



PathHunter® Cell-Based Assays for Kinase Receptors



Functional, Dimerization, and Translocation Cellular Assays for Receptor Tyrosine Kinases (RTKs)

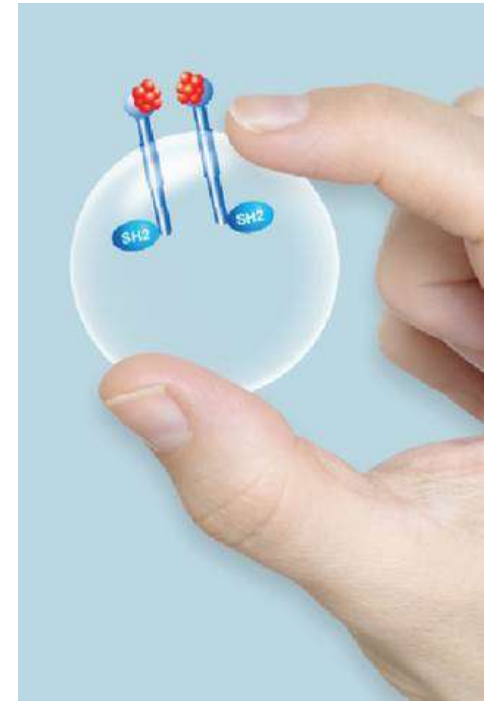
OUR EXPERTISE
IN YOUR HANDS.
DISCOVER
CONFIDENTLY.

Receptor Tyrosine Kinases

Coordinating a wide variety of cellular functions

Receptor Tyrosine Kinases (RTKs)

- ~20 Classes and ~60 known human RTKs
 - e.g. FGF, PDGF, Trk, Eph, IGF, INS
- Cell surface receptors that bind hormones, growth factors, and cytokines
- Regulators of cellular growth/proliferation, survival, development/differentiation, and disease progression
- Play a critical role the development and progression of several cancer types



Receptor Tyrosine Kinases

Understanding RTKs structure and signaling

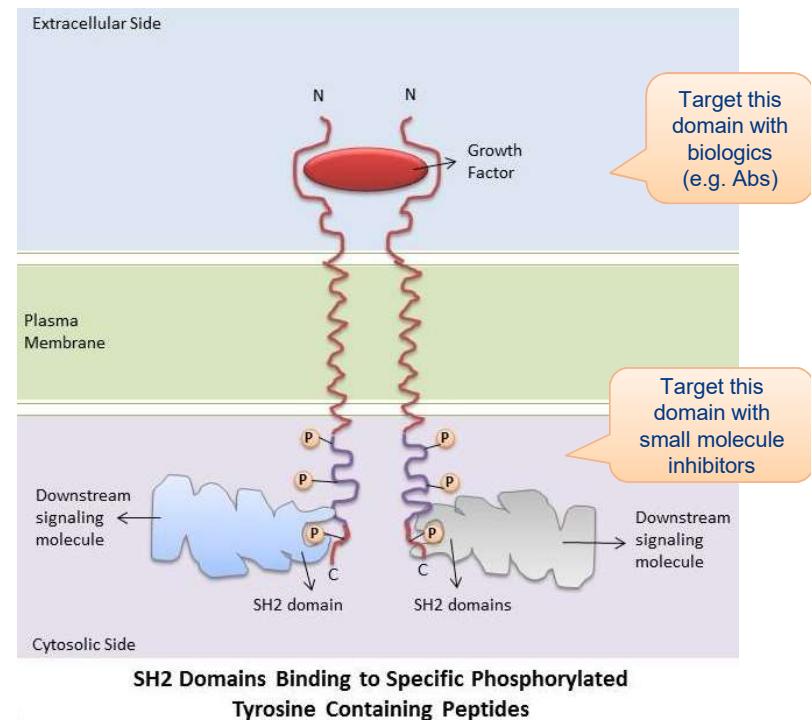
RTKs structure

- N-terminal extracellular ligand-binding domain
- C-terminal intracellular catalytic domain responsible for kinase activity
 - Autophosphorylation and tyrosine phosphorylation of RTK substrates

RTK cellular signaling

1. Ligand binding induces receptor dimerization (hetero- or homodimerization) and kinase domain activation
2. Autophosphorylation of RTKs
3. SH2-protein recruitment (at phosphotyrosine-containing motifs)
4. Promotion of further phosphorylation of downstream proteins (e.g. MAP kinases)
5. Ultimately stimulates enzymatic activity and associated with intended cellular response

Receptor Tyrosine Kinases Structure



Drug therapeutic relevance

RTKs are attractive targets for drug therapy due to their implications in a variety of cellular abnormalities and they are tractable targets for drug discovery medicinal chemistry

- Cancer, degenerative diseases, and cardiovascular diseases
- FDA has approved several anti-cancer drugs caused by activated RTKs

Example of FDA Approved Small Molecule Inhibitors and Antibodies Against RTKs* for Cancer Therapy

Drug	Target	Disease
Imatinib (Gleevec)	PDGFR, KIT, Abl, Arg	SML, GIST
Gefitinib (Iressa)	EGFR	Esophageal cancer, Glioma
Erlotinib (Tarceva)	EGFR	Esophageal cancer, Glioma
Sorafenib (Nexavar)	Raf, VEGFR, PDGFR, Flt3, KIT	Renal cell carcinoma
Sunitinib (Sutent)	KIT, VEGFR, PDGFR, Flt3	Renal cell carcinoma, GIST, Endocrine pancreatic cancer
Dasatinib (Sprycel)	Abl, Arg, KIT, PDGFR, Src	Gleevec-resistant CML
Nilotinib (Tasigna)	Abl, Arg, KIT, PDGFR	Gleevec-resistant CML
Lapatinib (Tykerb)	EGFR, ErbB2	Mammary carcinoma
Trastuzumab (Herceptin)	ErbB2	Mammary carcinoma
Cetuximab (Erbiximab)	EGFR	Colorectal cancer, Head and neck cancer
Bevacizumab (Avastin)	VEGF	Lung cancer, Colorectal cancer
Panitumumab (Vectibix)	EGFR	Colorectal cancer

Table adapted from "Cell signalling by receptor-tyrosine kinases," by Lemmon and Schlessinger's, 2010. Cell, 141, p. 1117-1134.

* Includes non RTK examples like Abl, Raf, ...

PathHunter® Cell-Based Assays
for Kinase Receptors

Functional, Dimerization, and Translocation Assays based on EFC technology

Investigate RTKs activity with high confidence

In general, it is difficult to create assays for cytokines and growth factors

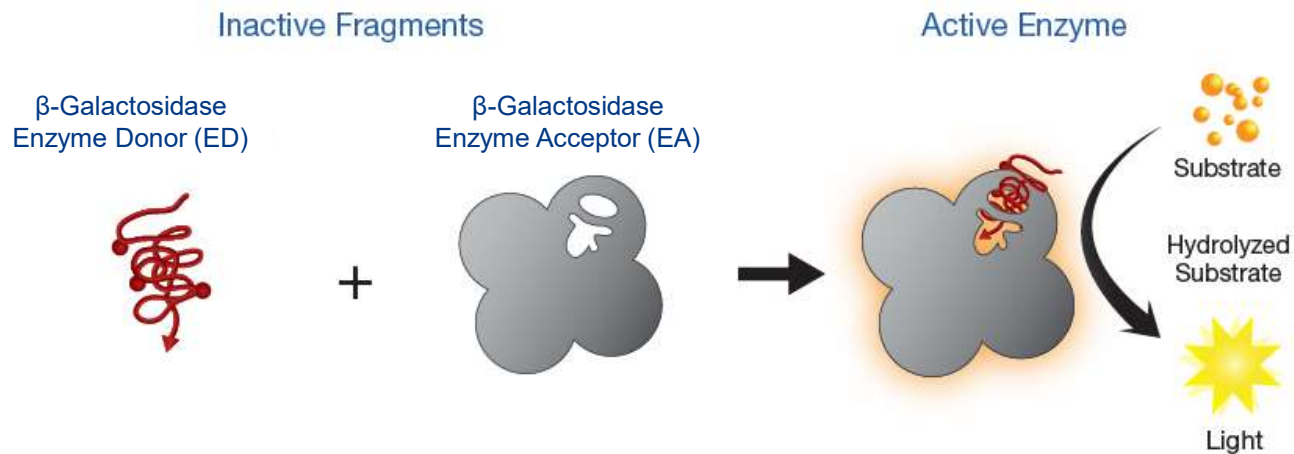
- Available assays have downstream readouts
- Lack of assay specificity
- Long complex protocols

Highlights of PathHunter cell-based assays for receptor kinases

- Provide solutions for analyzing functional activity, dimerization, translocation, and characterization
- Allow for screening, hit ID, and lead optimization of small molecules and biologics
- Give cellular context to activation and allow for the identification of novel inhibitors and therapeutic antibodies
 - E.g. Inhibitors of receptor dimerization and/or receptor function
- Based on proprietary DiscoverX enzyme complementation fragment technology platform

Enzyme Fragment Complementation (EFC)

PathHunter assays are based on the proprietary DiscoverX EFC Technology



EFC is a patented detection technology based on two recombinant β -galactosidase (β -gal) fragments

- Small peptide fragment (enzyme donor, ED)
- Large protein fragment (enzyme acceptor, EA)

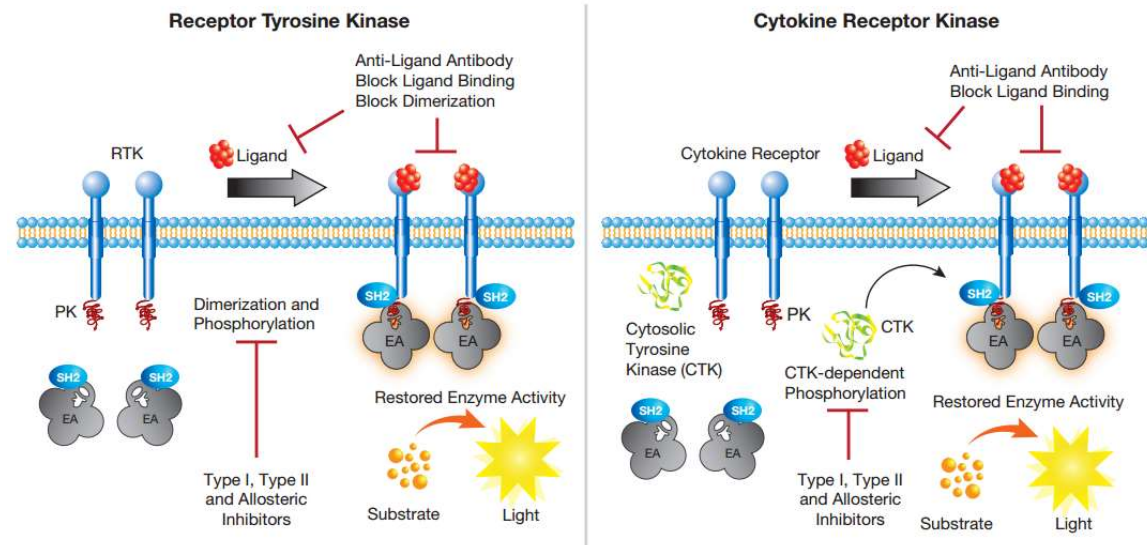
Separately, the fragments are inactive, but when combined, they form an active enzyme that hydrolyzes substrate to produce a chemiluminescent signal

Easily analyze kinase activation for multiple tyrosine kinases receptors

PathHunter cell-based functional assays for RTKs and cytokine receptors with CTKs

- Involves the target receptor (RTK or cytokine receptor kinase) tagged with PK and one of the many different partner proteins containing SH2 domains tagged with EA
- Upon ligand-induced activation, the receptors dimerize and cross-phosphorylate
- The SH2-EA fusion protein then binds the phosphorylated receptor-PK forcing the EFC interaction which is readout as a chemiluminescent signal

EFC Cell-Based Assays for RTKs, CTKs, & Cytokine Receptor Kinases



PathHunter® Cell-Based Dimerization Assays

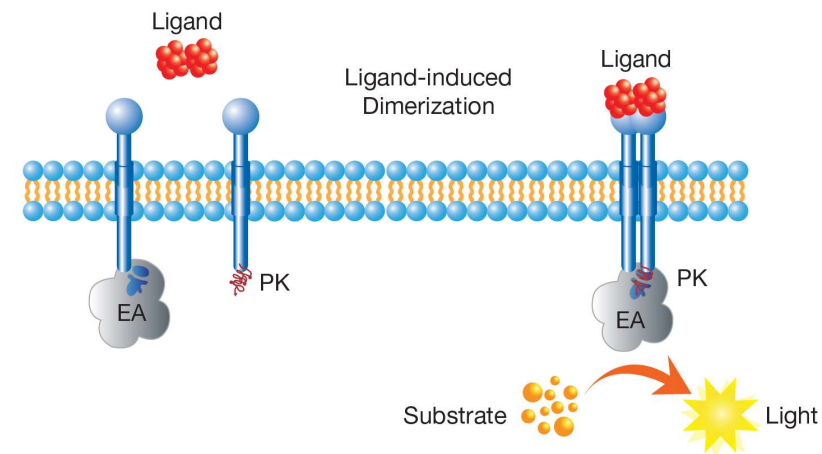
Measure heterodimerization, homodimerization, or co-receptor recruitment

PathHunter cell-based dimerization assays

- Involves two target receptors that are tagged with the EFC enzyme donor called ProLink™ (PK) or Enzyme Acceptor (EA)
- Upon ligand-induced activation (e.g. via biologics), the receptors dimerize forcing the two β -gal components to complement and create an active enzyme
- Active β -gal generates a chemiluminescent signal in the presence of substrate

Example heterodimer pairs and their therapeutic indications involvement

- EGFR-ErbB2 (Breast Cancer)
- ErbB2-ErbB3 (Cancer)



PathHunter® Cell-Based Translocation Assays

Track cellular movement of proteins to multiple membrane compartments

Monitor translocation of wild type, full length RTKs in an antibody-free, non-imaging, cell-based, homogeneous (no wash), and HTS-friendly assay format

Explore receptor movements from ER to cell membrane or cell membrane to endosome

- If these localizations are altered by a ligand, protein mutations, or aberrant signaling, undesirable effects may occur, often resulting in drug tolerance, unwanted side effects, and disease

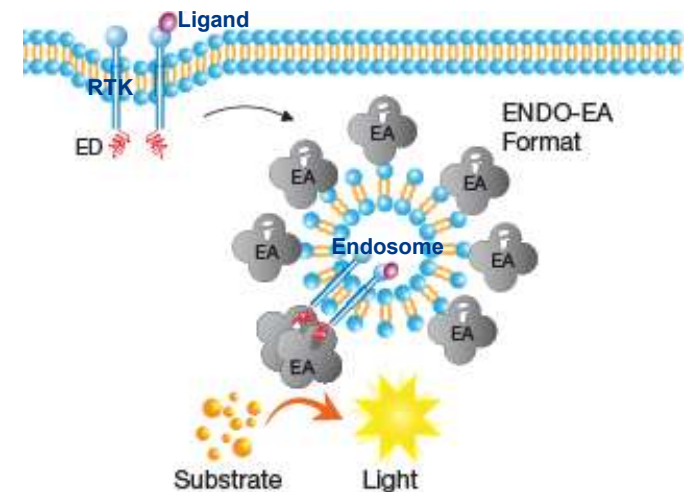
Study RTK recycling patterns

Discover compounds that have unique qualities

- e.g. agonists that behave as functional antagonists with respect to receptor internalization

Create your own quantitative cell-based assays to study translocation of any membrane protein





Tracking the movement of proteins between cellular compartments using EFC assays



Products and Services for Kinases

Technology access for analysis of RTKs, CTKs, and more

Products

-  • Stable cell lines
-   • Complete assay ready kits
 - eXpress and bioassay
-  • Engineered parental cell lines
 - For development of translocation and SH2-recruitment functional/dimerization assays

Services

- TKscan™ cell-based tyrosine kinase panel for screening and profiling
 - tkMAX, tkELECT, tkE/IC50ELECT

Related products and services

- InCELL cell-based compound-target engagement assays
- ADP accumulation fluorescent assays
- Ligands and inhibitors
- KINOMEscan®
- BioMAP® phenotypic screening

Custom assay development (CAD)



Products and Services for Kinases

Broad Range of assay targets available for analysis of RTKs, CTKs, and more

Dimerization Assays*	
ACVR1/ACVR2	CSF2RB/CSF2RA
ACVR1B/BMPR2	EGFR/EGFR
ACVR1C/ACVR2	EGFR/ErbB2
ACVR1C/ACVR2B	EGFR/ErbB3
ACVRL1/ACVR2	EpoR/EpoR
ACVRL1/ACVR2B	ErbB2/ErbB3
ACVRL1/BMPR2	ErbB2/ErbB4
BMPR1A/ACVR2	ErbB4/ErbB4
BMPR1A/ACVR2B	FGFR3(G380R)/FGFR3(G380R)
BMPR1A/BMPR2	FGFR3/FGFR3
BMPR1B/ACVR2A	KDR/KDR
BMPR1B/ACVR2B	TGFB1/ACVR2
BMPR1B/BMPR2	TGFB1/ACVR2B
c-MET/c-MET	TGFB1/TGFB2
CSF1R/CSF1R	TGFB1/TGFB2/ENG

RTK and CTK Functional Assays			
c-KIT	FGFR1	mINSRb	CTK
c-MET	FGFR1- α -Klotho	KDR	BLK
c-Ret-GFR α 1	FGFR1- β -Klotho	PDGFRa	CSF3R-JAK1
c-Ret-GFR α 2	FGFR2	PDGFRb	EpoR-JAK2
DDR1	FGFR4	TrkA	FGR
EphA4	FGFR4- α -Klotho	TrkA-P75	GHR-JAK1
EphA5	FGFR4- β -Klotho	rhTrkA-rhP75	GHR-JAK2
EphA7	Flt3	rTrkA-rP75	JAK3
EphB1	Flt4	Tie2	LCK
EphB2	IGF1R	TrkB	PRLR-JAK1
EphB3	INSRa	TrkB-P75	PRLR-JAK2
EphB4	INSRb	TrkC	SYK
ErbB1	clNSRa	TrkC-P75	TYK2
ErbB2-ErbB3	clNSRb	TYRO3	YES1
ErbB4	mINSRa		

Translocation and Functional Assays	
DIY/Toolbox* (Parental Cell Lines)	
ENDO-EA	MEM-EA
PLCG1(SH2)-EA	SHC1(SH2)-EA

Select assays shown. Please refer to the website for the current list of PathHunter [RTK/CTK](#) assays, [dimerization](#) assays, [translocation](#) assays, and [kinase profiling](#) services
 * Dimerization assays: interleukin receptors and other receptors are also available but not shown here.

Advantages of PathHunter[®] cell-based assays for kinase receptors

Accurate, Sensitive, and Reproducible - Superior quality, reproducible data with large assay windows, and robust performance

Broadly Applicable

- Identify various ligands including anti-receptor, anti-ligand, activating antibodies, non-ATP pocket binders (allosteric modulators or dimerization inhibitors), or ligand binding inhibitors (ATP-competitors)
- Study hetero- and homodimers

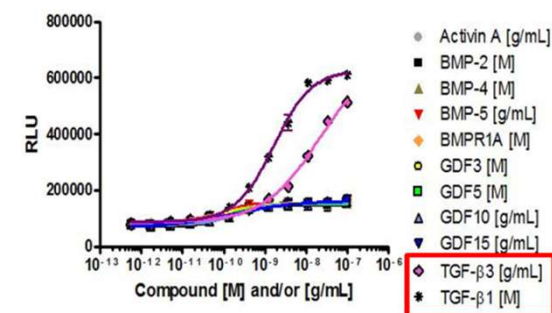
High Specificity - Tagged, full-length tyrosine kinase eliminates non-specificity due to background from endogenous tyrosine kinases

Easy-to-Use – Simple, one-step add and read protocol that is HTS-friendly

Universal, DIY Solution – Create your own quantitative cell-based assays to study translocation or function of RTK/CTKs

Example of Assay Specificity

TGF β R1-TGF β R2 Dimerization Assay



PathHunter® Cell-Based Assays for Kinase Receptors

Applications and Validation data for Functional, Dimerization, and Translocation (internalization) Assays

Inhibit the kinase domain of tyrosine kinases with small molecule inhibitors

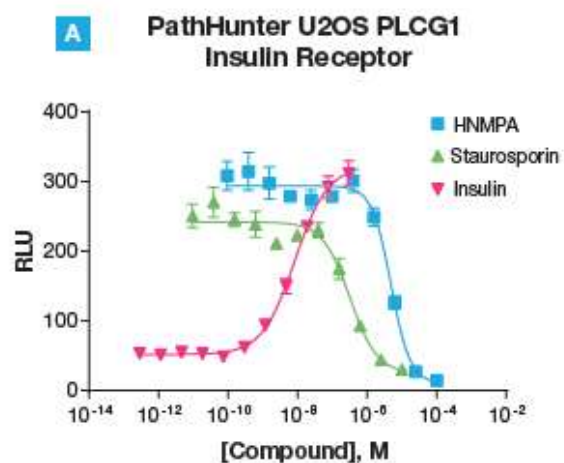
Select tyrosine kinase receptors and their small molecule inhibitors

Kinase ID	Compound ID	Historical IC ₅₀ (nM)
AXL	Crizotinib	195
BLK	EXEL-2880/XL-880/Foretinib	216
c-KIT	AG-013736 (Axitinib)	3.07
c-MET	EXEL-2880/XL-880/Foretinib	58.1
c-Ret-GFRa2	EXEL-2880/XL-880/Foretinib	7.88
DDR1	Dasatinib	0.679
ErbB1 (EGFR)	Gefitinib	98.8
ErbB2/ErbB3	Lapatinib	51.4
ErbB4	Lapatinib	39.4
FGFR2	AZD-4547	4.39
FGR	Dasatinib	0.997
FLT3 Activity	PKC-412	12.3
FLT3 Functional	EXEL-2880/XL-880/Foretinib	10.1
FLT4 Activity	EXEL-2880/XL-880/Foretinib	4
IGF1R	GSK 1838705A	14.2
INSR	GSK 1838705A	31.4

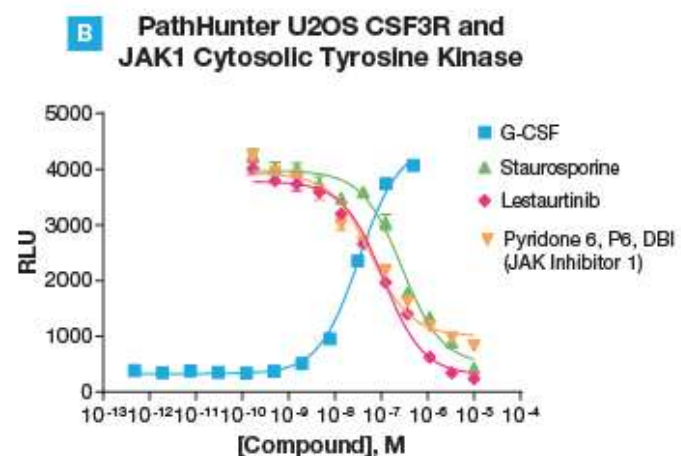
Kinase ID	Compound ID	Historical IC ₅₀ (nM)
JAK3	CP690550	1100
KDR	ABT-869	0.848
PDGFRa	AG-013736 (Axitinib)	10.2
PDGFRb	AG-013736 (Axitinib)	8.29
PRLR-JAK1	INCB018424	84.4
PRLR-JAK2	INCB018424	379
SYK	MNS	2030
TrkA	EXEL-2880/XL-880/Foretinib	9.51
TrkA-p75	EXEL-2880/XL-880/Foretinib	14.7
TrkB	EXEL-2880/XL-880/Foretinib	11.5
TrkB-p75	EXEL-2880/XL-880/Foretinib	15.4
TrkC	EXEL-2880/XL-880/Foretinib	6.56
TrkC-p75	EXEL-2880/XL-880/Foretinib	4.22
TYK2	INCB018424	1890

Evaluate multiple mechanisms of activation and small molecule inhibition

Insulin receptor (INSR) (SH2-adaptor protein = PLCG1)



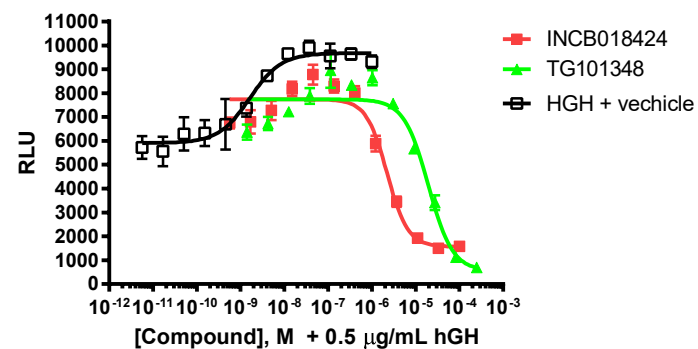
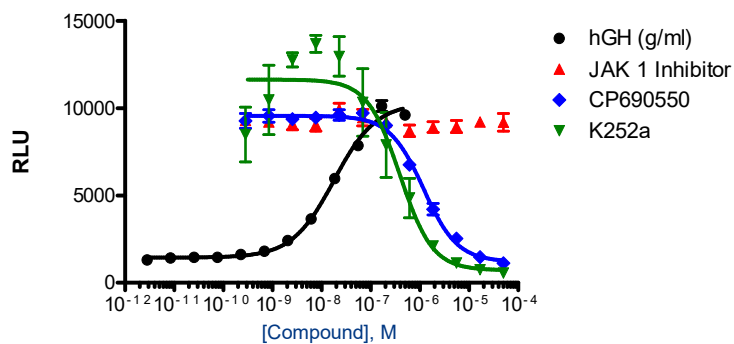
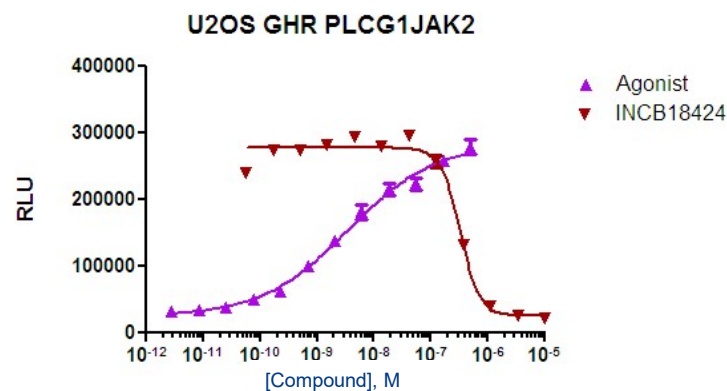
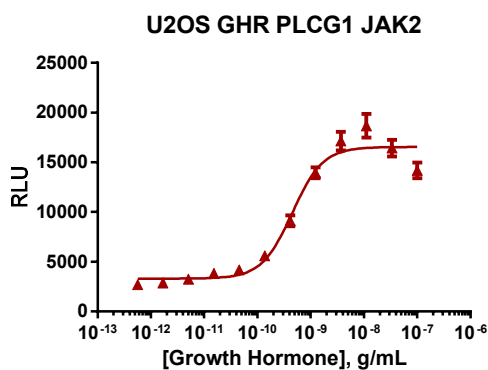
Colony stimulating factor 3 receptor (CSF3R) cytokine receptor (CTK = JAK1)



A. The insulin receptor is constitutively dimerized, but undergoes a conformational change to become active in the presence of insulin (agonist) and this can be inhibited by the small molecule antagonists staurosporine and HNMPA. B. The colony stimulating factor 3 receptor (CSF3R-JAK1) cell line was preincubated with antagonists staurosporine, lestaurtinib and pyridone 6, P6, DBI (a Jak inhibitor 1) and then challenged with EC₅₀ of agonist G-CSF. A dose dependent inhibition was observed with antagonist indicating that the assay can be used effectively to profile or screen inhibitors against the cytosolic tyrosine kinase JAK1.

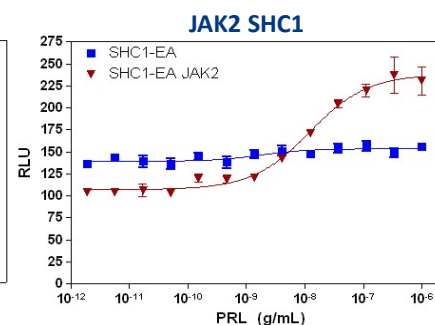
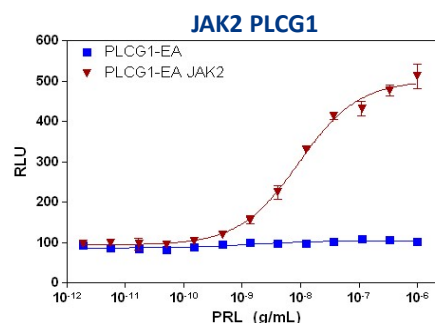
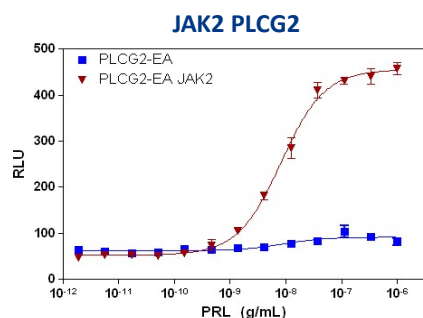
Characterize JAK-coupled cytokine receptor responses; identify JAK inhibitors

Growth Hormone Receptor examples

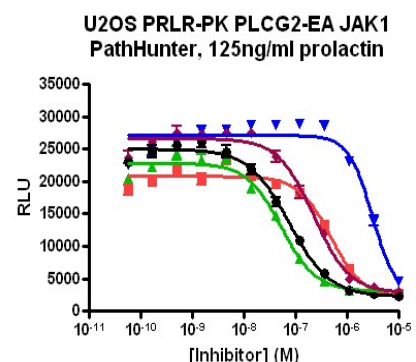
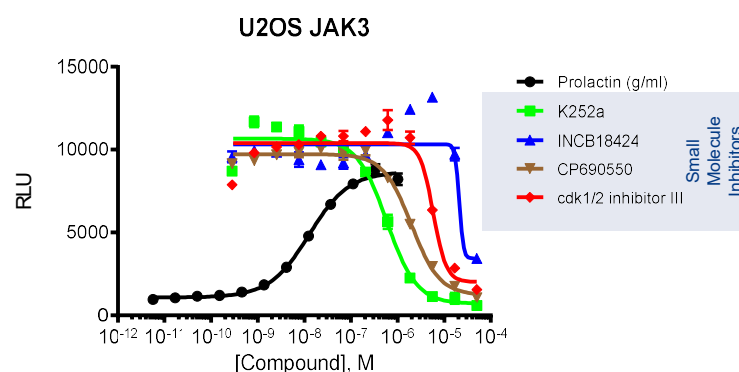
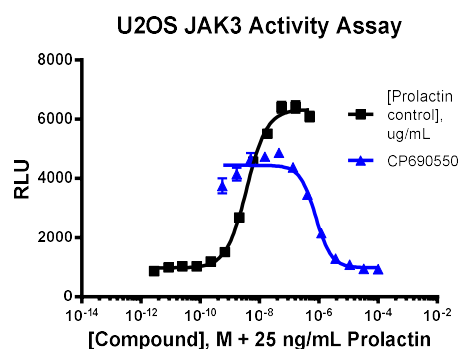


Characterize JAK-coupled cytokine receptor responses; identify JAK inhibitors

Prolactin Receptor examples



Red = + Jak2
 Blue = - Jak2

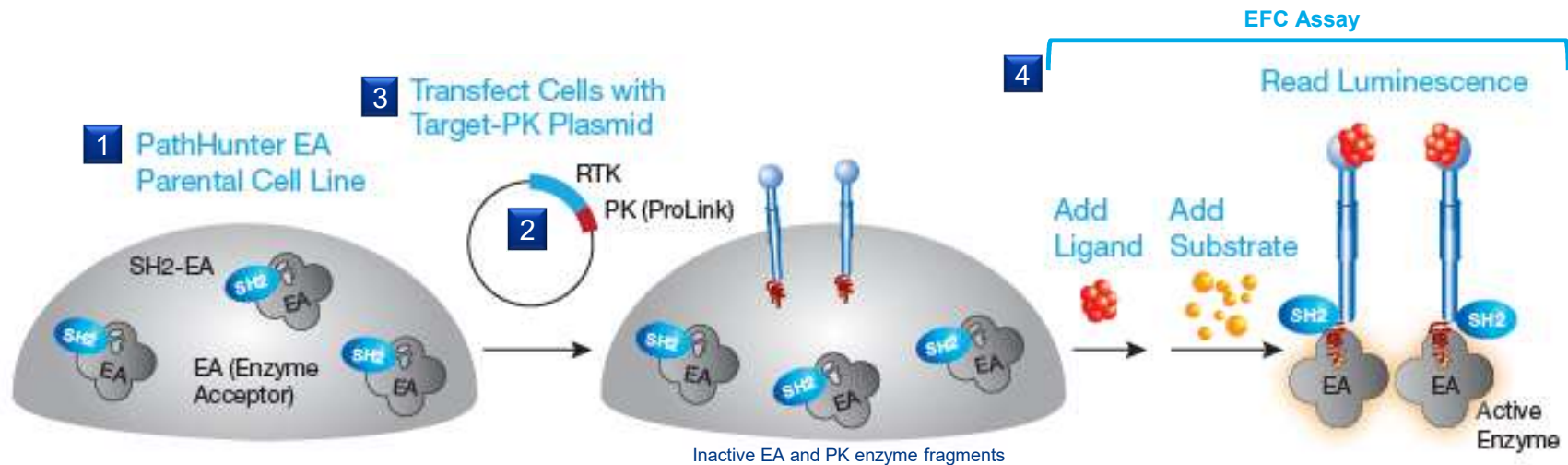


Inhibitor	Potency
AT-9283	+++
INCB18424	+++++
JNJ-28312141	+
CP-690550	++
AZD1480	++++
Agonist	ProLactin

+++++ = most potent
 + = least potent

Create Your Own PathHunter® Functional Assays

Workflow example for the development of an RTK cell-based functional assay



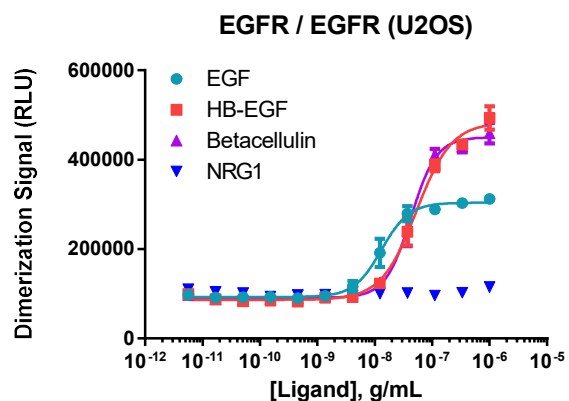
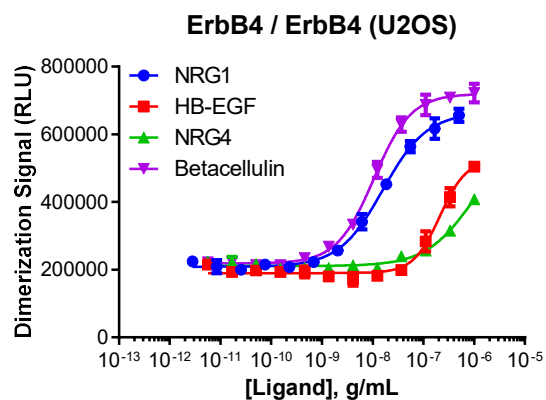
1. Start with an SH2-protein recruitment engineered parental cell line
2. Clone in protein of interest (e.g. RTK target) using PK (ED) cloning vector (plasmid)
 - Option to use a PK (ED) expression vector plasmid
3. Transfect plasmid into cells
4. Perform EFC assay (e.g. RTK functional assay) using ligand(s) and a PH Detection Kit (includes substrate)
 - Read chemiluminescent signal (obtain quantitative dose-response curve)

Similar workflow for translocation assays

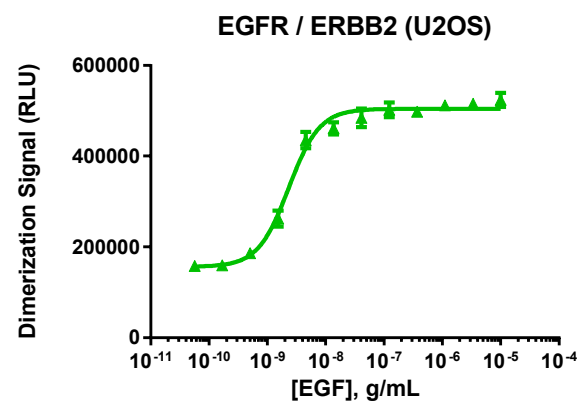
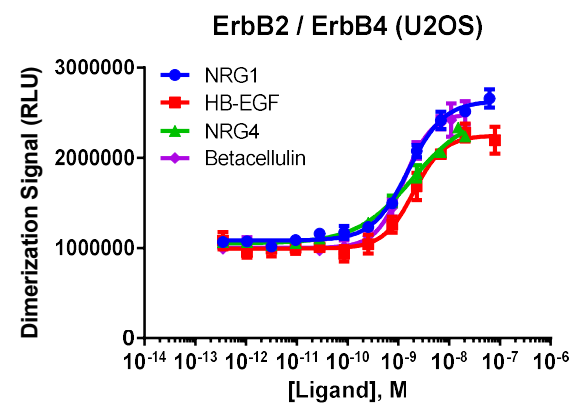
PathHunter® Cell-Based Dimerization Assays

Analyze ERB and EGF RTK dimerization

RTK Homodimers



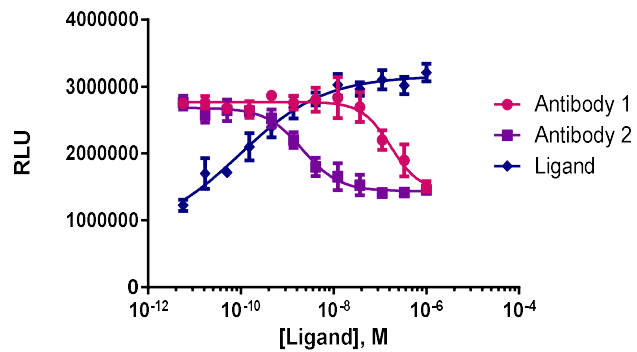
RTK Heterodimers



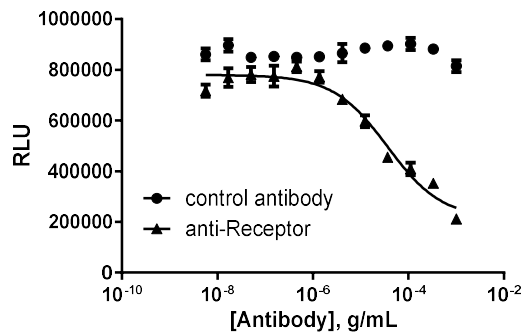
PathHunter® Cell-Based Dimerization Assays

Screen and profile anti-ligand and anti-receptor antibodies (biologics)

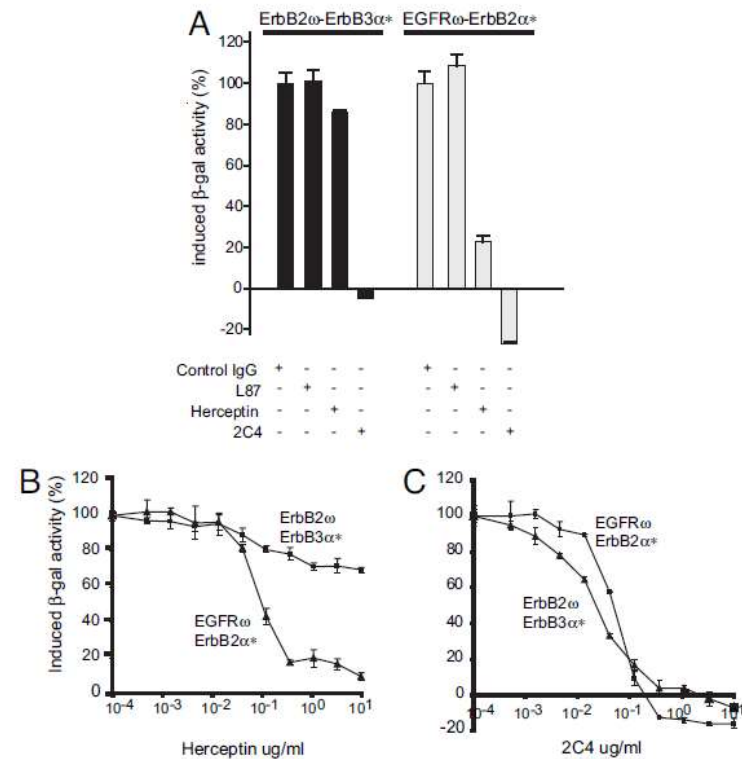
EGFR Homo-dimer Assay



TGFβR Hetero-dimer Assay



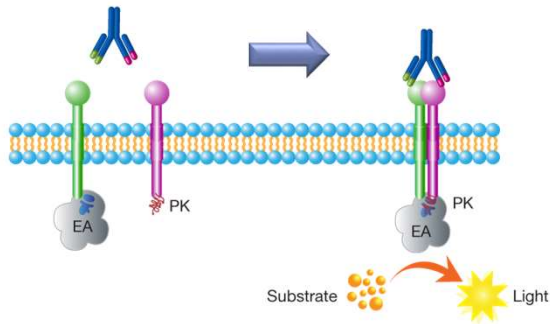
Dimerization assays demonstrate that Herceptin inhibits EGFR-ErbB2 not ErbB2-ErbB3*



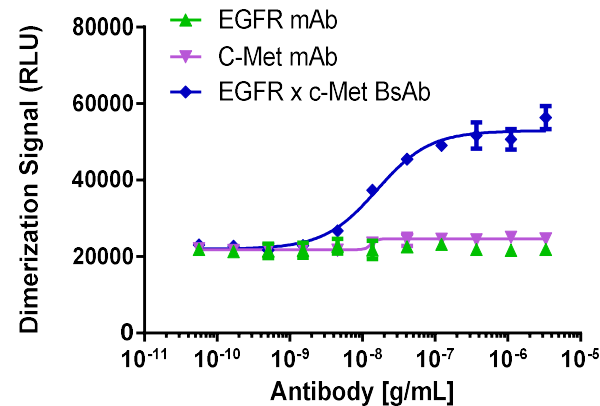
PathHunter[®] Cell-Based Dimerization Assays

Screen and optimize bispecific antibodies (biologics). Oncology relevant example.

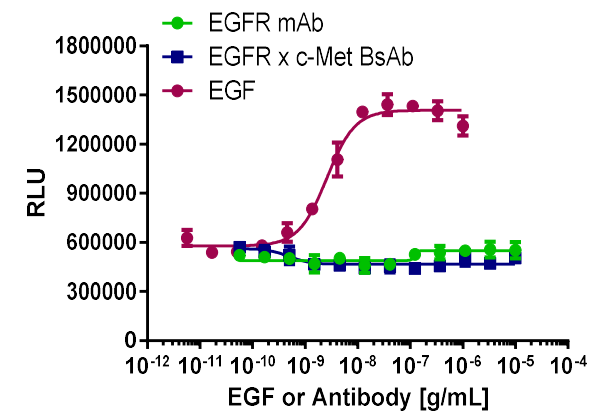
Bispecific Antibody-Induced Receptor Dimerization



EGFR/C-Met Bispecific Assay



EGFR Homodimer Assay



Assay developed to characterize bispecific antibodies

- Dimerization assay engineered to express two receptors (EGFR and c-Met) on same cell (e.g. mimic tumor model)
- Heterodimer model is specific for bispecific antibody (left panel); bispecific does not activate receptor homodimer assay (right panel)
- Published by S. W. Jarantow *et al.* in JBC (2015): Impact of Cell-surface Antigen Expression on Target Engagement and Function of an Epidermal Growth Factor Receptor × c-MET Bispecific Antibody

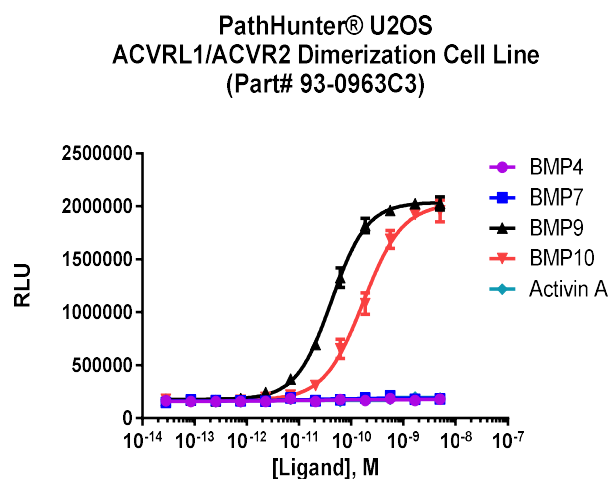
PathHunter® Cell-Based Dimerization Assays

Study other kinase receptors (example of serine/threonine kinase receptors)

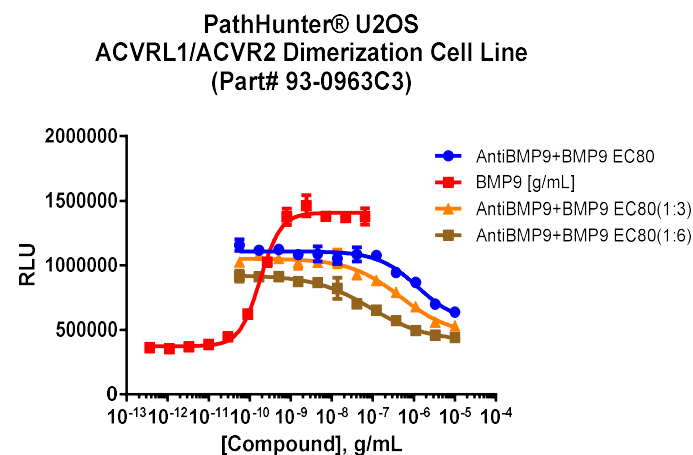
Serine/threonine-protein kinase receptors

- TGF β superfamily receptors (e.g. BMP and ACVRL1 receptors)
- Heterodimerize with activin-binding (e.g. ACVR2) receptors to form functional receptors
- Therapeutic targets in oncology and fibrosis

Specific Ligand Profiling



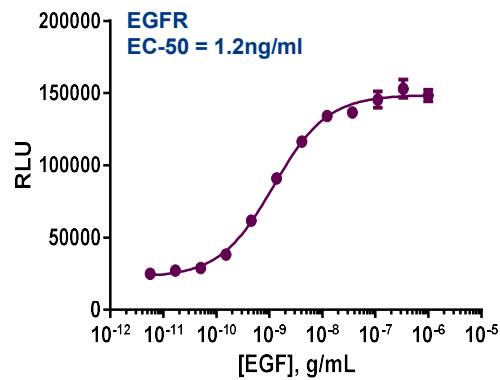
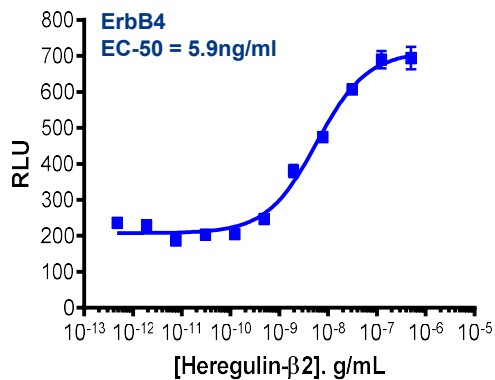
Antibody Profiling (anti-ligand, anti-receptor)



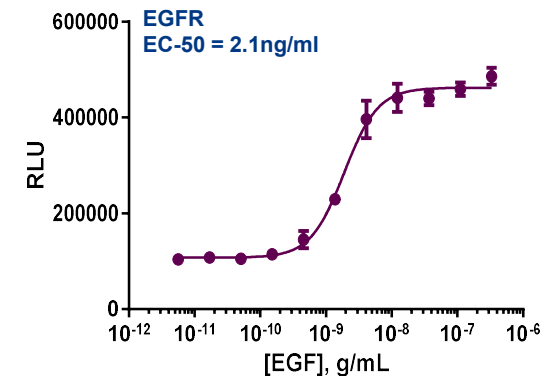
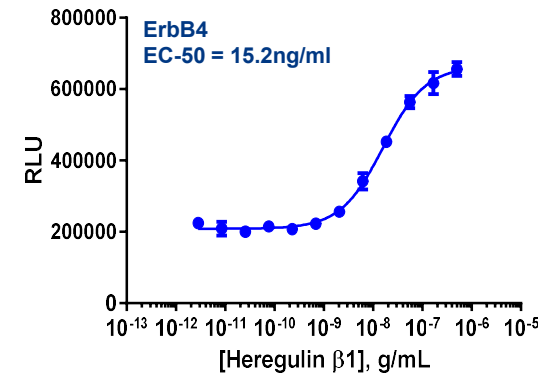
PathHunter® Cell-Based Assays for Kinase Receptors

Two assay formats provide receptor-proximal pathway responses at different levels (activation via adaptor recruitment vs. dimerization)

Functional Assays (Phosphorylation-dependent SH2-Recruitment Assays)



Dimerization Assays



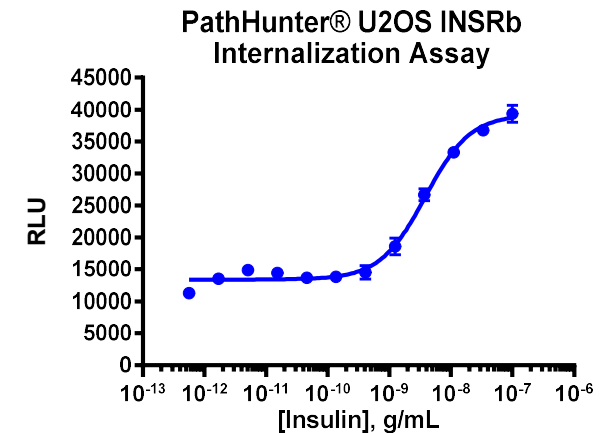
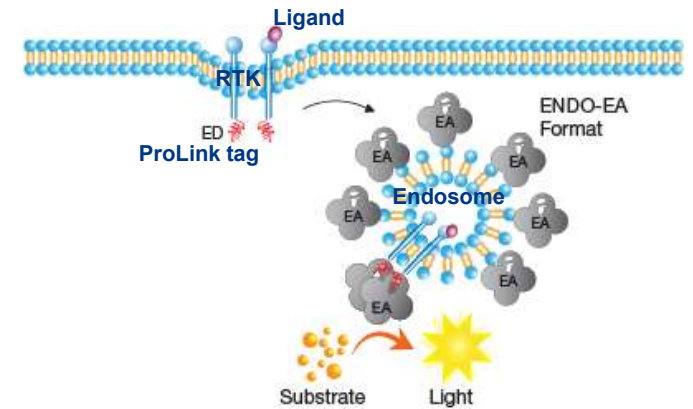
PathHunter® Cell-Based Translocation Assay

Measure internalization and trafficking of the RTKs

Provide a quantitative measurement of total endocytosed receptor protein using EFC

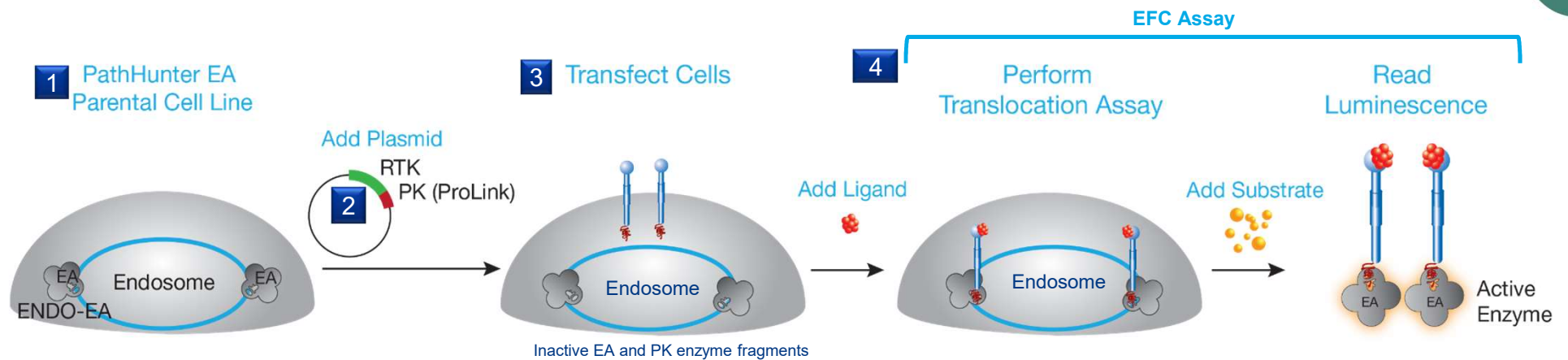
Example: Insulin receptor (e.g. INSRb)

- Cells are engineered to co-express the ProLink™ (PK)-tagged RTK and the Enzyme Acceptor (EA)-tagged endosome fusion proteins
- Activation of the RTK results (e.g. insulin) in internalization of the receptor and trafficking to cellular endosomes, where the two enzyme fragments complement and form a functional β -gal enzyme that is capable of hydrolyzing substrate and generating a chemiluminescent signal using the PathHunter Detection kit



Create Your Own PathHunter® Translocation Assays

Workflow example for the development of an RTK cell-based translocation assay



1. Start with an ENDO-EA engineered parental cell line
2. Clone in protein of interest (e.g. RTK target) using PK (ED) cloning vector (plasmid)
 - Option to use a PK (ED) expression vector plasmid
3. Transfect plasmid into cells
4. Perform EFC assay (e.g. translocation) using ligand(s) and a PH Detection Kit (includes substrate)
 - Read chemiluminescent signal (obtain quantitative dose-response curve)

Similar workflow as functional assays creation

PathHunter® cell-based assays for kinase receptors

Assay Highlights

- **Accurate, Sensitive and Reproducible** - Superior quality, reproducible data with large assay windows, and robust performance
- **Broadly Applicable** - Identify different ligand classes (small molecules and biologics, monospecific vs bispecific, agonist vs antagonist) and study hetero- and homodimers
- **High Specificity** - Tagged, full-length tyrosine kinase eliminates non-specificity from endogenous tyrosine kinases
- **Easy-to-Use** – Simple, one-step add and read protocol that is HTS-friendly
- **Universal, DIY Solution** – Create your own quantitative cell-based assays to study translocation or function of RTK/CTKs

Accessing kinase assays and solutions

[Kinase solutions page](#)

[Kinase targets](#)

[Kinase cell-based assays](#)

[Dimerization assays](#)

[Translocation assays](#)

[Kinase profiling services](#)

[Toolbox \(DIY\) solutions](#)