257B CSKS-AQT0257K	S/T	SignalChem (M01-10G)	S
MAP2K1 (MEK1), Cascade	Mitogen-activated protein kinase kinase 1	NM_002755	MAP2K1	CSKS-AQT0490B CSKS-AQT0490K	S/T/Y	Carna (07-141)	VS ⁵
MAP2K1 (MEK1),Direct	Mitogen-activated protein kinase kinase 1	NM_002755	MAP2K1	CSKS-AQT0972B CSKS-AQT0972K	S/T/Y	Carna (07-141)	VS
MAP2K2 (MEK2), Cascade	Mitogen-activated protein kinase kinase 2	NM_030662	MAP2K2	CSKS-AQT0490B CSKS-AQT0490K	S/T/Y	Carna (07-142)	VS ⁵
MAP2K2 (MEK2), Direct	Mitogen-activated protein kinase kinase 2	NM_030662	MAP2K2	CSKS-AQT0972B CSKS-AQT0972K	S/T/Y	Carna (07-142)	VS
MAP3K1 (MEKK1)	mitogen-activated protein kinase kinase kinase 1	NM_005921	MAP3K1	CSKS-AQT0739B CSKS-AQT0739K	S/T	Carna (07-103)	M
MAP3K2 (MEKK2)	mitogen-activated protein kinase kinase kinase 2	NM_006609	MAP3K2	CSKS-AQT0739B CSKS-AQT0739K	S/T	SignalChem (M10-10G)	M
MAP3K3 (MEKK3)	mitogen-activated protein kinase kinase kinase 3	NM_002401	MAP3K3	CSKS-AQT0743B CSKS-AQT0743K	S/T	Carna (07-105)	M
MAP3K5 (ASK1)	mitogen-activated protein kinase kinase kinase 5	NM_005923	MAP3K5	CSKS-AQT0557B CSKS-AQT0557K	S/T	SignalChem (M13-10G)	VS ⁶
MAP3K8 (Tpl2, COT)	mitogen-activated protein kinase kinase kinase 8	NM_005204	COT	CSKS-AQT0742B CSKS-AQT0742K	S/T	Carna (07-301)	VS
MAP3K9 (MLK1)	Mitogen-activated protein kinase kinase kinase 9	NM_033141	MLK1	CSKS-AQT0645B CSKS-AQT0645K	S/T	Carna (09-115)	VS
MAP3K11 (MLK3)	mitogen-activated protein kinase kinase kinase 11	NM_002419	MLK3	CSKS-AQT0646B CSKS-AQT0646K	S/T	Carna (09-017)	M
MAP3K19 (YSK4)	mitogen-activated protein kinase kinase kinase 19	NM_025052	YSK4	CSKS-AQT0742B CSKS-AQT0742K	S/T	SignalChem (Y06-11G)	VS
MAP3K20 (MLK7, ZAK)	mitogen-activated protein kinase kinase kinase 20	NM_016653	ZAK	CSKS-AQT0742B CSKS-AQT0742K	S/T	Carna (09-120)	M
MAP4K1 (HPK1)	Mitogen-activated protein kinase kinase kinase kinase 1	NM_007181	HPK1	CSKS-AQT0505B CSKS-AQT0505K	S/T	Thermo (PV6355)	VS
MAP4K2 (GCK)	Mitogen-activated protein kinase kinase kinase kinase 2	NM_004579	GCK	CSKS-AQT0505B CSKS-AQT0505K	S/T	SignalChem (M24-10G)	VS
MAP4K3 (GLK, KHS2)	Mitogen-activated protein kinase kinase kinase kinase 3	NM_003618	KHS2	CSKS-AQT0505B CSKS-AQT0505K	S/T	SignalChem (M25-11G)	S
MAP4K4 (HGK, ZC1)	Mitogen-activated protein kinase kinase kinase kinase 4	NM_004834	ZC1/HGK	CSKS-AQT0135B CSKS-AQT0135K	S/T	Carna (07-137)	VS
MAP4K5 (KHS1)	Mitogen-activated protein kinase kinase kinase kinase 5	NM_006575	KHS1	CSKS-AQT0178B CSKS-AQT0178K	S/T	SignalChem (M27-10G)	VS
MAPK1 (ERK2)	Mitogen-activated protein kinase 1	NM_002745	ERK2	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-143)	VS
MAPK3 (ERK1)	Mitogen-activated protein kinase 3	NM_002746	ERK1	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-142)	VS
MAPK7 (ERK5)	mitogen-activated protein kinase 7	NM_002750	ERK5	CSKS-AQT0373B CSKS-AQT0373K	S/T	Carna (04-146)	VS
MAPK8 (JNK1)	mitogen-activated protein kinase 8	NM_002750	JNK1	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-163)	VS
MAPK9 (JNK2)	mitogen-activated protein kinase 9	NM_002752	JNK2	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-164)	VS
MAPK10 (JNK3)	mitogen-activated protein kinase 10	NM_002753	JNK3	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-150)	VS
MAPK11 (p38 β)	mitogen-activated protein kinase 11	NM_002751	p38b	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-153)	VS
MAPK12 (p38 γ)	Mitogen-activated protein kinase 12	NM_002969	p38g	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-155)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
MAPK13 (p38δ)	Mitogen-activated protein kinase 13	NM_002754	p38d	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-154)	VS
MAPK14 (p38α)	mitogen-activated protein kinase 14	NM_001315	p38a	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-152)	VS
MAPKAPK2 (MK2)	Mitogen-activated protein kinase-activated protein kinase 2	NM_004759	MAPKAPK2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-142)	VS
MAPKAPK3 (MK3)	mitogen-activated protein kinase-activated protein kinase 3	NM_004635	MAPKAPK3	CSKS-AQT0425B CSKS-AQT0425K	S/T	Carna (02-143)	VS
MAPKAPK5 (MK5)	mitogen-activated protein kinase-activated protein kinase 5	NM_003668	MAPKAPK5	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-144)	VS
MARK1	microtubule affinity regulating kinase 1	NM_018650	MARK1	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-120)	VS
MARK2	Microtubule affinity regulating kinase 2	NM_004954	MARK2	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-121)	VS
MARK3	microtubule affinity regulating kinase 3	NM_002376	MARK3	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-122)	VS
MARK4	Microtubule affinity regulating kinase 4	NM_031417	MARK4	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-123)	VS
MASTL	microtubule associated serine/threonine kinase-like	NM_032844	MASTL	CSKS-AQT0693B CSKS-AQT0693K	S/T	SignalChem (M72-10G)	S
MELK	maternal embryonic leucine zipper kinase	NM_014791	MELK	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-124)	VS ¹
MERTK (MER)	MER proto-oncogene, tyrosine kinase	NM_006343	MER	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-108)	VS
MET (c-MET, HGFR)	MET proto-oncogene, receptor tyrosine kinase	NM_000245	MET	CSKS-AQT0702B CSKS-AQT0702K	Y	Carna (08-151)	VS
MINK1 (ZC3)	Misshapen-like kinase 1	NM_015716	ZC3/MINK	CSKS-AQT0183B CSKS-AQT0183K	S/T	SignalChem (M53-11G)	VS
MKNK1 (MNK1)	MAP kinase interacting serine/threonine kinase 1	NM_003684	MNK1	CSKS-AQT0426B CSKS-AQT0426K	S/T	SignalChem (M54-10G)	VS
MKNK2 (MNK2)	MAP kinase interacting serine/threonine kinase 2	NM_017572	MNK2	CSKS-AQT0426B CSKS-AQT0426K	S/T	SignalChem (M55-10G)	VS
MST1R (RON)	Macrophage stimulating 1 receptor	NM_016542	RON	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-152)	VS
MUSK	muscle associated receptor tyrosine kinase	NM_005592	MUSK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-153)	S
MYLK (MLCK)	Myosin light chain kinase	NM_053025	smMLCK	CSKS-AQT0151B CSKS-AQT0151K	S/T	SignalChem (M62-11G)	S ¹
MYLK2	Myosin light chain kinase 2	NM_033118	skMLCK	CSKS-AQT0517B CSKS-AQT0517K	S/T	SignalChem (M63-10H)	VS ¹
MYO3A	myosin IIIA	NM_017433	MYO3A	CSKS-AQT0505B CSKS-AQT0505K	S/T	SignalChem (M65-11G)	S
MYO3B	myosin IIIB	NM_138995	MYO3B	CSKS-AQT0135B CSKS-AQT0135K	S/T	SignalChem (M66-11G)	M
NEK1	NIMA related kinase 1	NM_012224	NEK1	CSKS-AQT0418B CSKS-AQT0418K	S/T	Carna (05-123)	VS
NEK2	NIMA related kinase 2	NM_002497	NEK2	CSKS-AQT0415B CSKS-AQT0415K	S/T	Carna (05-226)	VS
NEK3	NIMA related kinase 3	NM_002498	NEK3	CSKS-AQT0415B CSKS-AQT0415K	S/T	Carna (05-127)	M
NEK4	NIMA related kinase 4	NM_003157	NEK4	CSKS-AQT0418B CSKS-AQT0418K	S/T	Carna (05-128)	S
NEK5	NIMA related kinase 5	XM_292160	NEK5	CSKS-AQT0418B CSKS-AQT0418K	S/T	SignalChem (N07-11G)	VS
NEK7	NIMA related kinase 7	NM_133494	NEK7	CSKS-AQT0415B CSKS-AQT0415K	S/T	Carna (05-131)	M
NEK9	NIMA related kinase 9	NM_033116	NEK9	CSKS-AQT0415B CSKS-AQT0415K	S/T	Carna (05-133)	VS
NEK10	NIMA related kinase 10	NM_152534	NEK10	CSKS-AQT0415B CSKS-AQT0415K	S/T	SignalChem (N12-11G)	M
NIM1 (NIM1K)	NIM1 serine/threonine protein kinase	NM_153361	NIM1	CSKS-AQT0694B CSKS-AQT0694K	S/T	SignalChem (N14-10G)	S

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
NLK (LAK1)	nemo like kinase	NM_016231	NLK	CSKS-AQT0374B CSKS-AQT0374K	S/T	Carna (04-151)	S
NTRK1 (TRKA)	Neurotrophic receptor tyrosine kinase 1	NM_002529	TRKA	CSKS-AQT0104B CSKS-AQT0104K	Y	BPS (BS40280)	VS
NTRK2 (TRKB)	Neurotrophic receptor tyrosine kinase 2	NM_006180	TRKB	CSKS-AQT0104B CSKS-AQT0104K	Y	BPS (BS40281)	VS
NTRK3 (TRKC)	Neurotrophic receptor tyrosine kinase 3	NM_002530	TRKC	CSKS-AQT0104B CSKS-AQT0104K	Y	BPS (BS40282)	VS
NUAK1 (ARK5)	NUAK family kinase 1; (AMPK-related protein kinase 5)	NM_014840	NuaK1	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-126)	VS
NUAK2 (SNARK)	NUAK family kinase 2	NM_030952	NuaK2	CSKS-AQT0393B CSKS-AQT0393K	S/T	Carna (02-127)	VS
PAK1	p21 (RAC1) activated kinase 1	NM_002576	PAK1	CSKS-AQT0838B CSKS-AQT0838K	S/T	Carna (07-123)	S
PAK2	p21 (RAC1) activated kinase 2	NM_002577	PAK2	CSKS-AQT0838B CSKS-AQT0838K	S/T	Carna (07-124)	VS
PAK3	p21 (RAC1) activated kinase 3	NM_002578	PAK3	CSKS-AQT0838B CSKS-AQT0838K	S/T	Carna (07-125)	S
PAK4	p21 (RAC1) activated kinase 4	NM_005884	PAK4	CSKS-AQT0840B CSKS-AQT0840K	S/T	SignalChem (P05-10BG)	S
PAK5 (PAK7)	p21 (RAC1) activated kinase 5	NM_020341	PAK5	CSKS-AQT0840B CSKS-AQT0840K	S/T	SignalChem (P07-10G)	M
PAK6	p21 (RAC1) activated kinase 6	NM_020168	PAK6	CSKS-AQT0840B CSKS-AQT0840K	S/T	Carna (07-128)	M
PBK (TOPK)	PDZ binding kinase	NM_018492	PBK	CSKS-AQT0766B CSKS-AQT0766K	S/T	SignalChem (A05-11G)	S
PDGFRA (PDGFR α)	Platelet derived growth factor receptor alpha	NM_006206	PDGFR α	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-157)	VS
PDGFRB (PDGFR β)	Platelet derived growth factor receptor beta	NM_002609	PDGFR β	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-158)	VS
PDK1 (PDHK1)	pyruvate dehydrogenase kinase 1	NM_002610; NM_002613	PDHK1	CSKS-AQT0872B CSKS-AQT0872K	S/T	SignalChem (P57-10G)	M ⁷
PDK2 (PDHK2)	pyruvate dehydrogenase kinase 2	NM_002611	PDHK2	CSKS-AQT0896B CSKS-AQT0896K	S/T	SignalChem (P58-10G)	M ⁷
PDK3 (PDHK3)	pyruvate dehydrogenase kinase 3	NM_005391	PDHK3	CSKS-AQT0872B CSKS-AQT0872K	S/T	SignalChem (P59-10G)	M ⁷
PDK4 (PDHK4)	pyruvate dehydrogenase kinase 4	NM_002612	PDHK4	CSKS-AQT0872B CSKS-AQT0872K	S/T	SignalChem (P60-10G)	S ⁷
PDPK1 (PDK1)	3-phosphoinositide dependent protein kinase 1	NM_002610; NM_002613	PDHK1	CSKS-AQT0126B CSKS-AQT0126K	S/T	SignalChem (P14-10H)	VS
PHKG1	phosphorylase kinase catalytic subunit gamma 1	NM_006213	PHKg1	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-152)	S ¹
PHKG2	Phosphorylase kinase catalytic subunit gamma 2	NM_000294	PHKg2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-153)	S ¹
PIM1	Pim-1 proto-oncogene, serine/threonine kinase	NM_002648	PIM1	CSKS-AQT0474B CSKS-AQT0474K	S/T	Carna (02-054)	VS
PIM2	Pim-2 proto-oncogene, serine/threonine kinase	NM_006875	PIM2	CSKS-AQT0230B CSKS-AQT0230K	S/T	Carna (02-155)	VS
PIM3	Pim-3 proto-oncogene, serine/threonine kinase	NM_001001852.2	PIM3	CSKS-AQT0473B CSKS-AQT0473K	S/T	Carna (02-156)	VS
PINK1 (PINK)	PTEN induced putative kinase 1	NM_032409	PINK1	CSKS-AQT0686B CSKS-AQT0686K	S/T	BOSTON BIO(AP-182)	S
PKN1	protein kinase N1	NM_002741	PKN1	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-144)	VS
PKN2	protein kinase N2	NM_006256	PKN2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (01-145)	VS
PKN3	protein kinase N3	NM_013355	PKN3	CSKS-AQT0509B CSKS-AQT0509K	S/T	Carna (01-146)	S
PLK1	Polo like kinase 1	NM_005030	PLK1	CSKS-AQT0691B CSKS-AQT0691K	S/T	BPS (40033)	VS
PLK3	polo like kinase 3	NM_004073	PLK3	CSKS-AQT0691B CSKS-AQT0691K	S/T	SignalChem (P43-13G)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
PLK4	polo like kinase 4	NM_014264	PLK4	CSKS-AQT0686B CSKS-AQT0686K	S/T	Thermo (PV6394)	VS
PNCK (CAMK1β)	pregnancy up-regulated nonubiquitous CaM kinase	NM_001039582	CaMK1b	CSKS-AQT0420B CSKS-AQT0420K	S/T	Carna (02-105)	VS ¹
PRKAA1-B1-G1 (AMPKα1β1γ1)	Protein kinase AMP-activated catalytic subunit alpha 1	NM_006251	AMPKα1	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-113)	VS ⁸
PRKAA1-B2-G1 (AMPKα1β2γ1)	Protein kinase AMP-activated catalytic subunit alpha 1	NM_006251	AMPKα1	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-147)	VS ⁸
PRKAA2-B1-G1 (AMPKα2β1γ1)	protein kinase AMP-activated catalytic subunit alpha 2	NM_006252	AMPKα2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-114)	VS ⁸
PRKAA2-B2-G1 (AMPKα2β2γ1)	protein kinase AMP-activated catalytic subunit alpha 2	NM_006252	AMPKα2	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-148)	VS ⁸
PRKACA (PKACα)	Protein kinase cAMP-activated catalytic subunit alpha	NM_002730	PKACα	CSKS-AQT0458B CSKS-AQT0458K	S/T	Carna (01-127)	VS
PRKACB (PKACβ)	Protein kinase cAMP-activated catalytic subunit beta	NM_002731	PKACβ	CSKS-AQT0458B CSKS-AQT0458K	S/T	Carna (01-128)	VS
PRKACG (PKACγ, PKACg)	protein kinase cAMP-activated catalytic subunit gamma	NM_002732	PKACγ	CSKS-AQT0231B CSKS-AQT0231K	S/T	SignalChem (P53-10G)	VS
PRKCA (PKCα [alpha])	Protein kinase C alpha	NM_002737	PKCα	CSKS-AQT0316B CSKS-AQT0316K	S/T	Carna (01-133)	VS
PRKCB1 (PKCβ1 [beta1])	Protein kinase C beta	NM_002738	PKCβ	CSKS-AQT0316B CSKS-AQT0316K	S/T	Carna (01-134)	VS
PRKCB2 (PKCβ2 [beta2])	Protein kinase C beta	X07109	PKCβ	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-165)	VS
PRKCD (PKCδ [delta])	Protein kinase C delta	NM_006254	PKCδ	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-135)	VS
PRKCE (PKCε [epsilon])	Protein kinase C epsilon	NM_005400	PKCε	CSKS-AQT0465B CSKS-AQT0465K	S/T	Carna (01-136)	VS
PRKCG (PKCγ [gamma])	Protein kinase C gamma	NM_002739	PKCγ	CSKS-AQT0316B CSKS-AQT0316K	S/T	Carna (01-137)	VS
PRKCH (PKCη [eta])	Protein kinase C eta	NM_006255	PKCη	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-138)	VS
PRKCI (PKCι [iota])	Protein kinase C iota	NM_002740	PKCι	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-139)	VS
PRKCQ (PKCθ [theta])	Protein kinase C theta	NM_006257	PKCθ	CSKS-AQT0465B CSKS-AQT0465K	S/T	Carna (01-140)	VS
PRKCZ (PKCζ [zeta])	Protein kinase C zeta	NM_002744	PKCζ	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-141)	VS
PRKD1 (PKD1)	protein kinase D1	NM_002742	PKD1	CSKS-AQT0803B CSKS-AQT0803K	S/T	Carna (02-157)	VS
PRKD2 (PKD2)	protein kinase D2	NM_016457	PKD2	CSKS-AQT0803B CSKS-AQT0803K	S/T	Carna (02-158)	VS
PRKD3 (PKD3)	protein kinase D3	NM_005813	PKD3	CSKS-AQT0803B CSKS-AQT0803K	S/T	Carna (02-159)	VS
PRKDC (DNA-PK)	protein kinase, DNA-activated, catalytic polypeptide	NM_006904	DNAPK	CSKS-AQT0439B CSKS-AQT0439K	S/T	Thermo (PV5866)	VS
PRKG1 (PKG1, PGK)	protein kinase, cGMP-dependent, type I	NM_006258	DNAPK	CSKS-AQT0452B CSKS-AQT0452K	S/T	Carna (01-142)	VS ⁹
PRKG2 (PKG2, PGK2, CGK2)	protein kinase, cGMP-dependent, type II	NM_006259	PKG2	CSKS-AQT0452B CSKS-AQT0452K	S/T	Carna (01-143)	VS ⁹
PRKX	protein kinase, X-linked	NM_005044	PRKX	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (01-130)	S
PTK2 (FAK)	PTK2 protein tyrosine kinase 2	NM_153831	FAK	CSKS-AQT0061B CSKS-AQT0061K	Y	Carna (08-137)	S ²
PTK2B (PYK2)	Protein tyrosine kinase 2 beta	NM_004103	PYK2	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-138)	VS
PTK6 (BRK)	protein tyrosine kinase 6	NM_005975	BRK	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-165)	VS
Raf, b- (V600E), Cascade	B-Raf proto-oncogene, serine/threonine kinase	NM_002880	BRAF	CSKS-AQT0490B CSKS-AQT0490K	S/T	Thermo (PR7304A)	VS ¹⁰
Raf, b- (V600E), Direct	B-Raf proto-oncogene, serine/threonine kinase	NM_004333	BRAF	CSKS-AQT0682B CSKS-AQT0682K	S/T	Thermo (PR7304A)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
RAF1 (Y340E/Y341E), Cascade	Raf-1 proto-oncogene, serine/threonine kinase	NM_002880	RAF1	CSKS-AQT0490B CSKS-AQT0490K	S/T	Signal Chem (R01-13G-10)	VS ¹⁰
RAF1 (Y340E/Y341E), Direct	Raf-1 proto-oncogene, serine/threonine kinase	NM_002880	RAF1	CSKS-AQT0688B CSKS-AQT0688K	S/T	Signal Chem (R01-13G-10)	VS
RET	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-159)	VS
RET [G691S]	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-522)	VS
RET [M918T]	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-508)	VS
RET [S891A]	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-523)	VS
RET [Y791F]	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	Carna (08-521)	VS
RET:CCDC6 (CCDC6-RET)	Ret proto-oncogene	NM_020975	RET	CSKS-AQT0102B CSKS-AQT0102K	Y	SignalChem (R02-19BG)	VS
RIPK2	receptor interacting serine/threonine kinase 2	NM_003821	RIPK2	CSKS-AQT0209B CSKS-AQT0209K	S/T	Thermo (PV4213)	S
ROCK1	Rho associated coiled-coil containing protein kinase 1	NM_005406	ROCK1	CSKS-AQT0231B CSKS-AQT0231K	S/T	Carna (01-109)	VS
ROCK2	Rho associated coiled-coil containing protein kinase 2	NM_004850	ROCK2	CSKS-AQT0231B CSKS-AQT0231K	S/T	Carna (01-110)	VS
ROS1 (ROS)	ROS proto-oncogene 1, receptor tyrosine kinase	NM_002944	ROS	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-163)	VS
ROS1 [G2032R]	ROS proto-oncogene 1, receptor tyrosine kinase	NM_002944	ROS	CSKS-AQT0101B CSKS-AQT0101K	Y	SignalChem (R14-12BG)	VS
RPS6KA1 (RSK1)	Ribosomal protein S6 kinase A1	NM_002953	RSK3	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-149)	VS
RPS6KA2 (RSK3)	Ribosomal protein S6 kinase A2	NM_021135	RSK1	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-151)	VS
RPS6KA3 (RSK2)	Ribosomal protein S6 kinase A3	NM_004586	RSK2	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-150)	VS
RPS6KA4(MSK2)	ribosomal protein S6 kinase A4	NM_003942	MSK2	CSKS-AQT0535B CSKS-AQT0535K	S/T	Carna (01-148)	M
RPS6KA5 (MSK1)	ribosomal protein S6 kinase A5	NM_004755	MSK1	CSKS-AQT0535B CSKS-AQT0535K	S/T	Carna (01-147)	VS
RPS6KA6 (RSK4)	ribosomal protein S6 kinase A6	NM_014496	RSK4	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-152)	VS
RPS6KB1 (p70S6K)	Ribosomal protein S6 kinase, 70kDa, polypeptide 1	NM_003161	p70S6K	CSKS-AQT0231B CSKS-AQT0231K	S/T	Carna (01-158)	VS
RPS6KB2 (p70S6KB)	ribosomal protein S6 kinase B2	NM_003952	p70S6Kb	CSKS-AQT0231B CSKS-AQT0231K	S/T	SignalChem (R22-10G)	VS
SBK1 (SBK)	SH3 domain binding kinase 1	NM_001024401	SBK	CSKS-AQT0694B CSKS-AQT0694K	S/T	SignalChem (S56-10G)	M
SGK1 (SGK)	Serum/glucocorticoid regulated kinase 1	NM_005627	SGK	CSKS-AQT0536B CSKS-AQT0536K	S/T	Thermo (PV3818)	VS
SGK2	SGK2, serine/threonine kinase 2	NM_016276	SGK2	CSKS-AQT0536B CSKS-AQT0536K	S/T	Carna (01-159)	VS
SGK3	Serum/glucocorticoid regulated kinase family member 3	NM_013257	SGK3	CSKS-AQT0536B CSKS-AQT0536K	S/T	Carna (01-160)	VS
SIK1 (SIK)	salt inducible kinase 1	NM_173354	SIK	CSKS-AQT0508B CSKS-AQT0508K	S/T	Carna (02-131)	VS
SIK2 (QIK)	salt inducible kinase 2	NM_015191	QIK	CSKS-AQT0508B CSKS-AQT0508K	S/T	Carna (02-129)	VS
SIK3 (QSK)	salt inducible kinase 3	BC128510.1	QSK	CSKS-AQT0508B CSKS-AQT0508K	S/T	Dundee/MRC (DU 16623)	VS
SLK	STE20 like kinase	NM_014720	SLK	CSKS-AQT0505B CSKS-AQT0505K	S/T	Carna (07-129)	VS
SNRK	SNF related kinase	NM_017719	SNRK	CSKS-AQT0426B CSKS-AQT0426K	S/T	SignalChem (S16-10G)	S
SRC	SRC proto-oncogene, non-receptor tyrosine kinase	NM_005417	SRC	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-173)	VS

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
SRMS (SRM)	SRC-related kinase lacking C-terminal regulatory tyrosine and N-terminal myristylation sites;	NM_080823	SRM	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-174)	VS
SRPK1	SRSF protein kinase 1	NM_003137	SRPK1	CSKS-AQT0360B CSKS-AQT0360K	S/T	SignalChem (S21-10G)	VS
SRPK2	SRSF protein kinase 2	NM_182692	SRPK2	CSKS-AQT0360B CSKS-AQT0360K	S/T	Carna (04-161)	VS
SRPK3 (MSSK1)	SRSF protein kinase 3	NM_014370	MSSK1	CSKS-AQT0360B CSKS-AQT0360K	S/T	SignalChem (S41-10G)	S
STK3 (MST2)	Serine/threonine kinase 3	NM_006281	MST2	CSKS-AQT0135B CSKS-AQT0135K	S/T	Carna (07-117)	VS
STK4 (MST1)	Serine/threonine kinase 4	NM_006282	MST1	CSKS-AQT0135B CSKS-AQT0135K	S/T	SignalChem (S25-10G)	VS
STK10 (LOK)	serine/threonine kinase 10	NM_005990	LOK	CSKS-AQT0505B CSKS-AQT0505K	S/T	Carna (07-315)	VS
STK11 (LKB1, AMPKK)	serine/threonine kinase 11	NM_000455	LKB1	CSKS-AQT0728B CSKS-AQT0728K	S/T	Carna (02-119)	M
STK17A (DRAK1)	Serine/threonine kinase 17a	NM_004760	DRAK1	CSKS-AQT0173B CSKS-AQT0173K	S/T	Carna (02-137)	S
STK17B (DRAK2)	Serine/threonine kinase 17b	NM_004226	DRAK2	CSKS-AQT0174B CSKS-AQT0174K	S/T	SignalChem (S34-10G)	VS
STK24 (MST3)	Serine/threonine kinase 24	NM_003576	MST3	CSKS-AQT0505B CSKS-AQT0505K	S/T	Carna (07-118)	VS
STK25 (YSK1)	serine/threonine kinase 25	NM_006374	YSK1	CSKS-AQT0505B CSKS-AQT0505K	S/T	Carna(07-136)	VS
STK26 (MST4)	serine/threonine kinase 26	NM_016542	MST4	CSKS-AQT0505B CSKS-AQT0505K	S/T	Carna (07-119)	VS
STK38 (NDR1)	serine/threonine kinase 38	NM_007271	NDR1	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-125)	VS
STK38L (NDR2)	serine/threonine kinase 38 like	NM_015000	NDR2	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-126)	VS
SYK	Spleen associated tyrosine kinase	NM_003177	SYK	CSKS-AQT077B CSKS-AQT077K	Y	Carna (08-176)	VS
TAOK1 (TAO1)	TAO kinase 1	NM_020791	TAO1	CSKS-AQT0253B CSKS-AQT0253K	S/T	SignalChem (T24-11G)	M
TAOK2 (TAO2)	TAO kinase 2	NM_016151	TAO2	CSKS-AQT0253B CSKS-AQT0253K	S/T	Carna (07-133)	M
TAOK3 (TAO3)	TAO kinase 3	NM_016281	TAO3	CSKS-AQT0253B CSKS-AQT0253K	S/T	SignalChem (T26-11G)	M
TBK1	TANK-binding kinase 1	NM_013254	TBK1	CSKS-AQT0215B CSKS-AQT0215K	S/T	Carna (05-115)	VS
TEC	Tec protein tyrosine kinase	NM_003215	TEC	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-182)	VS
TEK (TIE2)	Tyrosine kinase with immunoglobulin like and EGF like domains 1	NM_000459	TIE2	CSKS-AQT0101B CSKS-AQT0101K	Y	Carna (08-185)	VS
TNIK (ZC2)	TRAF2 and NCK interacting kinase	NM_015028	ZC2/TNIK	CSKS-AQT0181B CSKS-AQT0181K	S/T	SignalChem (T27-11G)	VS
TNK2 (ACK)	tyrosine kinase non receptor 2	NM_005781	ACK	CSKS-AQT0066B CSKS-AQT0066K	Y	SignalChem (A05-11G)	M
TRPM7 (CHAK1)	transient receptor potential cation channel subfamily M member 7	NM_017672	ChaK1	CSKS-AQT0349B CSKS-AQT0349K	S/T	Carna (10-111)	VS
TSSK1B (TSSK1)	testis specific serine kinase 1B	NM_032028	TSSK1	CSKS-AQT0123B CSKS-AQT0123K	S/T	Carna (02-364)	S
TSSK2	testis specific serine kinase 2	NM_053006	TSSK2	CSKS-AQT0426B CSKS-AQT0426K	S/T	SignalChem (S38-10G)	VS
TSSK3	testis specific serine kinase 3	NM_052841	TSSK3	CSKS-AQT0426B CSKS-AQT0426K	S/T	Carna (02-166)	VS
TTK	TTK protein kinase	NM_003318	TTK	CSKS-AQT0897B CSKS-AQT0897K	S/T	Carna (05-169)	S
TTBK1	tau tubulin kinase 1	NM_032538	TTBK1	CSKS-AQT0686B CSKS-AQT0686K	S/T	SignalChem (T17-11G)	VS
TTBK2	tau tubulin kinase 2	NM_173500	TTBK2	CSKS-AQT0686B CSKS-AQT0686K	S/T	SignalChem (T18-11G)	S

HGNC Common Name	HGNC Approved Full Name	GenBank Acc.No.	Manning Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Enzyme Used in Sensor Peptide Development	Assay Strength*
TXK	TXK tyrosine kinase	NM_003328	TXK	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-183)	VS
TYK2 [JH1]	Tyrosine kinase 2	NM_003331	TYK2	CSKS-AQT0663B CSKS-AQT0663K	Y	Carna (08-147)	S
TYRO3	TYRO3 protein tyrosine kinase	NM_006293	TYRO3	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-109)	VS
ULK1	unc-51 like autophagy activating kinase 1	NM_003565	ULK1	CSKS-AQT0332B CSKS-AQT0332K	S/T	Academic Lab	VS
ULK2	unc-51 like autophagy activating kinase 2	NM_014683	ULK2	CSKS-AQT0332B CSKS-AQT0332K	S/T	Academic Lab	S
YES1 (YES)	YES proto-oncogene 1, Src family tyrosine kinase	NM_005433	YES	CSKS-AQT0104B CSKS-AQT0104K	Y	Carna (08-175)	VS
ZAP70	Zeta-chain (TCR) associated protein kinase 70kDa	NM_001079	ZAP70	CSKS-AQT0734B CSKS-AQT0734K	Y	Carna (08-177)	VS

Table 1.

Summary of PhosphoSens Kinase Target Assay Strength Based on Reaction Rates*					
Categories	Total	Very Strong (VS)	Strong (S)	Medium (M)	Weak (W)
Initial Rate (RFU/pmole/min)		>1000	300-999	100-299	<100
# Assays by Category	384	290	54	40	0
% Assays by Category		76%	14%	10%	0%
% Assays Very Strong or Strong		90%		Being improved	

Note: Substrate concentrations range from 10-15 μ M

Table 2.

Protocol Comments for Optimal Assay
1. 5 ng/ul CaM, 0.4 mM CaCl ₂
2. Activated (+0.5nM SRC)
3. Activated (+ poly(rI:C))
4. Activated (+ tRNA)
5. Uses inactive ERK2
6. Uses full-length ASK1
7. 10 mM KCl
8. 1 mM AMP
9. 1 mM cGMP
10. Uses inactive MEK1/ERK2