

KinSight Kinome Profiling Services

Continuous • Direct • Catalytic

A Powerful Approach for Understanding Kinase and Kinase Inhibitor Activity

400+ Unique Kinases for Comprehensive Profiling and Screening Results

Contact us at hello@assayquant.com to learn more about our Kinome Profiling Services powered by PhosphoSens Technology.

Add Mutants or Customize Your Panel!

Note: All kinases may not be included on every panel. Please reach out to our team for more information.

Recombinant Enzyme Name Complex, HGNC (Common) Name or [Form]

AAK1	CDK2/CycE1	DYRK4	HIPK3	MAP4K5 (KHS1)	NTRK2 (TRKB)	PRKCD (PKCδ [delta])	STK16 (MPSK1, PKL12, TSF1)
ABL1	CDK2/CycE2	EEF2K (eEF2K)	HIPK4	MAPK1 (ERK2)	NTRK3 (TRKC)	PRKCE (PKCε [epsilon])	STK17A (DRAK1)
ABL2 (ARG)	CDK2/CycO	EGFR (ERBB1)	HUNK	MAPK10 (JNK3)	NUAK1 (ARK5)	PRKCG (PKCγ [gamma])	STK17B (DRAK2)
AKT1 (PKB)	CDK3/cycE1	EIF2AK1 (HRI)	ICK	MAPK11 (p38b)	NUAK2 (SNARK)	PRKCH (PKCη [eta])	STK24 (MST3)
AKT2 (PKBb)	CDK3/CycE2	EIF2AK2 (PKR)	IGF1R	MAPK12 (p38y)	PAK1	PRKCI (PKCι [iota])	STK25 (YSK1)
AKT3 (PKBg)	CDK4 /CycD3	EIF2AK3 (PERK/PEK)	IKKB (IKKβ)	MAPK13 (p38δ)	PAK2	PRKCO (PKCθ [theta])	STK26 (MST4)
ALK	CDK4/CycD1	EIF2AK4 (GCN2)	IKKE (IKKε)	MAPK14 (p38a)	PAK3	PRKCZ (PKCζ [zeta])	STK3 (CLIK1, MST2)
AURKA (AURA)	CDK4/CycD2	EPHA1	INSR (IR)	MAPK15 (ERK7)	PAK4	PRKD1 (PKD1)	STK33
AURKB (AURB)	CDK5/p25	EPHA2	INSRR (IRR)	MAPK3 (ERK1)	PAK5 (PAK7)	PRKD2 (PKD2)	STK38 (NDR1)
AURKC (AURC)	CDK5/p35NCK	EPHA3	IRAK1	MAPK7 (BMK, ERK5)	PAK6	PRKD3 (PKD3)	STK38L (NDR2)
AXL	CDK6/CycD1	EPHA4	IRAK4	MAPK8 (JNK1)	PASK	PRKDC (DNA-PK)	STK4 (MST1)
BCKDK	CDK6/CycD2	EPHA5	ITK	MAPK9 (JNK2)	PBK (TOPK)	PRKG1 (PKG1, PGK)	SYK
BCR RET (fusion)	CDK6/CycD3	EPHA6	JAK1 [JH1]	MAPKAPK2 (MK2)	PDGFRA (PDGFRα)	PRKG2 (PKG2, CGK2)	TAOK1 (TAO1)
BLK	CDK7/CycH/MAT1	EPHA7	JAK1 [JH1-JH2]	MAPKAPK3 (MK3)	PDGFRB (PDGFRβ)	PRKX	TAOK2 (TAO2)
BMP2K (BIKE)	CDK7/CycH1	EPHA8	JAK2 [JH1]	MAPKAPK5 (MK5)	PDK1 (PDHK1)	PTK2 (FAK)	TAOK3 (TAO3)
BMX	CDK9 (TAK)/CycK	EPHB1	JAK2 [JH1-JH2]	MARK1	PDK2 (PDHK2)	PTK2B (PYK2)	TBK1
BRAF (BRAFA1, RAFA1)	CDK9 (TAK)/CycT1	EPHB2	JAK3 [JH1]	MARK2	PDK3 (PDHK3)	PTK6 (BRK)	TEC
BRSK1	CDK9 (TAK)/CycT2	EPHB3	KDR (VEGFR2)	MARK3	PDK4 (PDHK4)	RAF, b-[V600E]	TEK (TIE2)
BRSK2	CDKL5	EPHB4	KIT (SCFR, CD117)	MARK4	PDPK1 (PDK1)	RAF1 [Y340E/Y341E]	TESK2
BTK	CHEK1 (CHK1)	ERBB2 (HER2)	LATS1	MAST3	PHKG1	RET	TLK1
BUB1	CHEK2 (CHK2)	ERBB4 (HER4)	LATS2	MASTL	PHKG2	RIPK1	TLK2
BUB1B (BUBR1)	CHUK (IKKA, IKKα)	ERN1 (IRE1)	LCK	MATK (CTK)	PIM1	RIPK2	TNIK (ZC2)
CAMK1 (CAMK1a)	CIT (CRIK)	ERN2 (IRE2)	LIMK1	MELK	PIM2	RIPK3	TNK1
CAMK1D (CAMK1δ)	CLK1	FAM20C	LIMK2	MERTK (MER)	PIM3	ROCK1	TNK2 (ACK)
CAMK1G (CAMK1γ)	CLK2	FER	LRRK2	MET (c-MET, HGFR)	PINK1 (PINK)	ROCK2	TRPM7 (CHAK1)
CAMK2A (CAMK2α)	CLK3	FES	LTK	MINK1 (MAP4K6, MINK, ZC3)	PKN1	ROS1 (ROS)	TSSK1B (TSSK1)

CAMK2B (CAMK2β)	CLK4	FGFR1	LYNA (LYNa)	MLK4 (MAP3K21)	PKN2	RPS6KA1 (RSK1)	TSSK2
CAMK2D (CAMK2δ)	CSF1R (FMS)	FGFR2	LYNB (LYNb)	MNK1	PKN3	RPS6KA2 (RSK3)	TSSK3
CAMK2G (CAMK2γ)	CSK	FGFR3	MAK	MNK2	PLK1	RPS6KA3 (RSK2)	TSSK6 (SSTK)
CAMK4	CSNK1A1 (CK1A1, CK1α)	FGFR4	MAP2K1 (MEK1)	MST1R (RON)	PLK2	RPS6KA4 (MSK2)	TTBK1
CAMKK1	CSNK1A1L (CK1A1L, CK1a2)	FGR	MAP2K2 (MEK2)	mTOR (FRAP)	PLK3	RPS6KA5 (MSK1)	TTBK2
CAMKK2	CSNK1D (CK1D, CK1δ)	FLT1	MAP2K3 (MEK3)	MUSK	PLK4	RPS6KA6 (RSK4)	TTK
CDC42BP (MRCKA)	CSNK1E (CK1E, CK1ε)	FLT3	MAP3K1 (MEKK1)	MYLK (smMLCK)	PNCK (CAMK1β)	RPS6KB1 (p70S6K)	TXK
CDC42BPB (MRCKB)	CSNK1G1 (CK1G1, CK1γ1)	FLT4	MAP3K10 (MLK2)	MYLK2 (skMLCK)	PRKAA1-B1-G1 (AMPKα1β1γ1)	RPS6KB2 (p70S6KB)	TYK2 [JH1]
CDC42BPG (DMPK2)	CSNK1G2 (CK1G2, CK1γ2)	FRK	MAP3K11 (MLK3)	MYLK3 (caMLCK)	PRKAA1-B1-G2 (AMPKα1β1γ2)	SBK1 (SBK)	TYRO3
CDC7/DBF4 (CDC7)	CSNK1G3 (CK1G3, CK1γ3)	FYNa (FYN)	MAP3K12 (DLK)	MYLK4 (Sgk085)	PRKAA1-B1-G3 (AMPKα1β1γ3)	SGK1 (SGK)	ULK1
CDK1 (CDC2)/CycA1	CSNK2A1 (CK2A1, CK2α1)	FYNb (FYN)	MAP3K13 (LZK)	MYO3A	PRKAA1-B2-G1 (AMPKα1β2γ1)	SGK2	ULK2
CDK1 (CDC2)/CycE1	CSNK2A2 (CK2A2, CK2α2)	GAK	MAP3K14 (NIK)	MYO3B	PRKAA1-B2-G2 (AMPKα1β2γ2)	SGK3	ULK3
CDK1/CycA2 (CDC2)	DAPK1	GRK1 (RHOK)	MAP3K19 (YSK4)	NEK1	PRKAA1-B2-G3 (AMPKα1β2γ3)	SIK1 (SIK)	VRK1
CDK1/CycB1 (CDC2)	DAPK2 (DRP-1)	GRK2 (ADRBK1, BARK1)	MAP3K2 (MEKK2)	NEK10	PRKAA2-B1-G1 (AMPKα2β1γ1)	SIK2 (QIK)	VRK2
CDK12 (CRK7)/CycK	DAPK3 (ZIPK)	GRK3 (ADRBK2, BARK2)	MAP3K20 (MLK7, ZAK)	NEK11	PRKAA2-B1-G2 (AMPKα2β1γ2)	SIK3 (QSK)	WEE1
CDK13 (CHED)/CycK	DCLK1 (DCAMKL1)	GRK4 (GPRK4)	MAP3K3 (MEKK3)	NEK2	PRKAA2-B1-G3 (AMPKα2β1γ3)	SLK	YES1 (YES)
CDK14 (PFTAIRE1, PFTK1)/CycY	DCLK2 (DCAMKL2)	GRK5 (GPRK5)	MAP3K5 (ASK1)	NEK3	PRKAA2-B2-G1 (AMPKα2β2γ1)	SNRK	ZAP70
CDK15 (PFTAIRE2, PFTK2)/CycB1	DDR1	GRK6 (GPRK6)	MAP3K6 (ASK2)	NEK4	PRKAA2-B2-G2 (AMPKα2β2γ2)	SRC	New assays are continuously added! Please inquire if your target is not listed, or if you are interested in a custom panel.
CDK16/CycY (PCTAIRE1, PCTK1)	DDR2	GRK7 (GPRK7)	MAP3K7 (TAK1)	NEK5	PRKAA2-B2-G3 (AMPKα2β2γ3)	SRC N1	
CDK17 (PCTAIRE2, PCTK2)/CycY	DMPK (DMPK1)	GSK3A (GSK3α)	MAP3K8 (COT, Tpl2)	NEK6	PRKACA (PKACA, PKACα)	SRMS (SRM)	
CDK17 (PCTAIRE2, PCTK2)/p35NCK	DSTYK (RIPK5, SGK496)	GSK3B (GSK3β)	MAP3K9 (MLK1)	NEK7	PRKACB (PKACB, PKACβ)	SRPK1	
CDK18/CycY (PCTAIRE3, PCTK3)	DYRK1A	Haspin	MAP4K1 (HPK1)	NEK9	PRKACG (PKACG, PKACγ)	SRPK2	
CDK19 (CDK11)/CycC	DYRK1B	HCK	MAP4K2 (GCK)	NIM1 (NIM1K)	PRKCA (PKCα [alpha])	SRPK3 (MSSK1)	
CDK2/CycA1	DYRK2	HIPK1	MAP4K3 (GLK, KHS2)	NLK (LAK1)	PRKCB1 (PKCβ1 [beta1])	STK10 (LOK)	
CDK2/CycA2	DYRK3	HIPK2	MAP4K4 (H GK, ZC1)	NTRK1 (TRKA)	PRKCB2 (PKCβ2 [beta2])	STK11 (LKB1, AMPKK)	

For questions, please visit our website at www.assayquant.com or email hello@assayquant.com. To place an order, please email orders@assayquant.com.