



TARGET-SPECIFIC ASSAYS

- Kinase Assays
- Epigenetic Assays
- RAS Assays
- Targeted Protein Degradation Assays
- Protease Assays
- Phosphatase Assays
- Ion Channels
- and more





Reaction Biology has provided their service for numerous projects with us, especially on multiple kinase assays which have guided our way through drug development. I was very impressed by their level of expertise and professionalism. The insightful advice they provide has been instrumental to our successes. The reproducibility, quality, and reporting of the results have been outstanding throughout the many years we have been working together.

Dr Laurent Meijer

Chair & Chief Scientific Officer at
Perha Pharmaceuticals & ManRos Therapeutics



Target-Specific Assays

Reaction Biology offers a variety of target-specific assays for drug discovery. Most of our assays are high-throughput compatible and can be customized to your specific requirements.

Kinase Assays

Kinase Screening - FreeChoice	p. 15
Kinase Screening - Panels	p. 16
KinaseFinder	p. 18
Kinase SubstrateFinder	p. 19
NanoBRET Intracellular Kinase Assay	p. 20
Cellular Phosphorylation Assay	p. 22
BaF3 Cell Proliferation Assay	p. 24
In Vivo Kinase Activity Models	p. 26

Epigenetic Assays	
Reader Domain Assays p. 29
Methyltransferase Assays p. 33
Demethylase Assays p. 35
Histone Acetyltransferase Assays p. 36
Histone Deacetylase Assays p. 37
Cell-Based Epigenetic Assays p. 38
RAS-related Assays p. 39
Targeted Protein Degradation Assays p. 40
Ubiquitin-Proteasome Assays p. 41
PARP Assays p. 42
Protease Assays p. 43
Phosphatase Assays p. 44
Phosphodiesterase (PDE) Assays p. 45
Metabolic Pathway Assays p. 46
Ion Channel Assays p. 47
GPCR Assays p. 48
Nuclear Receptor Assays p. 51
Additional Assays p. 52
Customized Kinase Drug Discovery p. 53



Your trusted CRO for kinase drug discovery

Reaction Biology offers the most used and most valued kinase discovery platform in the industry comprising enzymatic, biophysical and cell-based assays as well as in vivo models that target kinase-related diseases.

Throughout the history of our company, our scientists were dedicated to use their expertise to advance kinase inhibitors which are now used in the clinic on a daily basis.

Our clients have identified us as their favorite CRO

HTStec Kinase Profiling Trends Survey has reported that the Reaction Biology kinase screening is the most used platform in industry. Almost half of those interviewed named Reaction Biology their favourite CRO.

Quality of assays

Reaction Biology uses state-of-the-art radioactivity-based assays for enzymatic kinase inhibitor screening. The assay directly measures the activity of the kinase with high reproducibility and avoids false negative and positives common with other kinase assay formats.

Customer service

We understand the needs of small and large research organizations alike. Our scientists are dedicated to providing immediate technical support and advice for every research project we undertake. Call us anytime.

Flexibility

We provide fee-for-service, customized projects as well as integrated drug discovery services. We can custom-tailor existing assays and develop new assays making us the CRO of choice for every project.

Expertise

Reaction Biology has been serving clients for over 20 years. Our staff includes expert scientists from a variety of research backgrounds who are available to assist you at each phase of the drug discovery process: hit identification, hit to lead and lead optimization.

Kinase assays at Reaction Biology screening facilities

With the largest portfolio of kinase assays available for drug discovery and 20 years of expertise in custom-tailored assay development, Reaction Biology is guaranteed to provide the kinase assay you need.

Explore our portfolio of 860+ kinase assays offered with a variety of assay formats:

Biochemical assays:

- Radiometric activity assays (HotSpot performed in US, ³³PanQinase performed in Germany)
- Luminescence-based activity assays (ADP-Glo performed in US and Germany)
- HTRF-based activity assay and Probe-Displacement assay (performed in US)

Cell-based assays:

- Kinase activity testing via Cell Phosphorylation assay format
- Target engagement via NanoBRET assay format

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
AAK1	NB, PD	ACVR1 (R206H)	HS, NB	ALK (F1174S)	HS, PQ, ADP-Glo
ABL1	HS, PQ, NB, ADP-Glo	ACVR1B	HS, PQ, NB	ALK (G1202R)	HS, PQ, ADP-Glo
ABL1 (E255K)	HS, PQ, NB, ADP-Glo	ACVR2A	PQ	ALK (G1269A)	HS
ABL1 (E255V)	HS	ACVR2B	PQ	ALK (G1269S)	HS
ABL1 (F317I)	HS, PQ, NB, ADP-Glo	ACVRL1	HS, PQ, NB	ALK (I1171N/D1203N)	HS
ABL1 (F317L)	HS, NB	ADK	NB	ALK (L1152R)	HS
ABL1 (G250E)	HS, PQ, ADP-Glo	AKT1	HS, PQ, NB, CPA, ADP-Glo	ALK (L1196M)	HS, PQ, ADP-Glo
ABL1 (H396P)	HS, PQ, NB, ADP-Glo	AKT1 (aa106-480)	PQ, ADP-Glo	ALK (L1196M/G1202R)	HS
ABL1 (M351T)	HS, PQ, NB, ADP-Glo	AKT1 (E17K)	HS, NB	ALK (R1275Q)	HS, PQ, ADP-Glo
ABL1 (Q252H)	HS, PQ, NB, ADP-Glo	AKT2	HS, PQ, NB, ADP-Glo	ALK (S1206R)	HS
ABL1 (T315I)	HS, PQ, NB, ADP-Glo	AKT2 (aa107-481)	PQ, ADP-Glo	ALK (T1151-L1152insT)	HS
ABL1 (V299L)	HS	AKT2 (E17K)	HS, NB	ALK (T1151M)	HS
ABL1 (Y245F)	HS	AKT3	HS, PQ, ADP-Glo	ALK-EML4	HS, PQ, ADP-Glo
ABL1 (Y245F/Y412F)	HS	AKT3 (aa106-479)	PQ, ADP-Glo	ALK-EML4 (Eex13Aex20)	HS
ABL1 (Y253F)	HS, PQ, NB, ADP-Glo	AKT3 (E17K)	HS, NB	ALK-EML4 (Eex14Aex20)	HS
ABL1 (Y253H)	HS	AKT3 (G171R)	HS, NB	ALK-EML4 (Eex20Aex20)	HS
ABL2	HS, PQ, NB, ADP-Glo	ALK	HS, PQ, CPA, ADP-Glo	ALK-EML4 (Eex6Aex20)	HS
ACVR1	HS, PQ, NB, ADP-Glo	ALK (C1156Y)	HS, PQ, ADP-Glo	ALK-KIF5B	HS
ACVR1 (G328V)	NB	ALK (F1174L)	HS, PQ, ADP-Glo	ALK-KLC1	HS
ACVR1 (G356D)	NB	ALK (F1174L)-EML4	HS, PQ, ADP-Glo	ALK-NPM1	HS, PQ, ADP-Glo
ACVR1 (Q207D)	HS, NB	ALK (F1174L)-NPM1	HS, PQ, ADP-Glo	ALK-STRN	HS

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
ALK-TFG	HS	BTK	HS, PQ, NB, ADP-Glo	CDK5-CDK5R1 (p25)	HS, PQ, ADP-Glo
ALK-TPM1	HS	BTK (C481S)	HS, NB	CDK5-CDK5R2	NB
ALK-TPM3	HS	BTK (E41K)	HS, NB	CDK6-CCND1	HS, PQ, NB, ADP-Glo
ARAF (Y301D/Y302D)	HS, PQ, ADP-Glo	BTK (P190K)	HS, NB	CDK6-CCND2	HS, PQ, ADP-Glo
ATM	HTRF, CPA	BTK (T474I)	HS	CDK6-CCND3	HS, PQ, NB, ADP-Glo
ATR	HTRF	BTK (Y485F)	HS	CDK7	NB
AURKA	HS, PQ, NB, ADP-Glo	BUB1B	PQ	CDK7-CCNH	NB
AURKB	HS, PQ, NB, CPA, ADP-Glo	CAMK1	HS, NB	CDK7-CCNH-MNAT1	HS, PQ, ADP-Glo
AURKB (G160L)	HS	CAMK1D	HS, PQ, NB	CDK8-CCNC	HS, PQ, NB, ADP-Glo
AURKC	HS, PQ, NB, ADP-Glo	CAMK1G	HS, NB	CDK9-CCNK	HS, PQ, NB, ADP-Glo
AXL	HS, PQ, NB, CPA, ADP-Glo	CAMK2A	HS, PQ, NB	CDK9-CCNT1	HS, PQ, NB, ADP-Glo
AXL (R199C)	HS	CAMK2B	HS, PQ	CDK9-CCNT2	HS, PQ, NB
BCR-ABL1	CPA	CAMK2D	HS, PQ, NB	CDK10-CCNL2	NB
BLK	HS, PQ, NB, ADP-Glo	CAMK2G	HS, PQ, NB	CDK10-CCNQ	PQ, ADP-Glo
BMP2K	NB	CAMK4	HS, PQ	CDK11A-CCNK	NB
BMPR1A	HS, PQ, NB, ADP-Glo	CAMKK1	HS, PQ	CDK11A-CCNL2	NB
BMPR1B	HS, PQ, ADP-Glo	CAMKK2	HS, PQ	CDK11B-CCNK	HS, PQ, ADP-Glo
BMPR2	HS, ADP-Glo, PD	CASK	PD	CDK12 (R722C)-CCNK	HS, PQ, ADP-Glo
BMX	HS, PQ, NB	CDC42BPA	HS, PQ	CDK12-CCNK	HS, PQ, NB, ADP-Glo
BRAF	HS, PQ, ADP-Glo	CDC42BPB	HS, PQ, ADP-Glo	CDK13-CCNK	HS, PQ, NB, ADP-Glo
BRAF (d485-489/P490Y)	HS	CDC42BPG	HS	CDK14-CCNY	HS, PQ, NB, ADP-Glo
BRAF (G464V)	HS	CDC7-DBF4	HS, PQ, ADP-Glo	CDK15-CCNA2	HS, PQ, ADP-Glo
BRAF (G469A)	HS	CDK1-CCNA1	PQ	CDK15-CCNB1	HS, PQ, ADP-Glo
BRAF (K601E)	HS	CDK1-CCNA2	HS, PQ, ADP-Glo	CDK15-CCNY	NB
BRAF (L597V)	HS	CDK1-CCNB1	HS, PQ, NB, ADP-Glo	CDK16-CCNY	HS, PQ, NB, ADP-Glo
BRAF (R506_K507 ^{ins} VLR)	HS	CDK1-CCNE1	HS, PQ, NB, ADP-Glo	CDK17-CCNY	HS, PQ, NB
BRAF (T599_V600 ^{ins} T)	HS	CDK2-CCNA1	HS, PQ, NB	CDK17-CDK5R1	PQ, ADP-Glo
BRAF (V600A)	HS	CDK2-CCNA2	HS, PQ, ADP-Glo	CDK18-CCNY	HS, PQ, NB, ADP-Glo
BRAF (V600D)	HS	CDK2-CCND1	PQ, ADP-Glo	CDK19-CCNC	HS, PQ, NB, ADP-Glo
BRAF (V600E)	HS, PQ, NB, CPA, ADP-Glo	CDK2-CCNE1	HS, PQ, NB, ADP-Glo	CDK20-CCNH	PQ, NB, ADP-Glo
BRAF (V600K)	HS	CDK2-CCNE2	HS, PQ	CDK20-CCNT1	PQ, ADP-Glo
BRAF-FAM131B	HS	CDK2-CCNO	HS	CDKL1	NB
BRAF-KIAA1549 (Kex15Bex9)	HS	CDK3-CCNC	HS, PQ, ADP-Glo	CDKL2	NB
BRAF-KIAA1549 (Kex16Bex9)	HS	CDK3-CCNE1	HS, PQ, NB, ADP-Glo	CDKL3	NB
BRAF-SRGAP3	HS	CDK3-CCNE2	HS, PQ	CDKL5	NB
BRSK1	HS, PQ, NB	CDK4-CCND1	HS, PQ, NB, ADP-Glo	CEP43-FGFR1	HS
BRSK2	HS, PQ, NB, ADP-Glo	CDK4-CCND2	HS, PQ, ADP-Glo	CHEK1	HS, PQ, NB, ADP-Glo
		CDK4-CCND3	HS, PQ, NB, ADP-Glo	CHEK2	HS, PQ, NB
		CDK5-CDK5R1	HS, PQ, NB, ADP-Glo	CHEK2 (I157T)	HS

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
CHUK	HS, PQ	DGKI	ADP-Glo (US)	EGFR (d747-749/A750P)	HS, PQ, CPA, ADP-Glo
CILK1	HS, PQ, NB, ADP-Glo	DGKK	ADP-Glo (US)	EGFR (d747-752/P753S)	HS, PQ, ADP-Glo
CIT	HS, PQ, ADP-Glo	DGKQ	ADP-Glo (US)	EGFR (d752-759)	HS, PQ, CPA, ADP-Glo
CK1α1 (E42C)	HS	DGKZ	ADP-Glo (US)	EGFR (D761Y)	HS
CK1α1 (I35C)	HS	DMPK	HS, PQ	EGFR (D770_N771 insG)	HS
CLK1	HS, PQ, NB, ADP-Glo	DSTYK	HS, PQ	EGFR (D770_N771 insNPG)	HS
CLK2	HS, PQ, NB	DYRK1A	HS, PQ, NB	EGFR (D770_N771 insNPG/ T790M)	HS
CLK3	HS, PQ, ADP-Glo	DYRK1B	HS, PQ, NB	EGFR (D770GY)	HS
CLK4	HS, PQ, NB	DYRK2	HS, PQ, NB	EGFR (G719C)	HS, PQ, ADP-Glo
CLK4 (I363V)	HS	DYRK3	HS, PQ, NB	EGFR (G719D)	HS
COQ8B	NB	DYRK4	HS, PQ	EGFR (G719S)	HS, PQ, CPA, ADP-Glo
CSF1R	HS, PQ, NB, ADP-Glo	EEF2K	HS, PQ	EGFR (K716A)	HS
CSK	HS, PQ, NB, ADP-Glo	EGFR	HS, PQ, CPA, BaF3, ADP-Glo	EGFR (K716A/C797S/L858R)	HS
CSNK1A1	HS, PQ, ADP-Glo	EGFR (A763_Y764insFHEA)	HS	EGFR (K716A/T790M/C797S/ L858R)	HS
CSNK1A1L	HS, NB	EGFR (A763_Y764insFQEA)	HS	EGFR (K716Q/L718Q)	HS
CSNK1D	HS, PQ, NB	EGFR (A767_S768insTLA)	HS	EGFR (K728A)	HS
CSNK1E	HS, PQ, NB	EGFR (C775S/T790M/L858R)	HS	EGFR (K728A/T790M/C797S/ L858R)	HS
CSNK1E (R178C)	HS	EGFR (C797A)	HS	EGFR (L718Q)	HS, PQ, ADP-Glo
CSNK1G1	HS, PQ	EGFR (C797G/L858R)	HS	EGFR (L747S)	HS
CSNK1G2	HS, PQ, NB	EGFR (C797S)	HS, PQ, ADP-Glo	EGFR (L792F)	HS
CSNK1G3	HS, PQ	EGFR (C797S/L858R)	HS, PQ, ADP-Glo	EGFR (L792F/L858R)	HS
CSNK2A1	HS, PQ, NB, ADP-Glo	EGFR (d746)	HS	EGFR (L792H)	HS
CSNK2A2	HS, PQ, NB, ADP-Glo	EGFR (d746-750)	HS, PQ, ADP-Glo	EGFR (L792H/C797S/L858R)	HS
DAPK1	HS, PQ, ADP-Glo	EGFR (d746-750/C775S/ T790M/L858R)	HS, PQ	EGFR (L792H/L858R)	HS
DAPK2	HS, PQ, NB	EGFR (d746-750/C797A)	HS	EGFR (L858R)	HS, PQ, CPA, ADP-Glo
DAPK3	HS, PQ	EGFR (d746-750/C797S)	HS, PQ, CPA, ADP-Glo	EGFR (L858R, T970M)	NB
DCLK1	HS	EGFR (d746-750/G724S)	HS	EGFR (L861Q)	HS, PQ, CPA, ADP-Glo
DCLK2	HS, PQ	EGFR (d746-750/S768I)	HS	EGFR (N771_P772insH)	HS
DCLK3	NB	EGFR (d746-750/T790M)	HS, CPA	EGFR (N771_P772insN)	HS
DDR1	HS, NB	EGFR (d746-750/T790M/ C797S)	HS, PQ, CPA, BaF3, ADP-Glo	EGFR (R999A)	HS
DDR2	HS, PQ, NB, ADP-Glo	EGFR (d746-750/T790M/ C797S/L858R)	HS, ADP-Glo	EGFR (S768I)	HS
DDR2 (N456S)	HS, PQ, NB, ADP-Glo	EGFR (d746-750/T790M/L792F)	HS	EGFR (T790M)	HS, PQ, CPA, ADP-Glo
DDR2 (T654M)	HS, PQ, ADP-Glo	EGFR (d746-750/T790M/ L792H)	HS	EGFR (T790M/C797G/L858R)	HS, ADP-Glo
DGKA	ADP-Glo (US)	EGFR (d746-750/T790M/L798I)	HS	EGFR (T790M/C797S)	HS
DGKB	ADP-Glo (US)	EGFR (d746-750/T790M/L858R)	HS	EGFR (T790M/C797S/L858R)	HS, PQ, CPA
DGKD	ADP-Glo (US)	EGFR (d747-749)	HS	EGFR (T790M/L792F/C797S/ L858R)	HS
DGKE	ADP-Glo (US)				
DGKG	ADP-Glo (US)				
DGKH	ADP-Glo (US)				

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
EGFR (T790M/L792F/L858R)	HS	ERN2	HS, NB	FLT3	HS, PQ, NB, CPA, ADP-Glo
EGFR (T790M/L792H/C797S/L858R)	HS	FER	HS, PQ, NB, ADP-Glo	FLT3 (D835H)	HS, NB
EGFR (T790M/L792H/L858R)	HS	FES	HS, PQ, NB	FLT3 (D835V)	HS, NB
EGFR (T790M/L858R)	HS, PQ, CPA, ADP-Glo	FGFR1	HS, PQ, NB, ADP-Glo	FLT3 (D835Y)	HS, PQ, NB, CPA, ADP-Glo
EGFR (V769_D770insGE)	HS	FGFR1 (V561M)	HS, PQ, ADP-Glo	FLT3 (F594_R595insR)	HS
EGFR (V948R)	HS	FGFR1 (W666R)	HS	FLT3 (F594_R595insREY)	HS
EIF2AK1	HS, PQ, ADP-Glo	FGFR2	HS, PQ, NB, CPA, ADP-Glo	FLT3 (F691L)	HS
EIF2AK2	HS, PQ	FGFR2 (C491A)	HS	FLT3 (F691L/D835Y)	HS
EIF2AK3	HS, PQ	FGFR2 (C491A/V564I)	HS	FLT3 (ITD)	HS, PQ, CPA, ADP-Glo
EIF2AK4	HS, PQ, NB, ADP-Glo	FGFR2 (C491A/V564L)	HS	FLT3 (ITD)-NPOS	HS
EPHA1	HS, PQ, NB, ADP-Glo	FGFR2 (C491F)	HS	FLT3 (ITD)-W51	HS
EPHA2	HS, PQ, NB, ADP-Glo	FGFR2 (C491S)	HS	FLT3 (K663Q)	NB
EPHA3	HS, PQ, NB, ADP-Glo	FGFR2 (C491S/V564L)	HS	FLT3 (N841I)	NB
EPHA4	HS, PQ, NB, ADP-Glo	FGFR2 (E565A)	HS	FLT3 (R595_E596insEY)	HS
EPHA5	HS, PQ, NB, ADP-Glo	FGFR2 (E565G)	HS	FLT3 (R834Q)	NB
EPHA6	HS, PQ, NB, ADP-Glo	FGFR2 (K526E)	HS	FLT3 (Y591_V592insVDFREYEYD)	HS
EPHA7	HS, PQ, NB, ADP-Glo	FGFR2 (K641R)	HS	FLT3 (Y591-V592insVDFREYEYD/D835Y)	HS
EPHA8	HS, PQ, NB	FGFR2 (K659N)	HS	FLT3 (Y591-V592insVDFREYEYD/F691L)	HS
EPHB1	HS, PQ, NB, ADP-Glo	FGFR2 (M420I)	HS	FLT4	HS, PQ, CPA, ADP-Glo
EPHB2	HS, PQ, NB, ADP-Glo	FGFR2 (N549H)	HS	FRK	HS, PQ, NB, ADP-Glo
EPHB3	HS, PQ, NB, ADP-Glo	FGFR2 (N549K)	HS	FYN	HS, PQ, NB, CPA, ADP-Glo
EPHB4	HS, PQ, NB, CPA, ADP-Glo	FGFR2 (R612T)	HS	FYN (Y531F)	HS, PQ, NB, ADP-Glo
ERBB2	HS, PQ, CPA, ADP-Glo	FGFR2 (V564F)	HS	FYN-A	HS
ERBB2 (775YVMA776)	PQ, ADP-Glo	FGFR2 (V564I)	HS	GAK	PQ, NB
ERBB2 (A775_G776insYVMA)	HS	FGFR2 (V564L)	HS	GRK1	HS
ERBB2 (D769H)	HS	FGFR3	HS, PQ, NB, ADP-Glo	GRK2	HS, PQ
ERBB2 (D769Y)	HS	FGFR3 (G697C)	HS, PQ, NB, ADP-Glo	GRK3	HS, PQ
ERBB2 (P1170A)	HS	FGFR3 (K650E)	HS, PQ, ADP-Glo	GRK4	HS, PQ
ERBB2 (P780-Y781insGSP)	HS	FGFR3 (K650M)	HS, PQ, ADP-Glo	GRK5	HS, PQ
ERBB2 (R896C)	HS	FGFR3 (K650Q)	HS	GRK6	HS, PQ
ERBB2 (V777_G778insCG)	HS	FGFR3 (V555M)	HS	GRK7	HS, PQ
ERBB2 (V777L)	HS	FGFR4	HS, PQ, NB, ADP-Glo	GSG2	HS, PQ, ADP-Glo
ERBB2 (V956R)	HS	FGFR4 (N535K)	HS, PQ, ADP-Glo	GSK3A	HS, PQ, NB
ERBB4	HS, PQ, CPA, ADP-Glo	FGFR4 (V550E)	HS, PQ, ADP-Glo	GSK3B	HS, PQ, NB, ADP-Glo
ERN1	HS, NB	FGFR4 (V550L)	HS	GSK3b (C199A)	HS
ERN1 (R727A)	HS	FGFR4 (V550M)	HS	HCK	HS, PQ, NB, ADP-Glo
ERN1 (R728A)	HS	FGR	HS, PQ, NB, ADP-Glo		
ERN1/IRE1 (R727A/R728A)	HS	FLT1	HS, PQ, NB, ADP-Glo		

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
HIPK1	HS, PQ	KIF20A	ADP-Glo	KSR2 (R676S)	HS
HIPK2	HS, PQ, NB	KIF22	ADP-Glo	LATS1	HS, PQ, NB
HIPK3	HS, PQ, NB, ADP-Glo	KIF23	ADP-Glo	LATS2	HS, PQ, NB
HIPK4	HS, PQ, NB	KIFC1	ADP-Glo	LCK	HS, PQ, NB, ADP-Glo
HUNK	HS, NB	KIFC3	ADP-Glo	LIMK1	HS, PQ, NB
IGF1R	HS, PQ, NB, CPA, ADP-Glo	KIT	HS, PQ, NB, CPA, ADP-Glo	LIMK2	HS, PQ, NB
IKKBK	HS, PQ, ADP-Glo	KIT (A829P)	HS, PQ, NB, ADP-Glo	LRRK2	HS, PQ, NB, CPA
IKBKE	HS, PQ, NB, ADP-Glo	KIT (d557-558)	HS, BaF3	LRRK2 (G2019S)	HS, PQ, NB
INSR	HS, PQ, NB, ADP-Glo	KIT (d557-558/V654A/A829P)	BaF3	LRRK2 (I2020T)	HS, PQ, NB
INSRR	HS, PQ	KIT (d557-558/V654A/D816A)	BaF3	LRRK2 (R1441C)	HS, PQ, NB
IRAK1	HS, PQ, NB	KIT (d557-558/V654A/D820A)	BaF3	LTK	HS, PQ, NB
IRAK2	HS	KIT (d557-558/V654A/D822K)	BaF3	LYN	HS, PQ, NB, ADP-Glo
IRAK3	NB, PD	KIT (d557-558/Y823D)	HS	LYN (d23-43)	HS
IRAK4	HS, PQ, NB, ADP-Glo	KIT (D816E)	HS, PQ	MAK	HS
IRAK4 (aa104-460) untagged	PQ, ADP-Glo	KIT (D816F)	HS	MAP2K1	HS, PQ, ADP-Glo
ITK	HS, PQ, NB, ADP-Glo	KIT (D816H)	HS, NB, ADP-Glo	MAP2K1 (F53L)	PQ, ADP-Glo
JAK1	PQ, ADP-Glo	KIT (D816I)	HS	MAP2K1 (P124L)	HS, PQ, ADP-Glo
JAK1 (aa850-1154)	PQ, ADP-Glo	KIT (D816V)	HS, PQ, NB, ADP-Glo	MAP2K1 (S218E/S222E)	PQ, ADP-Glo
JAK1 (aa866-1154)	HS	KIT (D816Y)	HS	MAP2K1/KRAS(G12C)	NB
JAK1 (S729C)	PQ, ADP-Glo	KIT (D820E)	HS	MAP2K1/MAP2K2	CPA
JAK1 JH2	PD	KIT (D820Y)	HS	MAP2K2	HS, PQ
JAK2	HS, PQ, NB, ADP-Glo	KIT (K642E)	HS	MAP2K2/KRAS(G12C)	NB
JAK2 (JH1&2)	HS, PQ	KIT (L576P)	NB	MAP2K3	HS, PQ, ADP-Glo
JAK2 (JH1)	NB	KIT (N822K)	HS	MAP2K4	HS, PQ, ADP-Glo
JAK2 (V617F)	HS, NB	KIT (T670I)	HS, PQ, ADP-Glo	MAP2K5	HS, PQ, ADP-Glo
JAK2 JH2	PD	KIT (V559A)	HS	MAP2K6	HS, NB
JAK3	HS, PQ, NB, ADP-Glo	KIT (V559D)	HS, PQ, NB, ADP-Glo	MAP2K6 (S207D/T211D)	PQ, ADP-Glo
JAK3 JH1 & JH2	HS	KIT (V559D/T670I)	HS, PQ, NB, ADP-Glo	MAP2K7	HS, PQ, ADP-Glo
KDR	HS, PQ, CPA, ADP-Glo	KIT (V559D/V654A)	HS, PQ, NB, ADP-Glo	MAP3K1	HS, PQ, ADP-Glo
KIF2C	ADP-Glo	KIT (V560G)	HS, PQ, ADP-Glo	MAP3K2	HS, PQ, NB, ADP-Glo
KIF3C	ADP-Glo	KIT (V560G/D816V)	HS	MAP3K3	HS, PQ, NB, ADP-Glo
KIF4A	ADP-Glo	KIT (V560G/N822K)	HS	MAP3K4	NB
KIF5B	ADP-Glo	KIT (V654A)	HS, PQ, ADP-Glo	MAP3K5	HS, PQ, ADP-Glo
KIF10/CENP-E	ADP-Glo	KIT (Y823D)	HS	MAP3K6	HS, PQ, ADP-Glo
KIF11/Eg5	ADP-Glo	KSR1	HS	MAP3K7	HS, PQ
KIF18A	ADP-Glo	KSR1 (A635F)	HS	MAP3K7-TAB1	PQ, ADP-Glo
KIF18B	ADP-Glo	KSR1 (L639F)	HS	MAP3K8	HS, PQ
KIF19	ADP-Glo	KSR2	HS	MAP3K9	HS, PQ, NB
				MAP3K10	HS, PQ, NB, ADP-Glo

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
MAP3K11	HS, PQ, NB, ADP-Glo	MAST4	NB	MET (Y1230H)	HS, PQ, NB, CPA, ADP-Glo
MAP3K12	HS, NB	MASTL	HS, PQ, ADP-Glo	MET (Y1230S)	HS
MAP3K13	NB	MATK	HS, PQ	MET (Y1235D)	HS, PQ, NB, ADP-Glo
MAP3K14	HS, PQ, ADP-Glo	MELK	HS, PQ, NB, ADP-Glo	MET-KIF5B	HS
MAP3K19	HS, NB	MELK (T460M)	HS, NB	MET-TFG	HS
MAP3K20	HS, PQ, NB	MERTK	HS, PQ, NB, ADP-Glo	MINK1	HS, PQ, ADP-Glo
MAP3K21	HS, PQ, NB, ADP-Glo	MERTK (A708S)	HS, NB	MKNK1	HS, PQ
MAP4K1	HS, PQ, NB, CPA, ADP-Glo	MET	HS, PQ, NB, CPA, ADP-Glo	MKNK2	HS, PQ, NB
MAP4K2	HS, PQ, NB	MET (D1213H)	HS	MOK	NB
MAP4K3	HS, PQ, NB	MET (D1228A)	HS	MPL	BaF3
MAP4K4	HS, PQ, ADP-Glo	MET (D1228G)	HS	MPL+CALR	BaF3
MAP4K5	HS, PQ, NB	MET (D1228H)	HS, PQ, NB, ADP-Glo	MPL-CALR (L367fs*46)	BaF3
MAPK1	HS, PQ, NB, ADP-Glo	MET (D1228N)	HS, PQ, NB, CPA, ADP-Glo	MST1R	HS, PQ, NB, CPA, ADP-Glo
MAPK1/KRAS(G12C)	NB	MET (D1228V)	HS	MTOR	HS, PQ
MAPK3	HS, PQ, NB, ADP-Glo	MET (D1228Y)	HS	MUSK	HS, PQ, NB, ADP-Glo
MAPK3/KRAS(G12C)	NB	MET (D1288H)	CPA	MYLK	HS, PQ
MAPK4	NB	MET (DelEx14)	HS	MYLK2	HS, PQ, NB, ADP-Glo
MAPK6	NB	MET (F1200I)	HS, PQ, NB, CPA, ADP-Glo	MYLK3	HS, PQ, NB
MAPK7	HS	MET (G1163R)	HS, PQ, ADP-Glo	MYLK4	HS, NB
MAPK7 (aa5-397)	PQ, ADP-Glo	MET (H1094L)	HS	MYO3A	HS
MAPK7 (CD)	HS	MET (H1094Y)	HS	MYO3B	HS
MAPK8	HS, PQ, NB, ADP-Glo	MET (K1244R)	HS	NEK1	HS, PQ, NB, ADP-Glo
MAPK9	HS, PQ, NB, ADP-Glo	MET (L1195F)	HS	NEK2	HS, PQ, NB, ADP-Glo
MAPK10	HS, PQ, NB, ADP-Glo	MET (L1195V)	HS, PQ, ADP-Glo	NEK3	HS, PQ, NB, ADP-Glo
MAPK11	HS, PQ, NB, ADP-Glo	MET (M1250I)	HS	NEK4	HS, PQ, NB
MAPK12	HS, PQ, ADP-Glo	MET (M1250T)	HS, PQ, NB, ADP-Glo	NEK5	HS, NB
MAPK13	HS, PQ, ADP-Glo	MET (P991S)	HS, NB	NEK6	HS, PQ, NB, ADP-Glo
MAPK14	HS, PQ, NB, ADP-Glo	MET (R1227K)	HS	NEK7	HS, PQ, ADP-Glo
MAPK14 (T106M)	HS, NB	MET (R970C)	HS	NEK8	HS, PQ, ADP-Glo
MAPK15	HS, PQ, ADP-Glo	MET (T1173I)	HS, NB	NEK9	HS, PQ, NB
MAPKAPK2	HS, PQ, ADP-Glo	MET (T992I)	HS, NB	NEK11	HS, PQ, NB
MAPKAPK3	HS, PQ, ADP-Glo	MET (V1092I)	HS, NB	NIM1K	HS, PQ, NB
MAPKAPK5	HS, PQ, ADP-Glo	MET (Y1230A)	HS, PQ, NB, CPA, ADP-Glo	NLK	HS, PQ, NB, ADP-Glo
MARK1	HS, PQ, ADP-Glo	MET (Y1230C)	HS, PQ, NB, ADP-Glo	NRK	NB
MARK2	HS, PQ, NB	MET (Y1230D)	HS, PQ, NB, CPA, ADP-Glo	NTRK1	HS, PQ, NB, ADP-Glo
MARK3	HS, PQ, NB, ADP-Glo			NTRK1 (A608D)	HS
MARK4	HS, PQ, NB			NTRK1 (F589L)	HS
MAST3	HS, NB				

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
NTRK1 (G595R)	HS	PDGFRB-TPM3	HS	PIK3CA (H1047R/Q859H)-PIK3R1	NB
NTRK1 (G595R/A608D)	HS	PDK1	HS, PQ, ADP-Glo	PIK3CA (H1047R/Q859K)-PIK3R1	NB
NTRK1 (G595R/G667A)	HS	PDK2	HS	PIK3CA (H1047R/R88Q)-PIK3R1	NB
NTRK1 (G595R/G667C)	HS	PDK3	HS	PIK3CA (H1047R/W780R)-PIK3R1	NB
NTRK1 (G595R/G667S)	HS	PDK4	HS	PIK3CA (H1047Y)-PIK3R1	NB
NTRK1 (G595R/L657M)	HS	PDPK1	HS	PIK3CA (I197K)/PIK3R1	NB
NTRK1 (G667A)	HS	PEAK1	HS	PIK3CA (I800L)-PIK3R1	NB
NTRK1 (G667C)	HS, PQ, NB, ADP-Glo	PHKG1	HS, PQ, NB	PIK3CA (M1004I)/PIK3R1	NB
NTRK1 (G667S)	HS	PHKG2	HS, PQ, NB	PIK3CA (M1043I)-PIK3R1	NB
NTRK1 (L657M)	HS	PI3KC2B	ADP-Glo	PIK3CA (Q546K)-PIK3R1	NB
NTRK1-TFG	HS	PI3KC2G/PIK3C2G	ADP-Glo	PIK3CA (Q859H)/PIK3R1	NB
NTRK1-TPM3	HS	PI4K2A	ADP-Glo	PIK3CA (Q859K)/PIK3R1	NB
NTRK1-TPR	HS	PI4K2B	ADP-Glo	PIK3CA (R88Q)/PIK3R1	NB
NTRK2	HS, PQ, NB, ADP-Glo	PI4KA	ADP-Glo	PIK3CA (S576C)/PIK3R1	NB
NTRK3	HS, PQ, ADP-Glo	PI4KB	ADP-Glo	PIK3CA (W780R)/PIK3R1	NB
NTRK3 (G623E)	HS	PIK3C2A	ADP-Glo	PIK3CA-PIK3R1	ADP-Glo, NB
NTRK3 (G623R)	HS	PIK3C2B	ADP-Glo	PIK3CA-PIK3R1/p65 α	ADP-Glo
NTRK3 (G623R/L686M)	HS	PIK3C2G	ADP-Glo	PIK3CB (D1067A)-PIK3R1	ADP-Glo
NTRK3 (G696A)	HS	PIK3C3	ADP-Glo, NB	PIK3CB (D1067V)-PIK3R1	ADP-Glo
NTRK3 (L686M)	HS	PIK3CA (C420R)-PIK3R1	NB	PIK3CB (D1067Y)-PIK3R1	ADP-Glo
NUAK1	HS, PQ, NB, ADP-Glo	PIK3CA (D549N)-PIK3R1	NB	PIK3CB (E1051K)-PIK3R1	ADP-Glo
NUAK2	HS, PQ, NB, ADP-Glo	PIK3CA (D725G)-PIK3R1	NB	PIK3CB (E633K)-PIK3R1	ADP-Glo
OXSR1	HS	PIK3CA (E172Q)-PIK3R1	NB	PIK3CB-PIK3R1	ADP-Glo, NB
PAK1	HS, PQ, ADP-Glo	PIK3CA (E418K)-PIK3R1	NB	PIK3CB-PIK3R2	ADP-Glo
PAK2	HS, PQ, ADP-Glo	PIK3CA (E542K)-PIK3R1	ADP-Glo, NB	PIK3CD-PIK3R1	ADP-Glo, NB
PAK2 (Y443N)	HS	PIK3CA (E545A)-PIK3R1	NB	PIK3CG	ADP-Glo
PAK3	HS, PQ, ADP-Glo	PIK3CA (E545K)-PIK3R1	ADP-Glo, NB	PIK3CG (L1049R)-PIK3R1	ADP-Glo
PAK4	HS, PQ, NB, ADP-Glo	PIK3CA (E545K/D549N)-PIK3R1	NB	PIKFYVE	NB, ADP-Glo
PAK5	HS, PQ, NB, ADP-Glo	PIK3CA (E545K/D725G)-PIK3R1	NB	PIM1	HS, PQ, CPA, ADP-Glo
PAK6	HS, PQ, NB, ADP-Glo	PIK3CA (E545K/E418K)-PIK3R1	NB	PIM2	HS, PQ, CPA, ADP-Glo
PASK	HS, PQ	PIK3CA (E545K/E453K)-PIK3R1	NB	PIM3	HS, PQ, NB, CPA, ADP-Glo
PBK	HS, PQ, ADP-Glo	PIK3CA (E545K/I197K)-PIK3R1	NB	PIP4K2A	ADP-Glo
PDGFRA	HS, PQ, ADP-Glo	PIK3CA (E545K/S576C)-PIK3R1	NB	PIP4K2B	ADP-Glo
PDGFRA (D842V)	HS, PQ, ADP-Glo	PIK3CA (H1047L)-PIK3R1	NB	PIP4K2C	ADP-Glo, NB
PDGFRA (G680R)	HS	PIK3CA (H1047R)-PIK3R1	ADP-Glo, NB	PIP5K1A	ADP-Glo
PDGFRA (T674I)	HS, PQ, ADP-Glo	PIK3CA (H1047R/M1004I)-PIK3R1	NB	PIP5K1B	ADP-Glo, HS, NB
PDGFRA (V561D)	HS, PQ, NB, ADP-Glo	PIK3CA (H1047R/Q859H)-PIK3R1	NB		
PDGFRA-FIP1L1	HS				
PDGFRB	HS, PQ, CPA, ADP-Glo				

Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
PIP5K1C	ADP-Glo	PRKCH	HS, PQ, ADP-Glo	RET (S904F)	HS
PKD2/PRKD2 (G848E)	HS	PRKCI	HS, PQ, ADP-Glo	RET (V738A)	HS
PKMYT1	HS, NB, ADP-Glo	PRKCCQ	HS, PQ, NB, ADP-Glo	RET (V778I)	HS
PKN1	HS, PQ, ADP-Glo	PRKCCZ	HS, PQ, ADP-Glo	RET (V804E)	HS, PQ, ADP-Glo
PKN1-ANXA4 (Aex2Pex13)	HS	PRKCCZ (aa184-592)	HS, PQ, ADP-Glo	RET (V804L)	HS, PQ, NB, ADP-Glo
PKN1-TECR	HS	PRKD1	HS, PQ, ADP-Glo	RET (V804L)-KIF5B	HS
PKN2	HS, PQ	PRKD2	HS, PQ	RET (V804M)	HS, PQ, NB, ADP-Glo
PKN3	HS, PQ, NB, ADP-Glo	PRKD2 (G870E)	HS	RET (V804M)-KIF5B	HS
PLK1	HS, PQ, NB, ADP-Glo	PRKD3	HS, PQ, ADP-Glo	RET (V804M/G810S)	HS
PLK2	HS, PQ, NB, ADP-Glo	PRKDC	HS, PQ	RET (Y791F)	HS, PQ, ADP-Glo
PLK3	HS, PQ, NB, ADP-Glo	PRKG1 (A)	HS, PQ	RET (Y806C)	HS
PLK4	HS, PQ, NB, ADP-Glo	PRKG1 (B)	HS	RET (Y806H)	HS, PQ, ADP-Glo
PNCK	HS	PRKG2	HS, PQ, NB	RET (Y806N)	HS
PRKAA1	PQ, NB, ADP-Glo	PRKX	HS, PQ, NB	RET-BCR	HS
PRKAA1 (aa1-312)	PQ, ADP-Glo	PTK2	HS, PQ, NB, CPA, ADP-Glo	RET-CCDC6	HS, PQ, ADP-Glo
PRKAA1-PRKAB1-PRKAG1	HS	PTK2 (aa411-686)	PQ, ADP-Glo	RET-KIF5B	HS
PRKAA1-PRKAB1-PRKAG2	HS	PTK2B	HS, PQ, NB, ADP-Glo	RET-NCOA4	HS
PRKAA1-PRKAB1-PRKAG3	HS	PTK6	HS, PQ, NB, ADP-Glo	RET-PRKARA1A	HS
PRKAA1-PRKAB2-PRKAG1	HS	RAF1 (R391W)	HS	RIOK2	NB
PRKAA1-PRKAB2-PRKAG2	HS	RAF1 (Y340D/Y341D)	HS, PQ, ADP-Glo	RIPK1	NB, ADP-Glo
PRKAA1-PRKAB2-PRKAG3	HS	RET	HS, PQ, NB, ADP-Glo	RIPK2	HS, PQ, NB, ADP-Glo
PRKAA2	NB	RET (A883F)	HS	RIPK3	HS
PRKAA2-PRKAB1-PRKAG1	HS	RET (E732K)	HS	RIPK4	HS, PQ, ADP-Glo
PRKAA2-PRKAB1-PRKAG2	HS	RET (E762Q)	HS, PQ, ADP-Glo	ROCK1	HS, PQ, NB, CPA
PRKAA2-PRKAB1-PRKAG3	HS	RET (G691S)	HS, PQ, ADP-Glo	ROCK2	HS, PQ, NB, CPA, ADP-Glo
PRKAA2-PRKAB2-PRKAG1	HS	RET (G810C)	HS, PQ, ADP-Glo	ROS1	HS, PQ, NB
PRKAA2-PRKAB2-PRKAG2	HS	RET (G810R)	HS, PQ, ADP-Glo	ROS1 (F2004C)	HS
PRKAA2-PRKAB2-PRKAG3	HS	RET (G810S)	HS, PQ, ADP-Glo	ROS1 (F2004I)	HS
PRKACA	HS, PQ, NB	RET (G810S/M918T)	HS	ROS1 (G2032R)	HS
PRKACA (L206R)	HS	RET (L730I)	HS, PQ, ADP-Glo	ROS1 (G2101A)	HS
PRKACA-DNAJB1	HS	RET (L730M)	HS, PQ, ADP-Glo	ROS1 (G2101C)	HS
PRKACB	HS, NB	RET (L790F)	HS	ROS1 (L2086F)	HS
PRKACG	HS	RET (M918T)	HS, PQ, NB, ADP-Glo	ROS1-GOPC	HS
PRKCA	HS, PQ, NB, ADP-Glo	RET (R749T)	HS, PQ, ADP-Glo	ROS1-TPM3	HS
PRKCB (1)	HS, PQ, NB, ADP-Glo	RET (R813Q)	HS, PQ, ADP-Glo	RPS6KA1	HS, PQ, NB
PRKCB (2)	HS, PQ, ADP-Glo	RET (R912P)	HS	RPS6KA2	HS, PQ, NB
PRKCD	HS, PQ, NB, ADP-Glo	RET (S891A)	HS, PQ, ADP-Glo	RPS6KA3	HS, PQ, NB
PRKCE	HS, PQ, NB, ADP-Glo	RET (S904A)	HS	RPS6KA3 (I416V)	HS, NB
PRKCG	HS, PQ, NB, ADP-Glo				

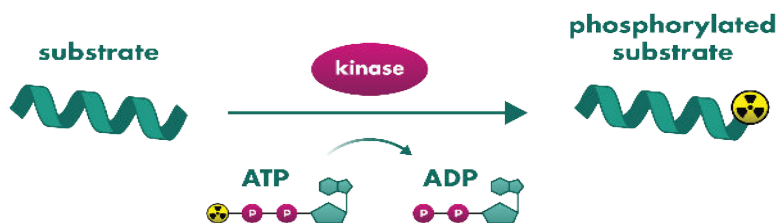
Kinase	Assay Format	Kinase	Assay Format	Kinase	Assay Format
RPS6KA3 (L608F)	HS, NB	STK32C	HS	TSSK4	HS
RPS6KA4	HS, PQ, NB	STK33	HS, PQ, NB	TSSK6	HS
RPS6KA5	HS, PQ	STK35	HS, NB	TTBK1	HS, PQ, ADP-Glo
RPS6KA6	HS, PQ, NB	STK36	NB	TTBK1 (aa1-480)	PQ, ADP-Glo
RPS6KB1	HS, PQ, ADP-Glo	STK38	HS, PQ, NB, ADP-Glo	TTBK2	HS, PQ
RPS6KB2	HS, PQ	STK38L	HS, PQ, NB, ADP-Glo	TTK	HS, PQ, NB, ADP-Glo
SBK1	HS, PQ	STK39	HS, PQ	TXK	HS, PQ, NB
SBK3	NB	SYK	HS, PQ, ADP-Glo	TYK2	HS, PQ, NB
SGK1	HS, PQ, NB, ADP-Glo	SYK (aa356-635)	PQ, ADP-Glo	TYK2 (JH1&2)	HS, PQ
SGK2	HS, PQ, NB	TAOK1	HS	TYK2 (JH1)	NB
SGK3	HS, PQ, ADP-Glo	TAOK2	HS, PQ	TYK2 (JH2)	NB, PD
SIK1	HS, PQ, NB, ADP-Glo	TAOK3	HS, PQ, ADP-Glo	TYRO3	HS, PQ, NB, ADP-Glo
SIK2	HS, PQ, NB, ADP-Glo	TBK1	HS, PQ, NB, ADP-Glo	ULK1	HS, PQ, NB, ADP-Glo
SIK3	HS, PQ, NB, ADP-Glo	TEC	HS, PQ, NB	ULK2	HS, PQ, NB, ADP-Glo
SLK	HS, PQ, NB	TEK	HS, PQ, NB, ADP-Glo	ULK3	HS, PQ, NB, ADP-Glo
SNRK	HS, NB	TEK (A1124V)	HS, NB	VRK1	HS, PQ, ADP-Glo
SPHK1	ADP-Glo	TEK (P883A)	HS, NB	VRK2	HS, PQ, ADP-Glo
SPHK2	ADP-Glo	TEK (R849W)	HS, PQ, NB, ADP-Glo	WEE1	HS, PQ, NB, ADP-Glo, PD
SRC	HS, PQ, NB, CPA, ADP-Glo	TEK (R915C)	HS	WEE2	HS, NB
SRC (T341M)	HS	TEK (Y1108F)	HS, PQ, NB, ADP-Glo	WNK1	HS, PQ, ADP-Glo
SRC (Y530F)	HS	TEK (Y897C)	HS, NB	WNK2	HS, PQ
SRC N1	HS	TEK (Y897H)	HS	WNK3	HS, PQ
SRMS	HS, PQ, NB	TEK (Y897H/R915C)	HS	YES1	HS, PQ, NB, ADP-Glo
SRPK1	HS, PQ, ADP-Glo	TEK (Y897S)	HS, PQ, NB, ADP-Glo	YES1 (T348I)	HS
SRPK2	HS, PQ, ADP-Glo	TESK1	HS, NB	ZAP70	HS, PQ, ADP-Glo
SRPK3	HS, PQ	TESK2	HS	ZAP70 (Y319F)	HS
STK3	HS, PQ, NB, ADP-Glo	TGFBR1	HS, PQ, ADP-Glo		
STK4	HS, PQ, NB, ADP-Glo	TGFBR2	HS, PQ, NB		
STK10	HS, NB	TIE1	NB		
STK11	HS, PQ, NB	TLK1	HS, PQ, NB, ADP-Glo		
STK16	HS, PQ, NB, ADP-Glo	TLK2	HS, PQ, NB, ADP-Glo		
STK17A	HS, PQ	TNIK	HS, PQ, ADP-Glo		
STK17B	PQ, NB, ADP-Glo	TNK1	HS, PQ, NB, ADP-Glo		
STK24	HS, PQ, NB	TNK2	HS, PQ, NB, ADP-Glo		
STK25	HS, PQ	TNNI3K	NB		
STK26	HS, PQ, NB, ADP-Glo	TRPM7	HS		
STK32A	HS, PQ, NB	TSSK1B	HS, PQ, NB		
STK32B	HS, PQ, NB	TSSK2	HS, PQ, ADP-Glo		
		TSSK3	HS, ADP-Glo		

HS : Radiometric HotSpot Kinase Activity Assay
HTRF : Homogeneous Time Resolved Fluorescence
NB : NanoBRET Intracellular Kinase Assay
PD : Probe-Displacement Assay
PQ : Radiometric ³³PanQinase Kinase Activity Assay
CPA : Cell Phosphorylation Assay
BaF3 : BaF3 Cell Proliferation Assay

Kinase Screening – Free Choice

With over 860+ kinases, Reaction Biology offers the largest selection of kinases available for screening and profiling services. It is the most used kinase screening service in the industry according to HTStec Kinase Profiling Trends Survey.

- Get the highest quality data possible with the gold standard radiometric assay format
- Any class of inhibitor can be tested including ATP competitive and non-competitive as well as allosteric inhibitors
- High-throughput compatible
- Customized assay development possible
- Deliverables: single concentration % inhibition; IC_{50} and/or K_i values
- A reference compound is included in every study for no additional cost



Assay formats

Compound screening on protein kinases is performed with highly sensitive radiometric assays. Phosphate from ^{33}P -labelled ATP is transferred onto a substrate and directly measured avoiding false positives and negatives common with other assay formats.

We offer two radiometric assay formats which differ only in the way of substrate retention via a filter membrane (HotSpot™ assay, used in the US facility) or on a ScintiPlate surface (^{33}P anQinase™ assay, used in the German facility).

Lipid kinases are screened with ADP-Glo Platform from Promega.

Kinase Panel Screening

Screen against the largest selection of kinases and the most widely used panel in industry. Our kinase panels are run once or twice per month allowing us to offer screening with a turnaround time of only two weeks.

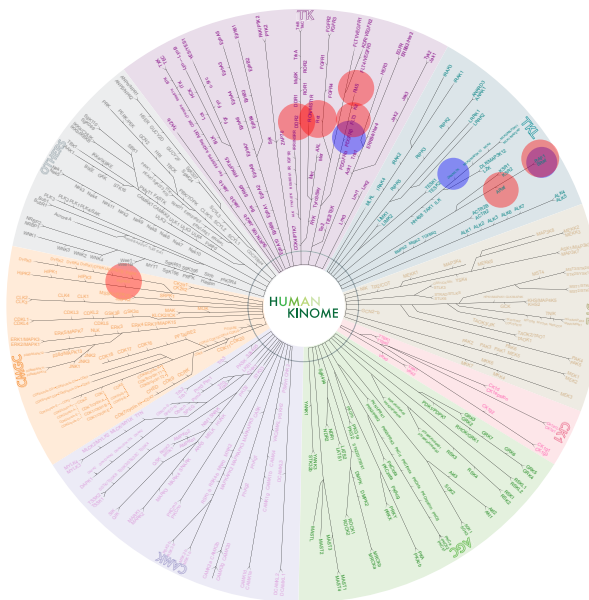
- Highest reproducibility
- Visualize your results with the kinase mapper
- Panels run with HotSpot™ assay include a free control compound's IC₅₀ for every assay
- Deliverable: % of inhibition (single point) or IC₅₀ value determination

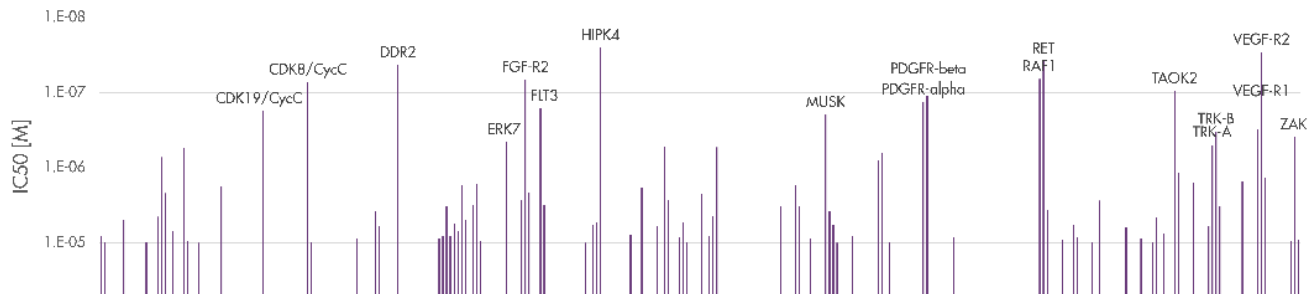
Panels run at US facility	# of kinases
Wild Type Kinase Panel	379
Mutant Kinase Panel	354
Tyrosine Kinase Panel	79
Diversify Panel	70
Lipid Kinase Panel	26
Atypical Kinase Panel	24
Diacylglycerol Kinase Panel	10

Panels run at German facility	# of kinases
Wild Type Kinase Panel	368
Mutant Kinase Panel	96
Diversify Panel	70
Lipid Kinase Panel	18

Kinase Mapper

A kinase mapper tool can be used by customers for graphic presentation of kinase screening results. Shown is an example with Sorafenib profiled with the Wild Type Kinase Panel. Kinases that were inhibited more than 90% are highlighted in red circles, those inhibited more than 75% are shown in blue.

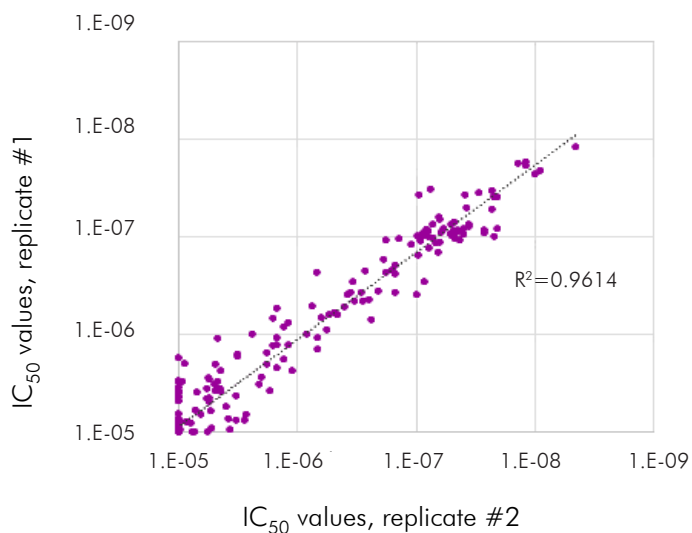




Example of IC₅₀ value determination for sorafenib with the Wild Type Kinase Panel by using the ³³PanQinase™ assay format

Sorafenib activity was determined with 6 concentrations on 320 wild type protein kinases for IC₅₀ value determination.

Using the IC₅₀ value determination setup yields a true value of the inhibition of the compound for every individual kinase. False positives or false negatives that may occur when testing with a suboptimal concentration will be avoided.



High Reproducibility

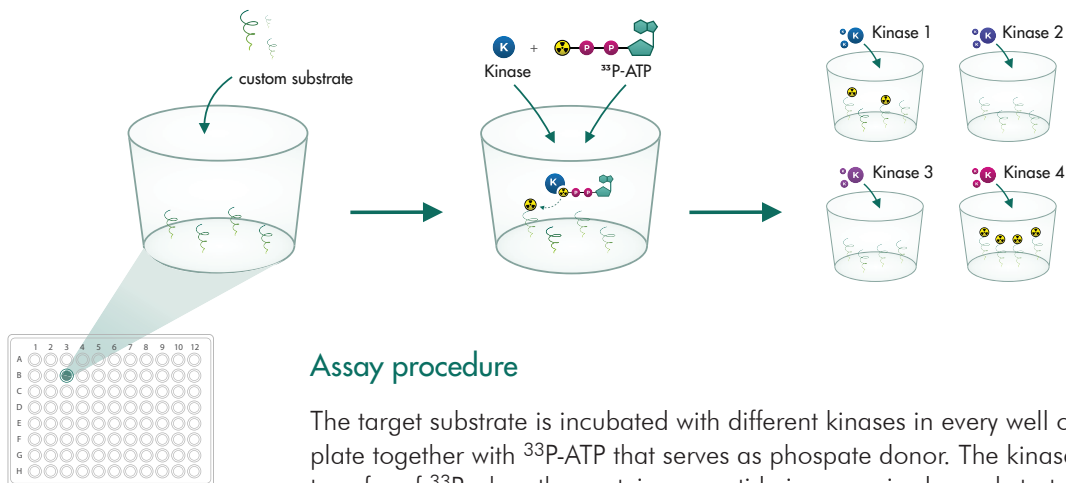
Selectivity profiles from two independent experiments with ponatinib on 320 wild type protein kinases.

KinaseFinder

The KinaseFinder can identify kinases that phosphorylate a substrate of interest. This service is ideal for the characterization of physiological pathways.

- Potential substrates can be peptides or proteins
- Option to compare peptides with phospho-site mutations
- Option to follow up with an SDS-PAGE and autoradiogram to visualize the phosphorylated substrate
- Deliverable: Absolute activity measurement of each kinase on your substrate

Type of KinaseFinder	# of kinases	Assay format
Tyr kinases	94	³³ PanQinase™
Ser/Thr kinases	245	³³ PanQinase™
Ser/Thr & Tyr kinases	339	³³ PanQinase™
Wild Type Kinase Panel	380	HotSpot™
custom panel	selected by customer	HotSpot™



Assay procedure

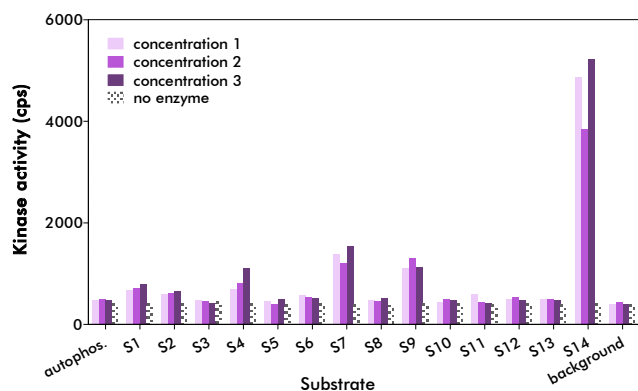
The target substrate is incubated with different kinases in every well of a multi-well plate together with ³³P-ATP that serves as phosphate donor. The kinase will catalyse the transfer of ³³P when the protein or peptide is recognized as substrate. Quantification of phosphorylated substrate is performed via scintillation counting

Kinase SubstrateFinder

The Kinase SubstrateFinder can identify suitable substrates for a specific kinase of interest. The generic substrate panels comprise various proteins, whereas the physiological substrate panels include biotinylated peptide libraries.

- Testing of generic substrates with ATP consumption assay ADP-Glo (Promega)
- Testing of physiologic substrates with radiometric assay using ^{33}P -ATP
- Deliverable: absolute activity measurement of kinase with each substrate

Type of Kinase SubstrateFinder	# of substrates	Assay format
Tyr Generic Substrate Panel	19	ADP-Glo assay
Ser/Thr Generic Substrate Panel	39	ADP-Glo assay
Ser/Thr & Tyr Generic Substrate Panel	58	ADP-Glo assay
Tyr Physiologic Substrate Panel	145	^{33}P PanQinase™
Ser/Thr & Tyr Physiologic Substrate Panel	720	^{33}P PanQinase™



Example of MELK assay development

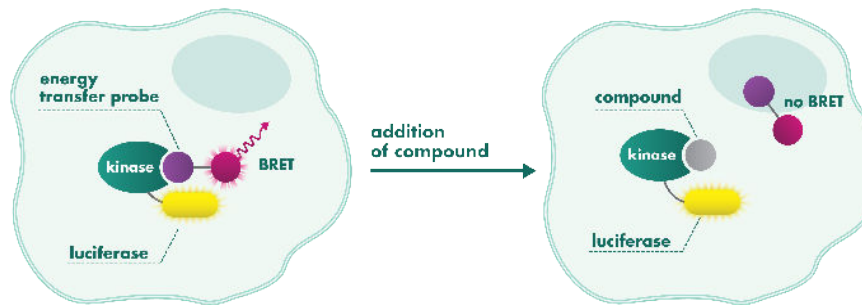
MELK kinase activity was measured with a variety of substrates. Controls are no-enzyme controls of each substrate and autophosphorylation of the MELK kinase without substrate.

Substrate 14 is to be most suitable for establishment of a MELK assay.

NanoBRET Intracellular Target Engagement Kinase Assay

Reaction Biology offers target engagement assays using Promega's NanoBRET technology that enables the quantitative determination of kinase inhibitor occupancy in live cells, without disruption of cellular membrane integrity.

- Intact cells with physiological ATP concentration, protein complex, co-factors and pH values.
- High-throughput compatible
- Deliverable: apparent binding affinity of inhibitor (IC_{50})

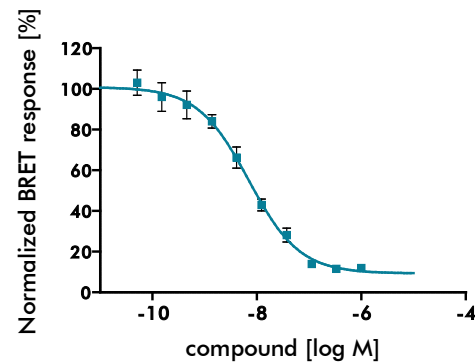


Assay principle

The assay is a compound competition assay that relies on bioluminescence resonance energy transfer (BRET) between a luciferase-tagged kinase and a fluorescent tracer. Quantitation and specificity are key attributes of the NanoBRET system.

DDR1 inhibition by Dasatinib

HEK293 cells transiently expressing NanoLuc®- DDR1 fusion vector were treated with the Tracer K-4 and reference compound Dasatinib for 1 hour. The BRET signal was measured on an EnVision 2104 multilabel microplate reader.



NanoBRET TE Intracellular CDK Panel Screening

Reaction Biology offers the NanoBRET Intracellular CDK Panel Screening service, which includes 20 clinically relevant and recognized CDK targets. This is an excellent platform for assessing test compound's binding affinity and selectivity across a diverse panel of CDK kinases in the physiological environment of intact cells, and it provides qualitative data on the test compound's permeability, affinity, selectivity, and residence time.

- 10-dose IC50 duplicate format
- Performed on a regular basis and economical
- Fast turnaround time: get your results in around 4 weeks
- Semi-automated processing for highly reproducible data

20 CDK Targets:

Target	Synonyms
CDK1+Cyclin B1	CDK1-CCNB1, CDC2, p34
CDK2+Cyclin E1	CDK2-CCNE1, p33
CDK3+Cyclin E1	CDK3-CCNE1, CDKN3
CDK4+Cyclin D1	CDK4-CCND1, CMM3, PSK-J3
CDK5+CDK5R1	CDK5-CDK5R1, PSSALRE
CDK6+Cyclin D1	CDK6-CCND1, PLSTIRE
CDK7+Cyclin H	CDK7-CCNH, CAK1, CDKN7, STK1, p39MO15
CDK8+Cyclin C	CDK8-CCNC, K35
CDK9+Cyclin K	CDK9-CCNK, C-2k, CDC2L4, PITALRE, TAK
CDK10+Cyclin L2	CDK10-CCNL2, PISSLRE

Target	Synonyms
CDK11A+Cyclin K	CDK11A-CCNK, PITSLRE
CDK12+Cyclin K	CDK12-CCNK, CRKRS, CRK7
CDK13+Cyclin K	CDK13-CCNK, CDC2L, CDC2L5, CHED
CDK14+Cyclin Y	CDK14-CCNY, PFTK1, PFTAIRE1
CDK15+Cyclin Y	CDK15-CCNY, ALS2CR7, PFTK2, PFTAIRE2
CDK16+Cyclin Y	CDK16-CCNY, PCTAIRE, PCTGAIRE, PCTK1, PCTAIRE1
CDK17+Cyclin Y	CDK17-CCNY, PCTK2, PCTAIRE2
CDK18+Cyclin Y	CDK18-CCNY, PCTK3, PCTAIRE, PCTAIRE3
CDK19+Cyclin C	CDK19-CCNC, CDC2L6, CDK11
CDK20+Cyclin H	CDK20-CCNH, CAK-kinase p42, CCRK, CDCH, p42, PNQALRE

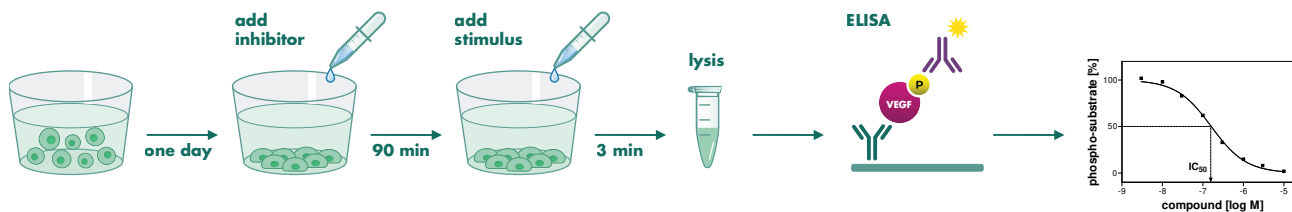
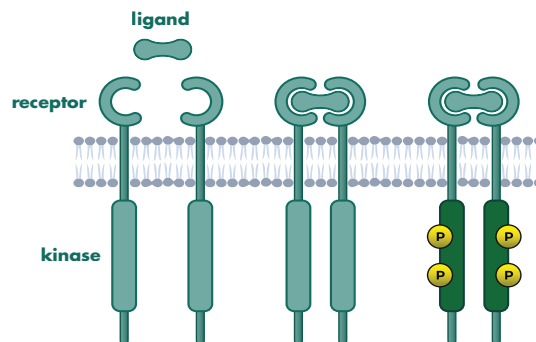
Cellular Phosphorylation Assay

The Cellular Phosphorylation Assay quantifies changes in the phosphorylation state of a substrate as a result of treatment with your inhibitor in intact cells. The assays have been designed to address compound activity in a physiological environment on a physiological substrate.

- Physiological kinase, substrate and ATP concentrations
- Assay can be performed with blood containing drug for plasma-inhibitory study
- Deliverable: % inhibition of kinase activity and IC_{50} determination

Assays based on endogenous kinases

Assays are performed with cells expressing the kinase of interest which is either overexpressed or constitutively active, or kinase activity is triggered by ligand administration.

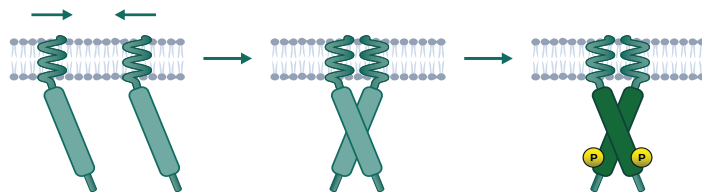


Example of VEGF-R2 signaling

Human endothelial cells are known to express VEGF-R2. The cells incubate with the test compound for 90 minutes to allow for target binding. After a 3-minute stimulation with ligand VEGF-A, cells are lysed and the substrate phosphorylation is quantified by ELISA with pan-phospho-tyrosine antibodies on captured VEGF-R2. The assays are performed with 8 compound concentrations in duplicate for IC_{50} value determination.

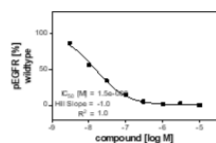
Assays based on exogenous kinases

Rat 1 fibroblasts were transfected to stably express the intracellular domain of EGF-R mutants fused to an artificial transmembrane domain. Dimerization of the receptors causes constitutive auto-phosphorylation that can be quantified via ELISA.

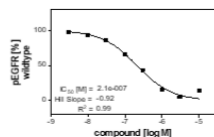


EGF-R Wild Type

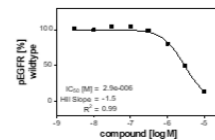
lapatinib



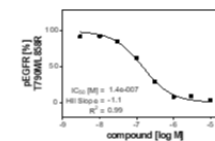
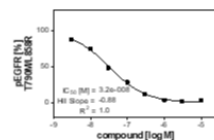
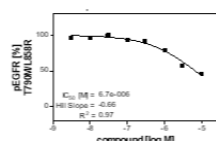
osimertinib



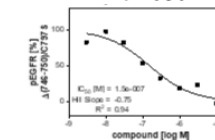
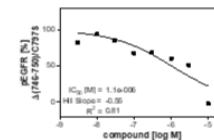
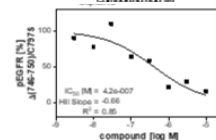
brigatinib



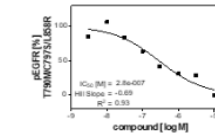
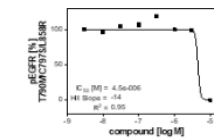
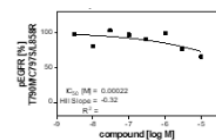
EGF-R T790M/L858R



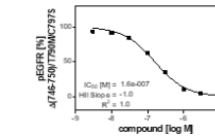
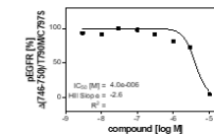
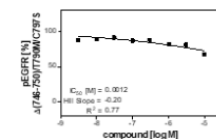
EGF-R Δ746750/C797S



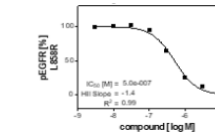
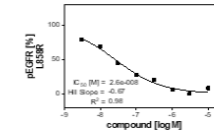
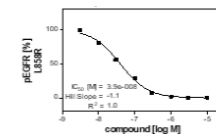
EGF-R T790M/C797S/L858R



EGF-R Δ746-750/T790M/C797S



EGF-R L858R



Example of EGF-R mutant analysis

Rat1 fibroblasts express the intracellular domain of EGF-R containing disease-relevant mutations and a transmembrane domain. The cells were incubated with three EGF-R-specific inhibitors and their potency was quantified via ELISA.

BaF3 Cell Proliferation Assay

The BaF3 Cell Proliferation Assay is performed with BaF3 cells, a pro-B-cell line that is dependent on interleukin 3 for its survival and proliferation. Transgenic overexpression of oncogenic kinases, in particular receptor tyrosine kinases, can transform the cell line to become independent of interleukin 3. This tool cell line is suitable to investigate the potency of kinase oncogenes and the downstream effects of kinase inhibition.

- Kinase inhibitor screening performed in the physiological environment of intact cells
- Determine the effects of compound treatment on the signaling activities of the target kinase
- Readout: Impact of kinase inhibition on cell proliferation

Available Assays

Target	Synonyms
EGFR wild type	ERBB, mENA, ERBB1
EGFR (d746-750/T790M/C797S)	ERBB, mENA, ERBB1
cKIT (d557-558/V654A/D816A)	KIT (d557-558/V654A/D816A)
cKIT (d557-558/V654A/D820A)	KIT (d557-558/V654A/D820A)
cKIT (d557-558/V654A/D822K)	KIT (d557-558/V654A/D822K)
cKIT (d557-558/V654A/A829P)	KIT (d557-558/V654A/A829P)
cKIT (d557-558)	KIT (d557-558)
MPL	Thrombopoietin receptor
MPL+CALR	Calreticulin
MPL-CALR (L367fs*46)	MPL-CALR (L367fs*46)
EpoR	Erythropoietin receptor

Assay Principle

The BaF3 cell line proliferates in the presence of interleukin 3.

The overexpression of receptor tyrosine kinases enables the BaF3 cell line to grow without the supplement of interleukin 3. The cell growth is driven by the signaling of the kinase.

The inhibition of the activity of the transforming kinase leads to the loss of growth stimuli resulting in cell apoptosis.

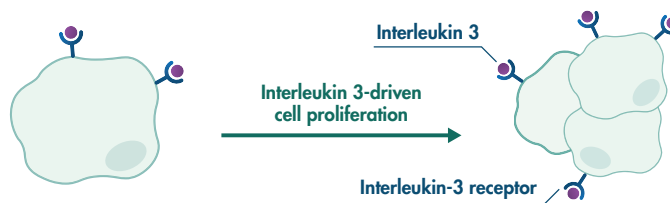
Assay Development

The BaF3 Cell Proliferation Assay can be performed with constructs of a large variety of receptor tyrosine kinases that can act as oncogenes driving cell survival, growth, and proliferation.

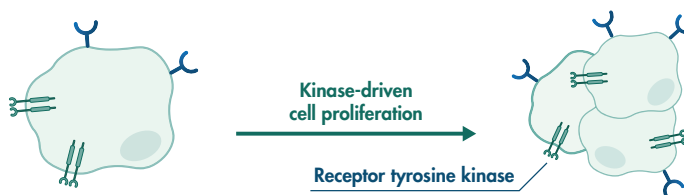
Study Example

BaF3 cells stably expressing EGFR wild type and EGFR (d746-750/T790M/C797S), respectively, were treated with kinase inhibitor Afatinib for 72 hours before quantification of live cells via Cell Titer Glo. The graph depicts the percentage of viable cells in relative to vehicle control (100 %) and staurosporine treatment (0 %).

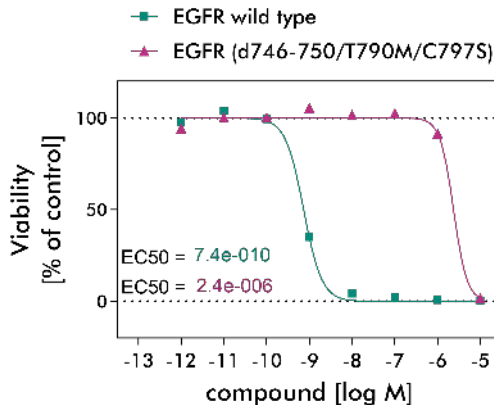
A. Untransfected BaF3 cells



B. BaF3 cells transfected with kinase oncogene



C. Kinase oncogene inhibition



In Vivo Kinase Tumor Models

Genetically engineered tumor models are well suited for investigation of a single driver of tumor growth such as an overexpressed or constitutively expressed kinase.

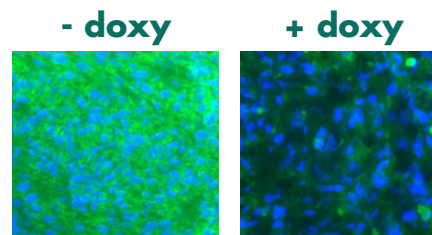
Reaction Biology offers two models based on fibroblast cells which were engineered to express an exogenous receptor kinase under the control of an inducible promotor. These models make excellent tools for the investigation of inhibitors in the in vivo setting.

- Target a human kinase in mice with intact immune system
- Implantation of engineered cells for comparable tumor growth
- Assess compound efficacy and evaluate mechanisms of drug resistance

Kinase	Cell line
human IGF receptor	MEF (mouse)
human ErbB2 receptor	NIH3T3 (mouse)

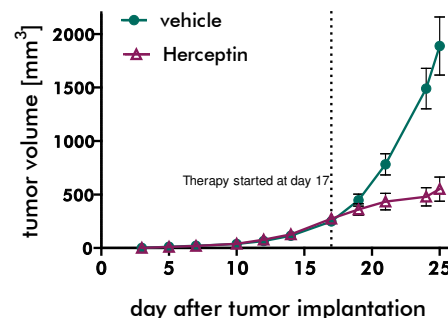
Example: ErbB2 Model

NIH3T3 cells were stably transfected to express human ErbB2 under the control of a Tet-inducible promoter. ErbB2 is expressed in the absence of doxycycline and expression is inhibited in the presence of doxycycline.



Example: ErbB2 Inhibition with Herceptin

NIH3T3-ErbB2-Rrep cells were implanted subcutaneously into mice. At an average tumor size of 400 mm³, mice were treated with the anti-ErbB2 antibody Herceptin resulting in tumor regression.



Customized Kinase Drug Discovery

High-Throughput Screening

Bring your own compound library or use one of our libraries for high-throughput screening with our well validated kinase assays. Contact us to talk about the best approach for a successful screening project.

Custom Assay Development

Our experience of establishing more than 730 kinase assays is the basis for the successful development of the custom-tailored assay for your drug discovery project. We will be happy to provide guidance in construct selection, protein production, substrate requirements and assay condition optimization.

ATP and Substrate Competition Assay

To determine whether a compound's mechanism of action is ATP competitive or substrate competitive, we determine the IC_{50} or K_i values at various ATP and substrate concentrations.

Mechanism of Action Analysis

Using a variety of biochemical and biophysical methods, we can determine the kinetic behavior of your compound including binding affinity, residence time, on- and off-rates, that are crucial to your compound's therapeutic efficacy.

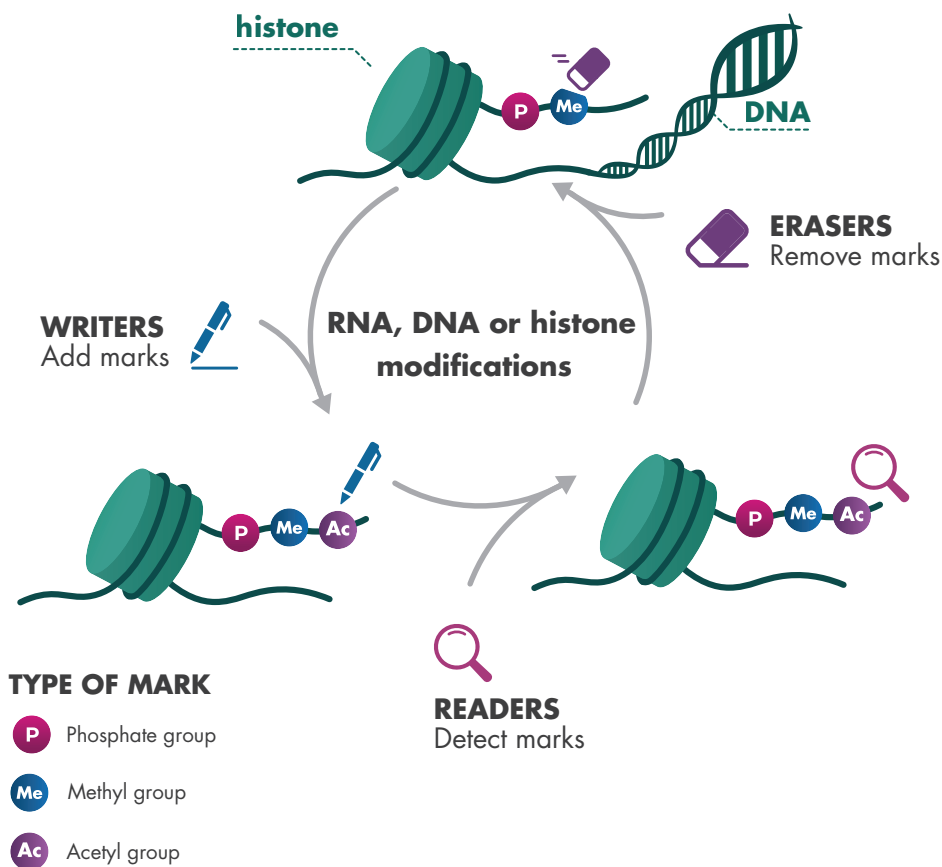
Kinetics, Binding affinity

SPR is commonly used to determine the kinetics of target-analyte binding kinetics. The assay detects changes in the molecular mass of a target after binding of the analyte. The target is immobilized to a sensor chip and the analyte flows to the target. Target binding is monitored in real-time for both: association and dissociation.

Kinase Activation Assay

The Kinase Activation Assay is suitable for the discovery of allosteric compounds that inhibit the activation of a target kinase by an upstream kinase in a so called cascade assay.

EPIGENETIC ASSAYS



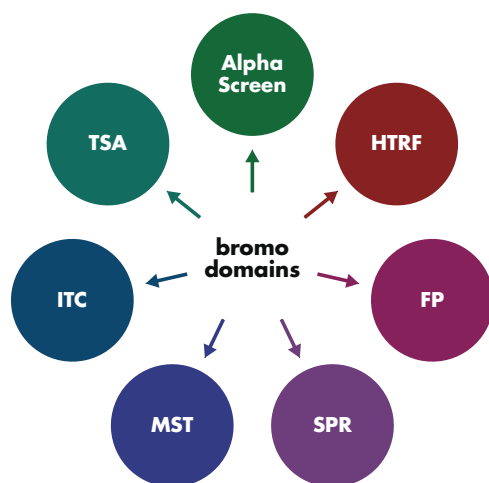
Reaction Biology offers extensive epigenetic drug discovery services including protein production, assay development, high-throughput screening, SAR support, mechanism of action analyses and cell-based assays.

The target families include proteins that regulate post-translational processes such as methylation, acetylation and phosphorylation.

Reader Domain Assays

Reaction Biology offers both biochemical and biophysical assays to study epigenetic reader domains. More than 100 assays have been established for screening, lead optimization or selectivity profiling for reader domain inhibitors.

- All reader domain proteins are produced at our facility and are available for purchase.
- Extensive coverage of the bromodomain family.
- Visualize your bromodomain profiling results with the mapper tool.



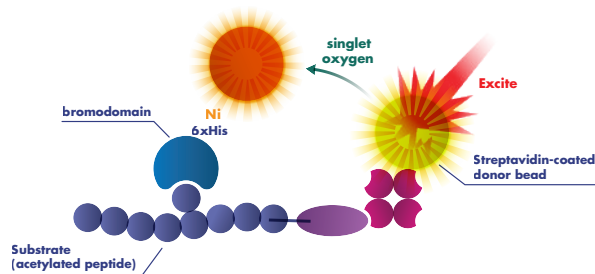
Assay formats available at Reaction Biology for bromodomain targets

Biochemical assay formats to quantify compound binding:

- AlphaScreen
- HTRF
- Fluorescence polarization (FP)

Biophysical assay formats for determination of binding affinity, on- and off-rates and parameters of agent-target interaction on the molecular level

- Surface plasmon resonance (SPR)
- Microscale thermophoresis (MST)
- Isothermal titration calorimetry (ITC)
- Thermal shift assay (TSA)



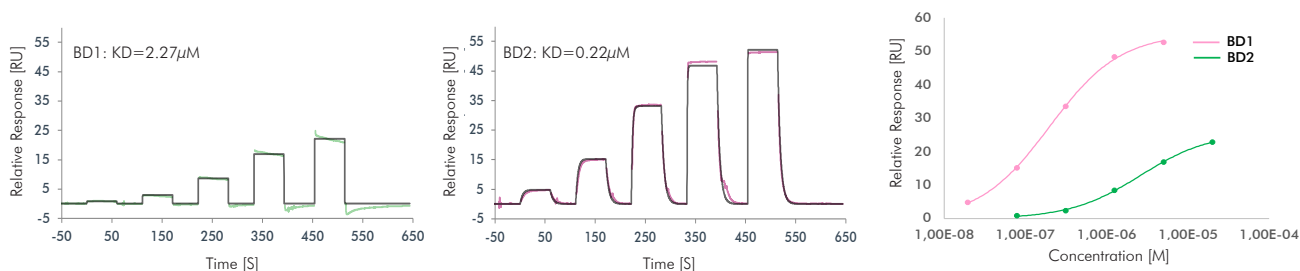
AlphaScreen assay

For bromodomain screening, an acetylated peptide substrate and a bromodomain protein are captured on AlphaScreen beads. A binding interaction between the substrate and protein brings the beads into close proximity. Laser excitation of the complex results in a sequence of chemical reactions resulting in an enhanced fluorescent emission. The presence of an inhibitor interferes with substrate/protein binding resulting in a reduced fluorescent signal. Reaction Biology offers AlphaScreen assays for many reader domains for screening and profiling.

Surface plasmon resonance (SPR) assay

SPR measures biomolecular interactions in real time for screening of targets that are enzymes as well as non-enzymatic targets such as bromodomains. The target proteins are immobilized on the surface of a sensor chip. The compound flows over the sensor chip and binds to the target increasing the molecular mass of the protein upon binding which is measured with an optical readout.

Using SPR we can describe several parameters of the inhibitor-bromodomain interaction: 1. Which of my compounds bind? 2. Is the compound specific to my target? 3. How strong is the binding? 4. What are the association and dissociation rates? 5. Where does it bind?



Example of BRD4 domain interaction with RVX-208 as detected by SPR

Human bromodomain BRD4 contains tandem bromodomains (BD1 and BD2) that have unique biological functions. Inhibitors that selectively bind to one of the domains may affect different biological outcomes. By using recombinant BRD4 with individual bromodomains, SPR revealed that RVX-208 is about 10-times more selective for BD2 over BD1.

Reader Domain	Assays
ASH1L-[BRD]-GST	TS
ATAD2-His	TS, AS
ATAD2B-GST	TS
ATAD2B-His	TS, AS
BAZ1A-GST	TS
BAZ1B-His	TS, AS
BAZ2A-GST	TS
BAZ2A-His	AS
BAZ2B-His	TS, AS
BPTF-[BRD]-His	TS, AS
BPTF-[PHD-BRD]-His	TS, AS
BRD1-GST	TS
BRD1-His	TS, AS
BRD2-1-GST	TS
BRD2-1-His	TS, AS
BRD2-2-GST	TS
BRD2-2-His	TS, AS
BRD2-Tndm-His	TS, AS
BRD3-1-GST	TS
BRD3-1-His	TS, AS
BRD3-2-GST	TS
BRD3-2-His	TS, AS
BRD3-Tndm-GST	TS
BRD3-Tndm-His	TS, AS
BRD4 Full length	TS, AS
BRD4-1	SPR
BRD4-1-GST	TS
BRD4-1-His	TS, AS
BRD4-2	SPR
BRD4-2-GST	TS
BRD4-2-His	TS, AS
BRD4-Tndm-GST	TS
BRD4-Tndm-His	TS, AS
BRD7-GST	TS

Reader Domain	Assays
BRD9-GST	TS
BRD9-His	TS, AS
BRDT-1-His	TS, AS
BRDT-2-His	TS
BRDT-Tndm-His	TS, AS
BRPF1 α	TS
BRPF1b-GST	TS
BRPF1b-His	TS, AS
BRPF3-GST	TS
BRPF3-His	TS, AS
BRWD1-2-GST	TS
BRWD1-2-His	TS, AS
BRWD3-2-GST	TS
CBX7-[CHR]-GST	TS
CDg1-[CHR]-GST	TS
CECR2-GST	TS
CECR2-His	TS, AS
CHD1-[CHR]-GST	TS
CHD1-[CHR]-His	TS
CHD2-[CHR]-GST	TS
CHD2-[CHR]-His	TS
CHD4-[CHR]-GST	TS
CHD4-[PHD-CHR]-GST	TS
CHD7-[CHR]-GST	TS
CREBBP-GST	TS
CREBBP-His	TS, AS
EED	HTRF
EP300-GST	TS
EP300-His	TS
HP1alpha-[CHR]-His	TS
HP1alpha-GST	TS
HP1beta-[CHR]-GST	TS
HP1beta-[CHR]-His	TS, AS

Reader Domain	Assays
HP1beta-GST	TS
HP1beta-His	TS, AS
HP1beta-Strep	TS
HP1gamma-GST	TS
HP1gamma-His	TS
KAT2A	TS
KAT2B	TS
KAT5-2-[CHR]-His	TS
KAT5-3-[CHR]-GST	TS
KAT5-3-[CHR]-His	TS
L3MBTL1	TS
L3MBTL1-His	AS
L3MBTL3-His	AS
MPP8-[CHR]-GST	TS
PB1-1	TS
PB1-2	TS
PB1-3	TS
PB1-4	TS
PB1-5	TS
PB1-6	TS
PHIP-2	TS
PHIP-Tndm	TS
SMARCA2 α bromodomain	HTRF
SMARCA2 α -His	TS
SMARCA2b-His	TS, AS
SMARCA4-His	TS, AS
SP100-GST	TS
SP100-His	TS
SP110c-GST	TS
SP140-GST	TS
SP140-His	TS, AS
SP140L-GST	TS
SP140L-His	TS, AS

Reader Domain	Assays
TAF1-1-GST	TS
TAF1-2-GST	TS
TAF1-2-His	AS
TAF1L-1	TS
TAF1L-2	TS
TAF1L-Tndm-GST	TS
TRIM24	TS
TRIM28	TS
TRIM33 α	TS
TRIM33b-His	TS
TRIM66	TS
UHRF1 Full length	TS
UHRF1-[PHD]	TS
UHRF1-[PHD]-His	AS
UHRF1-[SRA]	TS
UHRF1-[TDR-PHD]	TS
UHRF1-[TDR-PHD]-His	AS
UHRF1-[TDR]-His	TS, AS
UHRF1-His Full length	AS
YTHDC1	HTRF
YTHDC2	HTRF
YTHDF1	HTRF
YTHDF2	HTRF
YTHDF2-YTH	HTRF
YTHDF3	HTRF
YTHDF3-YTH	HTRF

AS...
AlphaScreen

TS....
Thermal Shift
Assay

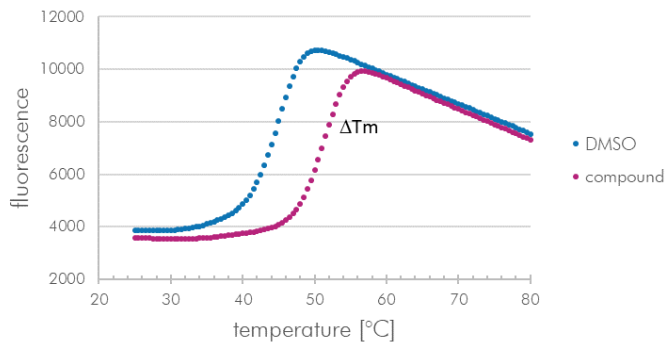
HTRF
Homogenous
Time-Resolved
Fluorescence

SPR
Surface
Plasmon
Resonance

BromoMELT

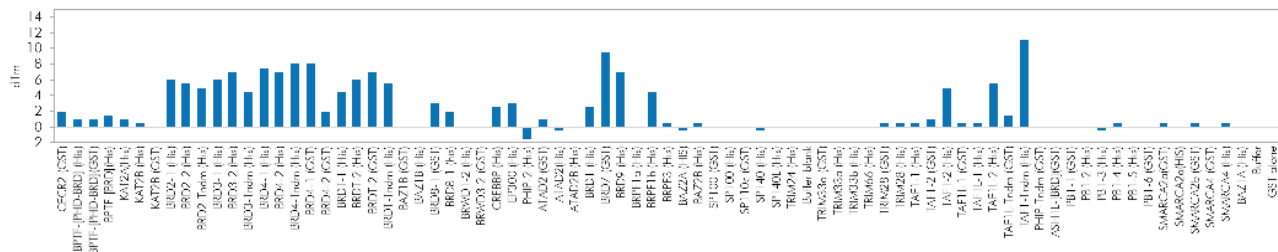
BromoMELT is a thermal shift assay for bromodomain targets that measures the change in protein melting temperature upon the binding of a ligand. Protein melting measurements are useful for identifying ligands, buffer conditions, co-factors and drugs affecting protein stability.

- Available as service or as kit to easily perform the assay in your own lab using a qPCR machine
- Includes 77 proteins representing 63 bromodomains
- Any inhibitor can be characterized within hours
- High-throughput compatible



Assay principle

The thermal shift assay determines the melting temperature at which there is 50% denaturation of the target protein. The difference between the melting temperatures of protein only (blue line) and protein plus ligand (purple line) is proportional to the binding affinity of the interaction.



Example of the selectivity profile of bromosporine

The binding of bromosporine to 77 bromodomain proteins was characterized using the BromoMELT assay kit. The difference in melting temperatures of target proteins bound to bromosporine versus DMSO control is proportional to the binding affinity of the protein/bromosporine interaction.

Methyltransferase Assays

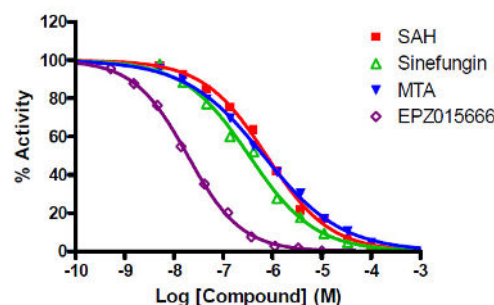
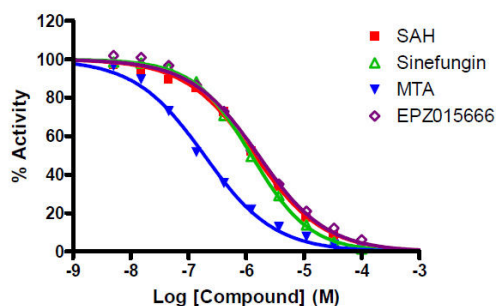
Reaction Biology offers radiometric activity assays and recombinant proteins for over 30 methyltransferases.

- Direct measurement of enzyme activity via radiometric assay
- Detection of inhibitors with varying binding modes
- Substrates can be nucleosomes, histones, peptides or other substrates
- Deliverable: % inhibition (single or multiple concentration) or IC_{50} values



Assay principle

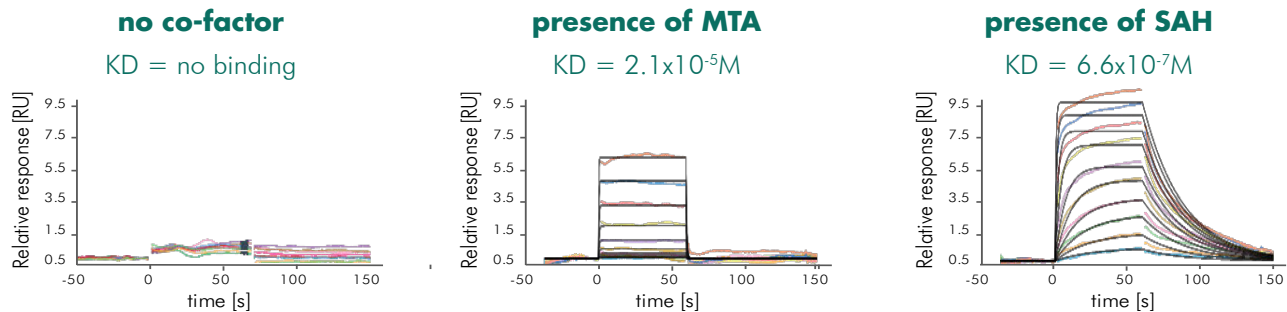
Methyltransferases use tritium-labeled S-adenosyl-L-methionine (SAM) as the methyl donor that is converted to S-adenosyl-L-homocysteine (SAH) during the transfer of the radioactive methyl group to the histone substrate.



Example: Inhibition of PRMT5/MEP50 activity

Concentration-dependent inhibition of PRMT5/MEP50 activity by inhibitors in comparison to SAH on either histone HA2 substrate (HotSpot assay, left) or H4-biotin substrate (FlashPlate assay, right).

Biophysical assays such as surface plasmon resonance (SPR) can be used to determine the binding affinities of inhibitors to epigenetic targets including enzymes and non-enzymatic proteins.



Example of a co-factor analysis by SPR

EPZ015666 is a substrate-competitive inhibitor that binds to its target PRMT5/MEP50 only in the presence of SAM or SAM analogues such as MTA and SAH.

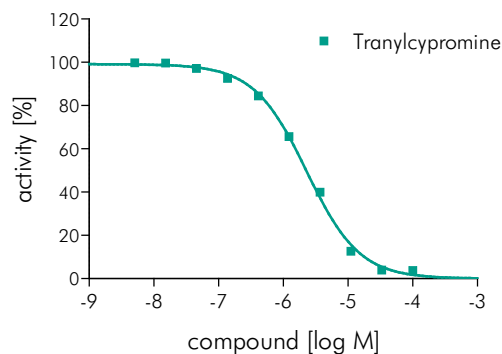
Methyltransferase	Protein available	Methyltransferase	Protein available	Methyltransferase	Protein available
ASH1L	√	MLL1 Complex	√	PRMT5/MEP50	√
COMT	√	MLL2 Complex	√	PRMT6	-
COMT (V108M)	√	MLL3 Complex	√	PRMT7	√
DNMT1	√	MLL4 Complex	√	PRMT8	√
DNMT3a	√	NRMT1	√	SET1B	√
DNMT3b	√	NSD1	√	SET7	√
DNMT3b/DNMT3L	√	NSD2	√	SET8	√
DOT1L	√	NSD2 (E1099K)	√	SETD2	√
EZH1 Complex	√	NSD2 (T1150A)	√	SETDB1	√
EZH2 (Y641F) Complex	√	NSD3	√	SMYD2	√
EZH2 Complex	√	PRDM9	√	SMYD3	√
G9a	√	PRMT1	√	SUV39H1	√
GLP	√	PRMT3	√	SUV39H2	√
METTL21A	-	PRMT4	√	SUV420H1-tv2	√
METTL3/METTL14	√	PRMT5 (C449S)/MEP50	√		

Demethylase Assays (KDM)

Reaction Biology offers assays for both Jumonji C-domain containing (JmjCs) and lysine-specific demethylases (LSD) histone demethylase subfamilies. LSDs are flavin-dependent monoamine oxidases that catalyze demethylation of Kme2 or Kme1 producing peroxide (H_2O_2) and formaldehyde (H_2CO) in the process. JmjC are Fe(II)/2-oxoglutarate-dependent dioxygenases that use a reactive Fe(IV)-oxoferryl species to catalyze hydroxylation reactions.

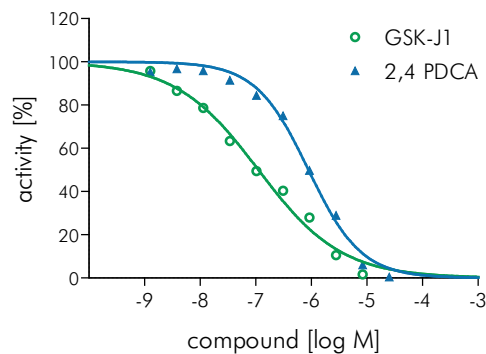
- Three assay formats are available to measure the activity of demethylases
- Customized conditions are available
- Deliverable: % inhibition (single or multiple test concentrations) or IC_{50} determination

Demethylase	Assay format	Protein available
KDM4A	HTRF	✓
KDM4C	HTRF	-
KDM4D	AlphaLISA	-
KDM5A	HTRF	✓
KDM5B	HTRF	✓
KDM5C	HTRF	✓
KDM6B	AlphaLISA	✓
LSD1	Amplex Red	✓



Example of LSD1 inhibition

LSDs activity was detected by quantification of H_2O_2 using Amplex Red reagent for IC_{50} value determination of a reference inhibitor.



Example of KDM5C inhibition

KDM5C activity was detected using HTRF (homogeneous time resolved fluorescence) technology with KDM5C and substrate-specific antibody for IC_{50} value determination of two reference inhibitors.

Histone Acetyltransferase Assays (HAT)

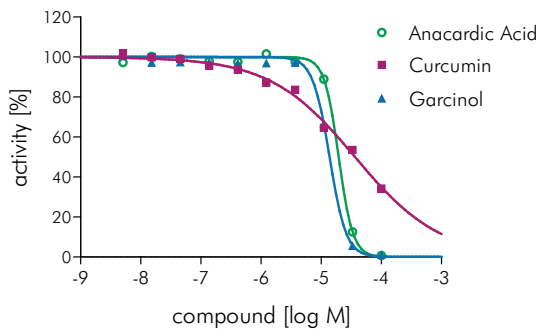
Reaction Biology offers radiometric activity assays for HAT enzymes using tritiated acetyl-Coenzyme A as cofactor.

- Direct measurement of enzyme activity via radiometric assay
- Customized conditions are available
- Deliverable: % inhibition (single or multiple test concentrations) or IC_{50} determination



Assay principle

Histone acetyltransferases acetylate lysines on histones and other proteins using tritium-labelled acetyl-Coenzyme A as the acetyl donor. The tritium-acetyl group is transferred onto histone substrate that is measured directly to reflect the enzyme activity.



Example of CBP inhibition

Full concentration-response of three reference inhibitors against CREP-binding protein CBP activity.

HAT	Protein available
CBP	✓
KAT2A	✓
KAT2B	✓
KAT5	✓
KAT6A	-
KAT6B	-
KAT7	-
KAT8	-
p300	✓

Histone Deacetylase (HDAC) and Sirtuin Assays

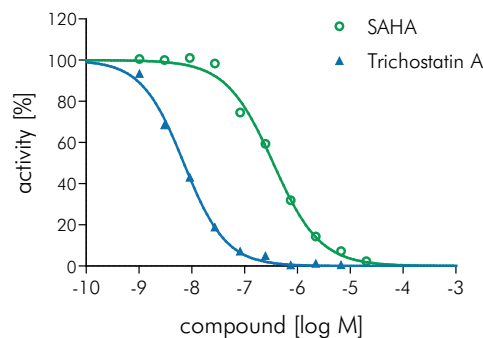
Reaction Biology offers fluorescence-based assays for both Zn^{2+} -dependent HDACs and NAD^{+} -dependent sirtuins.

Each assay is optimized based on its specific substrate:

- HDAC 1, 2, 3, 6, 10, and Sirt 1, 2, 3: p53 residues 379-382 (RHKKAc)
 - HDAC 8: p53 residues 379-382 (RHKAckAc)
 - HDAC 4, 5, 7, 9, 11: Trifluoroacetyl lysine
 - Sirt 5: Ac-Lys(Succinyl)-AMC
- Customized conditions and kinetic studies are available
 - Deliverable: % inhibition (single or multiple test concentrations) or IC_{50} determination

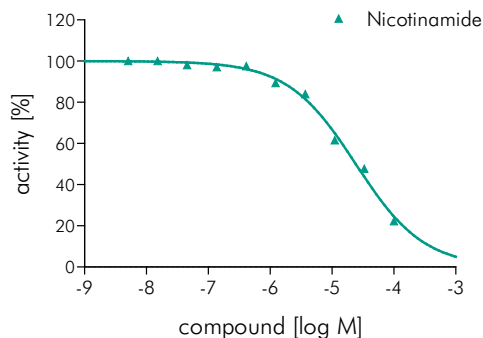
Assay principle

The assay is performed with purified human protein and a fluorogenic acetylated peptide substrate specifically designed for each enzyme. The deacetylated fluorogenic substrate is susceptible to cleavage by a protease to yield fluorescence.



Example of HDAC1 inhibition

Full concentration-response of two reference inhibitors against HDAC1 activity.



Example of SIRT5 inhibition

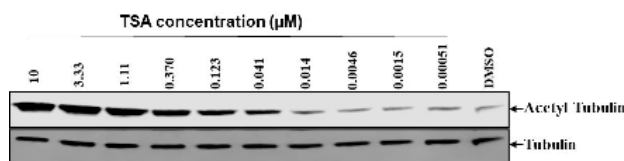
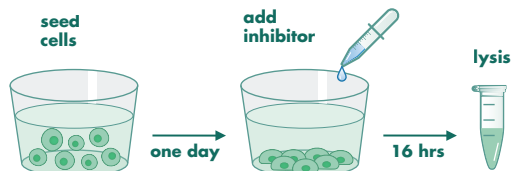
Full concentration-response of Nicotinamide, a pan-SIRT inhibitor, against SIRT5 activity.

Cell-based Epigenetic Assays

Cell-based assays are valuable tools for evaluating inhibitor potency to affect acetylation and/or methylation changes of substrates in a physiological environment using intact cells. The detection options include ELISA, Western Blot, NanoBRET and HDAC-Glo assays.

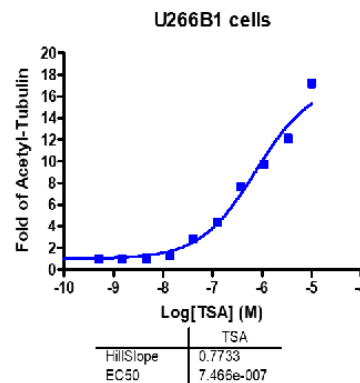
- Evaluate compound activity in intact cells.
- ELISA and Western Blot detect the endogenous substrates for direct activity measurement
- Deliverable: IC_{50} values of epigenetic enzyme inhibition

Readout	Assay format
Histone & Tubulin Deacetylation	Western Blot; ELISA
Histone Deacetylation (Class I/II)	HDAC-Glo
Histone Methylation	Western Blot
Histone Phosphorylation	ELISA



Example for detection of HDAC activity in cancer cells

Lysates of U266B1 cells treated with Trichostatin A (TSA) were subjected to Western Blotting and quantification via infrared imaging. The results are plotted in a curve for EC_{50} determination.



RAS PATHWAY ASSAYS

Reaction Biology provides a variety of services to discover new inhibitors targeting the RAS pathway. The small GTPase, RAS, is a known oncogene that is mutated in a large percentage of cancers and is associated with poor disease prognosis. Mutated RAS is locked in the activated GTP bound state and facilitates enhanced RAS signaling in cancer cells.

Most of our assays are available with wildtype and mutated RAS variants.

Available Assay Formats	Description
Nucleotide Exchange Assay	Measuring of SOS1/2 mediated exchange of fluorescently labeled GDP to GTP Alternative readout: Observation of an increase in HTRF upon binding of fluorescent GTP to K-RAS
Protein-Protein Interaction of RAS and SOS1	HTRF based assay for testing of compounds that disrupt SOS1 binding to RAS.
Protein-Protein Interaction of RAS and cRAF	HTRF-based assay for testing of compounds that disrupt cRAF binding to RAS. This assay is also suited for quantification of nucleotide exchange reaction.
Thermal Shift Direct Binding Assay	Compound binding affinity measurement suited for measurement of compound selectivity to RAS mutant panel.
SPR Direct Binding Assay	Surface Plasmon Resonance (SPR) determines the kinetics of compounds binding RAS and RAS mutants or SOS.
NanoBRET Target Engagement RAS Assay	Intracellular measurement of the binding affinity of compounds via competitive displacement of a switch I/II pocket tracer

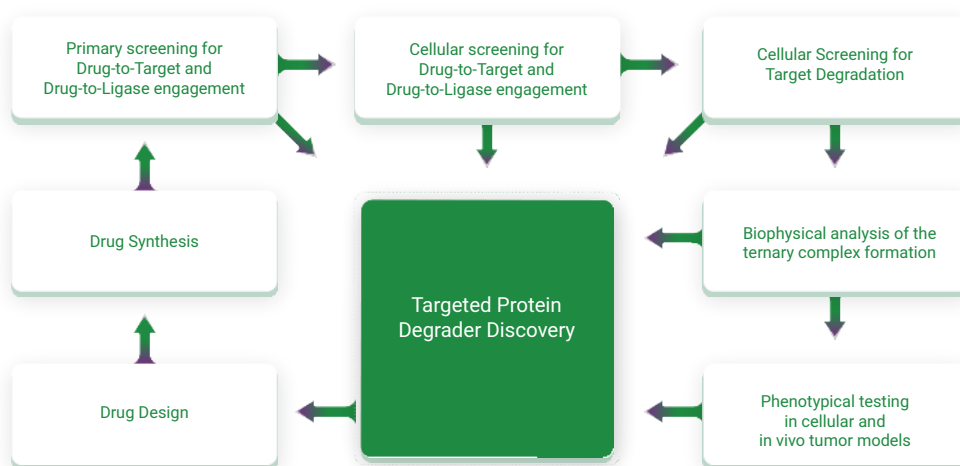
In addition to RAS specific assays we offer more than 80 RAS Pathway related assays for testing inhibitors against the ERK/MAPK and the PI3K signaling pathway as well as upstream pathways such as EGFR signaling including kinases, phosphatases and transcription factors.

TARGETED PROTEIN DEGRADATION ASSAYS

Rethinking PROTAC: An AI based platform to support your drug discovery project

Together with Medinoah, a medicinal chemistry provider, and PMRBioinfo, an AI computational company, we have created the Targeted Protein Degradation drug discovery platform. Medinoah has been one of the first CROs in synthesizing protein degradation molecules and advanced two PROTAC molecules into the IND stage. The AI powered Targeted Protein Degradation discovery platform will enable prediction of the binding mode of the ternary complex to significantly reduce the number of compounds and screening cycles needed for the generation of potent and optimized protein degradation molecules.

Workflow of discovery of new Targeted Protein Degradation molecules platform:

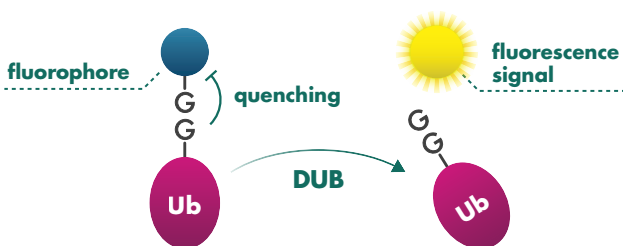


We are currently talking to biotech and pharma companies to perform fee-for-service, shared-cost collaboration or FTE-based work for the development of new Targeted Protein Degradation molecules. Please reach out to us to discuss options for supporting your Targeted Protein Degradation drug discovery program.

UBIQUITIN-PROTEASOME PATHWAY ASSAYS

Deubiquitinase enzymes (DUBs) are a group of proteases that cleave ubiquitin from proteins and other molecules. Ubiquitination of proteins affects protein degradation, cellular location, activities and protein-protein interactions. Reaction Biology offers fluorescent-based assays for screening of DUB inhibitors.

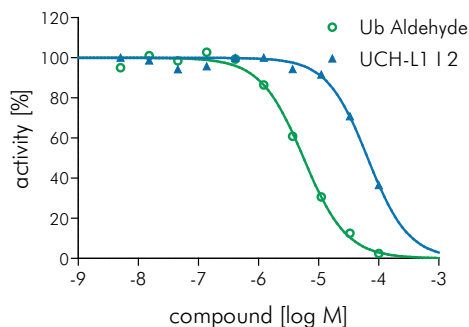
- Fluorescence-based quantification of deubiquitinase and 20S proteasome activity
- Custom assay development
- Deliverable: % inhibition (single point) or IC_{50} profiling, K_i determination



Assay procedure

Fluorescent Ubiquitin-AMC is a substrate containing the fluorophore, 7-amido-4-methylcoumarin (AMC), that is quenched when ubiquitinated. Upon incubation with a deubiquitinase, AMC is released and its fluorescence can be measured.

Available Targets	
20S Proteasome	USP5
A20	USP7
Ataxin3	USP8
BAP1	USP9X
MYSM1	USP10
NEDP1	USP11
SEN1	USP13
SEN2	USP14
UCHL1	USP15
UCHL3	USP20
UCHL5	USP25
USP2	USP28
USP4	USP30



Example of A20/TNFAIP3 inhibition

Concentration-dependent inhibition of deubiquitinase A20 by two inhibitors. IC_{50} value determination is based on 10 compound concentrations.

PARP ASSAYS

Poly (ADP-ribose) polymerase (PARP) is a family of proteins that transfer ADP-ribose units from NAD⁺ onto target nuclear proteins forming long branched Poly ADP-ribose chains. PARPs play a role in epigenetic regulation, for example, by poly ADP-ribosylation of histone substrates.

- Gold standard assay format: radiometric activity assay
- High-throughput compatible
- Custom assay development
- Deliverable: % inhibition (single point) or IC₅₀ profiling

PARP	Protein available
ARH3	✓
PARG	✓
PARP1	✓
PARP2	✓
PARP5B	-

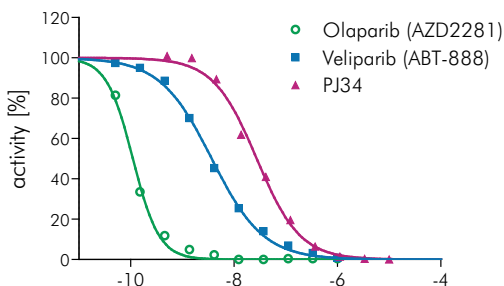


Assay procedure

Adenylate-NAD⁺ serves as co-factor for transfer of ³²P-labelled ADP-ribose units onto histones which will be quantified via scintillation counting.

NanoBRET TE PARP

PARP1
PARP2
PARP5A
PARP5B
PARP7
PARP11
PARP12



Example of PARP2 inhibition

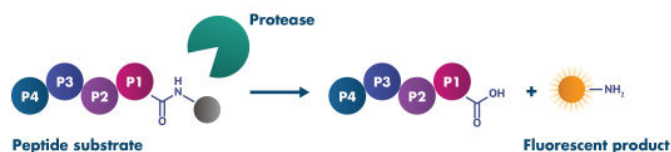
Dose-dependent inhibition of PARP2 by three reference inhibitors. IC₅₀ value determination is based on 10 compound concentrations.

PROTEASE ASSAYS

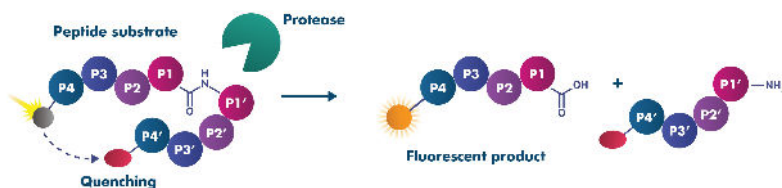
Reaction Biology offers a 65-member protease panel for routine profiling. Members include cysteine proteases, serine proteases, metalloproteases, aspartyl proteases, dipeptidases and others. Over 80 proteases are available for customized orders.

- Fluorescence-based activity assay
- High-throughput compatible
- Custom-assay development
- Deliverable: % inhibition (single point) or IC₅₀ value determination

Fluorogenic peptide substrate



FRET/quencher peptide substrate



Assay formats

Upper image: Fluorogenic peptide substrate is cleaved by target protease releasing a fluorescent product.

Lower image: The peptide substrates contain a fluorophore and a quencher suppressing fluorescence unless substrate is cleaved.

Proteases		
ACE1, 2	Furin	Plasmin
Activated Protein C	Granzyme B	Proteinase A, K
ADAM10, TS4, TS5	Hepsin	Renin
BACE1	HIV-1	SARS-CoV-2 Mpro, PLpro
Calpain1	HRV-3C	TACE
Caspase 1, 2, 3, 4, 5, 6, 7, 8, 9, 10, 11, 14	Kallikrein 1, 2, 3/PSA, 4, 5, 6, 7, 8, 11, 12, 13, 14	Thrombin
Cathepsin B, C, D, E, G, H, K, L, S, V, Z	MALT1	TMPRSS11D
Chymotrypsin	Matriptase	TMPRSS2
Chymase	Matriptase 2	tPA
DPP-III, IV, VIII, IX	MMP 1, 2, 3, 7, 8, 9, 10, 12, 13, 14	Trypsin
Elastase	Nepriylsin	Tryptase a/b1, g1
Factor VIIa, IXa, Xa, XIa, XIIa	Papain	Urokinase
FAP	Plasma Kallikrein	

PHOSPHATASE ASSAYS

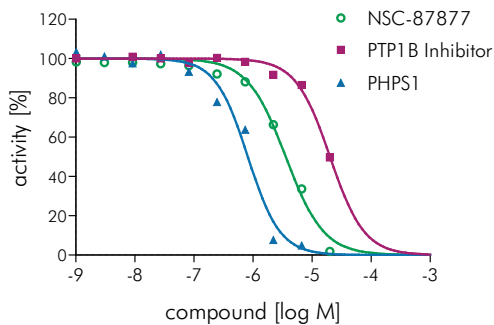
Protein phosphatases play important roles in cell signalling processes in interplay with kinases. Different than kinases, phosphatases are less specific for their substrates. All phosphatases catalyze the same basic hydrolysis reaction.

- Fluorescence-based activity assay
- High-throughput compatible
- Custom-assay development
- Deliverable: % inhibition (single point) or IC_{50} value determination



DiFMUP-based assay principle

The fluorinated MUP derivate is suitable as substrate for a large range of protein phosphatases. The reaction product of DiFMUP is fluorescent after dephosphorylation.



Example of PTPRC/CD45 inhibition

Full concentration-response of 3 reference inhibitors.

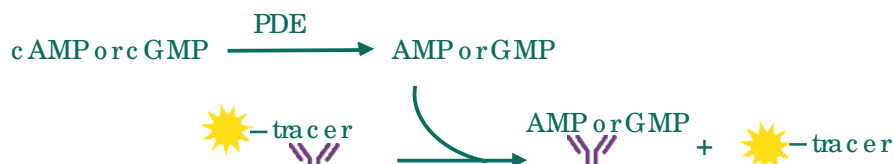
Phosphatases

DUSP22/MKPX
PP1A
PP1B
PP2A alpha/PPP2R1A Complex
PP2C alpha
PP2C gamma
PPAC (ACP1/LMW-PTP-A)
PPAC (ACP1/LMW-PTP-B)
PTEN
PTPN1/PTP1B-CD
PTPN1/PTP1B-FL
PTPN2/TC-PTP
PTPN6/SHP1
PTPN7/LC-PTP
PTPN11/SHP2 (E76K)-FL
PTPN11/SHP2-CD
PTPN11/SHP2-FL
PTPN12/PTP-PEST
PTPRB
PTPRC/CD45
PTPRF/LAR
PTPRJ/CD148

PHOSPHODIESTERASE (PDE) ASSAYS

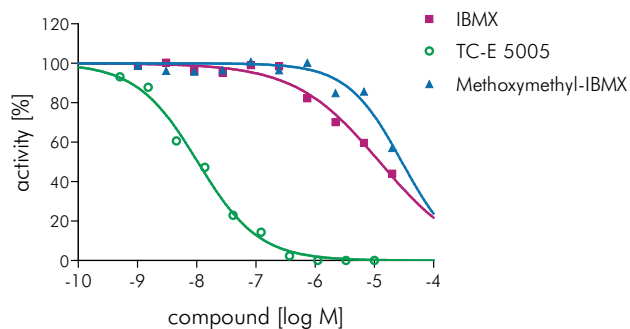
Phosphodiesterases (PDEs) catalyze the hydrolysis of cyclic AMP and cyclic GMP, thereby regulating these cyclic nucleotides' signaling pathways. PDE inhibitors have found utility in the treatment of a variety of conditions including pulmonary hypertension, acute refractory cardiac failure, erectile dysfunction, etc.

- Activity of enzymes is measured with the Transcreener AMP2/GMP2 FP PDE assay platform (BellBrook labs)
- High-throughput compatible
- Custom-assay development
- Deliverable: % inhibition (single point) or IC₅₀ values



Assay principle

PDE converts cAMP or cGMP to AMP or GMP which displace a fluorescent tracer from an antibody selective for AMP and GMP resulting in reduction of the fluorescence polarization signal.



Full concentration-response for three reference inhibitors of PDE 10A.

PDEs
PDE1A
PDE1B
PDE1C
PDE2A
PDE3A
PDE3B
PDE4A
PDE4B
PDE4C
PDE4D
PDE4D2
PDE5A
PDE7A
PDE7B
PDE8A
PDE8B
PDE9A
PDE10A

METABOLIC PATHWAY ASSAYS

Acetyl-CoA Carboxylase (ACC)

ACC is a biotin-dependent enzyme that catalyzes the ATP-dependent carboxylation of acetyl-CoA to malonyl-CoA. ACC is a crucial metabolic enzyme and attractive drug target. Reaction Biology provides compound screening against ACC by detecting the production of ADP.

Isocitrate dehydrogenase - IDH

Isocitrate dehydrogenases 1 and 2 (IDH1 and IDH2) are key metabolic enzymes that catalyze the conversion of isocitrate to α -ketoglutarate (α KG) and co-factor NADPH. Reaction Biology provides compound screening against IDH by measuring enzyme activity in a coupled system wherein NADPH produced in the initial reaction is a co-factor in the conversion of resazurin to fluorescent resorufin in a secondary reaction.

IDHs are also available for purchase.

NAD(P)H dehydrogenase [quinone] 1 - NQO

NQOs are involved in detoxification and biosynthetic pathways. Reaction Biology provides compound screening against NQOs by monitoring enzyme activity in an analogous coupled reaction as described above for IDHs.

Nucleotide Metabolism Pathway Assays

Nucleotide metabolism is the process in which nucleic acids (RNA, DNA, and cellular bioenergetics) are synthesized and degraded. DHODH (dihydroorotate dehydrogenase) synthesizes orotate from dihydroorotate (DHO) in the de novo pyrimidine synthesis pathway. The enzyme was shown to induce the differentiation of acute myeloid lymphomas and is therefore interesting as a drug target.

Our portfolio includes methylenetetrahydrofolate dehydrogenase (MTHFD) assays essential for purine and thymidylate biosynthesis. MTH1 (NUDT1) nucleotide pool sanitization assays enable discovery of compounds exploiting cancer's oxidative stress vulnerability by preventing removal of damaged nucleotides.

Carboxylase

ACC1

ACC2

IDH

IDH1 G97D IDH2 R140K

IDH1 R100A IDH2 R140Q

IDH1 R100Q IDH2 R172Q

IDH1 R132C IDH2 WT

IDH1 R132H

IDH1 Y139D

IDH1 WT

NQO

NQO1

NQO2

DHODH

DHODH

dDHODH (dog)

mDHODH (mouse)

rDHODH (rat)

MTH1

MTHFD1

MTH1FD2

ION CHANNEL ASSAYS

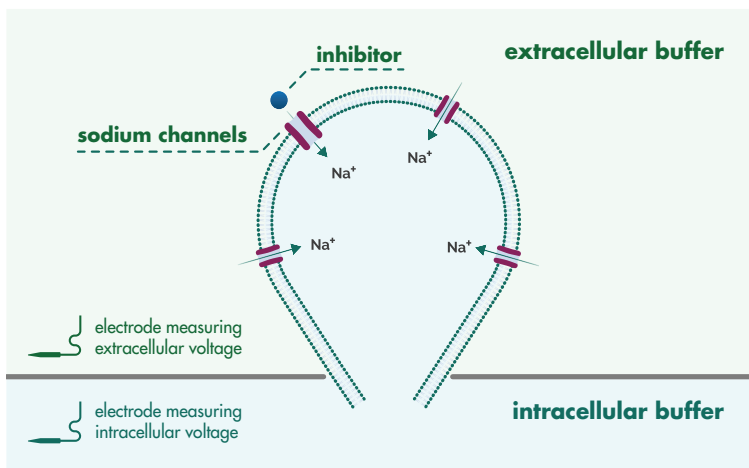
Reaction Biology offers cell-based ion channel testing for drug discovery and evaluation of drug safety.

- Cardiac-safety panel is available for pre-clinical evaluation of compound safety
- Investigation of voltage-gated and ligand-gated ion channels
- Single-concentration screen, vehicle and positive controls in every assay
- Deliverable: IC₅₀ value of inhibition of ion channel activity (6-point dose-response)

Manual patch clamp (MultiClamp 700B)

Manual patch clamp electrophysiology remains the regulatory gold standard for characterizing ion channel block. Gigaohm seal formation and whole-cell voltage clamp provide the highest fidelity measurements of compound effects on ionic currents, essential for IND-enabling cardiac safety packages.

Ion channels
hERG
Nav1.5
Cav1.2
Kv7.1/minK



Assay principle

Schematic of a cell expressing sodium channels in a planar patch-clamp setup with recording in the whole-cell format.

G-PROTEIN-COUPLED RECEPTOR (GPCR) ASSAYS

GPCRs represent the largest individual family of targets for currently approved medications. Recent advances in GPCR pharmacology, including biased signaling and allosteric modulators, have become increasingly important tools in drug discovery. Reaction Biology offers services to progress drug discovery research in the area of GPCR biology and pharmacology.

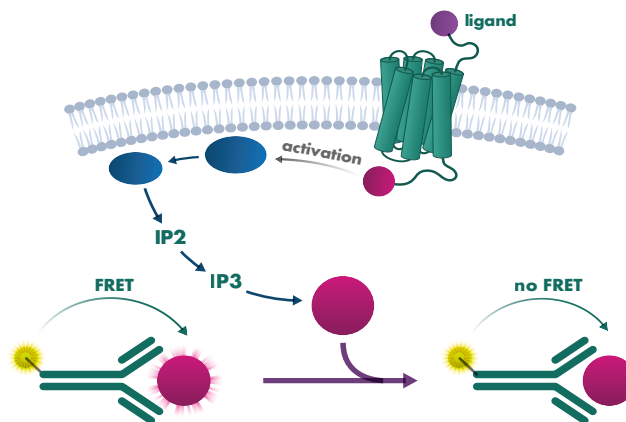
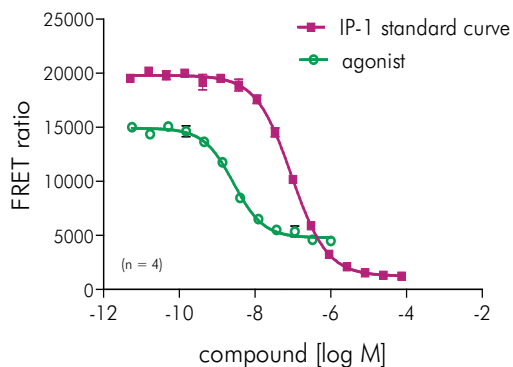
- We offer assay development, high-throughput screening, SAR support services.
- Transmembrane signaling assay formats can be readily established for your receptor of interest including calcium mobilization, β -arrestin translocation, cAMP generation and inositol 1-monophosphate (IP1) generation

Our dedicated team of GPCR experts will enable drug screening with assays tailored to your specific needs:

1. Define the needs and scope for the project together with our assay development team.
2. Make us familiar with the goals for your research project and define timelines to ensure goal-oriented work right from the start.
3. We will acquire or generate a cell line appropriate to the project needs.
4. The same high standards we use for our off-the-shelf assays will apply to newly developed assays for your project.
5. We guarantee fast turn-around times for data generation.
6. During every step of the process, you will be in close contact with your project manager for regular updates on the study progress.

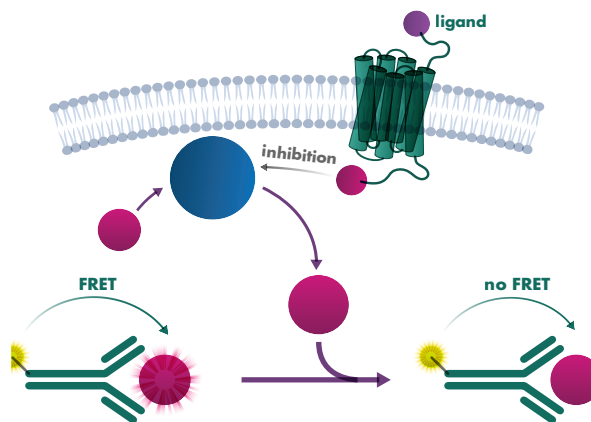
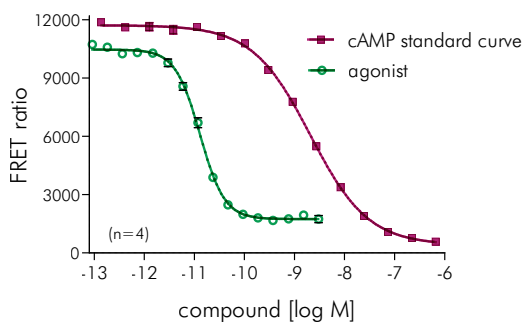
Agonist-induced IP1 generation

Cells engineered to express Gq-coupled GPCRs of interest are stimulated with an agonist for IP1 accumulation. The generated IP1 is measured using a competitive immunoassay wherein cellular IP1 competes with a labeled IP1 for binding to an anti-IP1-cryptate generating a FRET signal.



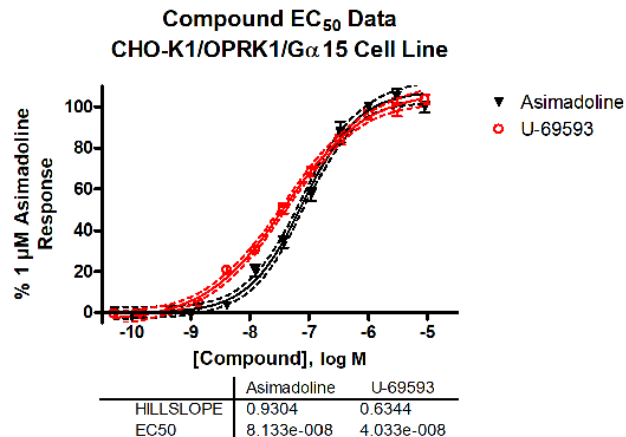
Agonist-induced cAMP generation

Cells engineered to express Gs-coupled GPCRs of interest are stimulated with an agonist to affect cellular activation. cAMP can be accurately measured by a variety of standard detection methods including a competitive immunoassay wherein cellular cAMP competes with a labeled cAMP probe to bind to an anti-cAMP-cryptate generating a FRET signal.



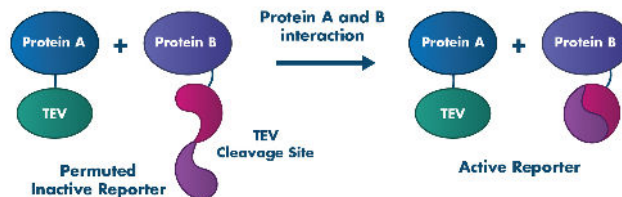
FLIPR calcium mobilization assay

The concentration-response of agonist stimulation for calcium mobilization is shown.

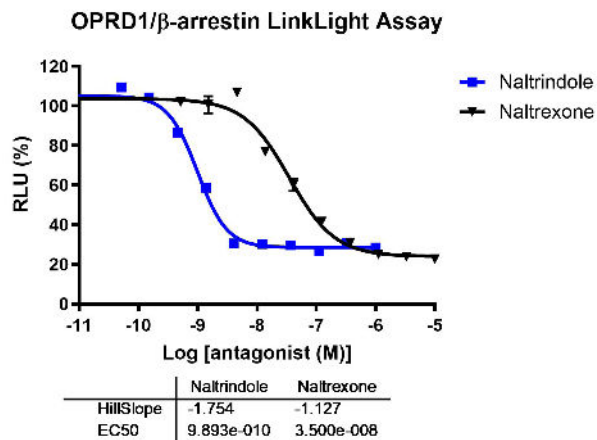


LinkLight™: Protein-Protein Interaction

Cells engineered to express GPCRs of interest are stimulated with test compounds to trigger receptor activation. Upon ligand binding, the GPCR undergoes conformational changes leading to G-protein signaling followed by GRK-mediated phosphorylation. The phosphorylated receptor recruits β-arrestin, bringing TEV protease (fused to the GPCR) into proximity with permuted luciferase (fused to β-arrestin). TEV cleavage activates the luciferase, generating a persistent luminescent signal that can be accurately measured by standard luminometry detection methods.



Levels of β-arrestin recruitment were measured after stimulation of a GPCR expressing cell line with an undisclosed agonist using LinkLight TEV/luciferase technology. The irreversible luminescent signal provides dose-dependent readout of regulatory pathway activation complementing cAMP/calcium data.



NUCLEAR RECEPTOR ASSAYS

Nuclear receptors are transcription factors that are regulated by small hydrophobic ligands such as hormones or vitamins. Upon ligand binding, nuclear receptors translocate into the nucleus to bind DNA and modulate the expression of their target genes to regulate a variety of cellular mechanisms such as growth, proliferation, metabolism, or homeostasis on a transcriptional level.

Reaction Biology offers a suite of assays to support the discovery of new drugs to modulate nuclear receptors with the following advantages:

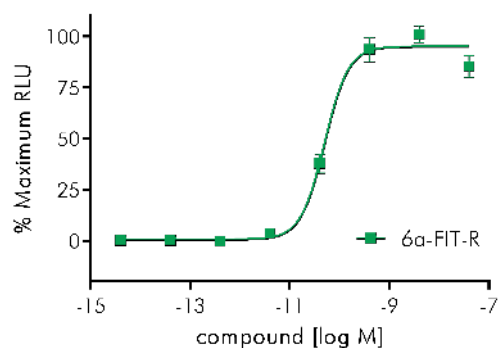
- Cell-based assays allow for drug testing in the physiological and complex environment of intact cells
- Any class of inhibitor can be tested such as modulators of translocation, activity, structural changes, dimerization, etc.
- Custom-assay development

Nuclear Receptors

Androgen receptor (AR)

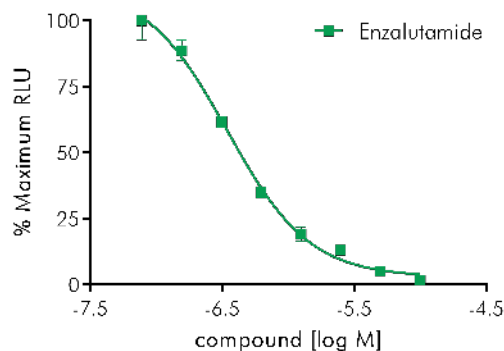
Farnesoid X receptor (FXR)

Pregnane X receptor (PXR)



AR Agonist Assay:

The androgen receptor reporter cell line was stimulated with various concentrations of agonist 6α-FIT-R for 24 hours before luminescence measurements were taken.



AR Antagonist Assay:

The androgen receptor reporter cell line was stimulated with a fixed concentration of agonist 6α-FIT-R and various concentrations of antagonist Enzalutamide for 24 hours before luminescence readout.

ADDITIONAL ASSAYS

Heat shock protein 90 – HSP90

HSP90 is an ATP-dependent molecular chaperone that stabilizes several regulatory molecules including many tyrosine kinases and transcription factors. Reaction Biology provides compound screening for both HSP90a and HSP90b isoforms.

HSP90	
HSP90a	
HSP90b	

Cytochrome P450 - CYP

CYPs are heme proteins that play key roles in the metabolism of drugs. Understanding a compound's inhibitory activity against key CYP proteins is essential for predicting drug-drug interactions. Reaction Biology provides compound profiling against the 14 most important CYP isoforms that affect drugs pharmacokinetics and responses.

CYP	
CYP 1A2	CYP 2D6
CYP 19A	CYP 2E1
CYP 2A6	CYP 2J2
CYP 2B6	CYP 3A4
CYP 2C8	CYP 3A5
CYP 2C9	CYP4A11
CYP 2C19	CYP4F3B

Kinesin ATPase

Kinesin superfamily proteins (KIFs) are a large family of molecular motor proteins, that share a highly conserved motor domain. Kinesins play important role in cell division and transport of vesicles and organelles within cells. Reaction Biology offers ATPase activity inhibitor screening assays for kinesin targets via Promega's ADP Glo™ detection system. Kinesin proteins available.

KIF	
KIF2C	KIF18B
KIF3C	KIF19
KIF4A	KIF20A
KIF5B	KIF22
KIF10/CENPE	KIF23
KIF11/Eg5	KIFC1
KIF18A	KIFC3

Apoptosis Pathway Assays

Apoptosis which is a highly regulated form of programmed cell death that is needed to remove damaged or aged cells. Modulating apoptosis has been exploited to fight malignancies for example by overcoming treatment resistance. The apoptosis assay technology measures the binding of the apoptotic protein such as BCL2 to a substrate peptide such as BAK. The interference of the binding by a test molecule can be detected via a fluorescence readout.

Apoptosis	
Bcl-2	MCL1
BCL-xL	p53/MDM2
clAP-1	XIAP
clAP-2	



CUSTOMIZED ASSAY DEVELOPMENT

Reaction Biology provides protein production and assay development based on customer requests. Targets include enzymes, protein-protein interaction, GPCRs, nuclear receptors, ion channels, and more.

Examples of assays developed for customers

Aldo-keto reductase	RNA polymerase
Dihydroorotate dehydrogenase	RNA epigenetic enzymes
DNA cytidine deaminase	PP2A/CIP2A
Ectonucleotide pyrophosphatase	RASGRP
GABA aminotransferase	Sentrin-specific proteases
RNAse H2	USP
Transcription factors	

Assay formats available for customized assay development

AlphaLisa	Radiometric assays using ^3H , ^{32}P or ^{33}P
AlphaScreen	Autoradiogram after SDS-PAGE
ELISA	HTFR
Thermal shift assay	ADP-Glo
Surface plasmon resonance	Fluorescent peptide screening
Microscale thermophoresis	NanoBRET
Isothermal titration calorimetry	Fluorescence Polarization
Flow cytometry	FRET
MSD	IncuCyte
Electrophysiology	

We have developed over 1,500 assays. Use our vast experience to develop an assay for the target of your interest.

NOTES



NOTES

LET'S DISCOVER TOGETHER.

Recombinant Proteins

- Kinase proteins
- Epigenetic proteins
- Substrates
- Custom-tailored protein production



Biochemical Assays

- Kinases, Epigenetic Enzymes
- Protein: Protein Interaction assays
- Metabolic & Pathway
- Receptors & Channels



Cell-Based Assays

- 2D and 3D proliferation assays
- Drug combination screening
- Migration assays
- Angiogenesis assay



Biophysical Assays

- Surface Plasmon Resonance
- Thermal Shift Assay
- Isothermal Titration Calorimetry
- Microscale Thermophoresis



In Vivo Pharmacology

- In Vivo Hollow Fiber Model
- Xenograft models
- Orthotopic models
- Metastasis models



Safety & Toxicology

- In Vitro Safety Panel
- Cardiac Safety Panel
- Cytochrome P450
- Maximum-Tolerated Dose
- GLP Regulatory Toxicology



Integrated Solutions

- Antibody Drug Conjugates
- RAS Drug Discovery
- Targeted Protein Degradation



Biomarker Discovery

- Genomic biomarkers
- Protein biomarkers
- Immunophenotyping



Immuno-Oncology

- In Vitro Killing Assays
- Syngeneic Mouse Models
- Proprietary Tumor Models
- Immunophenotyping



Reaction Biology | www.reactionbiology.com | requests@reactionbiology.com