



Accelerating Immunotherapy Drug Development Targeting Checkpoints with Robust MOA- Based Assays

Presented by Dr. Alexander Baumann

Head of European Business Development
Eurofins DiscoverX Products

Meet us at BOOTH #24

OUR EXPERTISE
IN YOUR HANDS.
DISCOVER
CONFIDENTLY.

Eurofins DiscoverX – Global Leader in Cell-Based Assays for Screening, Profiling, Potency, & Lot Release Programs

From Discovery to Development to Clinic to Post-Market

20+ Years of Enabling Drug Discovery and Development Programs



San Francisco Bay Area, California

R&D & Manufacturing

San Francisco Bay Area, California (HQ)
St. Charles, Missouri
Poitiers, France

10+

Druggable
target
classes

1500+

Stable cell line and
membrane preps

20+

Core patents

2000+

Publications across multiple
applications

55+

Qualified &
MOA- based
Bioassays

Validated

>30 Billion Data Points
screened in assay services with
same assays

**3 Certified
CRO Partners**

Scientific training to
enable global CROs

ICH Based Bioassay Qualification

Facilitate downstream
validation studies

Dedicated Scientific Support

Experienced team providing
scientific support

20+ Successful Assay Transfers

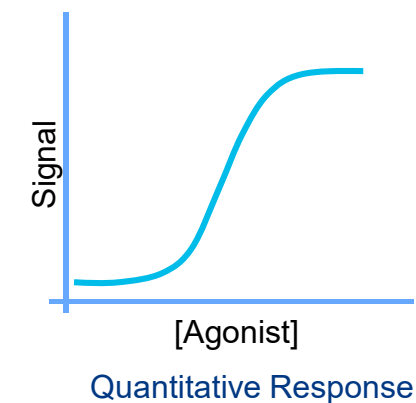
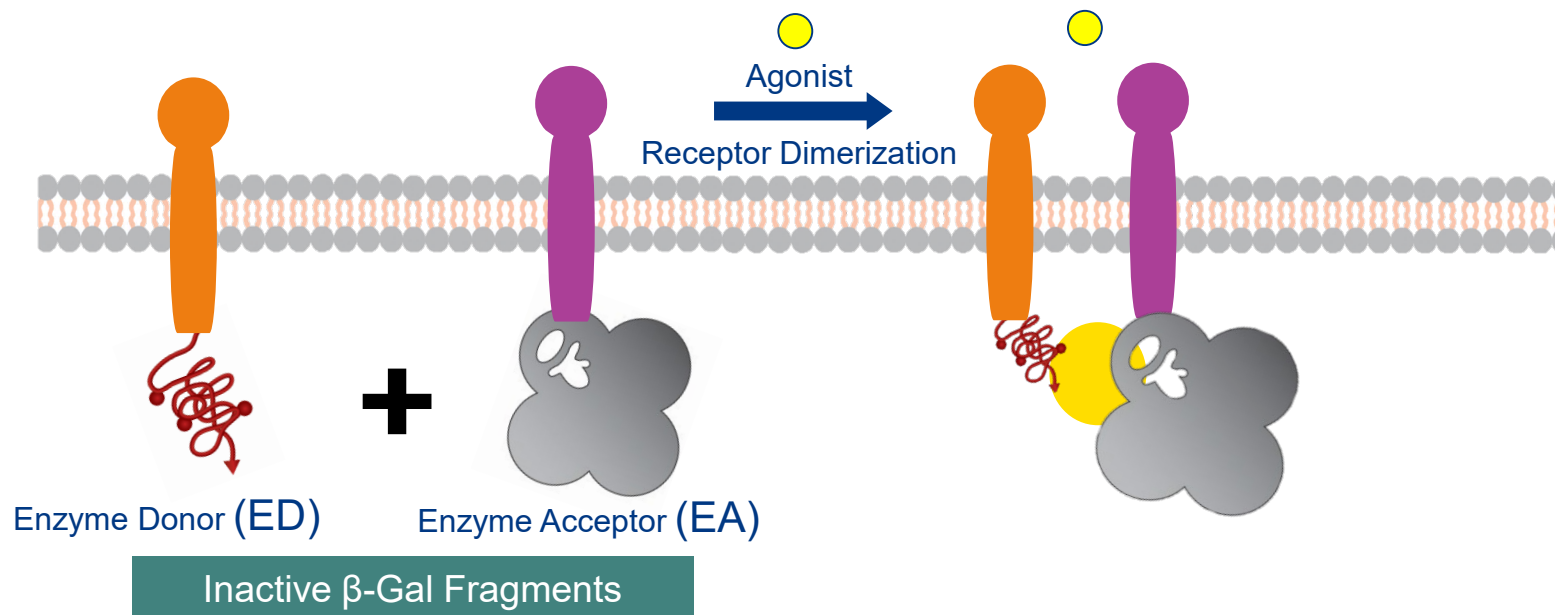
At clients/affiliated CRO sites

50+ Global Programs

For potency, stability and NAb
testing

Enzyme Fragment Complementation (EFC) Technology

Assay Design: e.g. Receptor Dimerization Assay (MOA for RTKs and Interleukins)



Homogenous

- Add-and-read assay format
- No washing steps
- No centrifugation step
- No shaking or filtration

Robust

- Enzymatically-amplified Assay
- Large signal-to-background
- High precision and reproducibility

Qualified and Validated

- Optimized for hundreds of targets
- Screened billions of data-points
- >2K peer-reviews publications

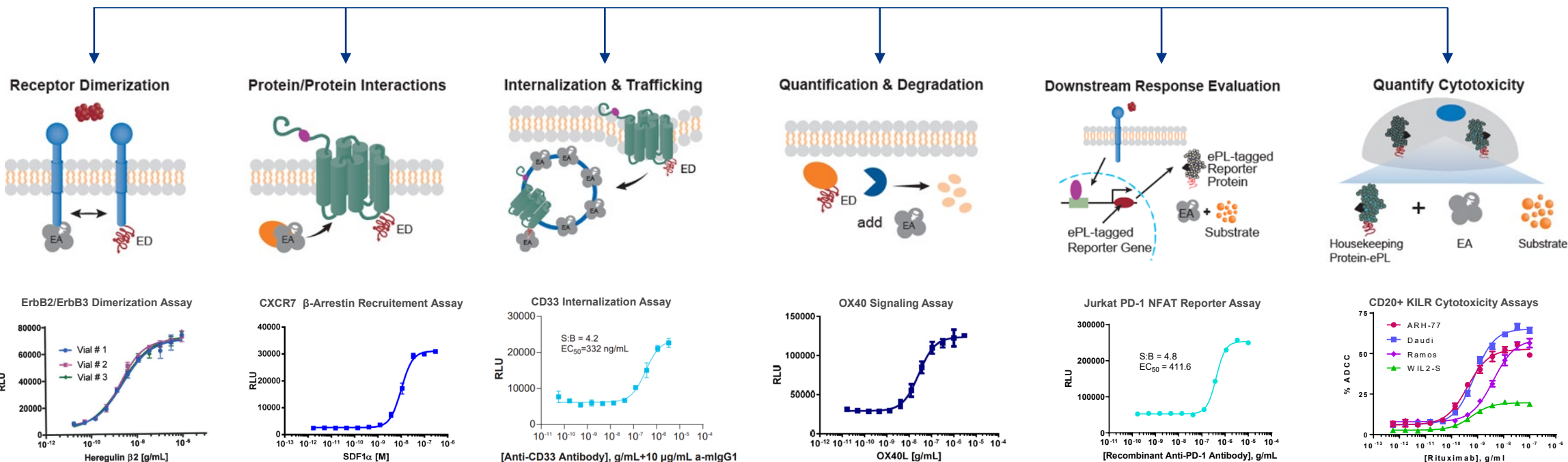
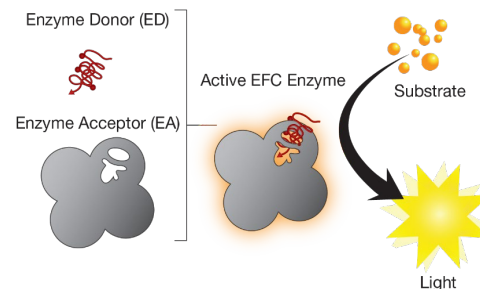
Transferrable

- Assay protocol is similar across platform
- Detailed user manuals
- Assays routinely transferred between multiple sites

Enzyme Fragment Complementation (EFC)

Lego®'s for Assay Developers!

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



PathHunter® Checkpoint Receptors Assays

Industry's Largest Menu of Stable Cell Lines and Bioassays

Functional Cell-Based Assays

Immunoglobulin Superfamily (IgSF)

SIRP α	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 γ R1a	Fc γ R1a	Fc γ R1b
-----------------	-----------------	-----------------

Early Access Cell Lines

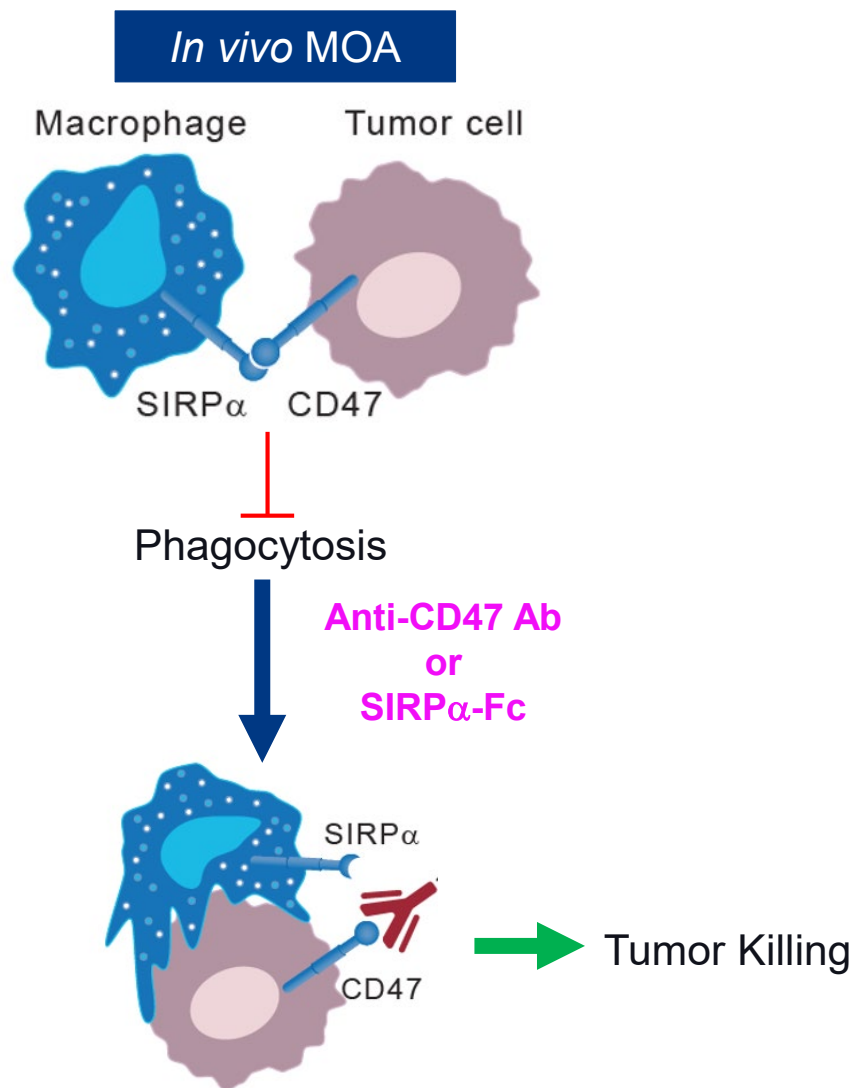
Binding Assays for Bi-specifics

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	

Case Study: PathHunter® SIRP α Signaling Bioassay Development

For Therapeutics Targeting CD47 | SIRP α Signaling Axis

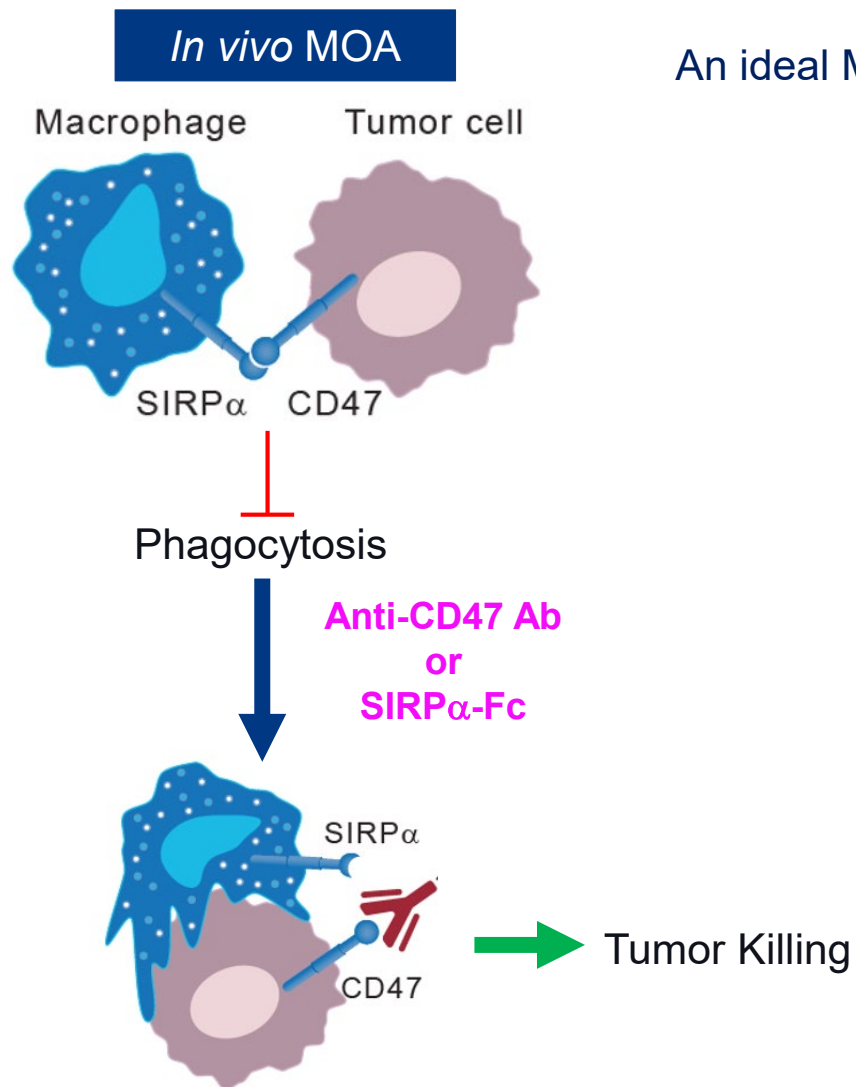
The SIRP α / CD47 Axis - Innate Immune Checkpoint



The SIRP α / CD47 Axis

- **SIRP α** : inhibitory receptor expressed on macrophages and dendritic cells that promotes phagocytosis of foreign objects
- **CD47**: ligand for SIRP α , expressed on nearly all cells, but significantly up-regulated in many tumor types
- **Blocking** the CD47 / SIRP α axis promotes tumor killing

The SIRP α / CD47 Axis - Innate Immune Checkpoint



An ideal MOA-reflective cell-based assay would quantify phagocytosis endpoint

Challenges with Phagocytosis Assays

- Typically evaluate co-localization of macrophage and target cells using flow cytometry or high content imaging
- Require use of human macrophages as effector cells
 - Macrophage cell lines not reliable as effectors
 - Differentiation from primary monocytes takes 7 days
 - Activated M1 macrophages difficult to lift (30' to 60')
- A more 'QC-friendly' MOA reflective assay is required

PathHunter® SIRPα Signaling Assay

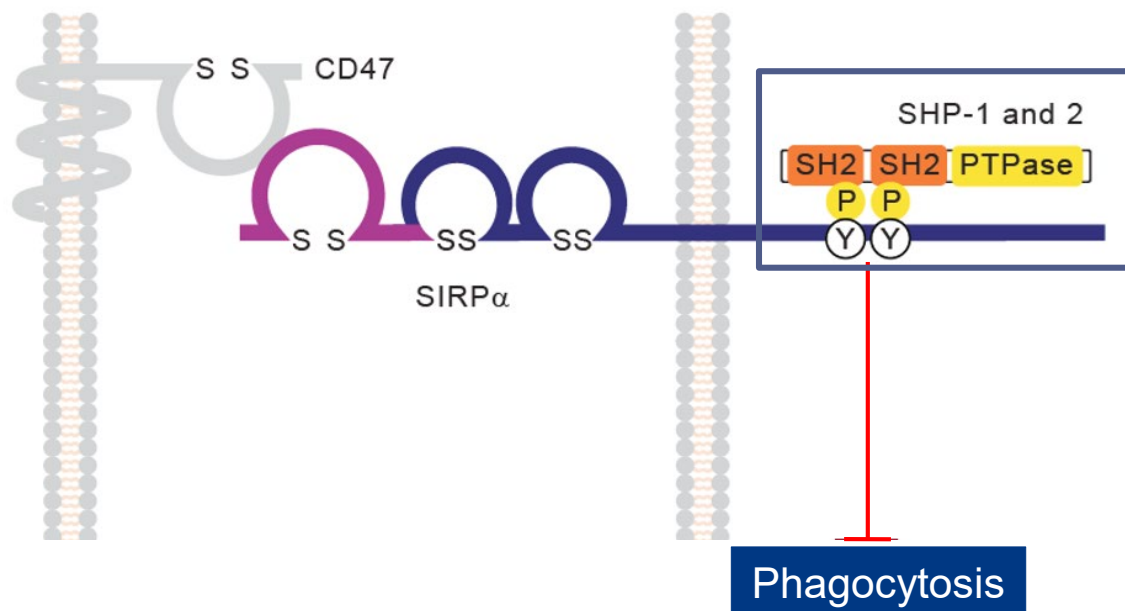
Assay Concept

Co-culture SHP recruitment model based on β -galactosidase enzyme fragment complementation

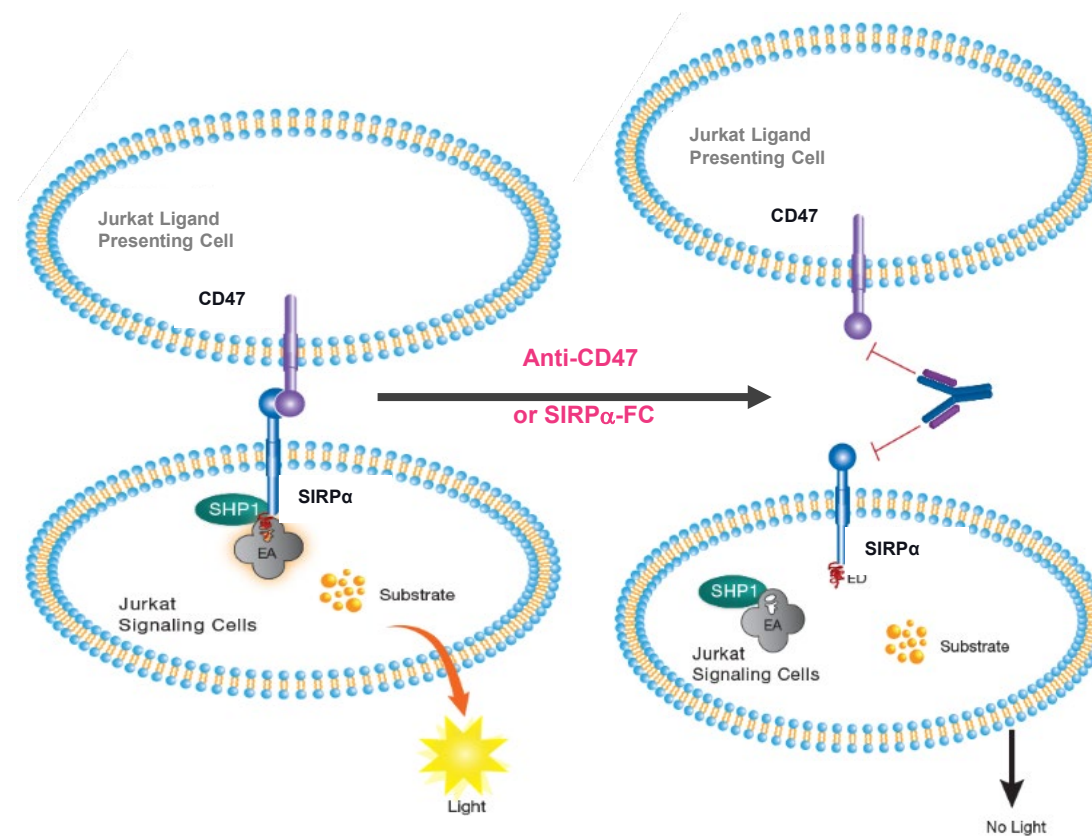
Molecular MOA

Tumor Cell

Macrophage



Co-Culture SHP Recruitment Model



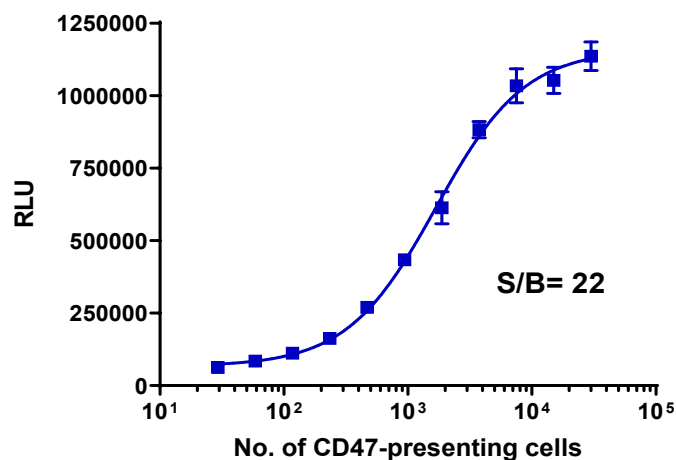
Adapted from Trends in Cell Biology, 2008. Vol 19, No. 2

PathHunter® SIRPα (CD47) Signaling Assay

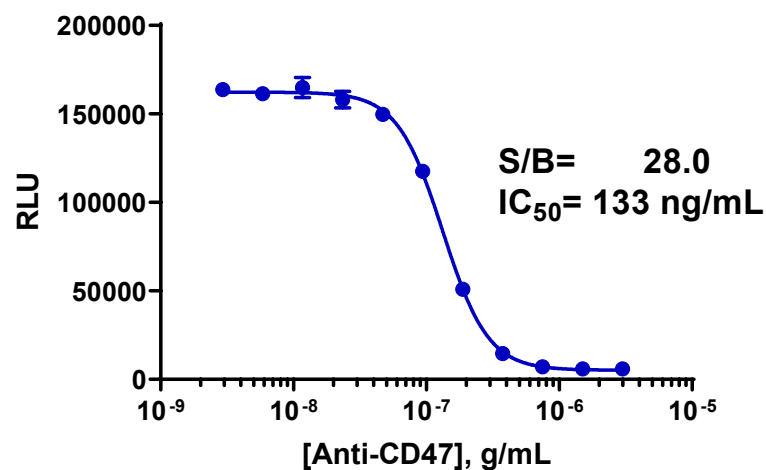
Co-culture model with stable surface expression of SIRPα and a stable functional response over 45+ passages

Cell-Presented Ligand

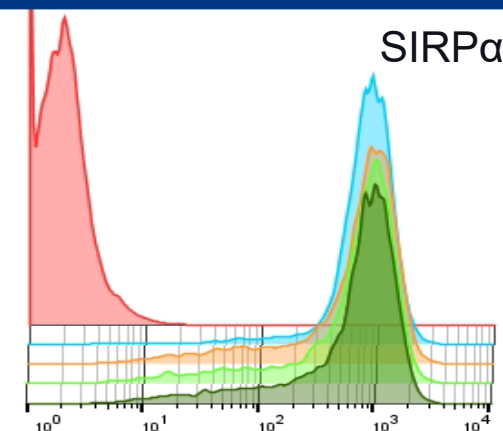
Agonist



Antagonist

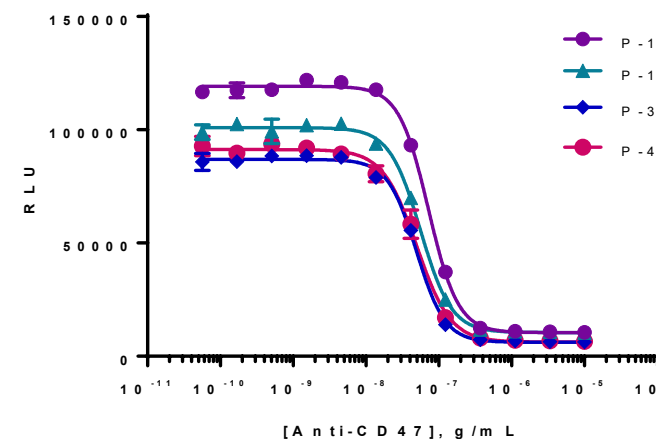


Stable SIRPα Surface Expression



SIRPα expression varies by <20% RSD over 45 passages

Stable Functional Response



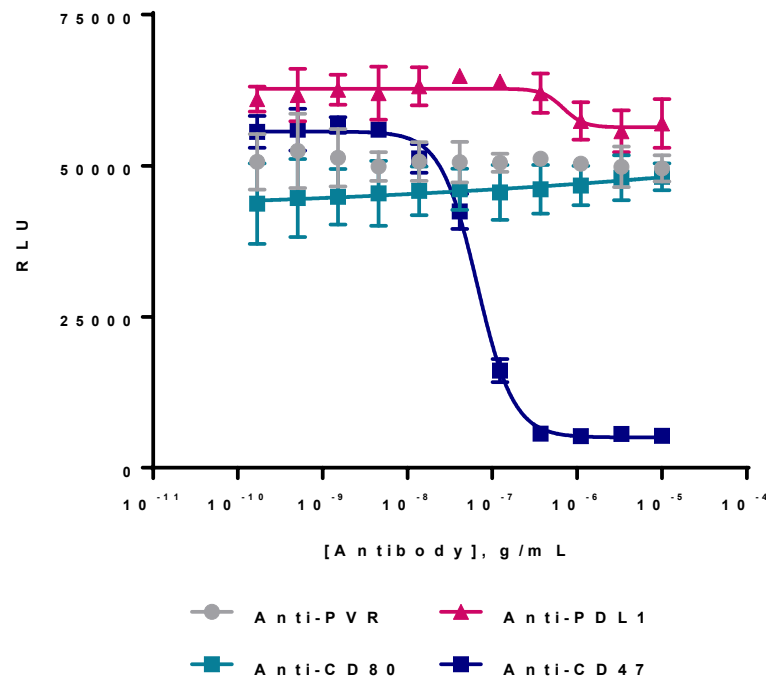
S/B : 17% RSD over 45 passages

IC_{50} : <19% RSD over 45 passages

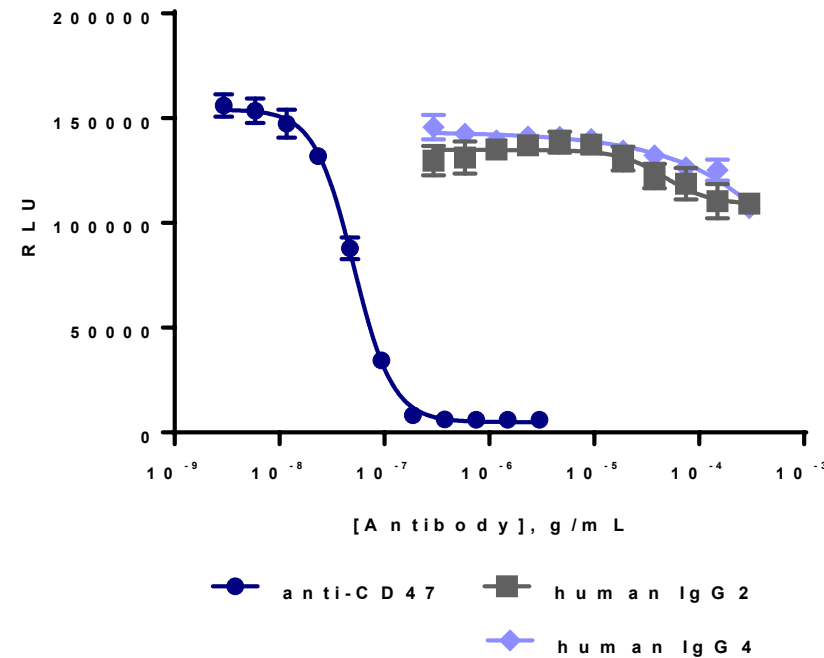
PathHunter® SIRP α (CD47) Signaling Assay

Excellent Specificity and Stability-Indicating Properties

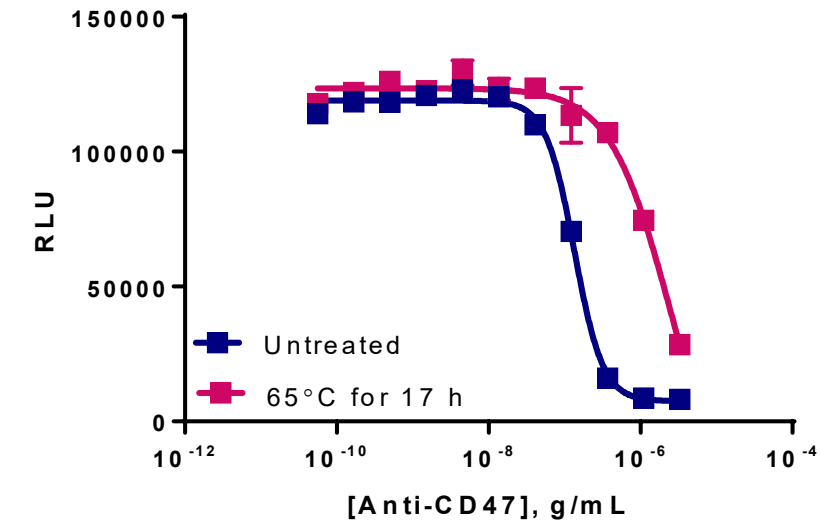
Assay Specificity



Assay Specificity



Stability-Indicating



RTU (Bioassay) Method

Thaw and add CD47 and SIRPα frozen cells (1 vial each) to assay plate

Prepare antibody dilutions and add to assay plate

24h

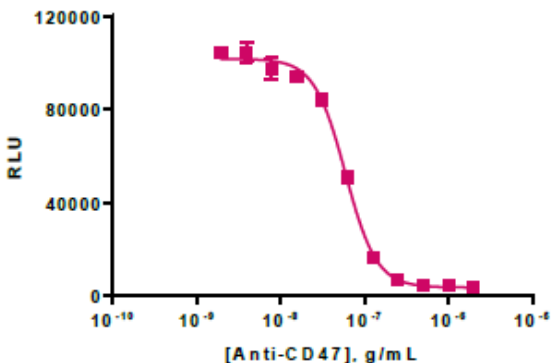
Add Detection Reagent to assay plate

1h

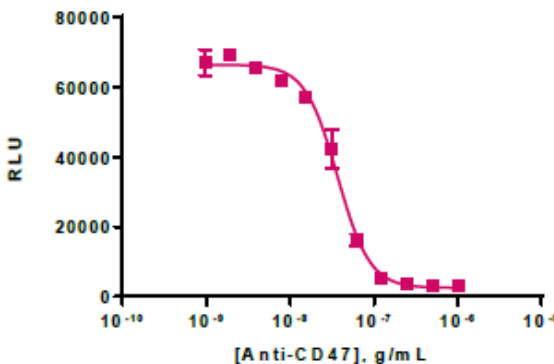
Read signal on plate reader

Total Assay Time:
~26 hours

Continuous Culture



Cryopreserved (RTU)

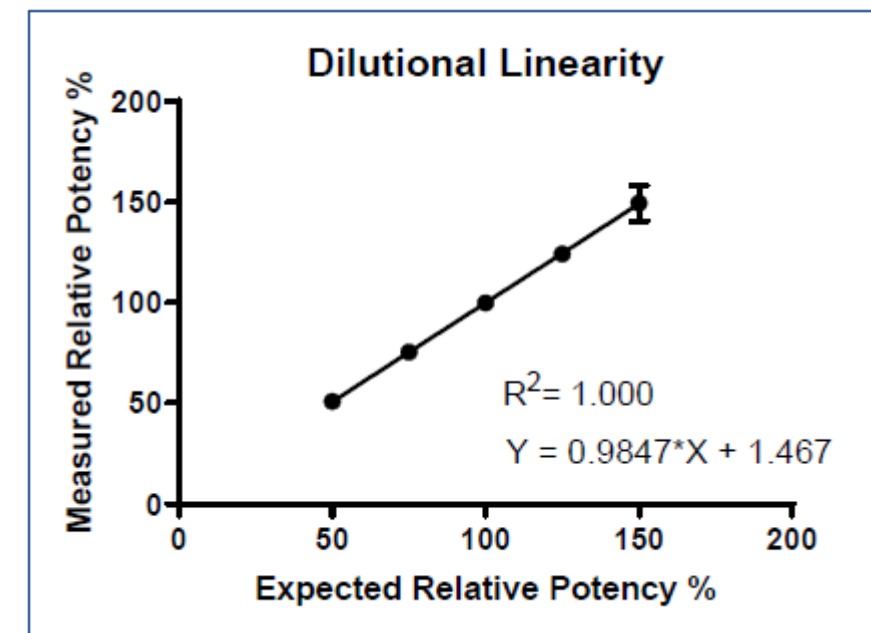


Comparable Performance to Continuous Culture Format

Format	HillSlope	IC ₅₀ (ng/mL)	S/B
Continuous Culture	-2.337	59.1	28
Cryopreserved (RTU)	-2.264	36.8	20

SIRPα Bioassay Qualification

Expected RP (%)	Exp #	Analyst #	Measured RP (%)	Average RP (%)	% RSD	% Accuracy	Relative Bias, %
150	1	1	164	149.5	5.96	99.7	-0.3
	2	1	144				
	3	1	145				
	4	2	140				
	5	2	148				
	6	2	156				
125	1	1	123	124.2	3.16	99.4	-0.6
	2	1	125				
	3	1	124				
	4	2	119				
	5	2	123				
	6	2	131				
100	1	1	102	99.8	3.66	99.8	0.2
	2	1	95				
	3	1	103				
	4	2	104				
	5	2	98				
	6	2	97				
75	1	1	75	75.3	5.15	100.4	0.4
	2	1	73				
	3	1	79				
	4	2	73				
	5	2	81				
	6	2	71				
50	1	1	55	50.8	6.51	101.6	1.6
	2	1	52				
	3	1	53				
	4	2	51				
	5	2	48				
	6	2	46				



Dilutional Linearity: $R^2 = 1.000$

Accuracy: 100.02%

Intermediate Precision: 6.5%

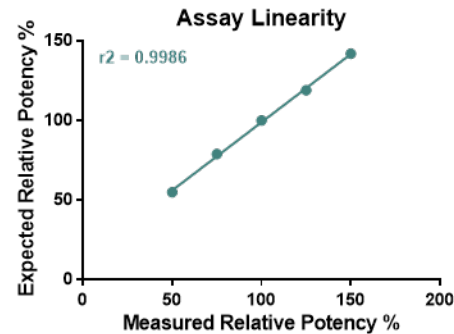
SIRPα Bioassay is highly accurate, precise and linear!

Examples of Qualified Bioassays for Precise Potency Determination for Therapeutic Drugs

Performance Qualification

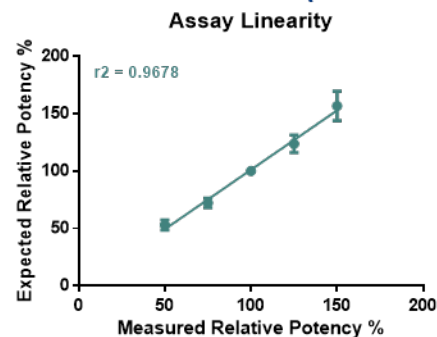
- ✓ Relative Accuracy
- ✓ Specificity
- ✓ Intermediate Precision
- ✓ Range

PD-1 Pembrolizumab (Keytruda)



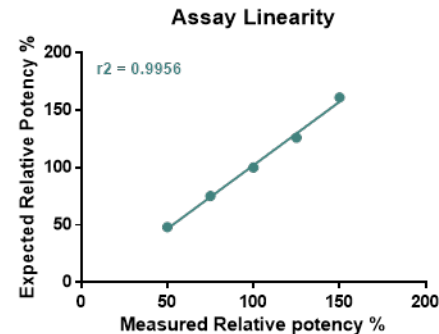
Pembrolizumab
Accuracy = 99.7%
Precision = 4.8%

KDR/KDR Ranibizumab (Lucentis)



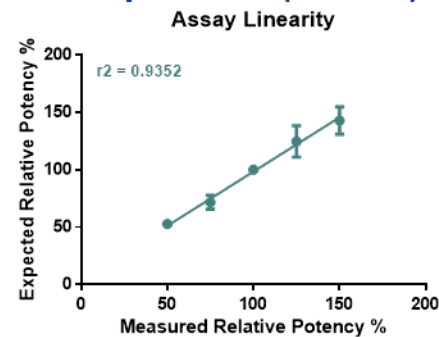
Ranibizumab
Accuracy = 100.5%
Precision = 5.4%

EGFR/ERBB2 Panitumumab (Vectibix)



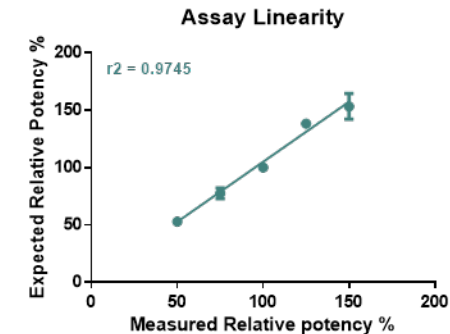
Panitumumab
Accuracy = 103.4%
Precision = 4.6%

PTHRI Teriparatide (Forteo)



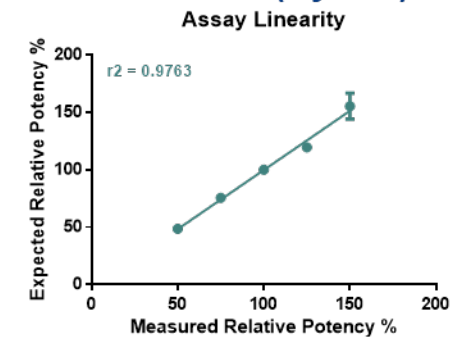
Teriparatide
Accuracy = 100.4%
Precision = 9.9%

IL6R/IL6ST Tocilizumab (Actemra)



Tocilizumab
Accuracy = 105.5%
Precision = 4.7%

GLP1R Exenatide (Byetta)



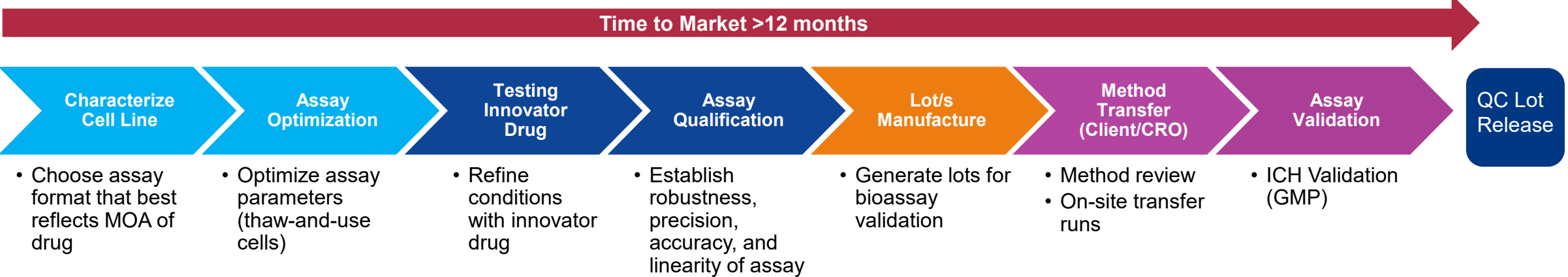
Exenatide
Accuracy = 100.8%
Precision = 5.2%

A central requirement for bioassays is the ability to accurately report drug potency using a functionally relevant readout

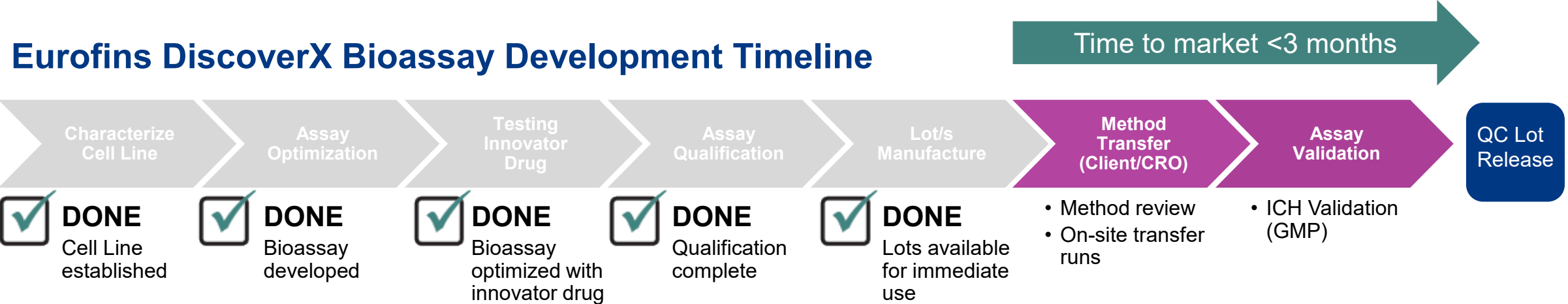
Eurofins DiscoverX bioassays are qualified with excellent linear potency as demonstrated with clinically relevant drugs.

Accelerating Implementation Phase for QC Lot Release with Qualified Bioassays

Typical Bioassay Development Timeline



Eurofins DiscoverX Bioassay Development Timeline



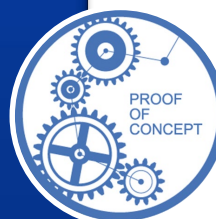
Eurofins DiscoverX Ready-To-Plate Bioassay Kits Save 9 Months of Assay Development Time

Thank You and Please Join Us at Booth #24

Options to suit your program needs



Purchase Membranes,
Cell Lines or Ready to
use Cells/Kits



Proof-of-Concept
(Feasibility Study) or
Custom Assay
Development



Cell Line Rental
(3-months Block)

Qualified Bioassays

Cell Lines

Custom Assay Development

Analytical Cell Banks

GPCRs

Checkpoint Receptors

Cytokine Receptors

Kinases

Signaling Pathways

TGF β Superfamily

ADCC Assays

ADCP Assays

CDC Assays

Target Cells

Effector Cells

For more information see discoverx.com/checkpoint or
email Alexanderbaumann@eurofins.com (EU contact) or heatherzanetakos@eurofins.com (US contact).