

A Powerful Approach for Understanding Kinase and Kinase Inhibitor Activity

Continuous • Direct • Catalytic

*For the end-point format, simply add Europium to the kinetic assay to yield a red-shifted, time-resolved assay to enable correction for compound auto-fluorescence and batch mode processing for HTS.

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
ABL1	ABL proto-oncogene 1, non-receptor tyrosine kinase	NM_005157	ABL1	CSKS-AQT0031B CSKS-AQT0031K	Y	Carna (08-001)
ABL2 (ARG)	ABL proto-oncogene 2, non-receptor tyrosine kinase	NM_005157	ARG	CSKS-AQT0030B CSKS-AQT0030K	Y	Carna (08-102)
AKT1 (PKB)	AKT Serine/Threonine Kinase 1	NM_005163	AKT1	CSKS-AQT0569B CSKS-AQT0569K	S/T	Carna (01-101)
AKT2 (PKBβ)	AKT Serine/Threonine Kinase 2	NM_001626	AKT2	CSKS-AQT0232B CSKS-AQT0232K	S/T	Carna (01-102)
AKT3 (PKBγ)	AKT Serine/Threonine Kinase 3	NM_005465	AKT3	CSKS-AQT0569B CSKS-AQT0569K	S/T	Carna (01-103)
AURKA (AURA)	Aurora kinase A	NM_003600	AurA	CSKS-AQT0455B CSKS-AQT0455K	S/T	Carna (05-101)
AURKB (AURB)	aurora kinase B	NM_004217	AurB	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (05-102)
AURKC (AURC)	aurora kinase C	NM_004217	AurB	CSKS-AQT0186B CSKS-AQT0186K	S/T	Carna (05-103)
BTK	Bruton tyrosine kinase	NM_000061	BTK	CSKS-AQT0663B CSKS-AQT0663K	Y	Carna (08-180)
CAMK4	calcium/calmodulin dependent protein kinase IV	NM_001744	CaMK4	CSKS-AQT0420B CSKS-AQT0420K	S/T	Carna (02-108)
CDC42BPG (DMPK2)	CDC42 binding protein kinase gamma	NM_017525	DMPK2	CSKS-AQT0837B CSKS-AQT0837K	S/T	SignalChem (D17-11G)
CDK1/CycE1	cyclin dependent kinase 1	NM_001786	CDC2	CSKS-AQT0296B CSKS-AQT0296K	S/T	Proqinase (0134-0055-1)
CDK2/CycE1	cyclin dependent kinase 2	NM_001798	CDK2	CSKS-AQT0255B CSKS-AQT0255K	S/T	Carna (04-165)
CDK4 /CycD3	cyclin dependent kinase 4	NM_000075	CDK4	CSKS-AQT0730B CSKS-AQT0730K	S/T	Carna (04-105)
CDK5/p25	cyclin dependent kinase 5	NM_004935	CDK5	CSKS-AQT0730B CSKS-AQT0730K	S/T	Carna (04-106)
CDK6/CycD3	cyclin dependent kinase 6	NM_001259	CDK6	CSKS-AQT0730B CSKS-AQT0730K	S/T	Carna (04-107)
CDK7/CycH/MAT1	cyclin dependent kinase 7	NM_001799	CDK7	CSKS-AQT0733B CSKS-AQT0733K	S/T	Carna (04-108)
CDK9 (TAK) w/CycT1	cyclin dependent kinase 9	NM_001261.2	CDK9	CSKS-AQT0449B CSKS-AQT0449K	S/T	Carna (04-110)
CDK14 (PFTAIRE1)/CycY	cyclin dependent kinase 14	NM_012395	PFTAIRE1	CSKS-AQT0730B CSKS-AQT0730K	S/T	SignalChem (P15-10G)
CDK16/CycY	cyclin dependent kinase 16	NM_006201	PCTAIRE1	CSKS-AQT0730B CSKS-AQT0730K	S/T	Carna (04-116)
CDK17 (PCTAIRE2)/CycY	cyclin dependent kinase 17	NM_002595	PCTAIRE2	CSKS-AQT0730B CSKS-AQT0730K	S/T	SignalChem (P10-10G)

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
CDK18/CycY (PCTAIRE3)	cyclin dependent kinase 18	NM_002596	PCTAIRE3	CSKS-AQT0730B CSKS-AQT0730K	S/T	Thermo (A30492)
CHEK1 (CHK1)	Check point kinase 1	NM_001274	CHK1	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-117)
CHEK2 (CHK2)	Check point kinase 2	NM_007194	CHK2	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-162)
EGFR (ERBB1)	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0663B CSKS-AQT0663K	Y	BPS (40187)
EGFR [C797S/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0663B CSKS-AQT0663K	Y	SignalChem (E10-122ZG)
EGFR [T790M/C797S/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0663B CSKS-AQT0663K	Y	SignalChem (E10-122VG)
EGFR [T790M/L858R]	Epidermal growth factor receptor	NM_005228	EGFR	CSKS-AQT0663B CSKS-AQT0663K	Y	Carna (08-510)
EIF2AK1 (HRI)	eukaryotic translation initiation factor 2 alpha kinase 1	NM_014413	HRI	CSKS-AQT0621B CSKS-AQT0621K	S/T	Carna (05-154)
EIF2AK4 (GCN2)	eukaryotic translation initiation factor 2 alpha kinase 4	NM_001013703	GCN2	CSKS-AQT0911B CSKS-AQT0911K	S/T	Carna (05-153)
ERN1 (IRE1)	endoplasmic reticulum to nucleus signaling 1	NM_001433	IRE1	CSKS-AQT0715B CSKS-AQT0715K	S/T	Thermo (A30971)
FGFR1	Fibroblast growth factor receptor 1	NM_000604	FGFR1	CSKS-AQT0515B CSKS-AQT0515K	Y	Carna (08-133)
FGFR2	Fibroblast growth factor receptor 2	NM_022970.3	FGFR2	CSKS-AQT0515B CSKS-AQT0515K	Y	Carna (08-134)
FGFR3	Fibroblast growth factor receptor 3	NM_000142	FGFR3	CSKS-AQT0515B CSKS-AQT0515K	Y	Carna (08-135)
FGFR4	Fibroblast growth factor receptor 4	NM_002011	FGFR4	CSKS-AQT0026B CSKS-AQT0026K	Y	Carna (08-136)
GRK2 (ADRBK1, BARK1)	G protein-coupled receptor kinase 2	NM_001619	BARK1	CSKS-AQT0641B CSKS-AQT0641K	S/T	Thermo (PV3361)
GSK3B (GSK3β)	glycogen synthase kinase 3 beta	NM_001146156	GSK3B	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-141)
JAK1 [JH1]	Janus kinase 1	NM_002227	JAK1	CSKS-AQT0658B CSKS-AQT0658K	Y	Thermo (PV4774)
LRRK2	leucine rich repeat kinase 2	NM_198578	LRRK2	CSKS-AQT0615B CSKS-AQT0615K	S/T	Thermo (PV48783)
LRRK2 [G2019S]	leucine rich repeat kinase 2	NM_198578	LRRK2	CSKS-AQT0615B CSKS-AQT0615K	S/T	Thermo (PV4881)
MAP3K5 (ASK1)	mitogen-activated protein kinase kinase kinase 5	NM_005923	MAP3K5	CSKS-AQT0557B CSKS-AQT0557K	S/T	SignalChem (M13-10G)
MAP3K8 (Tpl2, COT)	mitogen-activated protein kinase kinase kinase 8	NM_005204	COT	CSKS-AQT0716B CSKS-AQT0716K	S/T	Carna (07-301)
MAP4K1 (HPK1)	mitogen-activated protein kinase kinase kinase kinase 1	NM_007181	HPK1	CSKS-AQT0080B CSKS-AQT0080K	S/T	Thermo (PV6355)
MAP4K4 (HGK, ZC1)	mitogen-activated protein kinase kinase kinase kinase 4	NM_004834	ZC1/HGK	CSKS-AQT0135B CSKS-AQT0135K	S/T	Carna (07-137)
MAPK1 (ERK2)	mitogen-activated protein kinase 1	NM_002745	ERK2	CSKS-AQT0649B CSKS-AQT0649K	S/T	Carna (04-143)
MAPK3 (ERK1)	mitogen-activated protein kinase 3	NM_002746	ERK1	CSKS-AQT0650B CSKS-AQT0650K	S/T	Carna (04-142)
MAPK8 (JNK1)	mitogen-activated protein kinase 8	NM_002750	JNK1	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-163)
MAPK9 (JNK2)	mitogen-activated protein kinase 9	NM_002752	JNK2	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-164)
MAPK10 (JNK3)	mitogen-activated protein kinase 10	NM_002753	JNK3	CSKS-AQT0376B CSKS-AQT0376K	S/T	Carna (04-150)
MAPK11 (p38β)	mitogen-activated protein kinase 11	NM_002751	p38b	CSKS-AQT0490B CSKS-AQT0490K	S/T	Carna (04-153)
MAPK12 (p38γ)	mitogen-activated protein kinase 12	NM_002969	p38g	CSKS-AQT0649B CSKS-AQT0649K	S/T	Carna (04-155)

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
MAPK13 (p38δ)	mitogen-activated protein kinase 13	NM_002754	p38d	CSKS-AQT0649B CSKS-AQT0649K	S/T	Carna (04-154)
MAPK14 (p38α)	mitogen-activated protein kinase 14	NM_001315	p38a	CSKS-AQT0649B CSKS-AQT0649K	S/T	Carna (04-152)
MASTL	microtubule associated serine/threonine kinase-like	NM_032844	MASTL	CSKS-AQT0703B CSKS-AQT0703K	S/T	SignalChem (M72-10G)
MET (c-MET, HGFR)	MET proto-oncogene, receptor tyrosine kinase	NM_000245	MET	CSKS-AQT0658B CSKS-AQT0658K	Y	Carna (08-151)
PAK1	p21 (RAC1) activated kinase 1	NM_002576	PAK1	CSKS-AQT0838B CSKS-AQT0838K	S/T	Carna (07-123)
PAK2	p21 (RAC1) activated kinase 2	NM_002577	PAK2	CSKS-AQT0839B CSKS-AQT0839K	S/T	Carna (07-124)
PAK3	p21 (RAC1) activated kinase 3	NM_002578	PAK3	CSKS-AQT0845B CSKS-AQT0845K	S/T	Carna (07-125)
PAK4	p21 (RAC1) activated kinase 4	NM_005884	PAK4	CSKS-AQT0840B CSKS-AQT0840K	S/T	SignalChem (P05-10BG)
PAK5 (PAK7)	p21 (RAC1) activated kinase 5	NM_020341	PAK5	CSKS-AQT0840B CSKS-AQT0840K	S/T	SignalChem (P07-10G)
PAK6	p21 (RAC1) activated kinase 6	NM_020168	PAK6	CSKS-AQT0840B CSKS-AQT0840K	S/T	Carna (07-128)
PKN1	protein kinase N1	NM_002741	PKN1	CSKS-AQT0509B CSKS-AQT0509K	S/T	Carna (01-144)
PKN2	protein kinase N2	NM_006256	PKN2	CSKS-AQT0462B CSKS-AQT0462K	S/T	Carna (01-145)
PKN3	protein kinase N3	NM_013355	PKN3	CSKS-AQT0509B CSKS-AQT0509K	S/T	Carna (01-146)
PRKD1 (PKD1)	protein kinase D1	NM_002742	PKD1	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-157)
PRKD2 (PKD2)	protein kinase D2	NM_016457	PKD2	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-158)
PRKD3 (PKD3)	protein kinase D3	NM_005813	PKD3	CSKS-AQT0694B CSKS-AQT0694K	S/T	Carna (02-159)
PRKG1 (PKG1, PGK)	protein kinase, cGMP-dependent, type I	NM_006258	DNAPK	CSKS-AQT0324B CSKS-AQT0324K	S/T	Carna (01-142)
Raf, b- (V600E), Direct	B-Raf proto-oncogene, serine/threonine kinase	NM_004333	BRAF	CSKS-AQT0682B CSKS-AQT0682K	S/T	Thermo (PR7304A)
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)
ROCK1	Rho associated coiled-coil containing protein kinase 1	NM_005406	ROCK1	CSKS-AQT0713B CSKS-AQT0713K	S/T	Carna (01-109)
ROCK2	Rho associated coiled-coil containing protein kinase 2	NM_004850	ROCK2	CSKS-AQT0713B CSKS-AQT0713K	S/T	Carna (01-110)
RPS6KA1 (RSK1)	Ribosomal protein S6 kinase A1	NM_002953	RSK3	CSKS-AQT0535B CSKS-AQT0535K	S/T	Carna (01-149)
RPS6KA2 (RSK3)	Ribosomal protein S6 kinase A2	NM_021135	RSK1	CSKS-AQT0535B CSKS-AQT0535K	S/T	Carna (01-151)
RPS6KA3 (RSK2)	Ribosomal protein S6 kinase A3	NM_004586	RSK2	CSKS-AQT0233B CSKS-AQT0233K	S/T	Carna (01-150)
RPS6KB1 (p70S6K)	Ribosomal protein S6 kinase, 70kDa, polypeptide 1	NM_003161	p70S6K	CSKS-AQT0535B CSKS-AQT0535K	S/T	Carna (01-158)
SGK1 (SGK)	Serum/glucocorticoid regulated kinase 1	NM_005627	SGK	CSKS-AQT0566B CSKS-AQT0566K	S/T	Thermo (PV3818)
SGK2	SGK2, serine/threonine kinase 2	NM_016276	SGK2	CSKS-AQT0566B CSKS-AQT0566K	S/T	Carna (01-159)
SGK3	Serum/glucocorticoid regulated kinase family member 3	NM_013257	SGK3	CSKS-AQT0566B CSKS-AQT0566K	S/T	Carna (01-160)
SIK3 (QSK)	salt inducible kinase 3	BC128510.1	QSK	CSKS-AQT0508B CSKS-AQT0508K	S/T	Dundee/MRC (DU 16623)
YES1 (YES)	YES proto-oncogene 1, Src family tyrosine kinase	NM_005433	YES	CSKS-AQT0515B CSKS-AQT0515K	Y	Carna (08-175)