



# Accelerating Immunotherapy Drug Development with Simple Cell-Based Assays for Immune Checkpoint Receptors

Jennifer Lin-Jones, Ph.D.

Senior Group Leader, Assay Development

OUR EXPERTISE  
IN YOUR HANDS.  
DISCOVER  
CONFIDENTLY.

October 2021

# Eurofins DiscoverX

## Strong Foundation | Technical Expertise | End-to-End Solutions



### Dedicated Operations

Supporting programs from Research, Discovery to Lot Release

- **Products Division headquartered in Fremont, CA**
- Additional sites in Missouri, USA and Poitiers, France
- **800+ off-the-shelf assays for in-house development**
- Over 10,000 customers in NA, APAC and EMEA

### Deep Domain Expertise

Over 45 years of cumulative technical experience in

- **Cell line engineering & characterization**
- **Bioassay development, optimization & qualification**
- Analytical Cell Banks
- Membrane Preps and Frozen Assay Ready Cells
- Bulk Enzyme Production

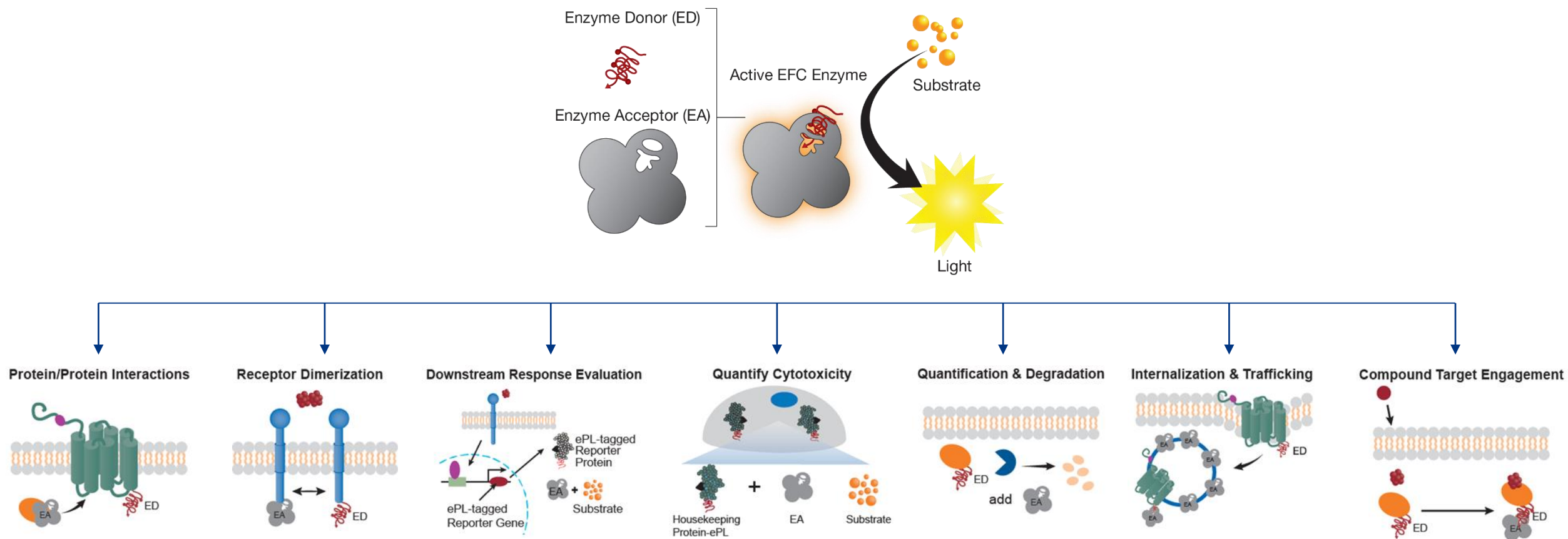
### Established Brand

Successfully implemented at global Pharma, Biotech & CRO

- **Products implemented in discovery & development**
- Over 50 billion data points screened
- **2,000+ publications**
- Several active Biotech/CRO-partnered programs
- **Implemented in lot release of several marketed biologics**

# Enzyme Fragment Complementation (EFC) Robust Platform for Cell-Based Assays

*Split  $\beta$ -galactosidase reporter system can be engineered to generate target-specific, homogeneous cell-based assays for immuno-oncology therapeutics*

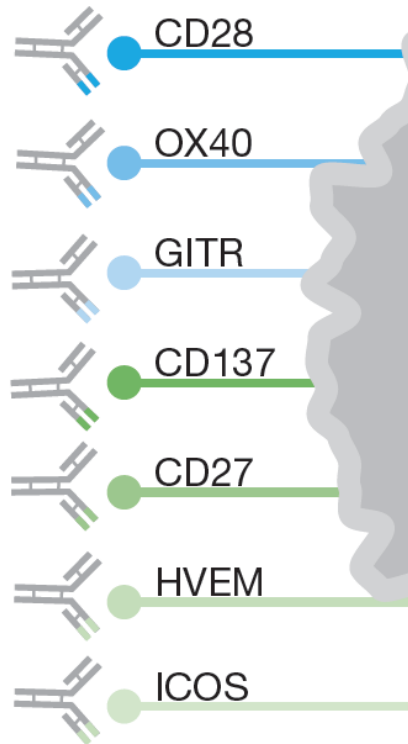


# Targeting T-Cell Co-Stimulatory and Inhibitory Checkpoint Receptors

*Tools are needed to screen for and develop new therapeutics*

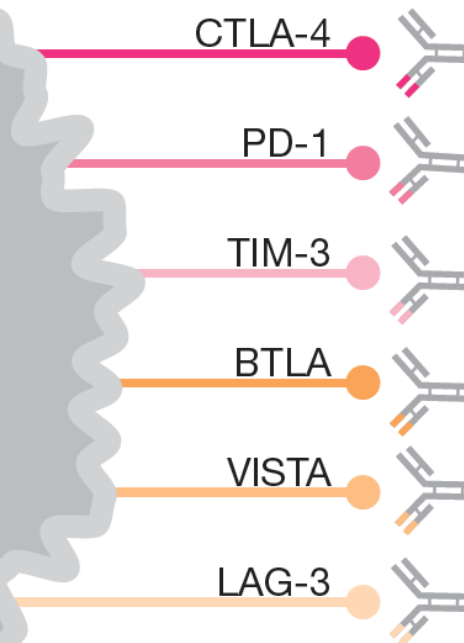
**Push the gas pedal  
on T cell activation  
to stimulate the  
immune system**

## Co-Stimulatory Receptors



Agonistic antibodies

## Co-Inhibitory Receptors



Blocking antibodies

**Remove the brakes  
inhibiting T cell  
activation to  
stimulate the immune  
system**

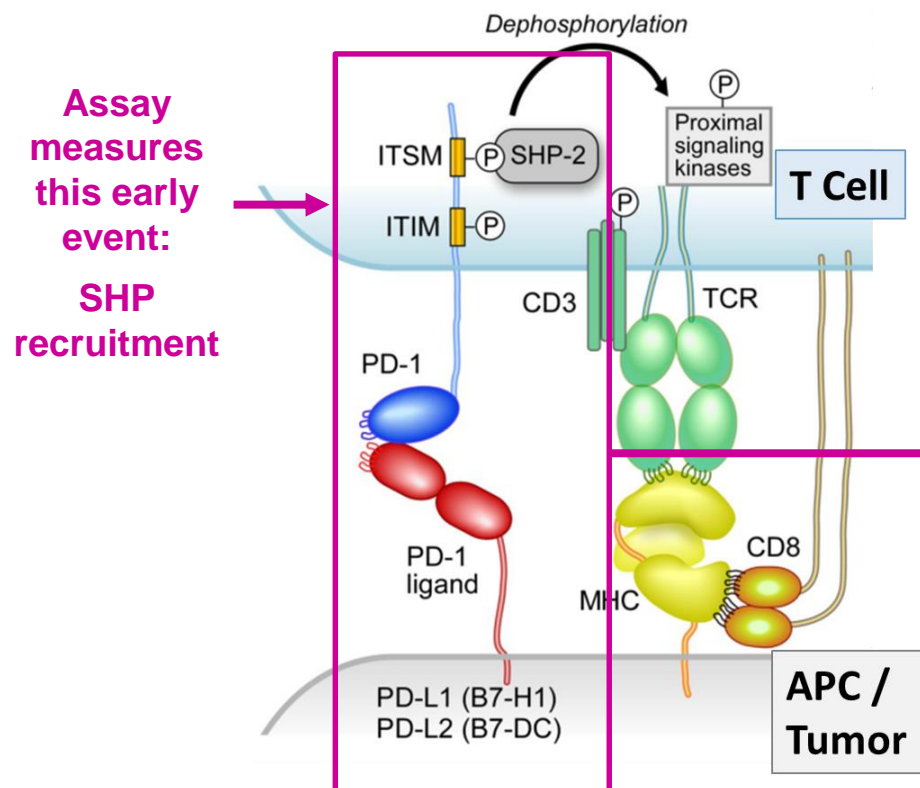
**Modulate immune response to destroy cancer cells**



# PD-1 Signaling Assay Concept

Quantify early step in PD-1 mediated inhibition of T cell activation: SHP recruitment

## Mechanism of Action



## Assay Design

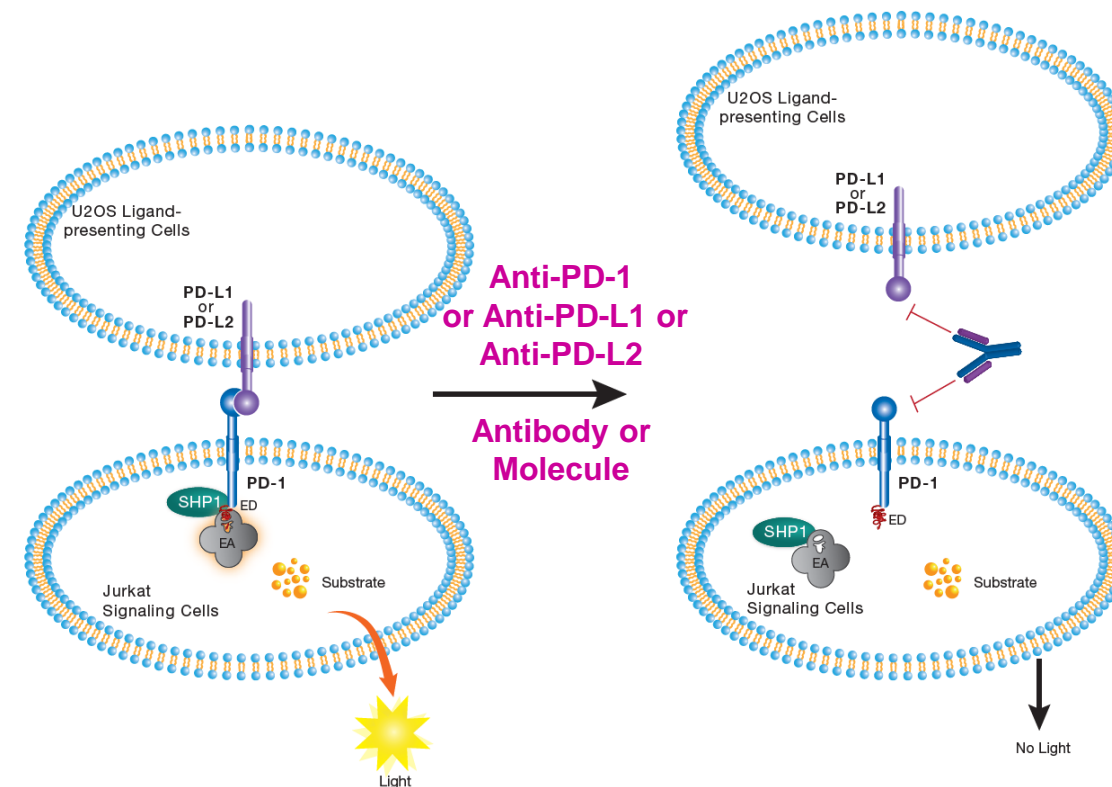
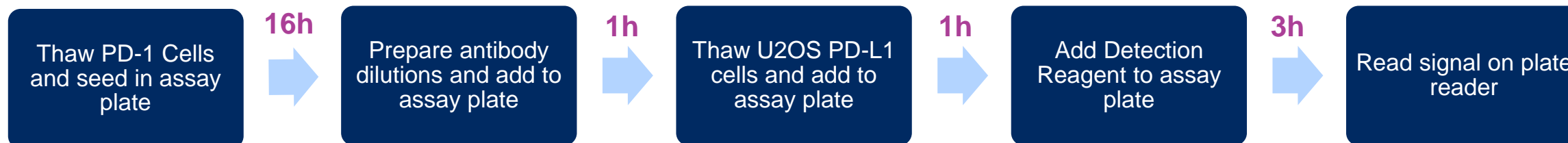


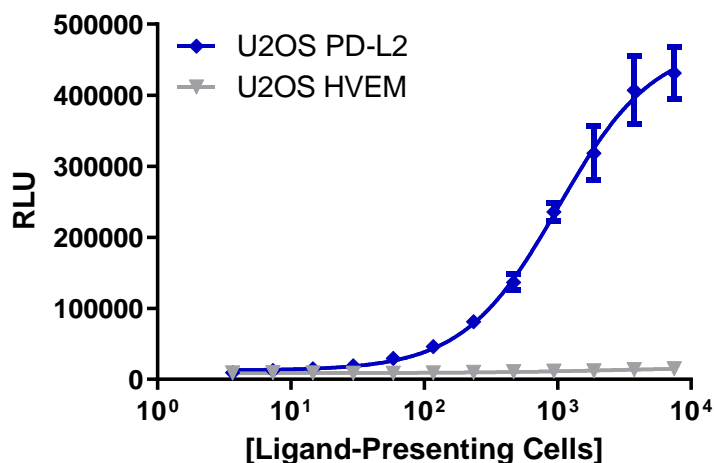
Figure from Science Webinar Series, Part 5: Gordon J. Freeman, Ph.D.

# PD-1 Signaling Assay is Highly Specific and Stability-Indicating

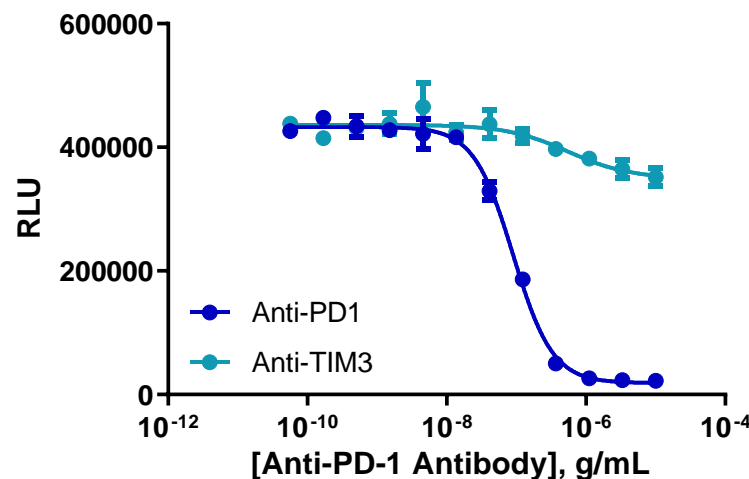
## Assay-Ready (RTU) Protocol



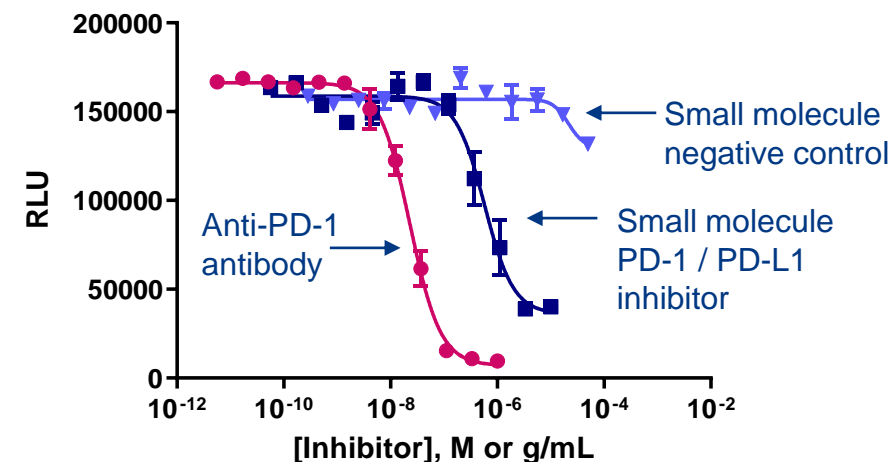
### Specificity (Agonist)



### Specificity (Antagonist)

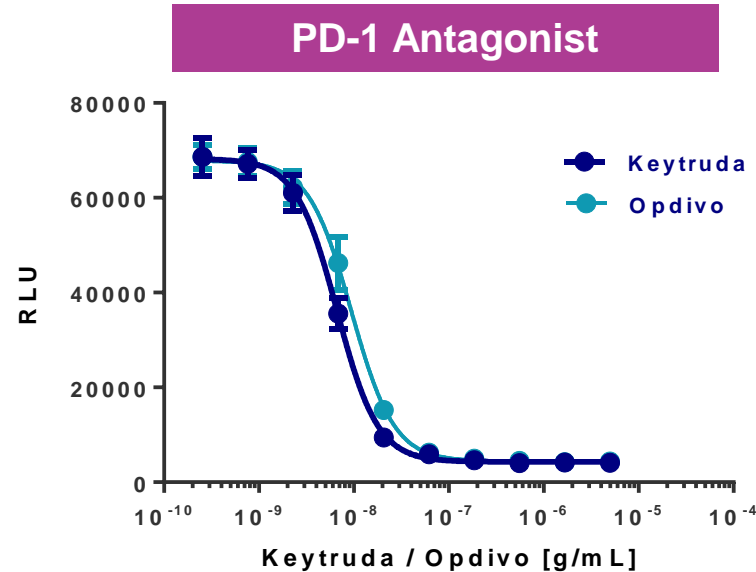


### Suitable for Small Molecules or Biologics

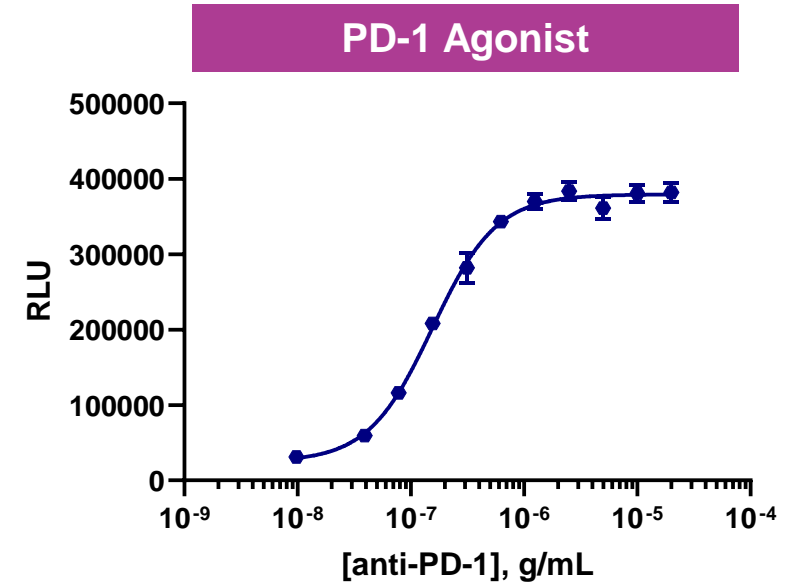


# PD-1 Signaling Assay Suitable for Therapeutic Development of Both PD-1 Agonists and Antagonists

## Immuno-oncology

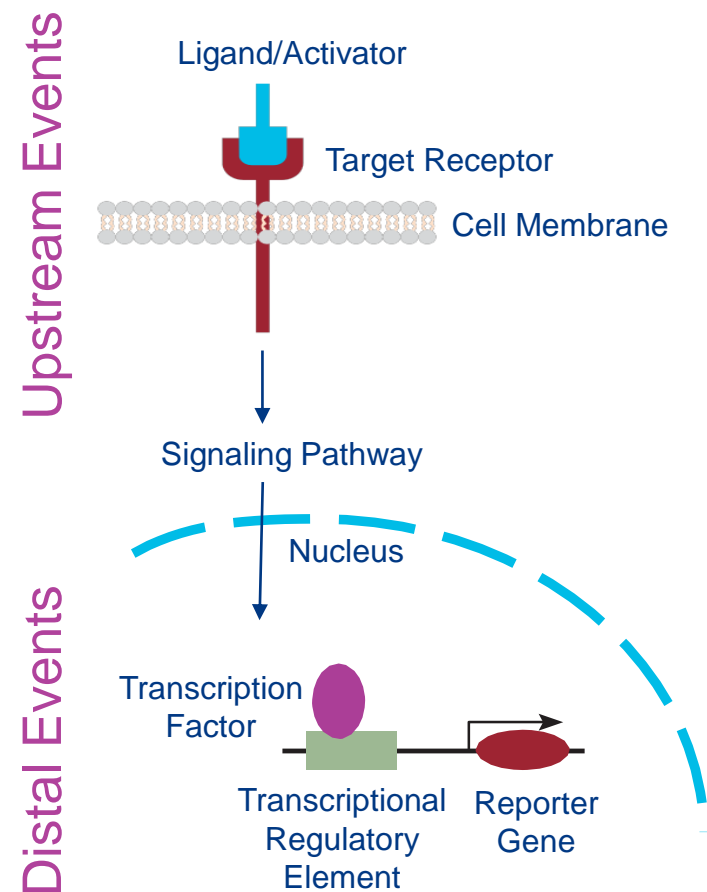


## Auto-immunity



*PD-1 Signaling assay is a versatile screening tool for PD-1 antagonists (I/O) or for PD-1 agonists (novel therapies for autoimmunity). The latter data were generated in co-culture with Fcγ receptor expressing cells.*

# Complementary Cell-Based Assays for Comprehensive Understanding of a Drug Molecule's MOA

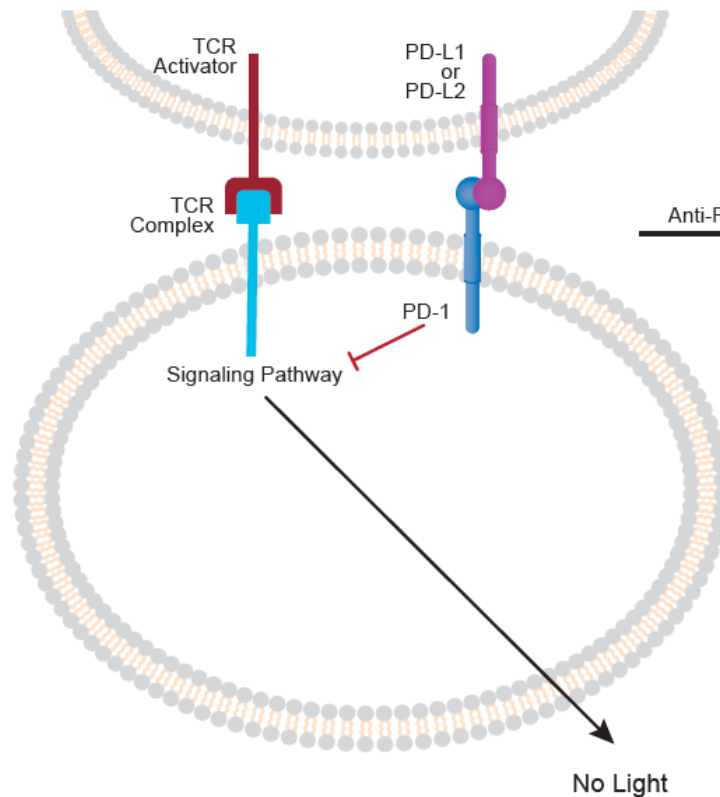


Cell-based Signaling Pathway Assays	
Upstream (Receptor-Proximal) Readout (Non-Reporter Gene Assays)	Downstream (Distal) Readout (Reporter Gene Pathway Assays)
Represent a readout more proximal to receptor activation and upstream of signal transduction and transcriptional changes.	Represent a distal pathway readout reflective of a phenotypic endpoint (e.g. proliferation)
Measure status of a specific functional event upstream of transcription	Measure changes in transcription that affect reporter protein expression
Results may more directly reflect drug MOA: measures a specific functional event	Results can integrate effects from multiple inputs on a signaling pathway: sum of downstream signaling events
Produce modest to large assay windows, with short assay times and low variability	



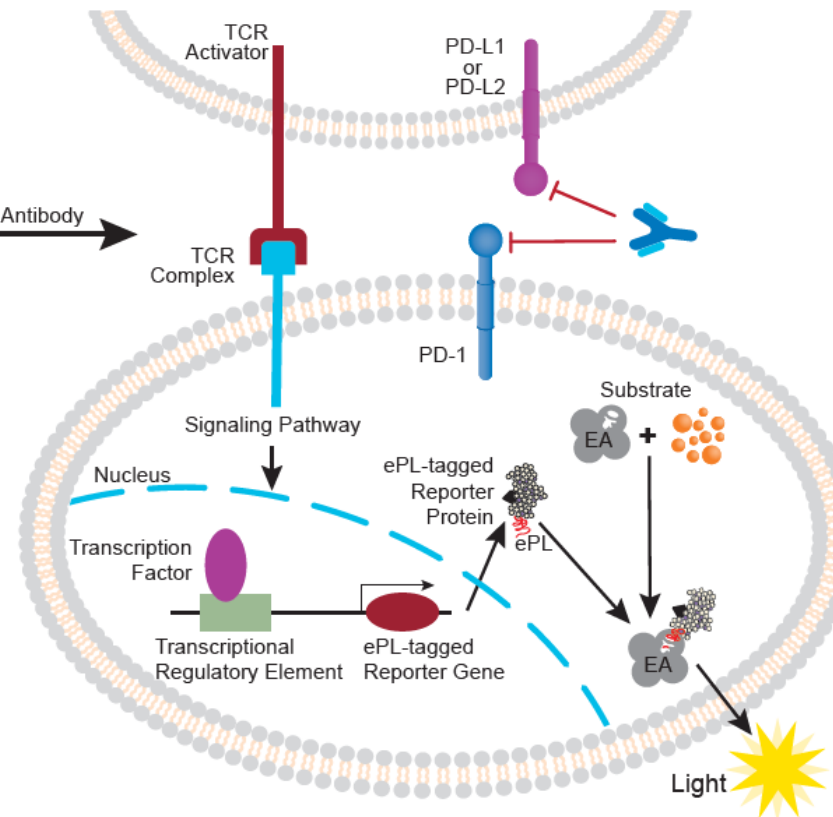
# Monitor PD-1 Inhibitor Activity Downstream of TCR Activation: Assay Principle

**U2OS PD-L1/TCR Activator cells +  
Jurkat PD-1-NFAT**



- TCR activation inhibited by PD-1 activation
- Reduction in reporter protein expression

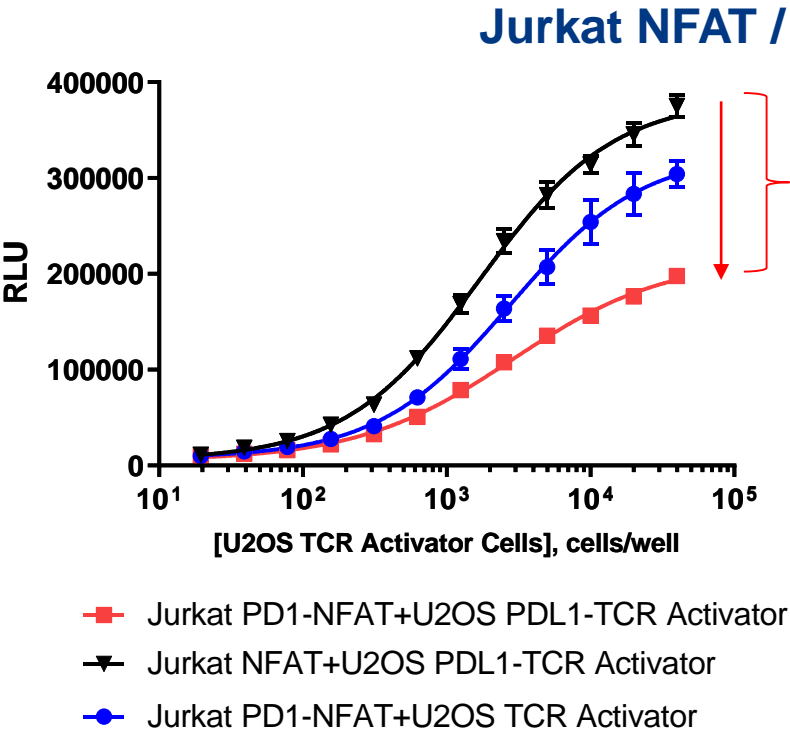
**U2OS PD-L1/TCR Activator cells +  
Jurkat PD-1-NFAT + anti-PD-1 Ab**



- Antibody inhibits PD-1 activation
- PD-1 inhibition of T cell activation released
- Increase in reporter protein expression

# PathHunter® Jurkat PD-1-NFAT Pathway Reporter Assay: Evaluate PD-1 Checkpoint Receptor Mediated Inhibition of T Cell Activation

- PD-1 inhibitory receptor introduced into PH Jurkat NFAT cells attenuates T cell activation
- PD-1 checkpoint receptor inhibition of T cell activation can be monitored by PD-1 receptor co-expression in Jurkat NFAT reporter cells in co-cultures with U2OS TCR activator cells co-expressing PD-L1 ligand



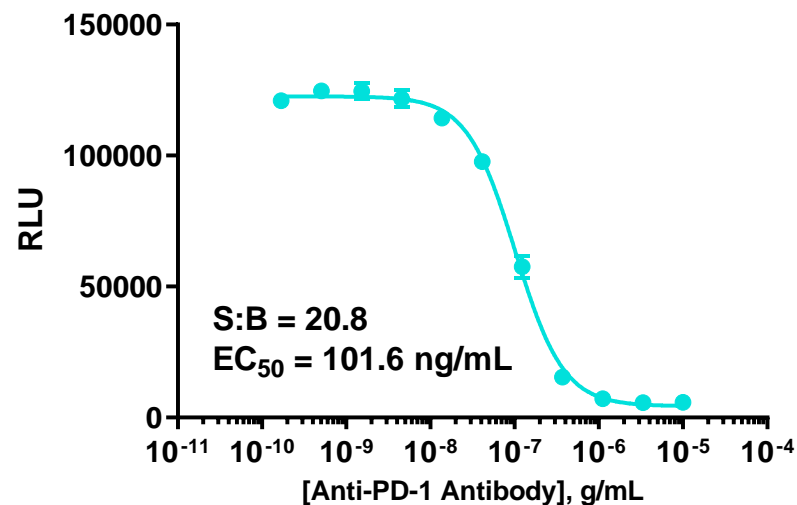
Decreased T cell activation when PD-1 is co-expressed in Jurkat NFAT cells and PD-L1 co-expressed with TCR activator in U2OS

	U2OS TCR activator	
Jurkat NFAT reporter	-PD-L1	+PD-L1
-PD1	(+++)	+++
+PD-1	++	+

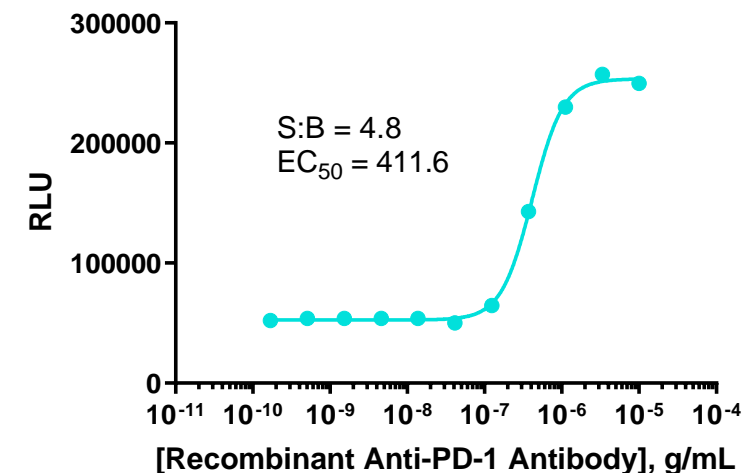
# Measure Proximal & Distal Events in the PD-1 Signaling Pathway

- Both PathHunter® PD-1 assays types provide a comprehensive understanding of the PD-1 pathway and the MOA of the anti-PD-1 Antibody
- Assays are robust and measure PD-1 inhibition with sensitive responses from an event either **proximal** (PD-1 SH2 recruitment, Jurkat PD-1 Signaling) to receptor activation, or **distal** (NFAT-regulated reporter, Jurkat PD-1-NFAT)

**PathHunter Jurkat PD-1 Signaling Assay**  
(93-1106C19)



**PathHunter Jurkat PD-1 NFAT Reporter Assay**  
(93-1141C19)

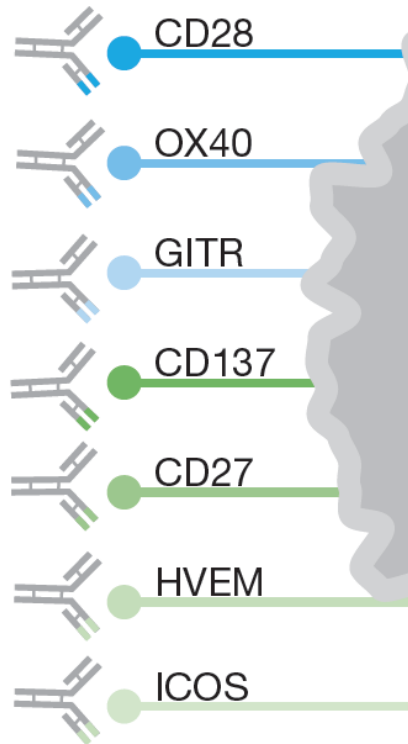


# Targeting T-Cell Co-Stimulatory and Inhibitory Checkpoint Receptors

*Tools are needed to screen for and develop new therapeutics*

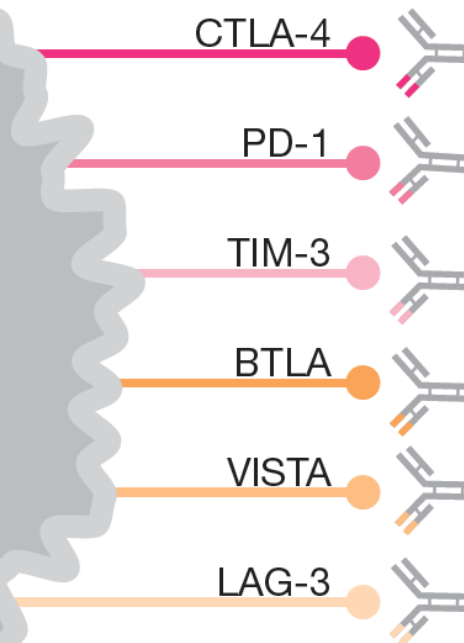
**Push the gas pedal  
on T cell activation  
to stimulate the  
immune system**

## Co-Stimulatory Receptors



Agonistic antibodies

## Co-Inhibitory Receptors



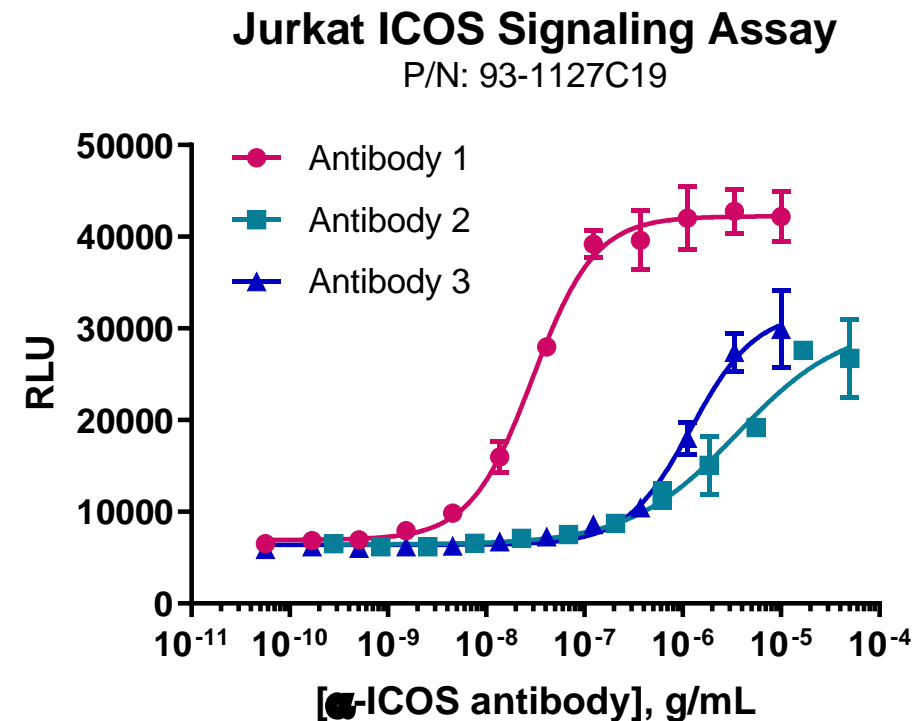
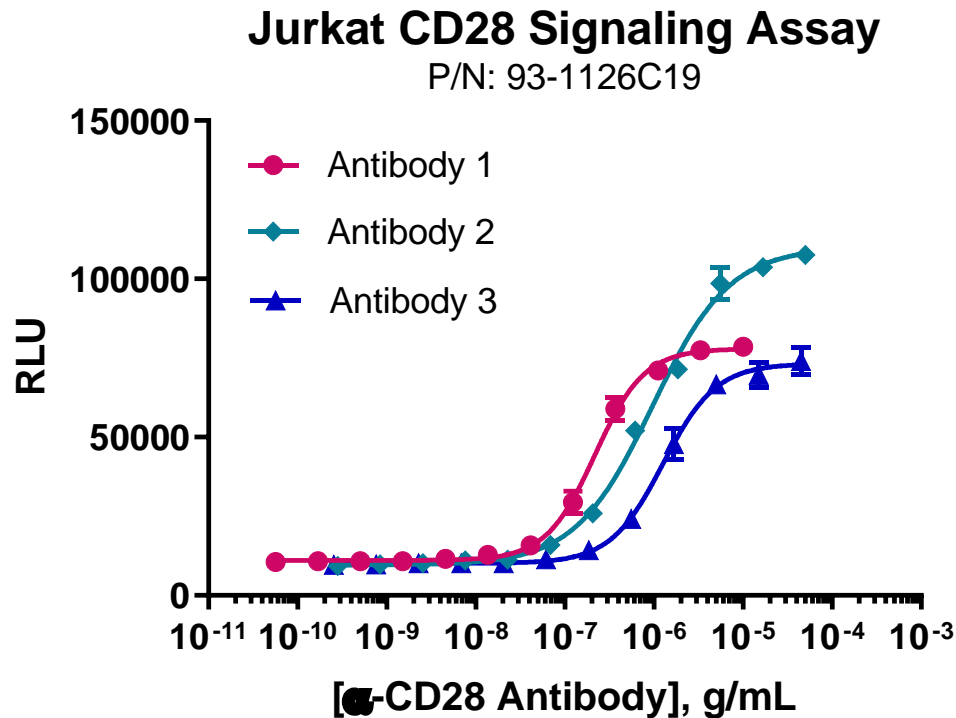
Blocking antibodies

**Remove the brakes  
inhibiting T cell  
activation to  
stimulate the immune  
system**

**Modulate immune response to destroy cancer cells**

# CD28 and ICOS Co-Stimulatory Receptors Also Signal Through SH2- Domain Containing Proteins

*Assays measure ligand-mediated recruitment of Grb2 (CD28) or p85 (ICOS)*

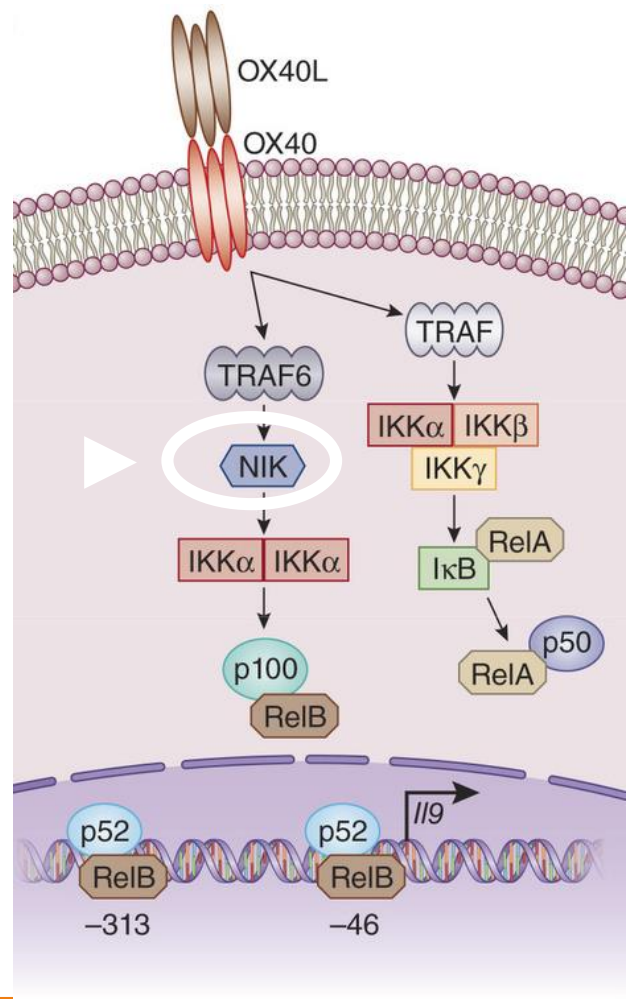


*Robust Assays for CD28 and ICOS Co-Stimulatory Receptors Enable Rank Ordering of Agonist Antibodies*

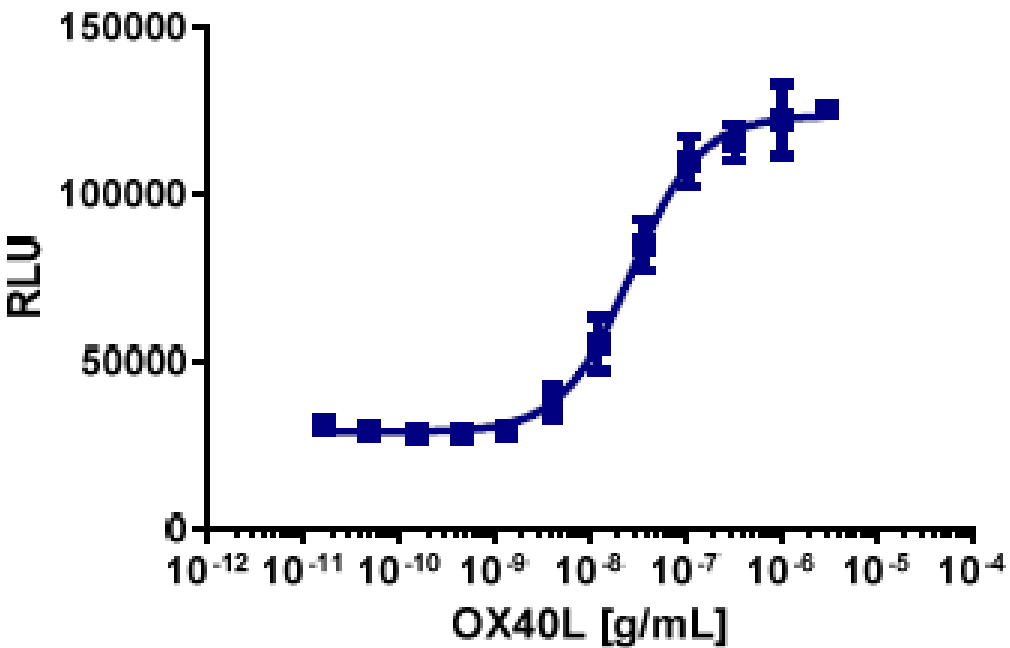


# PathHunter® OX40 Assay – Quantifies NIK Stabilization in Response to Pathway Activation

## Mechanism of Action



## OX40 Signaling Assay ([93-1080C3](#))

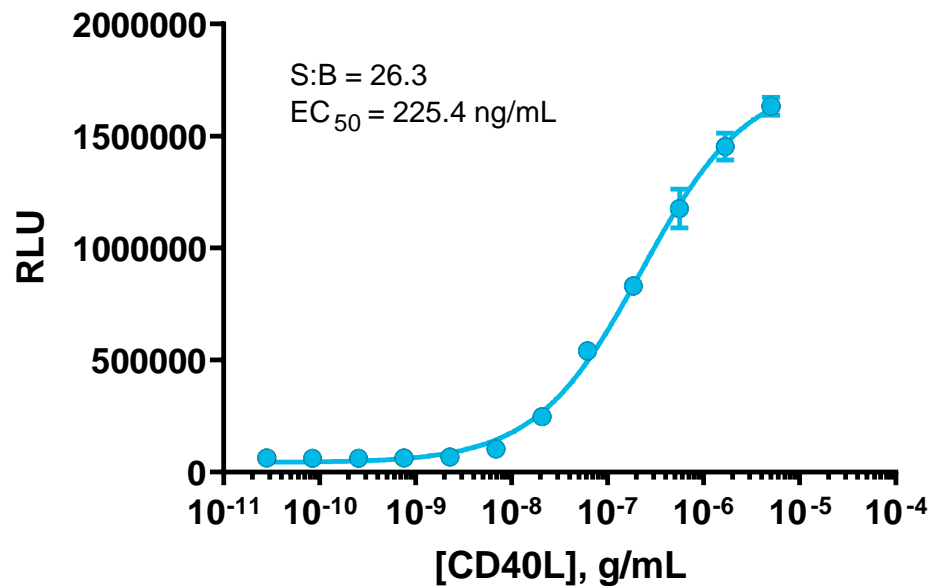


EC <sub>50</sub>	27.4 ng/mL
S:B	4
Assay Time	5 h

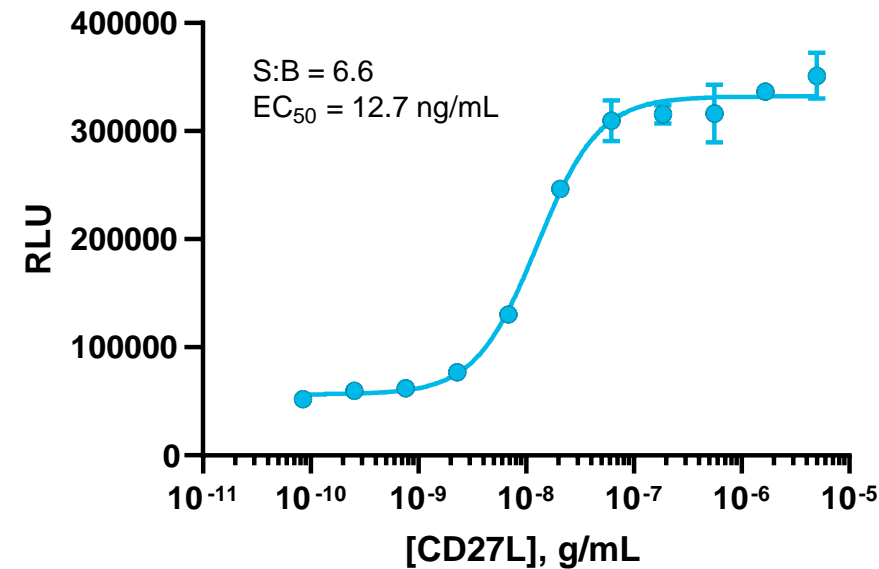
# PathHunter® Pathway Reporter Assays – Measure Pathway activation using Endogenous or Heterologously-Expressed Receptors

## Assays for Co-stimulatory TNFR superfamily receptors

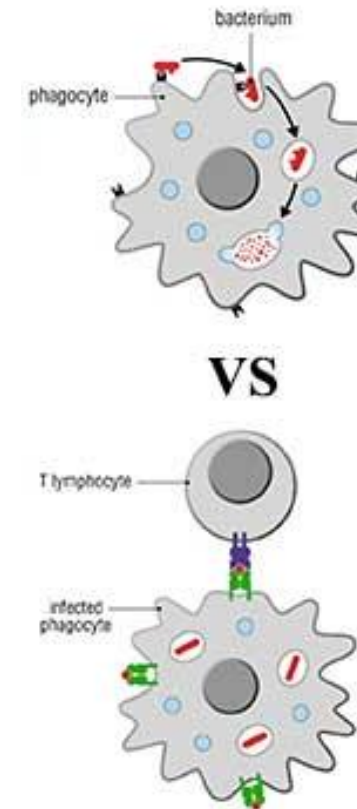
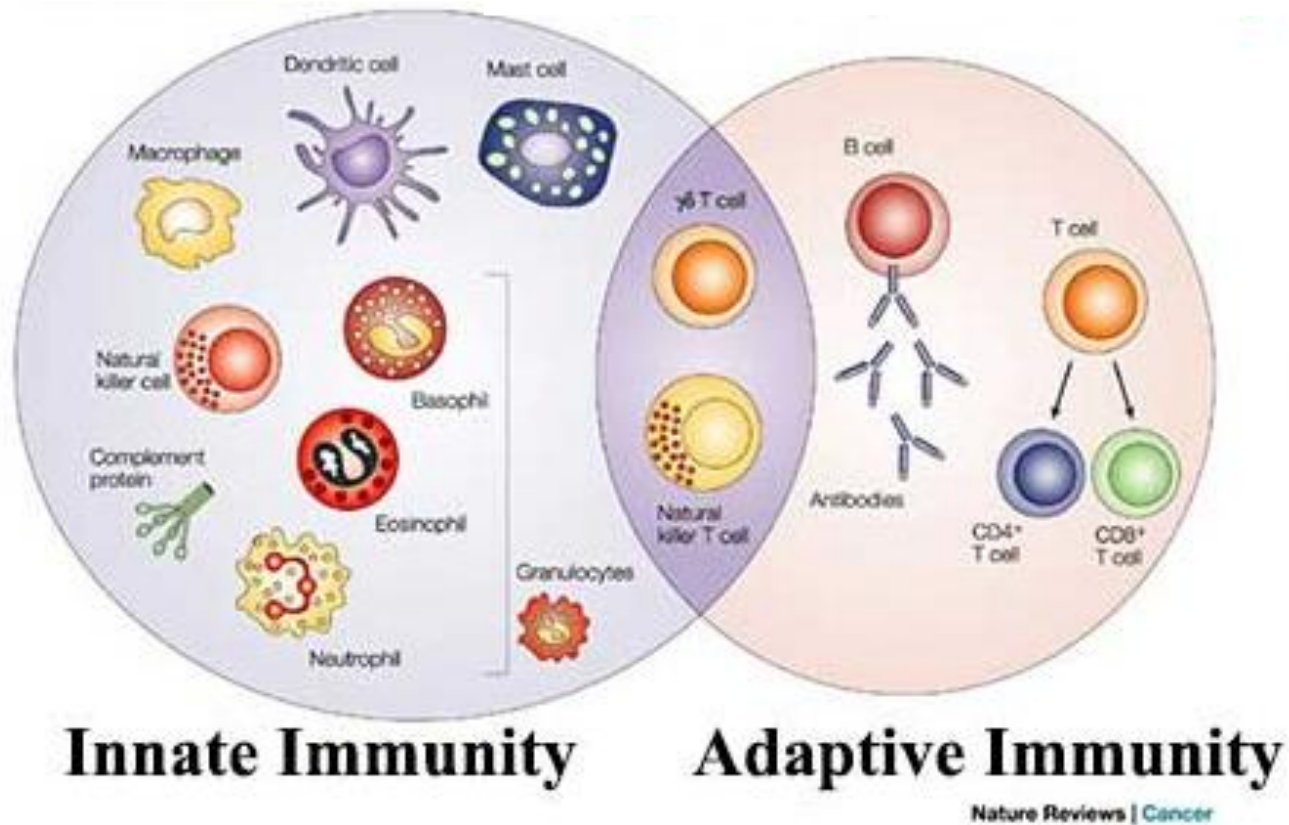
**PathHunter U2OS NF- $\kappa$ B  
Pathway Reporter Assay**



**PathHunter HEK293 CD27-NF- $\kappa$ B Pathway Reporter Assay**



# Immunotherapy Agents: Targeting Innate vs Adaptive Immunity



## Therapeutic Modalities

TLR Agonists  
STING Agonists  
**SIRPα / CD47**

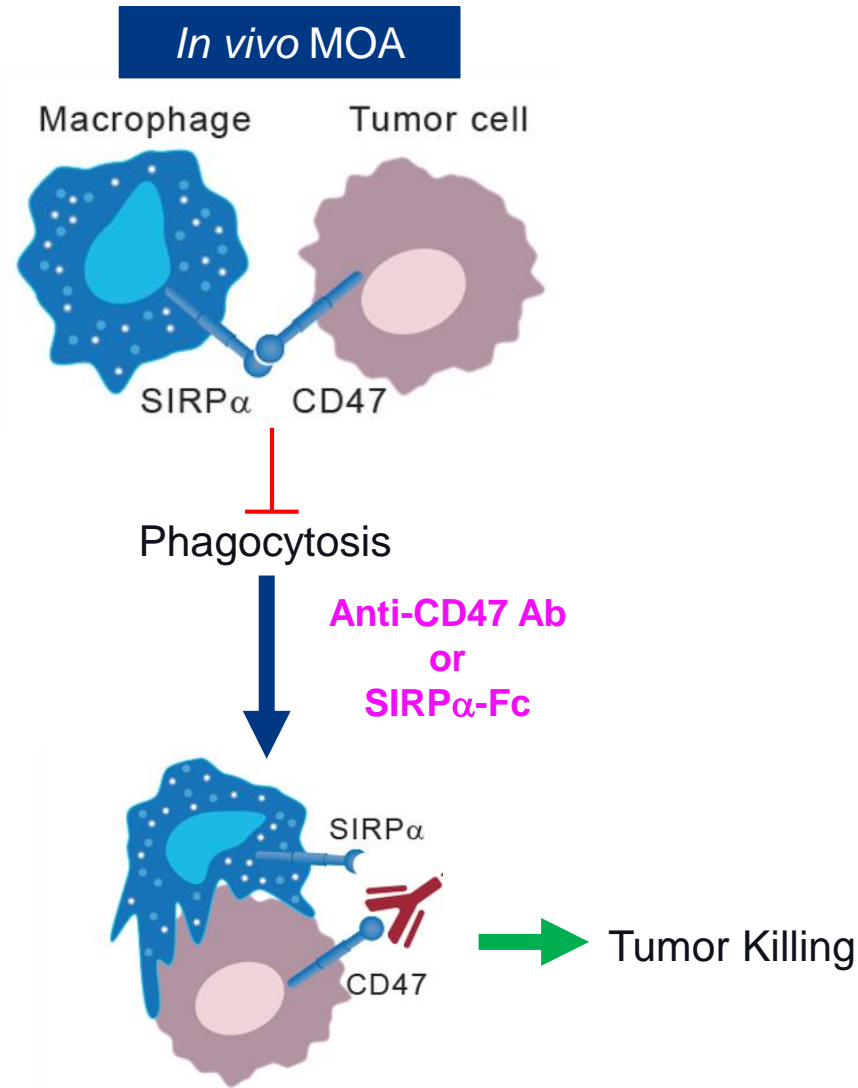
Checkpoint Inhibitors  
(**anti-PD-1/PD-L1**)

Checkpoint agonists  
(**OX40, CD137, ICOS**)

BiTEs, TRIKEs, etc

# The SIRP $\alpha$ / CD47 Axis

## An Innate Immune Checkpoint



- SIRP $\alpha$  is an inhibitory receptor expressed on macrophages and dendritic cells that promotes phagocytosis of foreign objects
- CD47, the ligand for SIRP $\alpha$ , is expressed on nearly all cells, but is significantly up-regulated in many tumor types, especially hematological malignancies such as AML and MDS
- ‘Don’t eat me’ signal that represses signaling via SIRP $\alpha$ , preventing myosin-IIA accumulation at the phagocytic synapse, leading to inhibition of phagocytosis
- Blocking the CD47 / SIRP $\alpha$  axis (e.g. with anti-CD47 antibodies, engineered receptor decoys, anti-SIRP $\alpha$  antibodies and bispecific agents) promotes tumor killing
  - phagocytosis of the tumor
  - Anti-CD47 blockade has also been shown to enhance adaptive immunity (e.g. prime an anti-tumor cytotoxic T cell response)

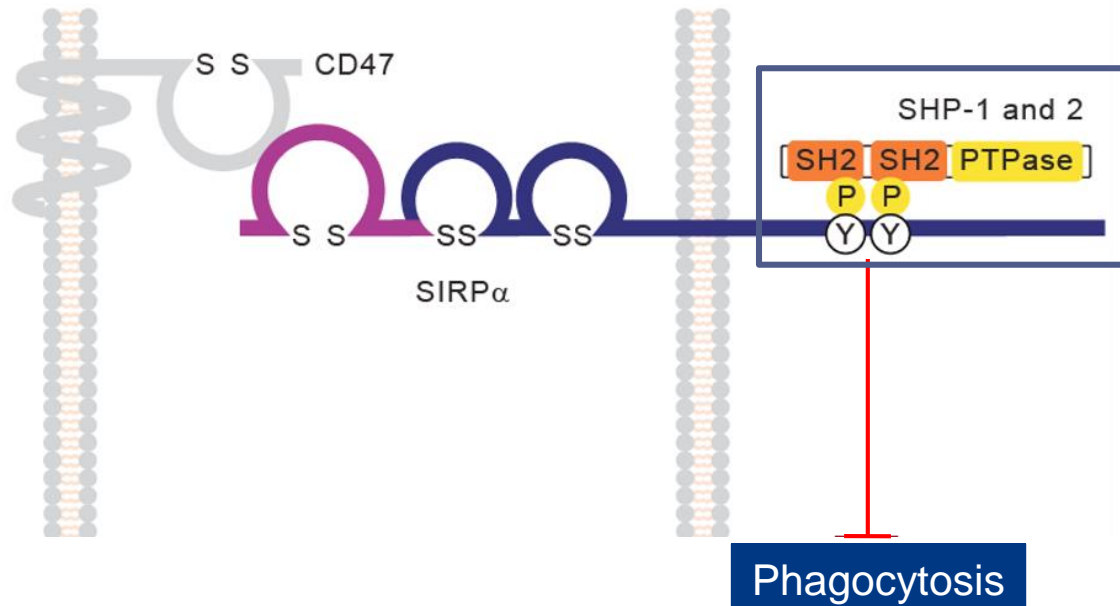
# PathHunter® SIRPα Signaling Assay: Assay Concept

*Co-culture SHP recruitment model based on  $\beta$ -galactosidase enzyme fragment complementation*

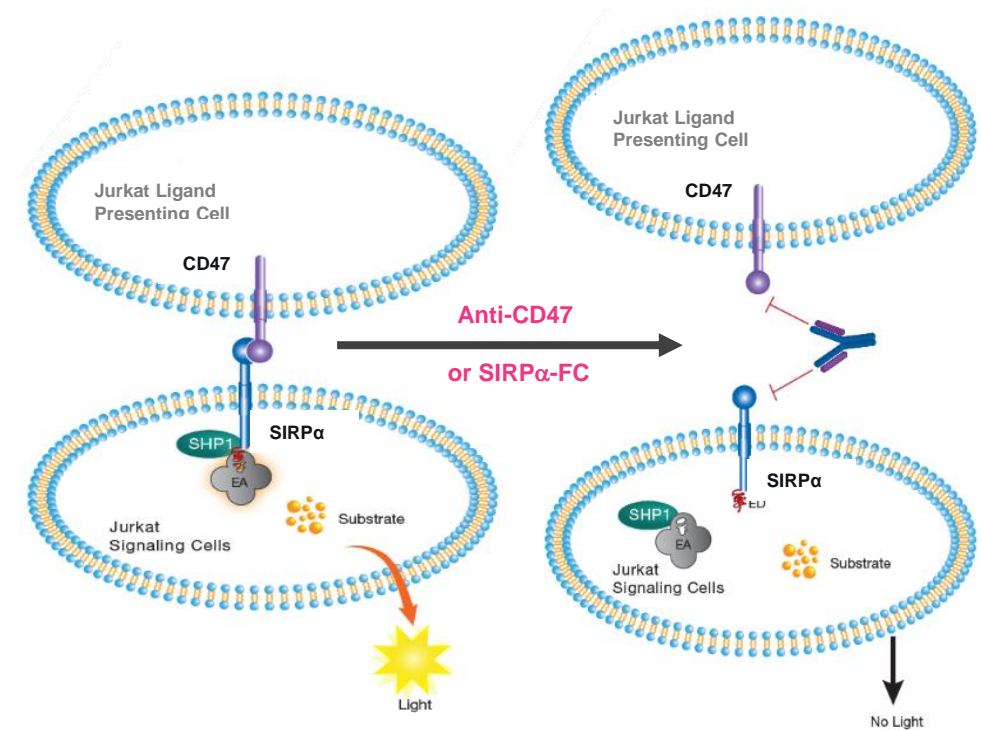
## Molecular MOA

### Tumor Cell

### Macrophage



## Co-Culture SHP Recruitment Model



Adapted from Trends in Cell Biology, 2008. Vol 19, No. 2

Assay quantifies ligand-induced recruitment of SHP-1 to ITIM motifs in C-terminal tail of SIRPα in response to phosphorylation

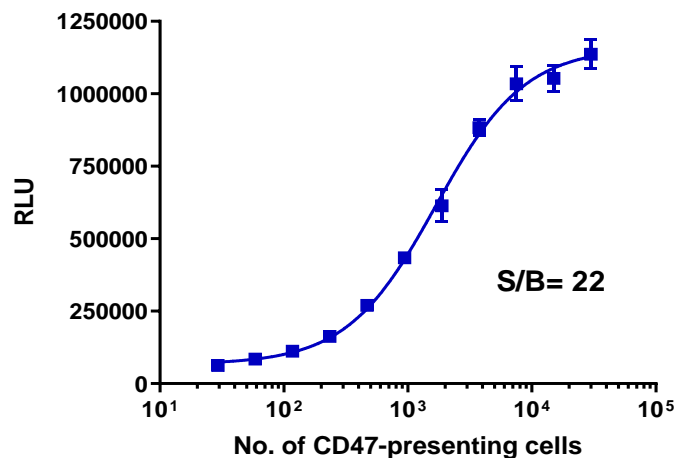


# PathHunter® SIRPα (CD47) Signaling Assay

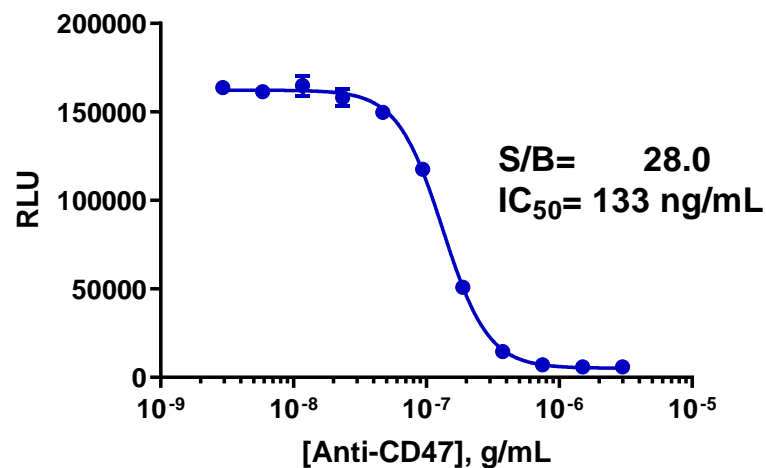
*Co-culture model with stable surface expression of SIRPα and a stable functional response over 45+ passages*

## Cell-Presented Ligand

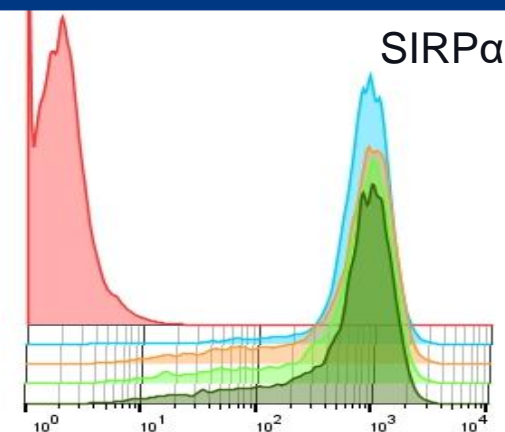
Agonist



Antagonist

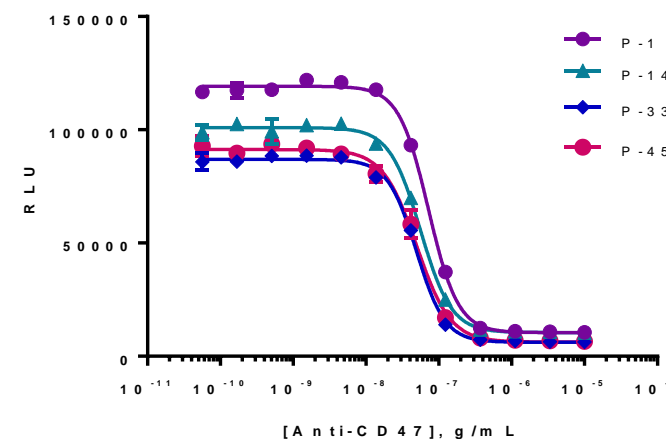


## Stable SIRPα Surface Expression



SIRPα expression varies by <20% RSD over 45 passages

## Stable Functional Response

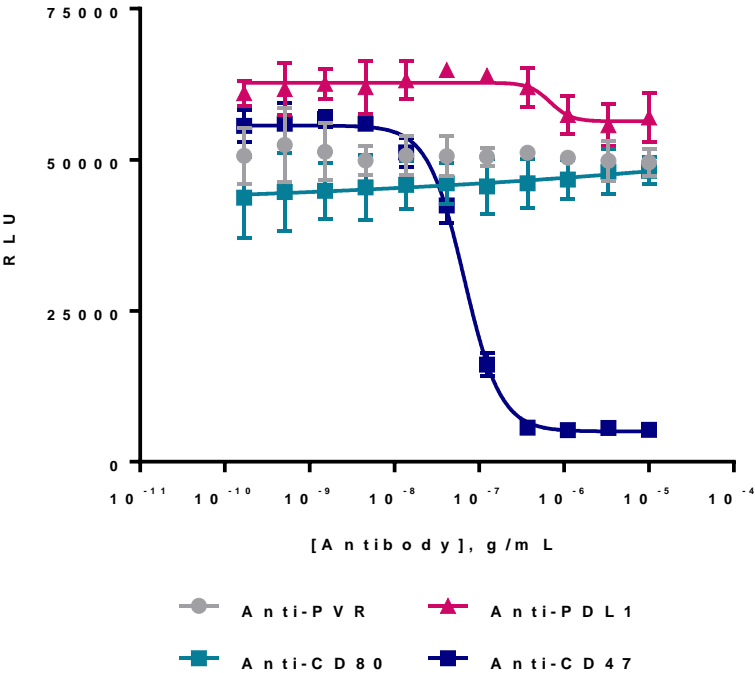


S/B : 17% RSD over 45 passages

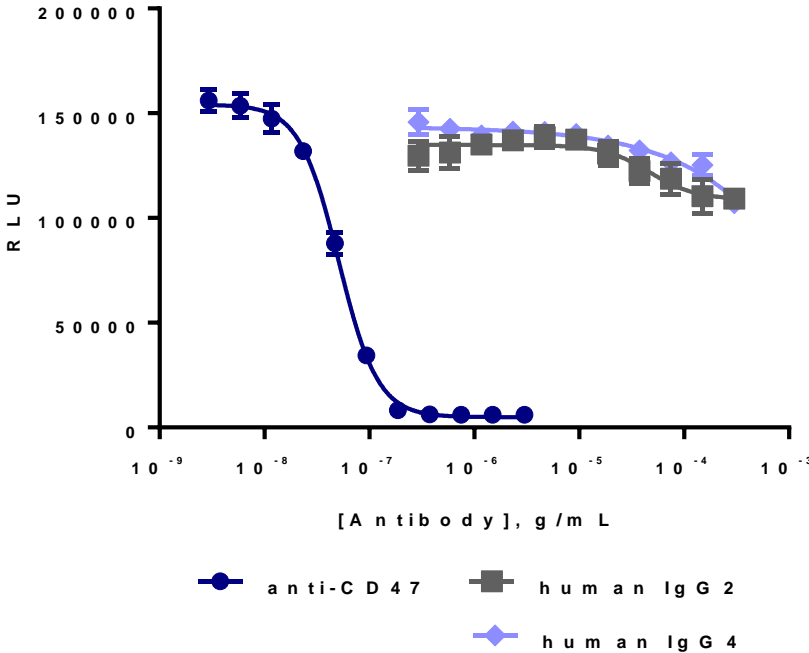
IC<sub>50</sub> : <19% RSD over 45 passages

# PathHunter SIRP $\alpha$ Signaling Assay: Excellent Specificity and Stability-Indicating Properties

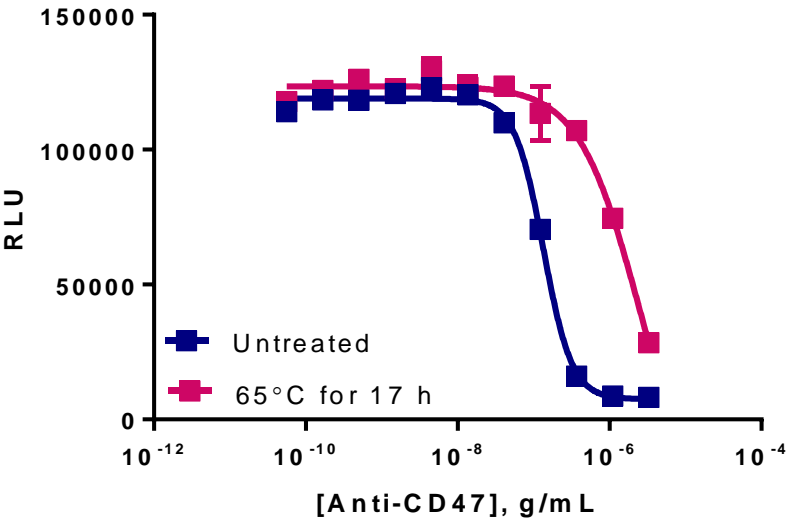
Assay Specificity



Assay Specificity

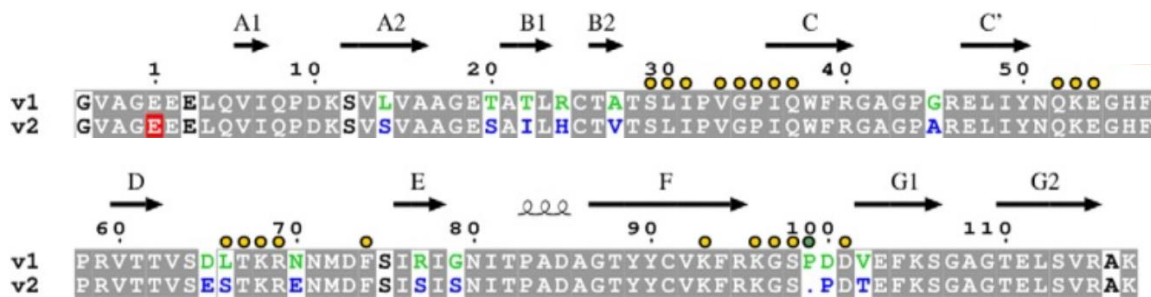


Stability-Indicating



# Signaling Assays for Most Common SIRP $\alpha$ Variants: V1 and V2

- At least 10 SIRP $\alpha$  variants identified
- SIRP $\alpha$  variants 1 and 2 (V1, V2) are most prevalent; differ by 15 amino acids



JOURNAL OF  
BIOLOGICAL  
CHEMISTRY

Published by the American Society for  
Biochemistry and Molecular Biology

*J Biol Chem*. 2014 Apr 4; 289(14): 10024–10028.

PMCID: PMC3974974

Published online 2014 Feb 18. doi: [10.1074/jbc.M114.550558](https://doi.org/10.1074/jbc.M114.550558)

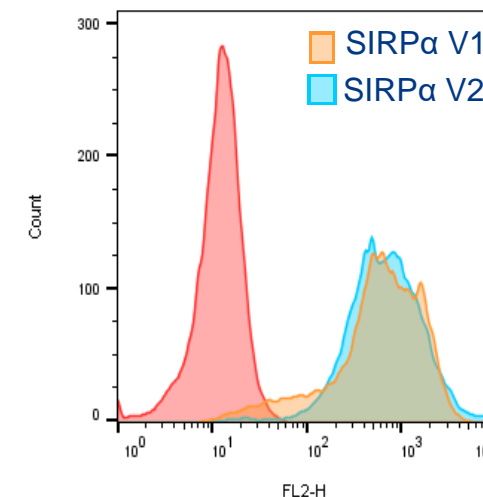
PMID: [24550402](https://pubmed.ncbi.nlm.nih.gov/24550402/)

Polymorphisms in the Human Inhibitory Signal-regulatory  
Protein  $\alpha$  Do Not Affect Binding to Its Ligand CD47\*

Deborah Hatherley, Susan M. Lea, Steven Johnson, and A. Neil Barclay<sup>1</sup>

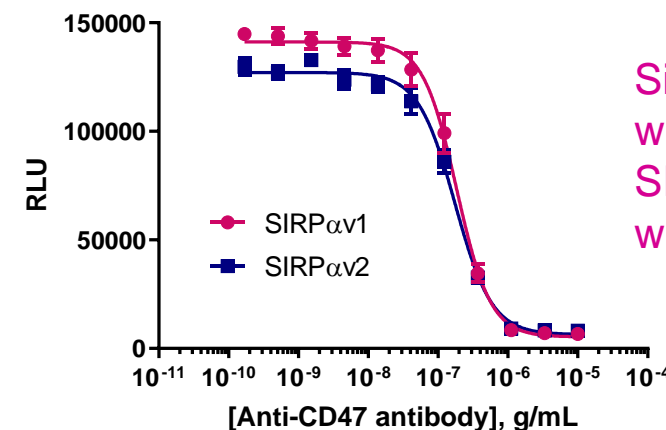
\* = This work was supported by Medical Research Council Grants G9826026 and G0900888 and by Wellcome Trust Senior Investigator Award 100298 (to the S. M. L. group).

## SIRP $\alpha$ Surface Expression



Comparable surface  
expression for SIRP $\alpha$   
V1 and V2

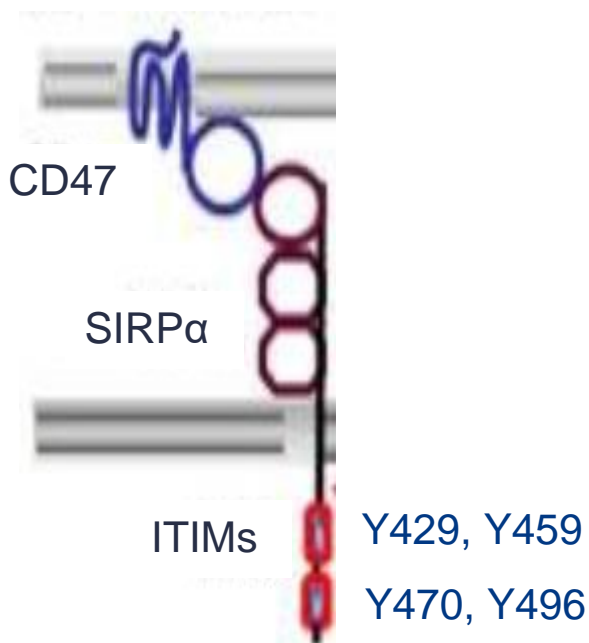
## SIRP $\alpha$ Functional Response



Similar assay  
window and IC<sub>50</sub> for  
SIRP $\alpha$  V1 and V2  
with anti-CD47

# Mutation of SIRP $\alpha$ ITIM Motifs Disrupts CD47-Mediated Signaling

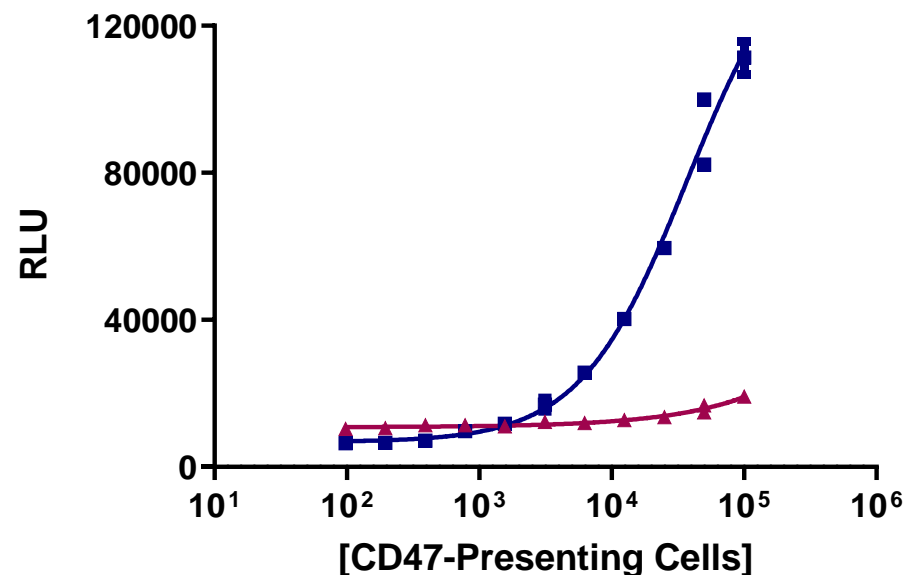
Evaluation of impact of mutations in the 4 tyrosine residues that are potential sites for phosphorylation: Y429, Y459, Y470 and Y496



Adapted from Zen, K. et al. Nat Commun 2013; 4: 2436

## Agonist

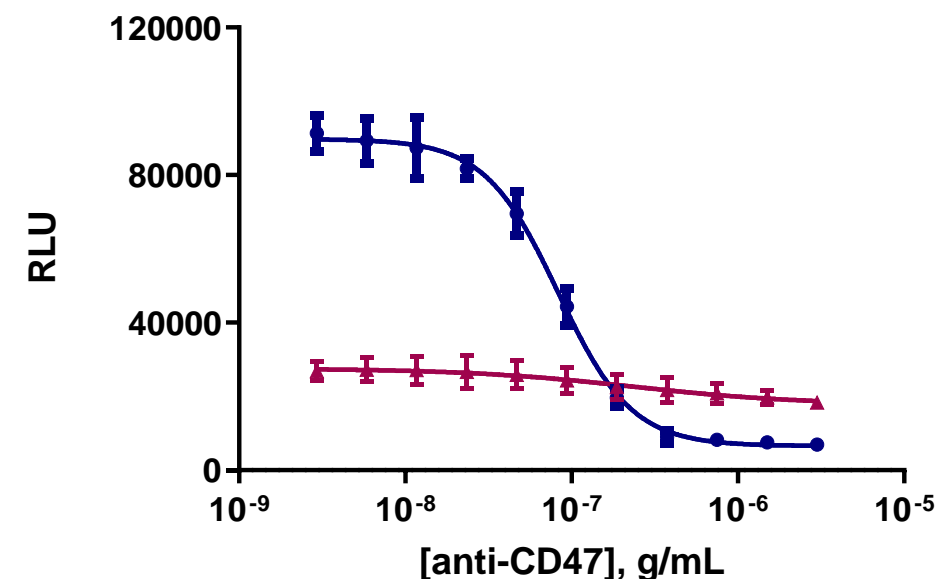
### Jurkat SIRP $\alpha$ Signaling Assay



—▲— SIRP $\alpha$  (Y429F, Y459F, Y470F, Y496F)  
—■— W/T SIRP $\alpha$

## Antagonist

### Jurkat SIRP $\alpha$ Signaling Assay



—▲— SIRP $\alpha$  (Y429F, Y459F, Y470F, Y496F)  
—■— W/T SIRP $\alpha$

*Single mutations disrupted signaling to different degrees (data not shown), but mutation of 3 or more tyrosine residues completely abrogated CD47-mediated SHP recruitment*

# PathHunter® Checkpoint Receptors Assays

## Industry's Largest Menu of Stable Cell Lines and Bioassays

### Functional Cell-based Assays

#### Immunoglobulin Superfamily (IgSF)

SIRP $\alpha$	CD47
ICOS	
CD28	
CD200R	CD200
PD1 (SHP1)	PD-L1
PD1 (SHP2)	PD-L2
BTLA	HVEM
CTLA4	CD86
mPD1	mPD-L1
mCTLA4	mCD80

Signaling Cell Lines

Ligand Cell Lines

#### TNFR Superfamily (TNFRSF)

CD137
CD40
OX40
CD27

### Tools for Testing Agonist Antibodies

#### Clustering Cell Lines

Fc $\gamma$ R1a	Fc $\gamma$ R1a	Fc $\gamma$ R1b
-----------------	-----------------	-----------------

#### Early Access Cell Lines

#### Binding Assays for Bi-specifics

PD-1/LAG3	PD-1/PD-L1
PD-1/CTLA4	PD-L1/CTLA4
PD-1/TIGIT	PD-L1/TIM3
PD-1/CEACAM1	TIM3/CEACAM1
PD-1/CD28	TIGIT/LAG3
mPD-1/mLAG3	mPD-1/mTIGIT
mPD-1/mCTLA4	



Thank You!

## Options to suit your program needs



Purchase Cell Line or  
Bioassay Kits



Quick Confirmation  
with an eXpress Kit



Proof-of-Concept  
Feasibility Study or  
Custom Assay  
Development



Cell Line Rental for  
in-house Testing

Stable Cell Lines

Qualified Bioassays

MOA-based Bioassays

Analytical Cell Banks

Custom Assay Development

GPCRs

Checkpoint Receptors

Cytokine Receptors

Kinases

Signaling Pathways

TGF $\beta$  Superfamily

ADCC Assays

ADCP Assays

CDC Assays

Target Cells

Effector Cells

Eurofins DiscoverX provides simple cell-based assays for immune checkpoint receptors to accelerate immunotherapy drug discovery and development

- **No primary cells** – Get biologically-relevant responses without primary cells
- **Easy-to-use protocol with fast results** – Increase efficiency with an “add-and-read” protocol and get results in 5-8 hours
- **Highly sensitive response** – Better sensitivity than competitor assays allows screening of early stage and dilute development samples
- **Multiple applications** – Drive development of biologic and small molecule drugs
- **Support broader drug program** – Cell-based assay for functional screening, lead optimization, and bioanalytical QC lot release applications

For further information: [www.discoverx.com/checkpoint](http://www.discoverx.com/checkpoint)