

KinaseSeeker™ Mixed Panel

Kinase	Group
AKT1	AGC
AKT2	AGC
AKT3	AGC
PDK1/PDPK1	AGC
PKA/PRKACA	AGC
PKC- ϵ /PRKCE	AGC
PKG1/PRKG1	AGC
PKX/PRKX	AGC
RPS6KA3/RSK2	AGC
RPS6KA4/MSK2	AGC
YANK2	AGC
AMPK- α 1/AMPK	CAMK
CAMK1D	CAMK
CAMK2D	CAMK
DAPK3	CAMK
MARK1	CAMK
MARK2	CAMK
PIM1	CAMK
SNF1LK/SIK1	CAMK
SNF1LK2/SIK2/QIK	CAMK
CLK2	CMGC
CDK5	CMGC
p38- α /MAPK14	CMGC
AAK1	Other
AURKA	Other

Kinase	Group
AURKB	Other
AURKC	Other
PLK4	Other
MST2	STE
SLK	STE
YSK1	STE
ABL1	TK
DDR1	TK
EPHA5	TK
EPHB2	TK
FLT1/VEGFR1	TK
FLT3	TK
HCK	TK
IGF1R	TK
ITK	TK
KIT	TK
MUSK	TK
PDGFRB	TK
PTK2B/PYK2	TK
SRC	TK
TNK1	TK
VEGFR2/KDR/FLK1	TK
ACVR2A/ACVR2	TKL
MLK2/MAP3K10	TKL
MLK3/MAP3K11	TKL