



## PRODUCTS TARGET LIST

Enabling Drug Discovery and Development Programs

- Checkpoint Receptors
- Cytokine Receptors/Interleukins
- Cytotoxicity
- Epigenetic Proteins
- GPCRs
- Ion Channels
- Kinases and Phosphatases
- Nuclear Hormone Receptors
- Signaling Pathways



Checkpoint Receptor Products		Product Type		
Target	Application	Cell Line	eXpress Kit	Bioassay Kit
BCMA	Internalization	●		
BTLA	Signaling	●		
CD28	Signaling	●		
CD33	Internalization	●		
CD47	Ligand-Presenting	●		
mCD80	Ligand-Presenting	●		
CD86	Ligand-Presenting	●		
CD137	Signaling	●		
CD200	Ligand-Presenting	● <i>New</i>		
CD200R	Signaling	● <i>New</i>		
CTLA4	Signaling	●		
mCTLA4	Signaling	●		
FcγR1a	Clustering	●		
FcγR11a	Clustering	●		
FcγR11b	Clustering	●		
mFcγR11b	Clustering	●		
HVEM	Ligand-Presenting	●		
ICOS	Signaling	●		
LAIR1	Signaling	●		
NIK	Signaling	●		
OX40	Signaling	●		
PD-1	Signaling	●		●
mPD-1	Signaling	●		
PD-L1	Ligand-Presenting	●		
mPD-L1	Ligand-Presenting	●		
PD-L2	Ligand-Presenting	●		
Siglec9	Internalization	●		
SIRPα	Signaling	●		●
TIM3/TIM3	Dimerization	●	●	
VISTA/VISTA	Dimerization	●	●	

Cytokine Receptor/Interleukin Products		Product Type		
Target	Application	Cell Line	eXpress Kit	Bioassay Kit
ACVR1/ACVR2	Dimerization	●		
ACVR1B/BMPR2	Dimerization	●	●	
ACVR1C/ACVR2	Dimerization	●	●	
ACVR1C/ACVR2B	Dimerization	●	●	
ACVRL1/ACVR2	Dimerization	●	●	
ACVRL1/ACVR2B	Dimerization	●	●	
ACVRL1/BMPR2	Dimerization	●	●	
BMPR1A/ACVR2	Dimerization		●	
BMPR1A/ACVR2A	Dimerization	●		
BMPR1A/ACVR2B	Dimerization	●		
BMPR1A/BMPR2	Dimerization	●	●	
BMPR1B/ACVR2A	Dimerization	●	●	
BMPR1B/ACVR2B	Dimerization	●		
BMPR1B/BMPR2	Dimerization	●		
c-MET/c-MET	Dimerization	●		
CSF1R/CSF1R	Dimerization	●	●	
CSF2RB/CSF2R	Dimerization			●
CSF2RB/CSF2RA	Dimerization	●	●	
EGFR/EGFR	Dimerization	●	●	
EGFR/ErbB2	Dimerization	●	●	
EGFR/ErbB3	Dimerization	●		
EpoR/EpoR	Dimerization	●	●	●
ErbB2/ErbB3	Dimerization	●		
ErbB2/ErbB4	Dimerization	●	●	
ErbB4/ErbB4	Dimerization	●	●	
FGFR3(G380R)/FGFR3(G380R)	Dimerization	●		
FGFR3/FGFR3	Dimerization	●		
IFNAR1/IFNAR2	Dimerization	●		
IFNGR1/IFNGR2	Dimerization	●	●	
IFNRL1/IL10RB	Dimerization	●	●	
IL1R/IL1R	Dimerization			●
IL1R1/IL1RAP	Dimerization		●	
IL1R1/ILRAP	Dimerization	●		
IL1RL1/IL1RAP	Dimerization	●	●	●
IL2	Signaling Reporter	● <i>New</i>		
IL2/IL2	Dimerization			●
IL2RB/IL2RG	Dimerization	●	●	
IL2RB/IL2RG/IL2RA	Dimerization	●	●	●
IL2	Signaling Reporter	● <i>New</i>		
IL3RA/CD131	Dimerization	●		
IL4R/IL13RA1	Dimerization	●	●	
IL4R/IL2RG	Dimerization	●	●	
IL5RA/CSF2RB	Dimerization	●	●	
IL6R/IL6ST	Dimerization			●
IL6RA/IL6ST	Dimerization	●	●	●
IL7R/CRLF2	Dimerization	●	●	
IL7R/IL2RG	Dimerization	●	●	●
IL9R/IL2RG	Dimerization	●		
IL10RA/IL10RB	Dimerization	●	●	●

Cytokine Receptor/Interleukin Products		Product Type		
Target	Application	Cell Line	eXpress Kit	Bioassay Kit
IL11R/IL6ST	Dimerization	●		
IL12RB1/IL12RB2	Dimerization	●	●	
IL-15 (IL2R $\beta$ /IL2R $\gamma$ )	Dimerization	● <i>New</i>		● <i>New</i>
IL17RA/IL17RB	Dimerization	●		
IL17RA/IL17RC	Dimerization	●	●	●
IL17RA/IL17RC (U2OS)	Dimerization		●	
IL17RD/TNFR2	Dimerization	●	●	
IL18	Signaling Reporter	●		
IL18R1/IL18RAP	Dimerization	●	●	
IL20R1/IL20R2	Dimerization	●		
IL20RA/IL10RB	Dimerization	●	●	
IL20RA/IL20RB	Dimerization		●	
IL21R/IL2RG	Dimerization	●	●	
IL22RA/IL10RB1	Dimerization	●		
IL22RA1/IL10RB	Dimerization		●	
IL23R/IL12RB1	Dimerization	●	●	
IL27RA/IL6ST	Dimerization	●		
IL31RA/OSMRb	Dimerization			●
IFN $\gamma$	Signaling Reporter	●		
KDR/KDR	Dimerization	●	●	
LEPR/LEPR	Dimerization	●		
OSMR/IL6ST	Dimerization	●	●	
OSMRb/IL31R	Dimerization	●		
OSMRb/IL31RA	Dimerization		●	
PD-1/PD-1	Dimerization	●		
PD-L1/PD-L1	Dimerization	●		
TGFBR1/ACVR2	Dimerization		●	
TGFBR1/ACVR2B	Dimerization	●	●	
TGFBR1/TGFBR2	Dimerization	●	●	
TGFBR1/TGFBR2/ENG	Dimerization	●		
TGFBR1A/ACVR2A	Dimerization	●		
TPOR/TPOR	Dimerization	●		

KILR® Cytotoxicity Products			Product Type			
Target	Cell Type	Application	Cell Line	Cell Pool	Primary Cell	Bioassay Kit
CD16	Effector Cell	Cytotoxicity			●	
EGFR	U2OS	Cytotoxicity	●			
LAG3	Jurkat	Cytotoxicity	●			
PD-1	Jurkat	Cytotoxicity	●			
PD-L1	U2OS	Cytotoxicity	●			
PD-L2	U2OS	Cytotoxicity	●			
TIM3	Jurkat	Cytotoxicity	●			
	4T1	Cytotoxicity		●		
	A498	Cytotoxicity		●		
	A549	Cytotoxicity		●		
	ARH-77	Cytotoxicity	●			
	BT-474	Cytotoxicity		●		
	CCRF-CEM	Cytotoxicity		●		
	COLO-205	Cytotoxicity		●		
	Daudi	Cytotoxicity		●		●
	DU-145	Cytotoxicity		●		
	EL4	Cytotoxicity		●		
	H322	Cytotoxicity		●		
	HCT-116	Cytotoxicity		●		
	HepG2	Cytotoxicity		●		
	HL-60	Cytotoxicity		●		
	HT-1080	Cytotoxicity		●		
	Hut78	Cytotoxicity		●		
	Jurkat	Cytotoxicity		●		
	K562	Cytotoxicity		●		
	MCF7	Cytotoxicity		●		
	MDA-MB-231	Cytotoxicity		●		
	MM-1R	Cytotoxicity		●		
	MOLT-4	Cytotoxicity		●		
	NCI-H292	Cytotoxicity		●		
	NCI-N87	Cytotoxicity		●		● <i>New</i>
	PANC-1	Cytotoxicity		●		
	Raji	Cytotoxicity		●		●
	Ramos	Cytotoxicity		●		
	RPMI 8226	Cytotoxicity		●		
	SKBR3	Cytotoxicity		●		
	SK-MEL-28	Cytotoxicity		●		
	SK-MEL-5	Cytotoxicity		●		
	SKOV3	Cytotoxicity		●		
	SR	Cytotoxicity		●		
	T2	Cytotoxicity		●		
	THP-1	Cytotoxicity		●		
	U118-MG	Cytotoxicity		●		
	WIL2-S	Cytotoxicity		●		

Epigenetic Protein Products		Product Type		
Target	Application	Cell Line	eXpress Kit	Protein
BAZ2A	Target Engagement	●		
BRD2(1)	Target Engagement	●	●	
BRD2(1,2)	Target Engagement	●		
BRD3(1)	Target Engagement	●		
BRD3(1,2)	Target Engagement	●	●	
BRD4(1)	Target Engagement	●	●	
BRD4	Target Protein Degradation	●		
BRDT(1)	Target Engagement	●		
BRDT(1,2)	Target Engagement	●	●	
DNMT1	Screening & Activity Analysis			●
DOT1L	Screening & Activity Analysis			●
EHMT1	Screening & Activity Analysis			●
EP300	Target Engagement	●		
G9a	Target Engagement	●	●	
G9a	Screening & Activity Analysis			●
GLP	Target Engagement	●	●	
HAT1	Screening & Activity Analysis			●
hPHF8	Screening & Activity Analysis			●
JMJD2A	Screening & Activity Analysis			●
JMJD2B	Screening & Activity Analysis			●
JMJD2C	Screening & Activity Analysis			●
JMJD2D	Screening & Activity Analysis			●
JMJD2E	Screening & Activity Analysis			●
JMJD3	Screening & Activity Analysis			●
KAT2A (GCN5L2)	Screening & Activity Analysis			●
NSD1	Screening & Activity Analysis			●
PRMT1	Screening & Activity Analysis			●
PRMT3	Target Engagement	●		
PRMT3	Screening & Activity Analysis			●
SETD2	Screening & Activity Analysis			●
SETD7	Screening & Activity Analysis			●
SETD8	Screening & Activity Analysis			●
Sirtuin 1	Screening & Activity Analysis			●
sirtuin 2	Screening & Activity Analysis			●
SMARCA4	Target Engagement	●	●	
SUV39H2	Screening & Activity Analysis			●
SUV4-20H2	Screening & Activity Analysis			●
TAF1L(2)	Target Engagement	●		
WHSC1 (NSD2)	Screening & Activity Analysis			●

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		ADCYAP1R1	cAMP & Calcium	●			
ADCYAP1R1	Internalization	●		●			
ADCYAP1R1	Binding & GTPγS Function						●
ADCYAP1R1 (PAC1)	cAMP & Calcium	●		●			
ADCYAP1R1 (PAC1)	β-Arrestin	●		●			
mADCYAP1R1 (PAC1)	β-Arrestin	●		●			
ADORA1	cAMP & Calcium	●					
ADORA1	β-Arrestin	●					
ADORA1	Binding & GTPγS Function						●
ADORA2A	cAMP & Calcium	●				●	
ADORA2B	cAMP & Calcium	●		●			
mADORA2B	β-Arrestin	●		●			
ADORA3	cAMP & Calcium	●					
ADORA3	β-Arrestin	●		●			
ADORA3	Binding & GTPγS Function						●
mADORA3	β-Arrestin	●					
ADRA1A	cAMP & Calcium	●				●	
ADRA1A	Binding & GTPγS Function						●
ADRA1B	cAMP & Calcium	●					
ADRA1B	β-Arrestin	●		●			
ADRA1B	cAMP & Calcium					●	
ADRA1B	Binding & GTPγS Function						●
mADRA1B	β-Arrestin		●				
ADRA1D	β-Arrestin	●					
ADRA1D (82-572)	cAMP & Calcium	●				●	
ADRA1D (82-572)	Binding & GTPγS Function						●
ADRA2A	cAMP & Calcium	●		●		●	
ADRA2A	β-Arrestin	●		●			
ADRA2A	Binding & GTPγS Function						●
ADRA2B	cAMP & Calcium	●				●	
ADRA2B	β-Arrestin	●		●			
ADRA2B	Binding & GTPγS Function						●
ADRA2C	cAMP & Calcium	●		●			
ADRA2C	Internalization	●					
ADRA2C	β-Arrestin	●					
ADRB1	cAMP & Calcium	●		●			
ADRB1	β-Arrestin	●		●			
ADRB1	Binding & GTPγS Function						●
ADRB2 (B2AR)	cAMP & Calcium	●		●		●	
ADRB2 (B2AR)	Internalization	●					
ADRB2 (B2AR)	β-Arrestin	●		●			
ADRB2 (B2AR)	Binding & GTPγS Function						●
ADRB2(W158A)	Pharmacotraficking	●					
ADRB3	cAMP & Calcium	●		●		●	
ADRB3	Binding & GTPγS Function						●
AGTR1	β-Arrestin	●		●	●		
AGTR1	cAMP & Calcium	●				●	
AGTR1	Binding & GTPγS Function						●

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		AGTR2	Binding & GTPγS Function				
AGTRL1	β-Arrestin	●		●	●		
AGTRL1	cAMP & Calcium	●		●		●	
AGTRL1	Internalization	●		●			
AGTRL1	Binding & GTPγS Function						●
mAGTRL1	β-Arrestin	●		●			
AVPR1A	cAMP & Calcium	●				●	
AVPR1A	Internalization	●					
AVPR1A	β-Arrestin	●		●			
AVPR1A	Binding & GTPγS Function						●
AVPR1B	cAMP & Calcium	●				●	
AVPR1B	Internalization	●		●			
AVPR1B	β-Arrestin	●		●			
AVPR1B	Binding & GTPγS Function						●
AVPR2	cAMP & Calcium	●		●		●	
AVPR2	Internalization	●					
AVPR2	β-Arrestin	●		●			
AVPR2	Binding & GTPγS Function						●
mAVPR2	β-Arrestin		●				
AVPR2(S167T)	Pharmacotraficking	●					
BAI1	β-Arrestin	●					
BAI2	β-Arrestin	●		●			
BAI3	β-Arrestin	●		●			
BDKRB1	cAMP & Calcium	●					
BDKRB1	β-Arrestin	●		●			
BDKRB2	β-Arrestin	●		●	●		
BDKRB2	cAMP & Calcium	●				●	
BDKRB2	Internalization	●					
BDKRB2	Binding & GTPγS Function						●
BRS3	cAMP & Calcium	●				●	
BRS3	β-Arrestin	●		●			
BRS3	Binding & GTPγS Function						●
mBRS3	β-Arrestin		●				
C3AR1	cAMP & Calcium	●				●	
C3AR1	Internalization	●					
C3AR1	β-Arrestin	●		●			
C3AR1	Binding & GTPγS Function						●
C5AR1	β-Arrestin	●		●	●		
C5AR1	cAMP & Calcium	●		●		●	
C5AR1	Internalization	●		●			
C5AR1	Binding & GTPγS Function						●
mC5AR1	β-Arrestin	●		●			
C5L2	Internalization	●					
C5L2	β-Arrestin	●		●			
C5R1	cAMP & Calcium	●					
CALCR	cAMP & Calcium	●					
CALCR	β-Arrestin	●		●			
CALCRL-RAMP1(CGRP1)	cAMP & Calcium	●		●			



GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		CALCRL-RAMP1(CGRP1)	Binding & GTPγS Function				
CALCRL-RAMP1(CGRP1)	β-Arrestin	●		●			
CALCRL-RAMP2(AM1)	β-Arrestin	●		●			
CALCRL-RAMP3(AM2)	cAMP & Calcium	●		●			
CALCRL-RAMP3(AM2)	β-Arrestin	●		●			
CALCR-RAMP1	cAMP & Calcium					●	
CALCR-RAMP1	Binding & GTPγS Function						●
CALCR-RAMP2(AMY2)	β-Arrestin	●		●			
CALCR-RAMP3(AMY3)	cAMP & Calcium	●					
CALCR-RAMP3(AMY3)	β-Arrestin	●		●			
CASR	cAMP & Calcium	●				●	
CCKAR	cAMP & Calcium	●				●	
CCKAR	Internalization	●		●			
CCKAR	β-Arrestin	●		●			
CCKAR	Binding & GTPγS Function						●
mCCKAR	β-Arrestin		●				
CCKBR	cAMP & Calcium	●				●	
CCKBR	Internalization	●					
CCKBR	β-Arrestin	●		●			
CCKBR	Binding & GTPγS Function						●
CCR1	cAMP & Calcium	●				●	
CCR1	Internalization	●		●			
CCR1	β-Arrestin	●					
CCR1	Binding & GTPγS Function						●
mCCR1	β-Arrestin	●					
CCR2	β-Arrestin	●		●	●		
CCR2	cAMP & Calcium	●				●	
CCR2	Internalization	●		●			
CCR2	Binding & GTPγS Function						●
mCCR2	β-Arrestin	●		●			
CCR3	cAMP & Calcium	●				●	
CCR3	β-Arrestin	●		●			
mCCR3	β-Arrestin	●		●			
CCR4	cAMP & Calcium	●				●	
CCR4	Internalization	●		●			
CCR4	β-Arrestin	●		●			
mCCR4	β-Arrestin	●		●			
CCR5	cAMP & Calcium	●		●		●	
CCR5	Internalization	●		●			
CCR5	β-Arrestin	●		●			
CCR5	Binding & GTPγS Function						●
mCCR5	β-Arrestin	●		●			
CCR6	cAMP & Calcium	●		●		●	
CCR6	Internalization	●		●			
CCR6	β-Arrestin	●		●			
CCR6	Binding & GTPγS Function						●
mCCR6	β-Arrestin	●		●			
CCR7	cAMP & Calcium	●		●		●	

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		CCR7	Internalization	●			
CCR7	β-Arrestin	●		●			
mCCR7	β-Arrestin	●		●			
CCR8	cAMP & Calcium	●					
CCR8	β-Arrestin	●		●			
CCR8	cAMP & Calcium					●	
CCR8	Binding & GTPγS Function						●
mCCR8	β-Arrestin	●					
CCR9	cAMP & Calcium	●				●	
CCR9	β-Arrestin	●		●			
CCR9	Binding & GTPγS Function						●
mCCR9	β-Arrestin	●		●			
CCR10	cAMP & Calcium	●				●	
CCR10	Internalization	●		●			
CCR10	β-Arrestin	●		●			
CCRL1	β-Arrestin	●					
CCRL2	β-Arrestin	●		●			
cEDG5	β-Arrestin	●					
CHRM1	cAMP & Calcium	●				●	
CHRM1	Internalization	●		●			
CHRM1	β-Arrestin	●		●			
CHRM1	Binding & GTPγS Function						●
CHRM2	cAMP & Calcium	●		●		●	
CHRM2	Internalization	●		●			
CHRM2	β-Arrestin	●		●			
CHRM2	Binding & GTPγS Function						●
CHRM3	cAMP & Calcium	●				●	
CHRM3	Internalization	●		●			
CHRM3	β-Arrestin	●		●			
CHRM3	Binding & GTPγS Function						●
CHRM4	cAMP & Calcium	●				●	
CHRM4	Internalization	●		●			
CHRM4	β-Arrestin	●		●			
CHRM4	Binding & GTPγS Function						●
rCHRM4	β-Arrestin	●		●			
CHRM5	cAMP & Calcium	●				●	
CHRM5	Internalization	●		●			
CHRM5	β-Arrestin	●		●			
CHRM5	Binding & GTPγS Function						●
CMKLR1	cAMP & Calcium	●		●		●	
CMKLR1	Internalization	●		●			
CMKLR1	β-Arrestin	●		●			
mCMKLR1	β-Arrestin	●		●			
CNR1 (CB1)	β-Arrestin	●		●	●		
CNR1 (CB1)	cAMP & Calcium	●		●		●	
CNR1 (CB1)	Binding & GTPγS Function						●
mCNR1 (CB1)	β-Arrestin	●		●			
CNR2 (CB2)	cAMP & Calcium	●		●		●	

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		CNR2 (CB2)	$\beta$ -Arrestin	●			
CNR2 (CB2)	Binding & GTP $\gamma$ S Function						●
mCNR2 (CB2)	$\beta$ -Arrestin	●		●			
rCNR2 (CB2)	$\beta$ -Arrestin	●					
CRHR1	cAMP & Calcium	●		●		●	
CRHR1	Internalization	●					
CRHR1	$\beta$ -Arrestin	●		●			
CRHR1	Binding & GTP $\gamma$ S Function						●
mCRHR1	$\beta$ -Arrestin	●		●			
CRHR2	cAMP & Calcium	●		●		●	
CRHR2	Internalization	●					
CRHR2	$\beta$ -Arrestin	●		●			
CRHR2	Binding & GTP $\gamma$ S Function						●
CRTH2	cAMP & Calcium	●					
CRTH2	$\beta$ -Arrestin	●		●			
mCRTH2	$\beta$ -Arrestin	●					
rCRTH2	$\beta$ -Arrestin	●		●			
CX3CR1	cAMP & Calcium	●				●	
CX3CR1	$\beta$ -Arrestin	●		●			
CX3CR1	Binding & GTP $\gamma$ S Function						●
mCX3CR1	$\beta$ -Arrestin	●		●			
CXCR1	cAMP & Calcium	●				●	
CXCR1	Binding & GTP $\gamma$ S Function						●
CXCR1 (IL8RA)	cAMP & Calcium	●		●			
CXCR1 (IL8RA)	Internalization	●					
CXCR1 (IL8RA)	$\beta$ -Arrestin	●		●			
CXCR2 (IL8RB)	cAMP & Calcium	●		●		●	
CXCR2 (IL8RB)	Internalization	●		●			
CXCR2 (IL8RB)	Binding & GTP $\gamma$ S Function						●
CXCR2 (IL8RB)	$\beta$ -Arrestin	●		●			
mCXCR2 (IL8RB)	$\beta$ -Arrestin	●		●			
CXCR3	cAMP & Calcium	●		●		●	
CXCR3	Internalization	●		●			
CXCR3	$\beta$ -Arrestin	●		●			
CXCR3	Binding & GTP $\gamma$ S Function						●
mCXCR3	$\beta$ -Arrestin	●		●			
CXCR4	cAMP & Calcium	●		●		●	
CXCR4	Internalization	●					
CXCR4	$\beta$ -Arrestin	●					
CXCR4	Binding & GTP $\gamma$ S Function						●
mCXCR4	$\beta$ -Arrestin	●		●			
CXCR5 (BLR1)	$\beta$ -Arrestin	●		●	●		
CXCR5 (BLR1)	cAMP & Calcium	●		●		●	
mCXCR5 (BLR1)	$\beta$ -Arrestin	●		●			
CXCR6	cAMP & Calcium	●				●	
CXCR6	Internalization	●		●			
CXCR6	$\beta$ -Arrestin	●		●			
CXCR6	Binding & GTP $\gamma$ S Function						●

GPCR Products	Application	Product Type					
		Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
mCXCR6	β-Arrestin	●		●			
CXCR7 (CMKOR1)	Internalization	●		●			
CXCR7 (CMKOR1)	Binding & GTPγS Function						●
CXCR7 (CMKOR1)	β-Arrestin	●		●			
mCXCR7 (CMKOR1)	β-Arrestin	●		●			
cyno CXCR1	β-Arrestin	●		●			
cyno CXCR2	β-Arrestin	●		●			
cyno GPR120	β-Arrestin	●					
CysLT1	cAMP & Calcium	●				●	
CYSLTR2	cAMP & Calcium	●				●	
DARC	β-Arrestin	●		●			
DRD1	cAMP & Calcium	●		●		●	
DRD1	Internalization	●					
DRD1	β-Arrestin	●		●			
DRD1	Binding & GTPγS Function						●
rDRD1	β-Arrestin			●			
rDRD1	cAMP & Calcium	●		●			
rDRD1	β-Arrestin	●					
DRD2	cAMP & Calcium	●				●	
DRD2	Binding & GTPγS Function						●
DRD2L	cAMP & Calcium	●					
DRD2L	β-Arrestin	●		●			
rDRD2L	cAMP & Calcium	●		●			
rDRD2L	β-Arrestin	●					
DRD2S	cAMP & Calcium	●					
DRD2S	β-Arrestin	●		●			
DRD2S	cAMP & Calcium			●			
rDRD2S	cAMP & Calcium	●					
rDRD2S	β-Arrestin	●					
DRD3	β-Arrestin	●		●			
DRD3	Binding & GTPγS Function						●
DRD4	cAMP & Calcium	●				●	
DRD4	β-Arrestin	●		●			
DRD5	cAMP & Calcium	●		●		●	
DRD5	β-Arrestin	●		●			
DRD5	Binding & GTPγS Function						●
mDRD5	β-Arrestin	●					
EBI2	cAMP & Calcium	●					
EBI2	Internalization	●					
EBI2	β-Arrestin	●		●			
mEBI2	β-Arrestin	●		●			
EDG1	cAMP & Calcium	●					
EDG1	Internalization	●		●			
EDG1	β-Arrestin	●		●			
mEDG1	β-Arrestin	●		●			
EDG2	cAMP & Calcium	●					
EDG2	Internalization	●					
EDG2	β-Arrestin	●					

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		EDG3	cAMP & Calcium	●			
EDG3	Internalization	●					
EDG3	β-Arrestin	●		●			
mEDG3	β-Arrestin	●					
EDG4	cAMP & Calcium	●					
EDG4	β-Arrestin	●		●			
EDG5	Internalization	●					
EDG5	β-Arrestin	●		●			
mEDG5	β-Arrestin	●					
rEDG5	β-Arrestin	●					
EDG6	β-Arrestin	●					
EDG7	Internalization	●					
EDG7	β-Arrestin	●		●			
EDNRA	cAMP & Calcium	●				●	
EDNRA	β-Arrestin	●		●			
EDNRA	Binding & GTPγS Function						●
mEDNRA	β-Arrestin	●		●			
EDNRB	cAMP & Calcium	●				●	
EDNRB	Internalization	●		●			
EDNRB	β-Arrestin	●		●			
EDNRB	Binding & GTPγS Function						●
mEDNRB	β-Arrestin	●		●			
F2R (PAR1)	β-Arrestin	●		●			
F2RL1	cAMP & Calcium	●				●	
F2RL1	Internalization	●		●			
F2RL1	Binding & GTPγS Function						●
F2RL1 (PAR2, GPR11)	β-Arrestin	●		●			
F2RL3 (PAR4)	Internalization	●					
F2RL3 (PAR4)	β-Arrestin	●		●			
FFAR1 (GPR40)	β-Arrestin	●					
FFAR1 (GPR40)	cAMP & Calcium	●				●	
mFFAR1 (GPR40)	β-Arrestin	●					
rFFAR1 (GPR40)	β-Arrestin	●					
FFAR2 (GPR43)	cAMP & Calcium	●				●	
FFAR2 (GPR43)	β-Arrestin	●					
FFAR3 (GPR41)	cAMP & Calcium	●		●		●	
FFAR4	cAMP & Calcium	●				●	
FPR1	cAMP & Calcium	●					
FPR1	Internalization	●		●			
FPR1	β-Arrestin	●		●			
mFPR1	β-Arrestin	●					
FPR2	cAMP & Calcium	●				●	
FPR2	Binding & GTPγS Function						●
mFPR2	β-Arrestin	●		●			
FPR3	cAMP & Calcium	●		●			
FPRL1	cAMP & Calcium	●		●			
FPRL1	Internalization	●		●			
FPRL1	β-Arrestin	●		●			

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		FSHR	cAMP & Calcium	●			●
FSHR	β-Arrestin	●					
GABBR1 - GABBR2	cAMP & Calcium	●				●	
GABBR1 - GABBR2	Binding & GTPγS Function						●
GALR1	cAMP & Calcium	●		●		●	
GALR1	Internalization	●		●			
GALR1	β-Arrestin	●		●			
GALR1	Binding & GTPγS Function						●
mGALR1	β-Arrestin		●				
GALR2	cAMP & Calcium	●				●	
GALR2	β-Arrestin	●		●			
GALR2	Binding & GTPγS Function						●
mGALR2	β-Arrestin	●		●			
GCGR	cAMP & Calcium	●		●		●	
GCGR	Internalization	●		●			
GCGR	β-Arrestin	●		●			
GCGR	Binding & GTPγS Function						●
mGCGR	β-Arrestin	●		●			
GHRHR	cAMP & Calcium	●		●			
GHSR	β-Arrestin				●		
GHSR	cAMP & Calcium	●				●	
GHSR	Binding & GTPγS Function						●
GHSR1a	β-Arrestin	●		●			
mGHSR1a	β-Arrestin	●		●			
GHSR1b	β-Arrestin	●		●			
GIPR	cAMP & Calcium	●		●		●	
GIPR	β-Arrestin	●		●			
GIPR	Binding & GTPγS Function						●
mGIPR	cAMP & Calcium	●		●			
mGIPR	β-Arrestin	●		●			
GLP1R	cAMP & Calcium	●		●	●	●	
GLP1R	Internalization	●		●			
GLP1R	β-Arrestin	●		●	●		
GLP1R	Binding & GTPγS Function						●
mGLP1R	cAMP & Calcium	●		●			
mGLP1R	β-Arrestin	●		●			
GLP2R	β-Arrestin	●		●	●		
GLP2R	cAMP & Calcium	●		●		●	
GLP2R	Internalization	●					
GNRHR	cAMP & Calcium	●				●	
GNRHR	Binding & GTPγS Function						●
GPBAR1	cAMP & Calcium	●		●	●	●	
GPBAR1	β-Arrestin	●					
GPR1	Internalization	●					
GPR1	β-Arrestin	●		●			
GPR3	β-Arrestin	●		●			
GPR4	β-Arrestin	●		●			
GPR6	β-Arrestin	●		●			

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		GPR12	β-Arrestin	●		●	
GPR15	β-Arrestin	●		●			
GPR17	cAMP & Calcium	●					
GPR17	β-Arrestin	●					
GPR18	β-Arrestin	●		●			
GPR20	β-Arrestin	●		●			
GPR22	β-Arrestin	●					
GPR23	β-Arrestin	●		●			
GPR25	β-Arrestin	●		●			
GPR26	β-Arrestin	●		●			
GPR27	β-Arrestin	●		●			
GPR30	β-Arrestin	●		●			
GPR31	β-Arrestin	●		●			
GPR32	β-Arrestin	●		●			
GPR34	β-Arrestin	●					
GPR35	cAMP & Calcium	●		●			
GPR35	Internalization	●		●			
GPR35	β-Arrestin	●					
rGPR35	β-Arrestin	●		●			
GPR37 (EDNRBL)	β-Arrestin	●		●			
GPR37L1	β-Arrestin	●		●			
GPR39	cAMP & Calcium	●					
GPR39	β-Arrestin	●		●			
GPR45	β-Arrestin	●		●			
GPR50	β-Arrestin	●		●			
GPR52	β-Arrestin	●		●			
GPR55	β-Arrestin	●		●			
GPR61	β-Arrestin	●		●			
GPR65	β-Arrestin	●		●			
GPR68	cAMP & Calcium	●				●	
GPR75	β-Arrestin	●		●			
GPR78	β-Arrestin	●		●			
GPR79	β-Arrestin	●		●			
GPR81	cAMP & Calcium	●					
GPR83	β-Arrestin	●		●			
GPR84	cAMP & Calcium	●					
GPR84	β-Arrestin	●					
mGPR84	β-Arrestin	●					
GPR85	β-Arrestin	●		●			
GPR88	β-Arrestin	●		●			
mGPR91 (SUCNR1)	β-Arrestin		●				
GPR92	β-Arrestin	●		●			
GPR97	β-Arrestin	●		●			
GPR101	β-Arrestin	●					
GPR103	β-Arrestin	●					
GPR107	β-Arrestin	●		●			
GPR109A (HM74A)	cAMP & Calcium	●		●			
GPR109A (HM74A)	β-Arrestin	●		●			

GPCR Products	Application	Product Type					
		Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
GPR109B	cAMP & Calcium	●		●			
GPR109B	β-Arrestin	●		●			
GPR119	β-Arrestin	●					
GPR120	β-Arrestin			●			
rGPR120	β-Arrestin	●		●			
GPR120L	Internalization	●					
GPR120L	β-Arrestin	●					
GPR120S	β-Arrestin	●		●			
GPR123	β-Arrestin	●		●			
GPR132	β-Arrestin	●		●			
GPR135 (HUMNP11Y20)	β-Arrestin	●		●			
GPR137 (TM7SF1)	β-Arrestin	●		●			
GPR139	β-Arrestin	●					
GPR141	β-Arrestin	●		●			
GPR142	β-Arrestin	●		●			
GPR143	β-Arrestin	●		●			
GPR146	β-Arrestin	●		●			
GPR148	β-Arrestin	●		●			
GPR149	β-Arrestin	●		●			
GPR150 (AXOR70)	β-Arrestin	●		●			
GPR151	β-Arrestin	●					
GPR152	β-Arrestin	●					
GPR157	β-Arrestin	●		●			
GPR161	β-Arrestin	●					
GPR162	β-Arrestin	●		●			
GPR171	β-Arrestin	●		●			
GPR173	β-Arrestin	●		●			
GPR176	β-Arrestin	●		●			
GPR182	β-Arrestin	●		●			
GRM1	cAMP & Calcium	●					
GRM2	cAMP & Calcium	●		●		●	
GRM2	Binding & GTPγS Function						●
GRM3	cAMP & Calcium	●					
GRM4	cAMP & Calcium	●		●			
GRM5	cAMP & Calcium	●					
GRM6	cAMP & Calcium	●		●			
GRPR	cAMP & Calcium	●				●	
GRPR	Internalization	●		●			
GRPR	β-Arrestin	●		●			
GRPR	Binding & GTPγS Function						●
HCR2	cAMP & Calcium	●				●	
HCRTR1	cAMP & Calcium	●				●	
HCRTR1	Internalization	●		●			
HCRTR1	β-Arrestin	●		●			
mHCRTR1	β-Arrestin		●				
HCRTR2	cAMP & Calcium	●				●	
HCRTR2	Internalization	●		●			
HCRTR2	β-Arrestin	●		●			



GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		HRH1	$\beta$ -Arrestin	●		●	
HRH1	Binding & GTP $\gamma$ S Function						●
HRH2	cAMP & Calcium	●		●			
HRH2	$\beta$ -Arrestin	●		●			
HRH2	Binding & GTP $\gamma$ S Function						●
mHTR2A	$\beta$ -Arrestin	●					
mHTR2C	$\beta$ -Arrestin	●					
HRH3	cAMP & Calcium	●		●			
HRH3	$\beta$ -Arrestin	●		●			
HRH4	$\beta$ -Arrestin	●		●			
HTR1A	cAMP & Calcium	●					
HTR1A	$\beta$ -Arrestin	●		●			
HTR1B	cAMP & Calcium	●					
HTR1B	$\beta$ -Arrestin	●		●			
HTR1B	Binding & GTP $\gamma$ S Function						●
HTR1E	$\beta$ -Arrestin	●					
HTR1F	cAMP & Calcium	●					
HTR1F	$\beta$ -Arrestin	●		●			
HTR2A	cAMP & Calcium	●				●	
HTR2A	Internalization	●					
HTR2A	$\beta$ -Arrestin	●		●			
HTR2B	cAMP & Calcium	●				●	
HTR2C	cAMP & Calcium	●				●	
HTR2C	Internalization	●		●			
HTR2C	$\beta$ -Arrestin	●		●			
HTR2C	Binding & GTP $\gamma$ S Function						●
rHTR2C	$\beta$ -Arrestin	●		●			
HTR2C (VGV)	$\beta$ -Arrestin	●		●			
HTR2C (VNV)	$\beta$ -Arrestin	●		●			
HTR2C (VSV)	$\beta$ -Arrestin	●		●			
HTR4B	cAMP & Calcium	●				●	
HTR4B	Binding & GTP $\gamma$ S Function						●
HTR5A	cAMP & Calcium	●					
HTR5A	$\beta$ -Arrestin	●		●			
HTR6	cAMP & Calcium	●				●	
HTR6	$\beta$ -Arrestin	●					
HTR6	Binding & GTP $\gamma$ S Function						●
HTR7A	cAMP & Calcium	●					
HTR7B	cAMP & Calcium	●					
HTR7D	cAMP & Calcium	●					
HTS059RTA	cAMP & Calcium					●	
KISS1R	cAMP & Calcium	●				●	
KISS1R	$\beta$ -Arrestin	●		●			
LGR4 (GPR48)	$\beta$ -Arrestin	●		●			
LGR5 (GPR49)	$\beta$ -Arrestin	●		●			
LGR6	$\beta$ -Arrestin	●		●			
LHCGR	cAMP & Calcium	●		●	●	●	
LHCGR	$\beta$ -Arrestin	●		●			

GPCR Products		Product Type						
		Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell
LPAR1	cAMP & Calcium	●					●	
LPAR1	Binding & GTPγS Function							●
LPAR2	cAMP & Calcium	●					●	
LPAR3	cAMP & Calcium	●					●	
LPAR5	cAMP & Calcium	●					●	
LTB4R	cAMP & Calcium	●		●			●	
LTB4R	Internalization	●						
LTB4R	β-Arrestin	●		●				
LTB4R	Binding & GTPγS Function							●
mLTB4R1	β-Arrestin	●						
MC1R	cAMP & Calcium	●						
MC1R	β-Arrestin	●		●				
MC1R	Binding & GTPγS Function							●
MC2R	cAMP & Calcium	●					●	
MC3R	cAMP & Calcium	●		●				
MC3R	Internalization	●		●				
MC3R	β-Arrestin	●		●				
MC3R	Binding & GTPγS Function							●
mMC3R	β-Arrestin		●					
MC4R	β-Arrestin	●		●	●			
MC4R	cAMP & Calcium	●		●				
mMC4R	β-Arrestin		●					
MC4R(T162I)	Pharmacotraficking	●						
MC5R	cAMP & Calcium	●					●	
MC5R	β-Arrestin	●		●				
MC5R	Binding & GTPγS Function							●
mMC5R	β-Arrestin		●					
MCHR1	β-Arrestin	●		●	●			
MCHR1	cAMP & Calcium	●		●			●	
MCHR1	Internalization	●		●				
MCHR1	Binding & GTPγS Function							●
mMCHR1	β-Arrestin	●						
MCHR2	cAMP & Calcium	●						
MCHR2	β-Arrestin	●		●				
MCHR2	cAMP & Calcium						●	
MLNR	cAMP & Calcium	●					●	
MLNR	Internalization	●						
MLNR	β-Arrestin	●		●				
MLNR	Binding & GTPγS Function							●
MRGPRD	cAMP & Calcium	●					●	
MRGPRD	β-Arrestin	●		●				
MRGPRD	Binding & GTPγS Function							●
MRGPRE	β-Arrestin	●		●				
MRGPRF	β-Arrestin	●		●				
MRGPRX1	cAMP & Calcium	●					●	
MRGPRX1	β-Arrestin	●						
MRGPRX2	β-Arrestin	●		●	●			
MRGPRX2	cAMP & Calcium	●					●	

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		cMRGPRX2	β-Arrestin	●			
MRGPRX4	β-Arrestin	●		●			
MTNR1A	cAMP & Calcium	●		●			
MTNR1A	Internalization	●					
MTNR1A	β-Arrestin	●		●			
MTNR1A	Binding & GTPγS Function						●
MTNR1B	cAMP & Calcium	●		●			
MTNR1B	Internalization	●					
MTNR1B	β-Arrestin	●		●			
NMBR	cAMP & Calcium	●				●	
NMBR	Internalization	●		●			
NMBR	β-Arrestin	●		●			
NMBR	Binding & GTPγS Function						●
NMU1R	cAMP & Calcium	●				●	
NMU1R	Internalization	●					
NMU1R	β-Arrestin	●		●			
NMU1R	Binding & GTPγS Function						●
NMU2	cAMP & Calcium	●				●	
NPBWR1	cAMP & Calcium	●				●	
NPBWR1	β-Arrestin	●		●			
NPBWR1	Binding & GTPγS Function						●
NPBWR2	cAMP & Calcium	●					
NPBWR2	β-Arrestin	●		●			
NPBWR2	Binding & GTPγS Function						●
NPFFR1	cAMP & Calcium	●		●			
NPFFR1	β-Arrestin	●					
NPSR1	β-Arrestin	●					
NPSR1b	β-Arrestin			●			
NPY1R	β-Arrestin	●		●			
mNPY1R	β-Arrestin	●					
NPY2R	cAMP & Calcium	●		●	●	●	
NPY2R	β-Arrestin	●		●	●		
NPY2R	Internalization	●					
NPY2R	Binding & GTPγS Function						●
mNPY2R	β-Arrestin	●					
NPY4R	cAMP & Calcium	●					
NPY4R	Binding & GTPγS Function						●
NTSR1	cAMP & Calcium	●				●	
NTSR1	Internalization	●		●			
NTSR1	β-Arrestin	●		●			
NTSR1	Binding & GTPγS Function						●
OPN5	β-Arrestin	●		●			
OPRD1	cAMP & Calcium	●		●		●	
OPRD1	Internalization	●		●			
OPRD1	β-Arrestin	●		●			
OPRD1	Binding & GTPγS Function						●
mOPRD1	β-Arrestin	●		●			
OPRK1	cAMP & Calcium	●		●		●	

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		OPRK1	Internalization	●		●	
OPRK1	β-Arrestin	●		●			
OPRK1	Binding & GTPγS Function						●
mOPRK1	β-Arrestin	●		●			
OPRL1	cAMP & Calcium	●		●		●	
OPRL1	Internalization	●					
OPRL1	β-Arrestin	●		●			
OPRL1	Binding & GTPγS Function						●
OPRM1	cAMP & Calcium	●		●		●	
OPRM1	Internalization	●		●			
OPRM1	β-Arrestin	●		●			
OPRM1	Binding & GTPγS Function						●
mOPRM1	β-Arrestin	●		●			
rOPRM1	β-Arrestin	●		●			
OXER1	cAMP & Calcium	●		●			
OXER1	Internalization	●					
OXER1	β-Arrestin	●					
OXGR1 (GPR80, GPR99)	cAMP & Calcium	●				●	
OXGR1 (GPR80, GPR99)	β-Arrestin	●		●			
OXTR	cAMP & Calcium	●				●	
OXTR	Internalization	●		●			
OXTR	β-Arrestin	●		●			
OXTR	Binding & GTPγS Function						●
mOXTR	β-Arrestin	●					
P2RY1	cAMP & Calcium	●				●	
P2RY1	Internalization	●		●			
P2RY1	β-Arrestin	●		●			
P2RY2	cAMP & Calcium	●				●	
P2RY2	Internalization	●		●			
P2RY2	β-Arrestin	●		●			
P2RY4	cAMP & Calcium	●				●	
P2RY4	Internalization	●		●			
P2RY4	β-Arrestin	●		●			
P2RY6	β-Arrestin	●		●			
mP2RY6	β-Arrestin	●					
P2RY8	β-Arrestin	●		●			
P2RY11	cAMP & Calcium	●					
P2RY11	β-Arrestin	●		●			
P2RY11	cAMP & Calcium					●	
P2RY12	Internalization	●					
P2RY12	β-Arrestin	●		●			
mP2RY12	β-Arrestin	●		●			
PPYR1 (NPY4R)	cAMP & Calcium	●		●			
PPYR1 (NPY4R)	β-Arrestin	●		●			
mPPYR1 (NPY4R)	β-Arrestin	●					
PRLHR	cAMP & Calcium	●				●	
PRLHR	Internalization	●					
PRLHR	β-Arrestin	●		●			

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		PRLHR	Binding & GTPγS Function				
PROKR1	cAMP & Calcium	●		●		●	
PROKR1	Internalization	●		●			
PROKR1	β-Arrestin	●		●			
PROKR1	Binding & GTPγS Function						●
rPROKR1	β-Arrestin	●		●			
PROKR2	cAMP & Calcium	●				●	
PROKR2	Internalization	●					
PROKR2	β-Arrestin	●		●			
PROKR2	Binding & GTPγS Function						●
rPROKR2	β-Arrestin	●		●			
PTAFR	cAMP & Calcium	●				●	
PTAFR	Internalization	●					
PTAFR	β-Arrestin	●		●			
mPTAFR	β-Arrestin	●		●			
PTGDR	cAMP & Calcium	●		●		●	
PTGDR	Binding & GTPγS Function						●
PTGDR2	Binding & GTPγS Function						●
PTGER1	cAMP & Calcium	●				●	
PTGER1	β-Arrestin	●					
PTGER1	Binding & GTPγS Function						●
PTGER2	cAMP & Calcium	●		●		●	
PTGER2	β-Arrestin	●		●			
PTGER2	Binding & GTPγS Function						●
PTGER3	cAMP & Calcium	●		●		●	
PTGER3	β-Arrestin	●		●			
PTGER3	Binding & GTPγS Function						●
PTGER4	cAMP & Calcium	●				●	
PTGER4	Internalization	●					
PTGER4	β-Arrestin	●		●			
PTGER4	Binding & GTPγS Function						●
mPTGER4	β-Arrestin	●		●			
rPTGER4	β-Arrestin	●					
PTGFR	cAMP & Calcium	●				●	
PTGFR	Internalization	●		●			
PTGFR	β-Arrestin	●		●			
PTGIR	cAMP & Calcium	●		●		●	
PTGIR	Internalization	●		●			
mPTGIR	β-Arrestin		●				
PTGIR	β-Arrestin	●					
PTH1R	cAMP & Calcium	●				●	
PTH1R	Binding & GTPγS Function						●
PTH2R	cAMP & Calcium	●				●	
PTHR	cAMP & Calcium				●		
PTHR	β-Arrestin				●		
PTHR1	cAMP & Calcium	●		●			
PTHR1	Internalization	●		●			
PTHR1	β-Arrestin	●		●			

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		PTHR2	cAMP & Calcium	●		●	
PTHR2	Internalization	●		●			
PTHR2	β-Arrestin	●		●			
QRFR	cAMP & Calcium	●				●	
mRHO(P23H)	Pharmacotraficking	●					
RXFP1 (LGR7)	cAMP & Calcium	●		●			
mRXFP1	cAMP & Calcium	●					
RXFP2 (LGR8)	cAMP & Calcium	●					
RXFP3 (SALPR)	Internalization	●					
RXFP3 (SALPR)	β-Arrestin	●		●			
RXFP3 (SALPR)	cAMP & Calcium	●		●			
mRXFP3	β-Arrestin	●		●			
RXFP4	cAMP & Calcium	●					
RXFP4	β-Arrestin	●		●			
mRXFP4	β-Arrestin	●		●			
S1PR1	cAMP & Calcium	●				●	
S1PR1	Binding & GTPγS Function						●
S1PR2	cAMP & Calcium	●				●	
S1PR2	Binding & GTPγS Function						●
S1PR3	cAMP & Calcium	●				●	
S1PR3	Binding & GTPγS Function						●
S1PR4	cAMP & Calcium	●				●	
S1PR5	cAMP & Calcium	●				●	
S1PR5	Binding & GTPγS Function						●
SCTR	cAMP & Calcium	●		●		●	
SCTR	Internalization	●					
SCTR	β-Arrestin	●		●			
sEDG5	β-Arrestin	●					
SMO(W535L)	Pharmacotraficking	●					
SSTR1	β-Arrestin	●		●			
SSTR1	Binding & GTPγS Function						●
SSTR2	β-Arrestin	●		●	●	●	
SSTR2	cAMP & Calcium	●		●			
SSTR2	Internalization	●		●			
SSTR2	Binding & GTPγS Function						●
mSSTR2	β-Arrestin	●		●			
SSTR3	cAMP & Calcium	●		●		●	
SSTR3	β-Arrestin	●		●			
SSTR3	Binding & GTPγS Function						●
SSTR4	cAMP & Calcium	●		●		●	
SSTR4	β-Arrestin	●		●			
SSTR4	Binding & GTPγS Function						●
SSTR5	cAMP & Calcium	●				●	
SSTR5	β-Arrestin	●		●			
SSTR5	Binding & GTPγS Function						●
mSSTR5	β-Arrestin	●					
SUCNR1	cAMP & Calcium	●				●	
TAAR1	cAMP & Calcium	●					

GPCR Products		Product Type					
Target	Application	Cell Line	Cell Pool	eXpress Kit	Bioassay Kit	Frozen Cell	Membrane Prep
		mTAAR1	cAMP & Calcium	●			
rTAAR1	cAMP & Calcium	●					
TAAR5	β-Arrestin	●		●			
TACR1	cAMP & Calcium	●				●	
TACR1	Internalization	●					
TACR1	β-Arrestin	●		●			
TACR1	Binding & GTPγS Function						●
mTACR1	β-Arrestin		●				
TACR2	Internalization	●		●			
TACR2	β-Arrestin	●		●			
TACR3	cAMP & Calcium	●		●		●	
TACR3	Internalization	●					
TACR3	β-Arrestin	●		●			
TBXA2R	cAMP & Calcium	●				●	
TBXA2R	Internalization	●					
TBXA2R	β-Arrestin	●		●			
TBXA2R	Binding & GTPγS Function						●
TRHR	cAMP & Calcium	●				●	
TRHR	Internalization	●		●			
TRHR	β-Arrestin	●		●			
TSHR(L)	cAMP & Calcium	●				●	
TSHR(L)	β-Arrestin	●					
TSHRL	cAMP & Calcium	●					
UTR2	Internalization	●		●			
UTR2	β-Arrestin	●		●			
mUTR2	β-Arrestin	●		●			
UTS2R	Binding & GTPγS Function						●
VIPR1	cAMP & Calcium	●		●		●	
VIPR1	Internalization	●		●			
VIPR1	β-Arrestin	●		●			
VIPR1	Binding & GTPγS Function						●
mVIPR1	β-Arrestin	●		●			
rVIPR1	β-Arrestin	●		●			
VIPR2	β-Arrestin	●		●	●		
VIPR2	cAMP & Calcium	●		●		●	
VIPR2	Internalization	●		●			
VIPR2	Binding & GTPγS Function						●
XCR1	cAMP & Calcium	●				●	

Ion Channel Products		Product Type	
Target	Application	Cell Line	Membrane Prep
ASIC3	Electrophysiological studies	●	
Cav2.2	Electrophysiological studies	●	
Cav3.2	Electrophysiological studies	●	
CFTR	Electrophysiological studies	●	
CFTR-ΔF508	Pharmacotraficking	●	
GABAA α1/β3/γ2	Electrophysiological studies	●	
GABAA α3/β3/γ2	Electrophysiological studies	●	
GABAA α4/β3/γ2	Electrophysiological studies	●	
GABAA α5/β3/γ2	Electrophysiological studies	●	
GABAA α6/β3/γ2	Electrophysiological studies	●	
GluR6	Electrophysiological studies	●	
GlyRA1	Electrophysiological studies	●	
HCN1	Electrophysiological studies	●	
HCN2	Electrophysiological studies	●	
HCN3	Electrophysiological studies	●	
HCN4	Electrophysiological studies	●	
hERG	Electrophysiological studies	●	
hERG	Pharmacotraficking	●	
hERG	Binding Studies		●
KCNH2(G601S)	Pharmacotraficking	●	
KCNQ1/hminK	Electrophysiological studies	●	
Kir2.1	Electrophysiological studies	●	
Kir6.2/SUR2A	Electrophysiological studies	●	
Kv1.1	Electrophysiological studies	●	
Kv1.2	Electrophysiological studies	●	
Kv1.3	Electrophysiological studies	●	
Kv1.4	Electrophysiological studies	●	
Kv1.5	Electrophysiological studies	●	
Kv1.6	Electrophysiological studies	●	
Kv1.7	Electrophysiological studies	●	
Kv1.8	Electrophysiological studies	●	
Kv12.2	Electrophysiological studies	●	
Kv2.1	Electrophysiological studies	●	
Kv2.1/Kv9.2	Electrophysiological studies	●	
Kv3.1	Electrophysiological studies	●	
Kv3.2	Electrophysiological studies	●	
Kv3.3	Electrophysiological studies	●	
Kv4.2/KChIP2	Electrophysiological studies	●	
Kv4.3/KChIP1	Electrophysiological studies	●	
Kv4.3/KChIP2	Electrophysiological studies	●	
Kv7.2/Kv7.3	Electrophysiological studies	●	
Kv7.3/Kv7.5	Electrophysiological studies	●	
Kv7.4	Electrophysiological studies	●	
Kv7.4/Kv7.5	Electrophysiological studies	●	
nAChR α1/β1/δ/ε	Electrophysiological studies	●	
nAChR α3/β4	Electrophysiological studies	●	
nAChR α4/β2	Electrophysiological studies	●	
nAChR α7/ric3	Electrophysiological studies	●	
Nav1.1	Electrophysiological studies	●	



Ion Channel Products		Product Type	
Target	Application	Cell Line	Membrane Prep
Nav1.2	Electrophysiological studies	●	
Nav1.3	Electrophysiological studies	●	
Nav1.4	Electrophysiological studies	●	
Nav1.5	Electrophysiological studies	●	
Nav1.6	Electrophysiological studies	●	
Nav1.7	Electrophysiological studies	●	
Nav1.8/ $\beta$ 1	Electrophysiological studies	●	
rNav1.8	Electrophysiological studies	●	
TRPA1	Electrophysiological studies	●	
TRPV1	Electrophysiological studies	●	
TRPV3	Electrophysiological studies	●	
TRPV4	Electrophysiological studies	●	

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	AAK1	Screening & Activity Analysis				●
Kinase	Abl	Screening & Activity Analysis				●
Kinase	ACK1	Screening & Activity Analysis				●
Kinase	ACTR2	Screening & Activity Analysis				●
Kinase	ActRIIB	Screening & Activity Analysis				●
Kinase	Akt1/PKB $\alpha$	Screening & Activity Analysis				●
Kinase	Akt2/PKB $\beta$	Screening & Activity Analysis				●
Kinase	Akt3/PKB $\gamma$	Screening & Activity Analysis				●
Kinase	ALK	Screening & Activity Analysis				●
Kinase	ALK1	Screening & Activity Analysis				●
Kinase	ALK2	Screening & Activity Analysis				●
Kinase	ALK4	Screening & Activity Analysis				●
Kinase	ALK6	Screening & Activity Analysis				●
Kinase	AMPK ( $\alpha$ 1, $\beta$ 1, $\gamma$ 1)	Screening & Activity Analysis				●
Kinase	A-Raf	Screening & Activity Analysis				●
Kinase	Arg	Screening & Activity Analysis				●
Kinase	Ask1	Screening & Activity Analysis				●
Kinase	ATM	Screening & Activity Analysis				●
Kinase	ATR/ATRIP	Screening & Activity Analysis				●
Kinase	Aurora A	Screening & Activity Analysis				●
Kinase	Aurora B	Screening & Activity Analysis				●
Kinase	Aurora C	Screening & Activity Analysis				●
Kinase	Axl	Screening & Activity Analysis				●
Kinase RTK	Axl	Functional	●			
Kinase	BIKe	Screening & Activity Analysis				●
Kinase	Blk	Target Engagement	●	●		
Kinase CTK	Blk	Activity	●	●		
Kinase	Blk	Screening & Activity Analysis				●
Kinase	BMPR2	Screening & Activity Analysis				●
Kinase	Bmx	Screening & Activity Analysis				●
Kinase	B-Raf	Screening & Activity Analysis				●
Kinase	BRK	Screening & Activity Analysis				●
Kinase	BrSK1	Screening & Activity Analysis				●
Kinase	BrSK2	Screening & Activity Analysis				●
Kinase	BTK	Target Engagement	●	●		
Kinase	BTK	Target Protein Degradation	●			
Kinase	BTK	Screening & Activity Analysis				●
Kinase	CaM	Screening & Activity Analysis				●
Kinase	CaMKI	Screening & Activity Analysis				●
Kinase	CaMKI $\gamma$	Screening & Activity Analysis				●
Kinase	CaMKI $\delta$	Screening & Activity Analysis				●
Kinase	CaMKII $\alpha$	Screening & Activity Analysis				●
Kinase	CaMKII $\beta$	Screening & Activity Analysis				●
Kinase	CaMKII $\delta$	Screening & Activity Analysis				●
Kinase	CaMKII $\gamma$	Screening & Activity Analysis				●
Kinase	CaMKIV	Screening & Activity Analysis				●
Kinase	CaMKK1	Screening & Activity Analysis				●
Kinase	CaMKK2	Screening & Activity Analysis				●
Phosphatase	CD45	Screening & Activity Analysis				●

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	Cdc7/cyclin B1	Screening & Activity Analysis				●
Kinase	Cdc25c	Screening & Activity Analysis				●
Kinase	CDK1/cyclin B	Screening & Activity Analysis				●
Kinase	CDK2/cyclin A	Screening & Activity Analysis				●
Kinase	CDK2/cyclin E	Screening & Activity Analysis				●
Kinase	CDK3/cyclin E	Screening & Activity Analysis				●
Kinase	CDK4/cyclin D3	Screening & Activity Analysis				●
Kinase	CDK5/p25	Screening & Activity Analysis				●
Kinase	CDK5/p35	Screening & Activity Analysis				●
Kinase	CDK6/cyclin D3	Screening & Activity Analysis				●
Kinase	CDK7/cyclin H/MAT1	Screening & Activity Analysis				●
Kinase	CDK9/cyclin T1	Screening & Activity Analysis				●
Kinase	CDK14/cyclin Y	Screening & Activity Analysis				●
Kinase	CDK16	Screening & Activity Analysis				●
Kinase	CDK17	Screening & Activity Analysis				●
Kinase	CDK18	Screening & Activity Analysis				●
Kinase	CDKL1	Screening & Activity Analysis				●
Kinase	CDKL2	Screening & Activity Analysis				●
Kinase	CDKL3	Screening & Activity Analysis				●
Kinase	CDKL4	Screening & Activity Analysis				●
Kinase	CDKN1A (p21)	Target Protein Degradation	●			
Kinase	ChaK1	Screening & Activity Analysis				●
Kinase	CHK1	Screening & Activity Analysis				●
Kinase	CHK2	Screening & Activity Analysis				●
Kinase RTK	cINSRa	Functional	●	●		
Kinase RTK	cINSRb	Functional	●	●		
Kinase	CK1α	Screening & Activity Analysis				●
Kinase	CK1ε	Screening & Activity Analysis				●
Kinase	CK1γ1	Screening & Activity Analysis				●
Kinase	CK1γ2	Screening & Activity Analysis				●
Kinase	CK1γ3	Screening & Activity Analysis				●
Kinase	CK1γδ	Screening & Activity Analysis				●
Kinase	CK2	Screening & Activity Analysis				●
Kinase	CK2α	Screening & Activity Analysis				●
Kinase	CK2α2	Screening & Activity Analysis				●
Kinase	c-Kit	Screening & Activity Analysis				●
Kinase RTK	c-Kit	Functional	●	●		
Kinase	CLIK1	Screening & Activity Analysis				●
Kinase	CLK1	Target Engagement	●	●		
Kinase	CLK1	Screening & Activity Analysis				●
Kinase	CLK2	Target Engagement	●	●		
Kinase	CLK2	Screening & Activity Analysis				●
Kinase	CLK3	Screening & Activity Analysis				●
Kinase	CLK4	Screening & Activity Analysis				●
Kinase	c-RAF	Screening & Activity Analysis				●
Kinase RTK	c-Ret-GFRα1	Functional	●	●		
Kinase RTK	c-Ret-GFRα2	Functional	●	●		
Kinase RTK	c-Ret-GFRα3	Functional	●			
Kinase	CRIK	Screening & Activity Analysis				●

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	CSF1R	Target Engagement	●	●		
Kinase CTK	CSF3R-JAK1	Functional	●	●		
Kinase	CSK	Screening & Activity Analysis				●
Kinase	CSNK1D	Target Engagement	●	●		
Kinase	DAPK1	Screening & Activity Analysis				●
Kinase	DAPK2	Screening & Activity Analysis				●
Kinase	DCAMKL1	Screening & Activity Analysis				●
Kinase	DCAMKL2	Screening & Activity Analysis				●
Kinase	DCAMKL3	Screening & Activity Analysis				●
Kinase	DDR1	Screening & Activity Analysis				●
Kinase RTK	DDR1	Functional	●	●		
Kinase	DDR2	Screening & Activity Analysis				●
Kinase	DMPK	Screening & Activity Analysis				●
Kinase	DNA-PK	Screening & Activity Analysis				●
Kinase	DRAK1	Screening & Activity Analysis				●
Kinase	DRAK2	Screening & Activity Analysis				●
Phosphatase	DUSP22	Screening & Activity Analysis				●
Kinase	DYRK1A	Screening & Activity Analysis				●
Kinase	DYRK1B	Screening & Activity Analysis				●
Kinase	DYRK2	Screening & Activity Analysis				●
Kinase	DYRK3	Screening & Activity Analysis				●
Kinase	eEF-2K	Screening & Activity Analysis				●
Kinase	EGFR	Screening & Activity Analysis				●
Kinase RTK	EGFR/ErbB2	Dimerization			●	
Kinase	EphA1	Screening & Activity Analysis				●
Kinase	EphA2	Screening & Activity Analysis				●
Kinase	EphA3	Screening & Activity Analysis				●
Kinase	EphA4	Screening & Activity Analysis				●
Kinase RTK	EphA4	Functional	●			
Kinase	EphA5	Screening & Activity Analysis				●
Kinase RTK	EphA5	Functional	●			
Kinase	EphA7	Screening & Activity Analysis				●
Kinase RTK	EphA7	Functional	●	●		
Kinase	EphA8	Screening & Activity Analysis				●
Kinase	EphB1	Screening & Activity Analysis				●
Kinase RTK	EphB1	Functional	●	●		
Kinase	EphB2	Screening & Activity Analysis				●
Kinase RTK	EphB2	Functional	●	●		
Kinase	EphB3	Screening & Activity Analysis				●
Kinase RTK	EphB3	Functional	●	●		
Kinase	EphB4	Screening & Activity Analysis				●
Kinase RTK	EphB4	Functional	●	●		
Kinase CTK	EpoR-JAK2	Functional	●	●		
Kinase RTK	ErbB1	Functional	●			
Kinase	ErbB2	Screening & Activity Analysis				●
Kinase RTK	ErbB2/ErbB3 (Her2)	Dimerization			●	
Kinase RTK	ErbB2/ErbB3 (Her2)	Functional	●	●		
Kinase	ErbB4	Screening & Activity Analysis				●
Kinase RTK	ErbB4	Functional	●			

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	FAK	Screening & Activity Analysis				●
Kinase	Fer	Screening & Activity Analysis				●
Kinase	Fes/Fps	Screening & Activity Analysis				●
Kinase	FGFR1	Screening & Activity Analysis				●
Kinase RTK	FGFR1	Functional	●	●		
Kinase RTK	FGFR1- $\alpha$ -Klotho	Functional	●			
Kinase RTK	FGFR1- $\beta$ -Klotho	Functional	●			
Kinase RTK	FGFR1v3B	Functional	●			
Kinase	FGFR2	Screening & Activity Analysis				●
Kinase RTK	FGFR2	Functional	●			
Kinase	FGFR3	Screening & Activity Analysis				●
Kinase RTK	FGFR3	Functional	●			
Kinase	FGFR4	Screening & Activity Analysis				●
Kinase RTK	FGFR4	Functional	●	●		
Kinase RTK	FGFR4- $\alpha$ -Klotho	Functional	●			
Kinase RTK	FGFR4- $\beta$ -Klotho	Functional	●	●		
Kinase RTK	FGFR4- $\beta$ -Klotho	SH2 Recruitment			●	
Kinase	FGR	Target Engagement	●			
Kinase CTK	FGR	Activity	●	●		
Kinase	Fgr	Screening & Activity Analysis				●
Kinase	Flt1	Screening & Activity Analysis				●
Kinase	Flt3	Screening & Activity Analysis				●
Kinase RTK	Flt3	Activity	●	●		
Kinase	Flt4	Screening & Activity Analysis				●
Kinase RTK	Flt4	Activity	●	●		
Kinase	Fms	Screening & Activity Analysis				●
Kinase	Fyn	Screening & Activity Analysis				●
Kinase	GAK	Screening & Activity Analysis				●
Kinase	GCK	Screening & Activity Analysis				●
Kinase	GCN2	Screening & Activity Analysis				●
Kinase RTK	GHR	SH2 Recruitment			●	
Kinase CTK	GHR-JAK1	Functional	●	●		
Kinase CTK	GHR-JAK2	Functional	●	●		
Kinase	GRK1	Screening & Activity Analysis				●
Kinase	GRK3	Screening & Activity Analysis				●
Kinase	GRK4	Screening & Activity Analysis				●
Kinase	GRK5	Screening & Activity Analysis				●
Kinase	GRK6	Screening & Activity Analysis				●
Kinase	GRK7	Screening & Activity Analysis				●
Kinase	GSK3 $\alpha$	Screening & Activity Analysis				●
Kinase	GSK3 $\beta$	Screening & Activity Analysis				●
Kinase	haspin	Screening & Activity Analysis				●
Kinase	HCK	Target Engagement	●	●		
Kinase	Hck	Screening & Activity Analysis				●
Phosphatase	HePTP	Screening & Activity Analysis				●
Kinase	HIPK1	Screening & Activity Analysis				●
Kinase	HIPK2	Screening & Activity Analysis				●
Kinase	HIPK3	Screening & Activity Analysis				●
Kinase	HIPK4	Screening & Activity Analysis				●

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	Histone	Screening & Activity Analysis				●
Kinase	HPK1	Screening & Activity Analysis				●
Kinase	HRI	Screening & Activity Analysis				●
Kinase	ICK	Screening & Activity Analysis				●
Kinase RTK	IGF1R	SH2 Recruitment			●	
Kinase RTK	IGF1R	Functional	●	●		
Kinase	IGF1R	Screening & Activity Analysis				●
Kinase	IKK $\alpha$	Screening & Activity Analysis				●
Kinase	IKK $\beta$	Screening & Activity Analysis				●
Kinase	IKK $\epsilon$	Screening & Activity Analysis				●
Kinase RTK	INSRa	Functional	●			
Kinase RTK	INSRb	SH2 Recruitment			●	
Kinase RTK	INSRb	Functional	●	●		
Kinase RTK	INSRb	Internalization	●			
Kinase	IR	Screening & Activity Analysis				●
Kinase	IRAK1	Screening & Activity Analysis				●
Kinase	IRAK4	Screening & Activity Analysis				●
Kinase	IRE1	Screening & Activity Analysis				●
Kinase	IRE2	Screening & Activity Analysis				●
Kinase	IRR	Screening & Activity Analysis				●
Kinase	Itk	Screening & Activity Analysis				●
Kinase	JAK1	Screening & Activity Analysis				●
Kinase	JAK2	Screening & Activity Analysis				●
Kinase CTK	JAK3	Activity	●	●		
Kinase	JAK3	Screening & Activity Analysis				●
Kinase	JNK1 $\alpha$ 1/SAPK1c	Screening & Activity Analysis				●
Kinase	JNK2 $\alpha$ 2/SAPK1a	Screening & Activity Analysis				●
Kinase	JNK3/SAPK1b	Screening & Activity Analysis				●
Kinase	KDR	Screening & Activity Analysis				●
Kinase RTK	KDR	Activity	●	●		
Kinase RTK	KDR/KDR	Dimerization			●	
Phosphatase	Lambda PP	Screening & Activity Analysis				●
Kinase	LATS1	Screening & Activity Analysis				●
Kinase	LATS2	Screening & Activity Analysis				●
Kinase CTK	LCK	Activity	●			
Kinase	Lck	Screening & Activity Analysis				●
Kinase	LIMK1	Screening & Activity Analysis				●
Kinase	LIMK2	Screening & Activity Analysis				●
Kinase	LKB1	Screening & Activity Analysis				●
Phosphatase	LMPTP-A	Screening & Activity Analysis				●
Phosphatase	LMPTP-B	Screening & Activity Analysis				●
Kinase	LOK	Screening & Activity Analysis				●
Kinase	LRRK2	Screening & Activity Analysis				●
Kinase	LTK	Screening & Activity Analysis				●
Kinase	Lyn	Screening & Activity Analysis				●
Kinase	MAK	Screening & Activity Analysis				●
Kinase	MAP4K3	Screening & Activity Analysis				●
Kinase	MAP4K4	Screening & Activity Analysis				●
Kinase	MAP4K5	Screening & Activity Analysis				●

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	MAPK1	Screening & Activity Analysis				●
Kinase	MAPK2	Screening & Activity Analysis				●
Kinase	MARK3	Screening & Activity Analysis				●
Kinase	MARK4	Screening & Activity Analysis				●
Kinase	MAPKAP-K2	Screening & Activity Analysis				●
Kinase	MAPKAP-K3	Screening & Activity Analysis				●
Kinase	MARK1	Screening & Activity Analysis				●
Kinase	MATK	Screening & Activity Analysis				●
Kinase	MEK1	Target Engagement	●	●		
Kinase	MEK1	Screening & Activity Analysis				●
Kinase	MEK2	Screening & Activity Analysis				●
Kinase	MEKK2	Screening & Activity Analysis				●
Kinase	MEKK3	Screening & Activity Analysis				●
Kinase	MELK	Screening & Activity Analysis				●
Kinase	Mer	Screening & Activity Analysis				●
Kinase RTK	MERTK	Functional	●			
Kinase	Met	Screening & Activity Analysis				●
Kinase	Met (D1246H)	Screening & Activity Analysis				●
Kinase	MINK	Screening & Activity Analysis				●
Kinase RTK	mINSRa	Functional	●	●		
Kinase RTK	mINSRb	Functional	●	●		
Kinase	MKK3	Screening & Activity Analysis				●
Kinase	MKK4	Screening & Activity Analysis				●
Kinase	MKK6	Screening & Activity Analysis				●
Phosphatase	MKP5	Screening & Activity Analysis				●
Kinase	MLCK	Screening & Activity Analysis				●
Kinase	MLK1	Screening & Activity Analysis				●
Kinase	MLK2	Screening & Activity Analysis				●
Kinase	MLK3	Screening & Activity Analysis				●
Kinase	MLK4	Screening & Activity Analysis				●
Kinase	Mnk2	Screening & Activity Analysis				●
Kinase	MOK	Screening & Activity Analysis				●
Kinase	MRCK $\alpha$	Screening & Activity Analysis				●
Kinase	MRCK $\beta$	Screening & Activity Analysis				●
Kinase	MRCK $\gamma$	Screening & Activity Analysis				●
Kinase	MSK1	Screening & Activity Analysis				●
Kinase	MSK2	Screening & Activity Analysis				●
Kinase	MSSK1	Screening & Activity Analysis				●
Kinase	MST1	Screening & Activity Analysis				●
Kinase	MST2	Screening & Activity Analysis				●
Kinase	MST3	Screening & Activity Analysis				●
Kinase	MST4	Screening & Activity Analysis				●
Kinase	MuSK	Screening & Activity Analysis				●
Kinase	MYLK2	Screening & Activity Analysis				●
Kinase	MYO3B	Screening & Activity Analysis				●
Kinase	NDR1	Screening & Activity Analysis				●
Kinase	NDR2	Screening & Activity Analysis				●
Kinase	NEK1	Screening & Activity Analysis				●
Kinase	NEK2	Screening & Activity Analysis				●



Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	NEK3	Screening & Activity Analysis				●
Kinase	NEK4	Screening & Activity Analysis				●
Kinase	NEK5	Screening & Activity Analysis				●
Kinase	NEK6	Screening & Activity Analysis				●
Kinase	NEK7	Screening & Activity Analysis				●
Kinase	NEK9	Screening & Activity Analysis				●
Kinase	NEK11	Screening & Activity Analysis				●
Kinase	NIM1	Screening & Activity Analysis				●
Kinase	NLK	Screening & Activity Analysis				●
Kinase	NuaK1 (ARK5)	Screening & Activity Analysis				●
Kinase	NUAK2	Screening & Activity Analysis				●
Kinase	OSR1	Screening & Activity Analysis				●
Kinase	p38 $\alpha$ /SAPK2a	Screening & Activity Analysis				●
Kinase	p38 $\beta$ 2/SAPK2b2	Screening & Activity Analysis				●
Kinase	p38 $\gamma$ /SAPK3	Screening & Activity Analysis				●
Kinase	p38 $\delta$ /SAPK4	Screening & Activity Analysis				●
Kinase	p53	Screening & Activity Analysis				●
Kinase	p70S6K	Screening & Activity Analysis				●
Kinase	PAK1	Screening & Activity Analysis				●
Kinase	PAK2	Screening & Activity Analysis				●
Kinase	PAK3	Screening & Activity Analysis				●
Kinase	PAK4	Screening & Activity Analysis				●
Kinase	PAK5	Screening & Activity Analysis				●
Kinase	PAK6	Screening & Activity Analysis				●
Kinase	PAR-1Ba	Screening & Activity Analysis				●
Kinase	PASK	Screening & Activity Analysis				●
Kinase	PDGFR $\alpha$	Screening & Activity Analysis				●
Kinase RTK	PDGFR $\alpha$	SH2 Recruitment			●	
Kinase RTK	PDGFR $\alpha$	Functional	●	●		
Kinase	PDGFR $\beta$	Screening & Activity Analysis				●
Kinase RTK	PDGFR $\beta$	SH2 Recruitment			●	
Kinase RTK	PDGFR $\beta$	Functional	●	●		
Kinase	PDHK2	Screening & Activity Analysis				●
Kinase	PDHK4	Screening & Activity Analysis				●
Kinase	PDK1	Screening & Activity Analysis				●
Kinase	PEK	Screening & Activity Analysis				●
Kinase	PhKg1	Screening & Activity Analysis				●
Kinase	PhKg2	Screening & Activity Analysis				●
Kinase	PI3K-C2 $\alpha$	Screening & Activity Analysis				●
Kinase	PI3K-C2 $\beta$	Screening & Activity Analysis				●
Kinase	PI3K-C2 $\gamma$	Screening & Activity Analysis				●
Kinase	PI3K $\alpha$ /p65 $\alpha$	Screening & Activity Analysis				●
Kinase	PI3K $\alpha$ /p85 $\alpha$	Screening & Activity Analysis				●
Kinase	PI3K $\beta$ /p85 $\alpha$	Screening & Activity Analysis				●
Kinase	PI3K $\beta$ /p85 $\beta$	Screening & Activity Analysis				●
Kinase	PI3K $\gamma$	Screening & Activity Analysis				●
Kinase	PI3K $\delta$	Target Engagement	●			
Kinase	PI3K $\delta$ /p85 $\alpha$	Screening & Activity Analysis				●
Kinase	PI4KIII $\alpha$	Screening & Activity Analysis				●



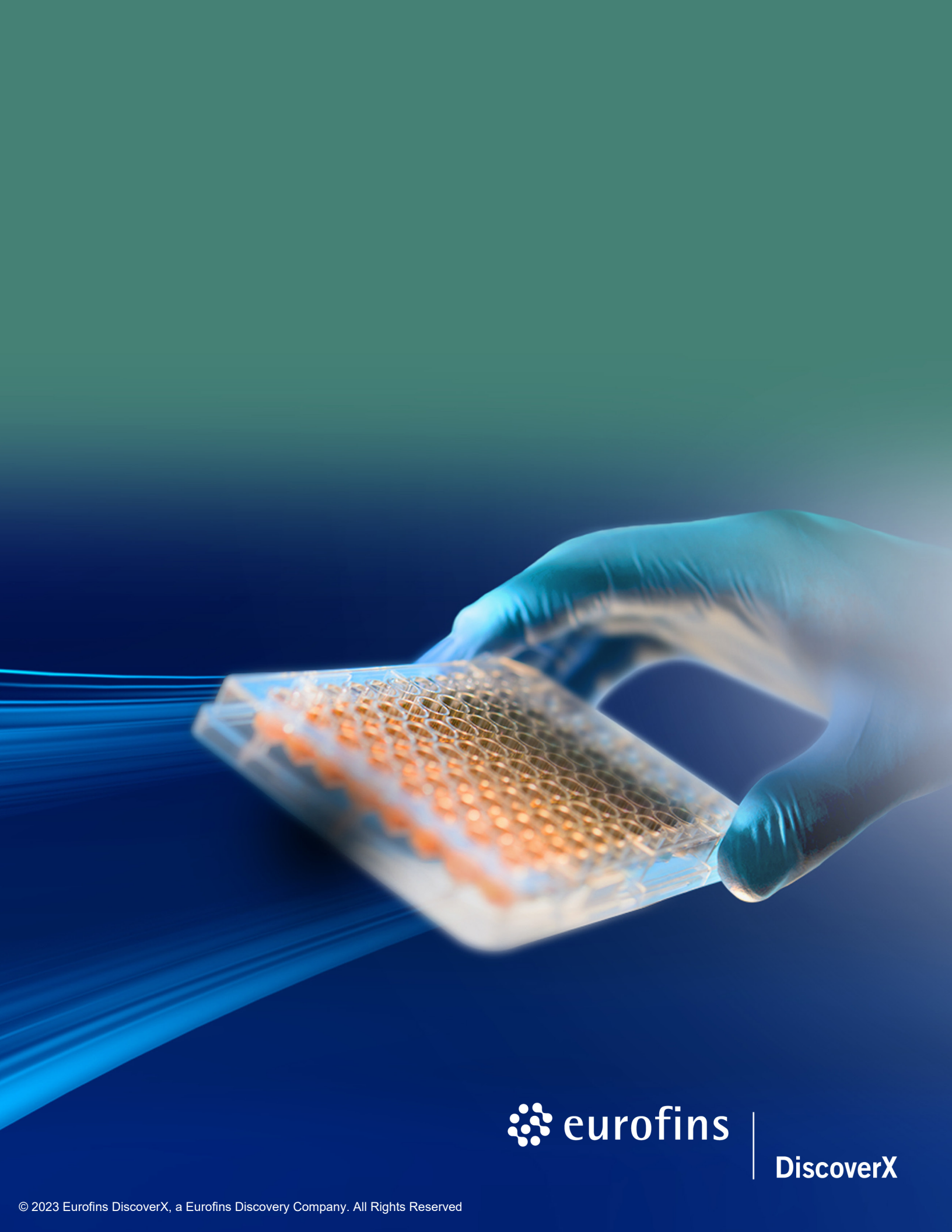
Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	PIK3C3	Screening & Activity Analysis				●
Kinase	PIKFYVE	Target Engagement	●			
Kinase	PIM1	Target Engagement	●	●		
Kinase	Pim-1	Screening & Activity Analysis				●
Kinase	Pim-2	Screening & Activity Analysis				●
Kinase	Pim-3	Screening & Activity Analysis				●
Kinase	PIP4K2α	Screening & Activity Analysis				●
Kinase	PIP5K1α	Screening & Activity Analysis				●
Kinase	PIP5K1γ	Screening & Activity Analysis				●
Kinase	PKA	Screening & Activity Analysis				●
Kinase	PKAcβ	Screening & Activity Analysis				●
Kinase	PKCμ	Screening & Activity Analysis				●
Kinase	PKCα	Screening & Activity Analysis				●
Kinase	PKCβI	Screening & Activity Analysis				●
Kinase	PKCβII	Screening & Activity Analysis				●
Kinase	PKCγ	Screening & Activity Analysis				●
Kinase	PKCδ	Screening & Activity Analysis				●
Kinase	PKCε	Screening & Activity Analysis				●
Kinase	PKCζ	Screening & Activity Analysis				●
Kinase	PKCη	Screening & Activity Analysis				●
Kinase	PKCθ	Screening & Activity Analysis				●
Kinase	PKCθ	Target Engagement	● <i>New</i>			●
Kinase	PKCι	Screening & Activity Analysis				●
Kinase	PKD2	Screening & Activity Analysis				●
Kinase	PKG1α	Screening & Activity Analysis				●
Kinase	PKG1β	Screening & Activity Analysis				●
Kinase	PKR	Screening & Activity Analysis				●
Kinase	Plk1	Screening & Activity Analysis				●
Kinase	PLK3	Screening & Activity Analysis				●
Phosphatase	PP1α	Screening & Activity Analysis				●
Phosphatase	PP5	Screening & Activity Analysis				●
Kinase	PRAK	Screening & Activity Analysis				●
Kinase	PRK2	Screening & Activity Analysis				●
Kinase	PrKX	Screening & Activity Analysis				●
Kinase CTK	PRLR-JAK1	Functional	●	●		
Kinase CTK	PRLR-JAK2	Functional	●	●		
Kinase	PTK5	Screening & Activity Analysis				●
Phosphatase	PTP-1B	Screening & Activity Analysis				●
Phosphatase	PTPMEG-1	Screening & Activity Analysis				●
Phosphatase	PTPMEG-2	Screening & Activity Analysis				●
Phosphatase	PTPN22	Screening & Activity Analysis				●
Phosphatase	PTPβ	Screening & Activity Analysis				●
Kinase	PYK2	Target Engagement	●	●		
Kinase	PYK2	Screening & Activity Analysis				●
Kinase	Ret	Screening & Activity Analysis				●
Kinase RTK	rhTrkA-rhP75	Functional	●	●		
Kinase	RIPK1	Target Engagement	●			
Kinase	RIPK1	Screening & Activity Analysis				●
Kinase	RIPK2	Target Engagement	●	●		

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase	RIPK2	Screening & Activity Analysis				●
Kinase	RIPK3	Target Engagement	●			
Kinase	RK2	Screening & Activity Analysis				●
Kinase	ROK $\alpha$ /ROCK-II	Screening & Activity Analysis				●
Kinase	ROK $\beta$ /ROCK-I	Screening & Activity Analysis				●
Kinase	Ron	Screening & Activity Analysis				●
Kinase	Ros	Screening & Activity Analysis				●
Phosphatase	RPTP $\mu$	Screening & Activity Analysis				●
Kinase	Rse	Screening & Activity Analysis				●
Kinase	Rsk1	Screening & Activity Analysis				●
Kinase	Rsk2	Screening & Activity Analysis				●
Kinase	Rsk3	Screening & Activity Analysis				●
Kinase	Rsk4	Screening & Activity Analysis				●
Kinase RTK	rTrkA-rP75	Functional	●	●		
Kinase	SAPK4	Screening & Activity Analysis				●
Kinase	SGK1	Screening & Activity Analysis				●
Kinase	SGK2	Screening & Activity Analysis				●
Kinase	SGK3	Screening & Activity Analysis				●
Phosphatase	SHP-1	Screening & Activity Analysis				●
Phosphatase	SHP-2	Screening & Activity Analysis				●
Kinase	SIK	Screening & Activity Analysis				●
Kinase	SIK1	Target Engagement	●	●		
Kinase	SIK2	Screening & Activity Analysis				●
Kinase	SIK3	Screening & Activity Analysis				●
Kinase	SNRK	Screening & Activity Analysis				●
Kinase	Src	Screening & Activity Analysis				●
Kinase	SRPK1	Screening & Activity Analysis				●
Kinase	SRPK2	Screening & Activity Analysis				●
Kinase	STK25	Screening & Activity Analysis				●
Kinase	STK32A	Screening & Activity Analysis				●
Kinase	STK33	Screening & Activity Analysis				●
Kinase	STK39	Screening & Activity Analysis				●
Kinase CTK	Syk	Activity	●	●		
Kinase	Syk	Screening & Activity Analysis				●
Kinase	Tak1	Screening & Activity Analysis				●
Kinase	TAO1	Screening & Activity Analysis				●
Kinase	TAO2	Screening & Activity Analysis				●
Kinase	TAO3	Screening & Activity Analysis				●
Kinase	TBK1	Screening & Activity Analysis				●
Phosphatase	TCPTP	Screening & Activity Analysis				●
Kinase	Tec	Screening & Activity Analysis				●
Kinase	TGFBR1	Screening & Activity Analysis				●
Kinase	Tie2	Screening & Activity Analysis				●
Kinase RTK	Tie2	Functional	●			
Kinase	TLK1	Screening & Activity Analysis				●
Kinase	TLK2	Screening & Activity Analysis				●
Phosphatase	TMDP	Screening & Activity Analysis				●
Kinase	mTOR (FRAP1)	Screening & Activity Analysis				●
Kinase	TrkA	Screening & Activity Analysis				●

Kinase and Phosphatase Products			Product Type			
Family	Target	Application	Cell Line	eXpress Kit	Bioassay Kit	Protein
Kinase RTK	TrkA	Functional	●	●		
Kinase RTK	TrkA-p75	SH2 Recruitment			●	
Kinase RTK	TrkA-P75	Functional	●	●		
Kinase	TrkB	Screening & Activity Analysis				●
Kinase RTK	TrkB	Functional	●	●		
Kinase RTK	TrkB-P75	SH2 Recruitment			●	
Kinase RTK	TrkB-P75	Functional	●	●		
Kinase	TrkC	Screening & Activity Analysis				●
Kinase RTK	TrkC	SH2 Recruitment			●	
Kinase RTK	TrkC	Functional	●	●		
Kinase RTK	TrkC-P75	Functional	●	●		
Kinase	TSSK1	Screening & Activity Analysis				●
Kinase	TSSK2	Screening & Activity Analysis				●
Kinase	TSSK3	Screening & Activity Analysis				●
Kinase	TSSK4	Screening & Activity Analysis				●
Kinase	TTBK2	Screening & Activity Analysis				●
Kinase	TTK	Screening & Activity Analysis				●
Kinase	Txk	Screening & Activity Analysis				●
Kinase CTK	TYK2	Activity	●	●		
Kinase	TYK2	Screening & Activity Analysis				●
Kinase RTK	TYRO3	Functional	●			
Kinase	ULK1	Screening & Activity Analysis				●
Kinase	ULK2	Screening & Activity Analysis				●
Kinase	ULK3	Screening & Activity Analysis				●
Phosphatase	VHR	Screening & Activity Analysis				●
Kinase	VRK2	Screening & Activity Analysis				●
Kinase	WEE1	Screening & Activity Analysis				●
Kinase	WNK2	Screening & Activity Analysis				●
Kinase	WNK3	Screening & Activity Analysis				●
Kinase	Yes	Screening & Activity Analysis				●
Kinase CTK	YES1	Activity	●	●		
Phosphatase	YopH	Screening & Activity Analysis				●
Kinase	ZAK	Screening & Activity Analysis				●
Kinase	ZAP-70	Screening & Activity Analysis				●
Kinase	ZIPK	Screening & Activity Analysis				●

Nuclear Hormone Receptor Products		Product Type
Target	Application	Cell Line
AR	Nuclear Translocation	●
AR	Protein Interaction	●
ARH	Protein Interaction	●
ER $\alpha$	Protein Interaction	●
FXR	Protein Interaction	●
GR	Nuclear Translocation	●
GR	Protein Interaction	●
LXR $\alpha$	Nuclear Translocation	●
LXR $\alpha$	Protein Interaction	●
LXR $\beta$	Protein Interaction	●
LXR $\beta$ -NCoR1	Protein Interaction	●
MR	Nuclear Translocation	●
MR	Protein Interaction	●
PPAR $\alpha$	Protein Interaction	●
PPAR $\gamma$	Protein Interaction	●
PPAR $\delta$	Protein Interaction	●
PR $\alpha$	Protein Interaction	●
PR $\beta$	Nuclear Translocation	●
PR $\beta$	Protein Interaction	●
RAR $\alpha$	Protein Interaction	●
RAR $\beta$	Protein Interaction	●
RXR $\alpha$	Protein Interaction	●
RXR $\gamma$	Protein Interaction	●
THR $\alpha$	Protein Interaction	●
THR $\beta$	Protein Interaction	●

Signaling Pathway Products		Product Type		
Target	Application	Cell Line	eXpress Kit	Bioassay Kit
AMHR2/BMPRI1A	Dimerization	●		
Bcl2-Bax	Protein Binding, Target Engagement	●		
Bcl2-Bim	Protein Binding, Target Engagement	●		
Bcl2L1-Bax	Protein Binding, Target Engagement	●	●	
Bcl2L1-Bim	Protein Binding, Target Engagement	●		
CD27 NF-kB	Reporter	●		
CDC25A	Degradation	●	●	
c-Myc	Target Protein Degradation	●		
GM-CSF	Reporter	●		
HepG2 STAT3	Reporter	●		
IκB	Signaling			●
IKZF3v1	Degradation	●		
IL11R STAT3	Reporter	●		
Keap1-NRF2	Nuclear Translocation	●	●	
MCL1-BIM	Protein Binding, Target Engagement	●		
NFAT	Reporter	●		
NFκB	Degradation	●	●	
NFκB	Nuclear Translocation	●	●	
NFκB	Reporter	●		
NPR1	Functional	●		
p53	Nuclear Translocation	●		
PD1	Reporter	●		
PI3	Nuclear Translocation	●	●	
RANK-IκB	Functional	●		
RANK-NFκB	Reporter	●		
SHP2	Protein Binding, Target Engagement	●		
SREBP1c	Nuclear Translocation	●		
SREBP2	Nuclear Translocation	●	●	
TREM1 DAP12	Reporter	●		
TORC1 (CRTC1)	Nuclear Translocation	●	●	
TORC2 (CRTC2)	Nuclear Translocation	●	●	
TORC3 (CRTC3)	Nuclear Translocation	●		
Wnt-FzGSK3β	Nuclear Translocation	●	●	
XBP1	Nuclear Translocation	●	●	



 eurofins

DiscoverX