

Technology Spotlight and Applications

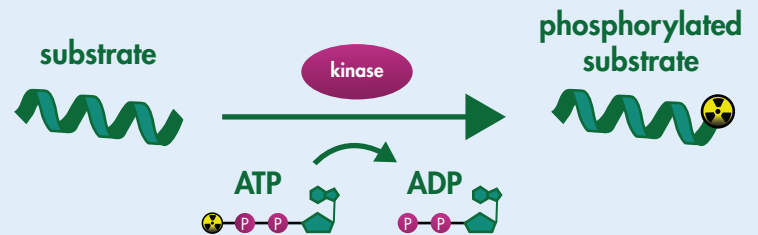
# Radiometric Filter Binding Assay

The gold standard for kinase profiling just got better

## Background Info

The radiometric filter binding assay is **the gold standard method** to study the activity of your target kinase by directly measuring its phosphorylated substrate products, without the use of modified substrates, coupling enzymes, or detection antibodies.

Radiometric assays are not kinase dependent and **can be applied to any kinase inhibitor** thus representing a truly universal assay that offers **consistent and robust results** with **low background signal**.



**1mM ATP now available on 355+ wild type kinase targets**

Reaction Biology's radiometric assay protocols are further optimised to deliver best-in-class quality data, minimise error rates and avoid false positive and negatives caused by other assay formats.

Our industry-leading radiometric assay can be run using **physiologically relevant** concentrations of ATP (1mM) in addition to previously available concentrations of 1µM, 10µM, or apparent ATP-K<sub>m</sub> up to 100µM.

## Why Choose a Radiometric Assay

	Measures Kinase Activity	Detects all types of inhibitors, including substrate-specific inhibitors	Accommodates both peptide and protein substrates	No modified substrates/Additional detection reagents	Universally applicable to all kinases
<b>Radiometric Filter Binding Assay</b>	✓	✓	✓	✓	✓
<b>FRET Peptide</b>	✓	✓	✗	✗	✓
<b>Luminescence</b>	✓	✓	✓	✗	✓
<b>Mobility Shift</b>	✓	✓	✗	✗	✗
<b>Competition Binding</b>	✗	Suboptimal	✗	✗	✗

## Why 1mM ATP?

### Relevant and meaningful results on cellular signaling and regulatory mechanisms

Selecting 1mM ATP concentration in radiometric kinase assays can provide a more accurate representation of physiological conditions, ensuring that the obtained results are more relevant and meaningful for understanding your compound's kinase inhibitory activity.

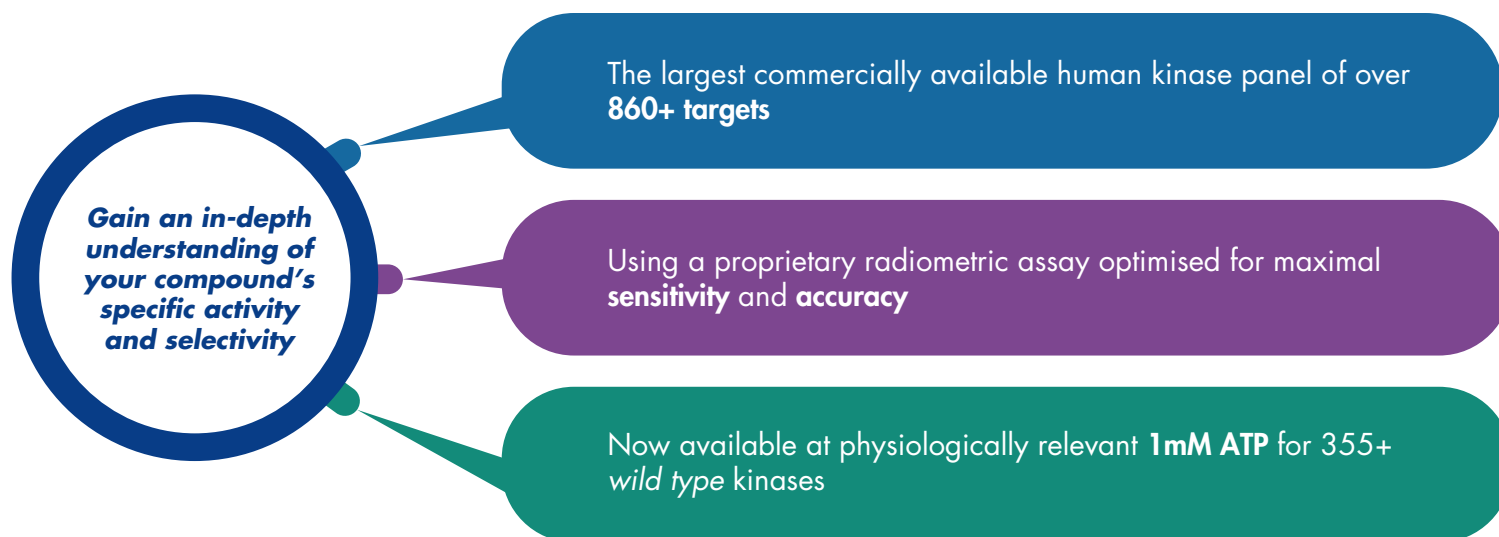
### Here's why our 1mM ATP kinase screening solution stands out:

- 1 True-to-Life Conditions:** our assays mimic typical intracellular ATP concentrations in mammalian cells, which are 1mM or higher
- 2 Maximize Your Kinase Inhibitor Potential:** our conditions guarantee an accurate estimation of your target's kinase inhibitory activity and kinetics.
- 3 Stability Guaranteed:** physiological ATP concentrations are required to maintain some kinase's stability, leading to accurate assay results.
- 4 Relevance to Cell and *In Vivo* Data:** obtain quick biochemical data applicable to follow up cell-based testing and *in vivo* studies

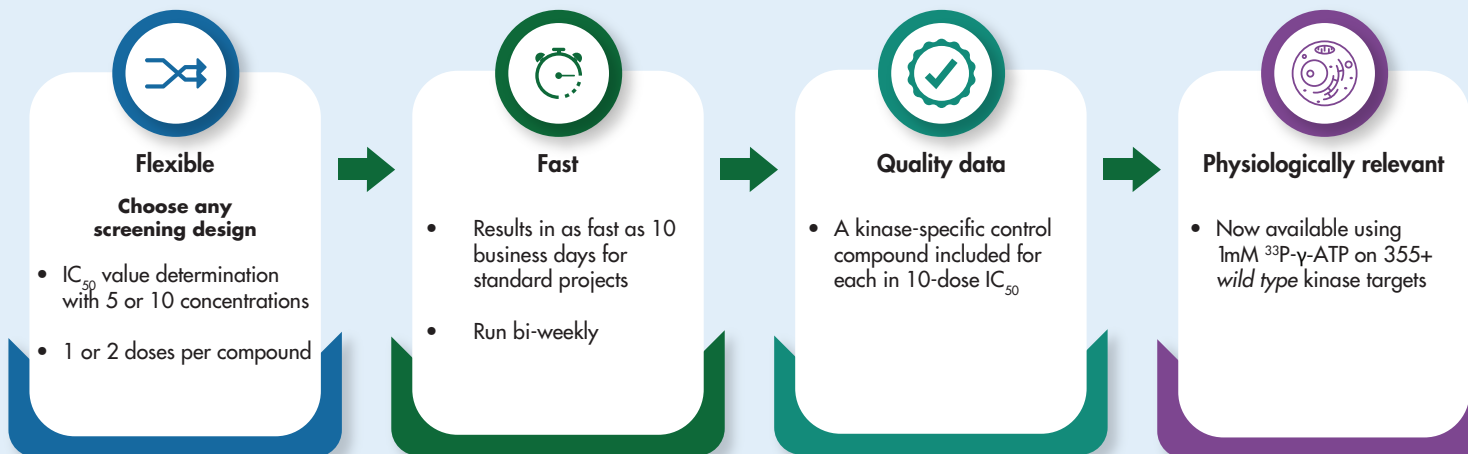
## Our Solution

### HotSpot™ Kinase Screening

*"The New Gold Standard, Only at Reaction Biology"*



## Discover the Benefits



# ATP-Max KinomeScreen

A full panel of 355+ 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 Kinome Targets
<b>AG</b>	MSK2/RPS6KA4
<b>AGC</b>	AKT1, AKT2, AKT3, DMPK2, GRK2, GRK3, GRK4, GRK5, GRK6, GRK7, LATS1, LATS2, LYN-B, MAST3, MASTL, MRCK $\alpha$ /CDC42BPA, MRCK $\beta$ /CDC42BPB, MSK1/RPS6KA5, P70S6K/RPS6KB1, P70S6KB/RPS6KB2, PDK1/PDPK1, PKA, PKACB, PKACG, PKC $\alpha$ , PKC $\beta$ 1, PKC $\beta$ 2, PKC $\delta$ , PKC $\epsilon$ , PKC $\eta$ , PKC $\gamma$ , PKC $\iota$ , PKC $\theta$ , PKC $\zeta$ , PKG1 $\alpha$ , PKG1 $\beta$ , PKG2/PRKG2, PKN1/PRK1, STK32B/YANK2, PRKX, ROCK1, ROCK2, RSK1, RSK2, RSK3, RSK4, SGK1, SGK2, SGK3/SGKL, STK21/CIT, STK32A/YANK1, STK38/NDR1, STK38L/NDR2
<b>CAMK</b>	ARK5/NUAK1, BRSK1, BRSK2, CAMK1A, CAMK1B, CAMK1D, CAMK1g, CAMK2 $\alpha$ , CAMK2b, CAMK2D, CAMK2G, CAMK4, CHK1, CHK2, DAPK1, DAPK2, DCAMKL1, DCAMKL2, DCAMKL3, HUNK, 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, TSSK4, ZIPK/DAPK3
<b>CK1</b>	CK1 $\alpha$ 1, CK1 $\alpha$ 1L, CK1 $\delta$ , CK1 $\epsilon$ , CK1 $\gamma$ 1, CK1 $\gamma$ 2, CK1 $\gamma$ 3, TTBK1, TTBK2, VRK2
<b>CMGC</b>	CDK1/CYCLIN A, CDK1/CYCLIN B, CDK1/CYCLIN E, CDK11B/CYCLIN K, CDK14/CYCLIN Y (PFTK1), CDK15/CYCLIN A2, CDK15/CYCLIN B1, CDK16/CYCLIN Y (PCTAIRE), CDK17/CYCLIN Y (PCTK2), CDK18/CYCLIN Y (PCTK3), CDK19/CYCLIN C, 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, CDK8/CYCLIN C, 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
<b>STE</b>	ASK1/MAP3K5, COT1/MAP3K8, GCK/MAP4K2, GLK/MAP4K3, HGK/MAP4K4, HPK1/MAP4K1, KHS/MAP4K5, LOK/STK10, NIK/MAP3K14, 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
<b>TK</b>	ABL1, ABL2/ARG, ACK1, ALK, AXL, BLK, BMX/ETK, BRK, BTK, C-KIT, C-MER, C-MET, CSK, C-SRC, CTK/MATK, DDR1, 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, LYNB, MUSK, PDGFRA, PDGFRB, PYK2, RET, RIPK1, RON/MST1R, ROS/ROS1, SRMS, SYK, TEC, TIE2/TEK, TNK1, TRKA, TRKB, TRKC, TXK, TYK1/LTK, TYK2, TYRO3/SKY, YES/YES1, ZAP70
<b>TKL</b>	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
<b>Other</b>	AUR-A, AUR-B, AUR-C, CAMKK2, DNA-PK, ERN1/IRE1, ERN2/IRE2, GAK, HASPIN, IKK $\beta$ /IKKBK, IKK $\epsilon$ /IKBKE, NEK1, NEK2, NEK3, NEK4, NEK5, NEK6, NEK7, NEK9, PKMYT1, PLK1, PLK2, PLK3, STK35/CLK1, TBK1, TLK1, TLK2, ULK1, ULK2, ULK3, WEE1, WEE2, 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
<b>AGC</b>	MAST3, ROCK1, PKA $\alpha$ , PKC $\alpha$ , PKC $\theta$ , Akt1/PKB $\alpha$ , RSK1, p70S6K, MSK2
<b>CAMK</b>	caMLCK, DAPK2, TSSK1, Pim3, Chk1, ARK5, MARK2, MNK2, CAMK1 $\alpha$ , DCAMKL1
<b>CK1</b>	CK1 $\alpha$ , CK2 $\alpha$ , CK1 $\gamma$ 1
<b>CMGC</b>	ERK1, p38 $\alpha$ , CDK4, JNK3, CDK7, CDK9, CDK1, GSK3 $\beta$ , CLK2, HIPK2, DYRK1A
<b>TKL</b>	Mlk1, ALK4, BRAF, TGF $\beta$ R2, LIMk1, IRAK4
<b>TK</b>	Abl1, BTK, TEC, Lyn, Src, EphA5, EphB2, IGF1R, TrkA, Met, Ret, FGFR2, FLT1/VEGFR1, Kit, FLT3, EGFR, Jak2
<b>STE</b>	MST1, LOK, PAK4, MEK1/MAP2K1, MEKK1
<b>Lipid Kinases</b>	PI3K $\alpha$ , PI3K $\beta$ , PI3K $\delta$ , PI3K $\gamma$
<b>Atypical Kinases</b>	AurA, Nek1, ULK1, IKK $\epsilon$ , PLK1

Discuss your project directly with our kinase experts



March 2026

1 Great Valley Parkway, Suite 2 Malvern, 19355, PA, USA  
requests@reactionbiology.com  
+1.610.722.0247  
www.reactionbiology.com  
Reaction Biology Corporation

Scan to learn more

