

ATP-Max HotSpot™ Kinase Panels

Discover the largest collection of kinase targets in the industry

Now available in radiometric filter binding assay format at 1mM ATP

ATP-Max KinomeScreen

A full panel of 340 kinase targets available at 1mM ATP, for comprehensive and deep interrogation of your compound's selectivity against the whole human kinome. Available as a full panel, or select targets to create your own custom panel.

Kinase Family	ATP-Max KinomeScreen Targets
AG	MSK2/RPS6KA4
AGC	AKT1, AKT2, AKT3, DMPK2, GRK2, GRK3, GRK4, GRK5, GRK6, GRK7, LATS1, LATS2, LYN-B, MAST3, MASTL, MRCK α /CDC42BPA, MRCK β /CDC42BPB, MSK1/RPS6KA5, P70S6K/RPS6KB1, P70S6KB/RPS6KB2, PDK1/PDPK1, PKA, PKACB, PKACG, PKC α , PKC β 1, PKC β 2, PKC δ , PKC ϵ , PKC η , PKC γ , PKC ι , PKC θ , PKC ζ , PKG1 α , PKG1 β , PKG2/PRKG2, PKN1/PRK1, STK32B/YANK2, PRKX, ROCK1, ROCK2, RSK1, RSK2, RSK3, RSK4, SGK1, SGK2, SGK3/SGKL, STK21/CIT, STK38/NDR1, STK38L/NDR2
CAMK	ARK5/NUAK1, BRK1, BRK2, CAMK1A, CAMK1B, CAMK1D, CAMK1g, CAMK2a, CAMK2b, CAMK2D, CAMK2G, CAMK4, CHK1, CHK2, DAPK1, DAPK2, DCAMKL1, DCAMKL2, LKB1, MAPKAPK2, MAPKAPK3, MAPKAPK5/PRAK, MARK1, MARK2/PAR-1BA, MARK3, MARK4, MELK, MLCK/MYLK, MLCK2/MYLK2, MNK1, MNK2, MYLK3, MYLK4, NIM1, PASK, PHKG1, PHKG2, PIM1, PIM2, PIM3, PKCMU/PRKD1, PKCNU/PRKD3, PKD2/PRKD2, SIK1, SIK2, SIK3, SNARK/NUAK2, SNRK, STK22D/TSSK1, TSSK2, TSSK3/STK22C, ZIPK/DAPK3
CK1	CK1 α 1, CK1 α 1L, CK1 δ , CK1 ϵ , CK1 γ 1, CK1 γ 2, CK1 γ 3, TTBK1, TTBK2, VRK2
CMGC	CDK1/CYCLIN A, CDK1/CYCLIN B, CDK1/CYCLIN E, CDK14/CYCLIN Y (PFTK1), CDK15/CYCLIN A2, CDK15/CYCLIN B1, CDK16/CYCLIN Y (PCTAIRE), CDK17/CYCLIN Y (PCTK2), CDK18/CYCLIN Y (PCTK3), CDK2/CYCLIN A, CDK2/CYCLIN A1, CDK2/CYCLIN E, CDK2/CYCLIN E2, CDK2/CYCLIN O, CDK3/CYCLIN C, CDK3/CYCLIN E, CDK3/CYCLIN E2, CDK4/CYCLIN D1, CDK4/CYCLIN D2, CDK4/CYCLIN D3, CDK5/P25, CDK5/P35, CDK6/CYCLIN D1, CDK6/CYCLIN D2, CDK6/CYCLIN D3, CDK7/CYCLIN H, CDK9/CYCLIN K, CDK9/CYCLIN T1, CDK9/CYCLIN T2, CK2A, CK2A2, CLK1, CLK2, CLK3, CLK4, DYRK1/DYRK1A, DYRK1B, DYRK2, DYRK3, ERK1, ERK2/MAPK1, ERK5/MAPK7, ERK7/MAPK15, GSK3A, GSK3B, HIPK1, HIPK2, HIPK3, HIPK4, JNK1, JNK2, JNK3, LCK2/ICK, MAK, MSSK1/STK23, P38A/MAPK14, P38B/MAPK11, P38d/MAPK13, P38G, SRPK1, SRPK2
STE	ASK1/MAP3K5, COT1/MAP3K8, GCK/MAP4K2, GLK/MAP4K3, HGK/MAP4K4, HPK1/MAP4K1, KHS/MAP4K5, LOK/STK10, MEK1, MEK3, MEK5, MEKK1, MEKK2, MEKK3, MINK/MINK1, MKK4, MKK6, MKK7, MST1/STK4, MST2/STK3, MST3/STK24, MYO3A, MYO3B, OSR1/OXSR1, PAK1, PAK2, PAK3, PAK4, PAK5, PAK6, SLK/STK2, STK25/YSK1, STK39/STLK3, TAOK1, TAOK2/TAO1, TAOK3/JIK, TNIK, YSK4/MAP3K19

Kinase Family	ATP-Max KinomeScreen Targets
TK	ABL1, ABL2/ARG, ACK1, ALK, AXL, BLK, BMX/ETK, BRK, BTK, C-KIT, C-MER, C-MET, CSK, C-SRC, CTK/MATK, DDR2, EGFR, EPHA1, EPHA2, EPHA3, EPHA4, EPHA5, EPHA6, EPHA7, EPHA8, EPHB1, EPHB2, EPHB3, EPHB4, ERBB2/HER2, ERBB4/HER4, FAK/PTK2, FER, FES/FPS, FGFR1, FGFR2, FGFR3, FGFR4, FGR, FLT1/VEGFR1, FLT3, FLT4/VEGFR3, FMS, FRK/PTK5, FYN, HCK, IGF1R, IR, IRR/INSRR, ITK, JAK1, JAK2, JAK3, KDR/VEGFR2, LCK, LYN, MUSK, PDGFRA, PDGFRB, PYK2, RET, RON/MST1R, ROS/ROS1, SRMS, SYK, TEC, TIE2/TEK, TRKA, TRKB, TRKC, TXK, TYK1/LTK, TYK2, TYRO3/SKY, YES/YES1, ZAP70
TKL	ALK4/ACVR1B, ALK5/TGFBR1, ALK6/BMPR1B, ARAF, BMPR2, BRAF, IRAK1, IRAK4, LIMK1, LIMK2, LRRK2, MLK1/MAP3K9, MLK2/MAP3K10, MLK3/MAP3K11, MLK4, PBK/TOPK, RAF1, RIPK2, RIPK4, RIPK5, TESK1, TESK2, TGFBR2, ZAK/MLTK
Other	AUR-A, AUR-B, AUR-C, CAMKK2, ERN1/IRE1, ERN2/IRE2, HASPIN, IKK β /IKKB, IKK ϵ /IKBKE, NEK1, NEK2, NEK3, NEK4, NEK5, NEK6, NEK7, NEK9, PLK1, PLK2, PLK3, TBK1, TLK1, TLK2, ULK1, ULK2, ULK3, WNK3

ATP-Max Diversify

A tailored panel of representative isoforms from each major kinase family available at 1 mM ATP, for quick understanding of your compound's selectivity

Kinase Family	ATP-Max Diversify Targets
AGC	MAST3, ROCK1, PKA α , PKC α , PKC θ , Akt1/PKB α , RSK1, p70S6K, MSK2
CAMK	caMLCK, DAPK2, TSSK1, Pim3, Chk1, ARK5, MARK2, MNK2, CAMK1 α , DCAMKL1
CK1	CK1 α , CK2 α , CK1 γ 1
CMGC	ERK1, p38 α , CDK4, JNK3, CDK7, CDK9, CDK1, GSK3 β , CLK2, HIPK2, DYRK1A
TKL	Mlk1, ALK4, BRAF, TGFBR2, LIMk1, IRAK4
TK	Abl1, BTK, TEC, Lyn, Src, EphA5, EphB2, IGF1R, TrkA, Met, Ret, FGFR2, FLT1/VEGFR1, Kit, FLT3, EGFR, Jak2
STE	MST1, LOK, PAK4, MEK1/MAP2K1, MEKK1
Lipid Kinases	PI3K α , PI3K β , PI3K δ , PI3K γ
Atypical Kinases	AURA, Nek1, ULK1, IKK ϵ , PLK1

Discuss your project directly with our kinase experts



v1.0 | Feb 2024

Reaction Biology Corp.

- 1 Great Valley Parkway, Suite 2 Malvern, 19355, PA, USA
- sales@reactionbiology.com
- +1.610.722.0247
- www.reactionbiology.com

