

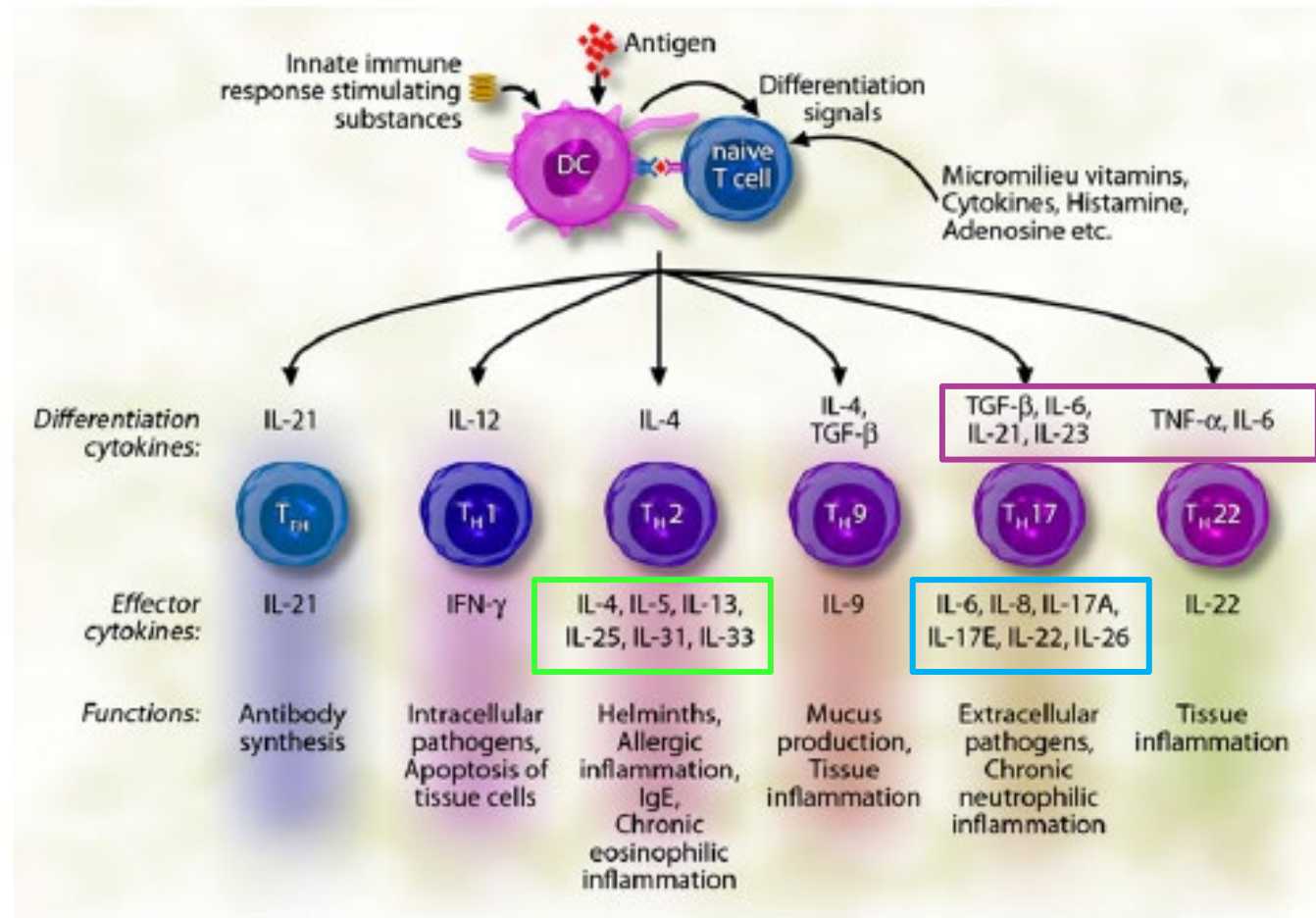
Advancing Therapeutics Targeting Cytokines with Functional Cell-based Assays

Jennifer Lin-Jones, Ph.D.

Associate Director, Assay Development

Cytokines Are Key Mediators of Inflammation

Cytokines produced in specific activated T cell populations are key mediators of inflammation and therapeutics



Approved Drugs Targeting Cytokines

Dupixent (Dupilumab)

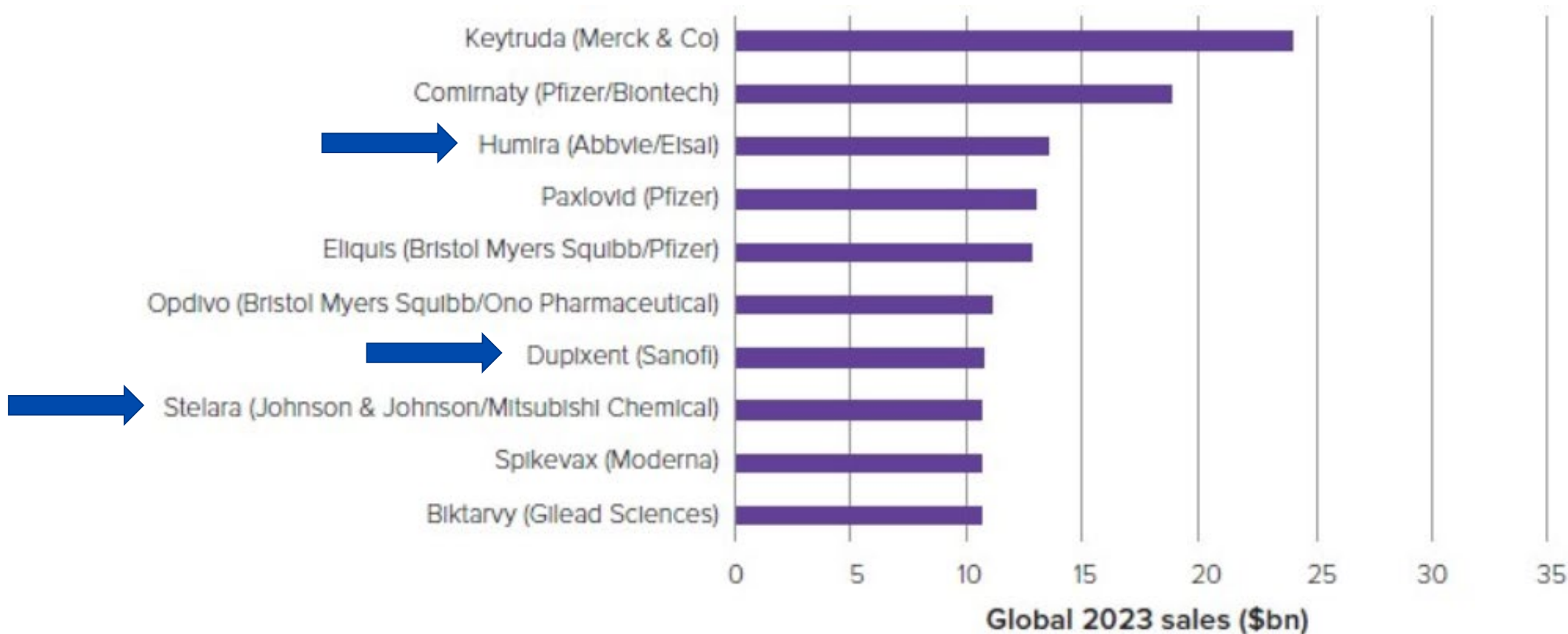
Actemra (Tocilizumab)
Cosentyx (Secukinumab)

Humira (Adalimumab)
Stelara (Ustekinumab)
Tremfya (Guselkumab)
Skyrizi (Rizankizumab)

And others in clinical development...

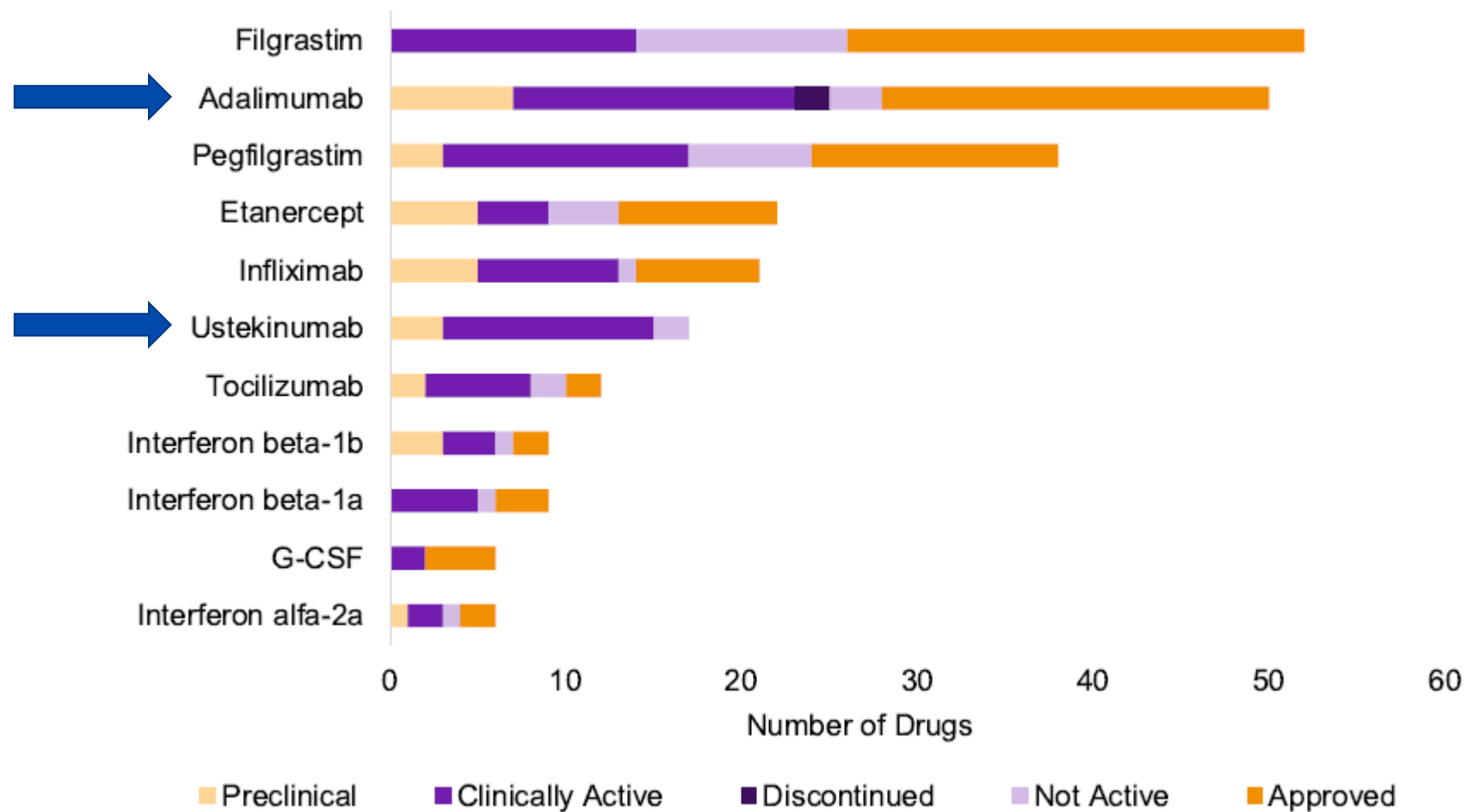
Source: Akdis *et al.* *J Allergy Clin Immunol*, Vol 127. All registered trademarks are the property of their respective owners. IL = interleukin.

Top Selling Therapeutics Targeted Cytokines



Source: Evaluate Vantage: The biggest-selling drugs of 2023. All registered trademarks are the property of their respective owners.

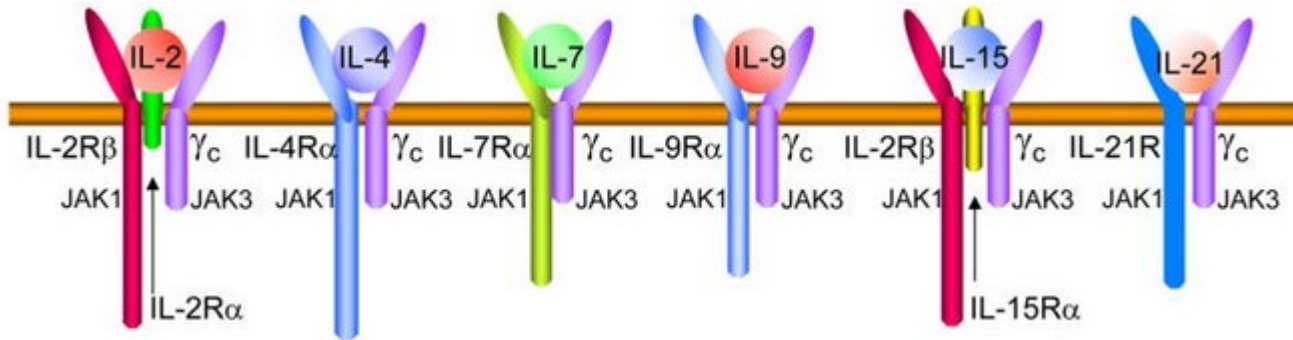
Cytokine Biosimilars Are Among the Top Selling Drugs



Source: Beacon Cytokine Webinar, 2023

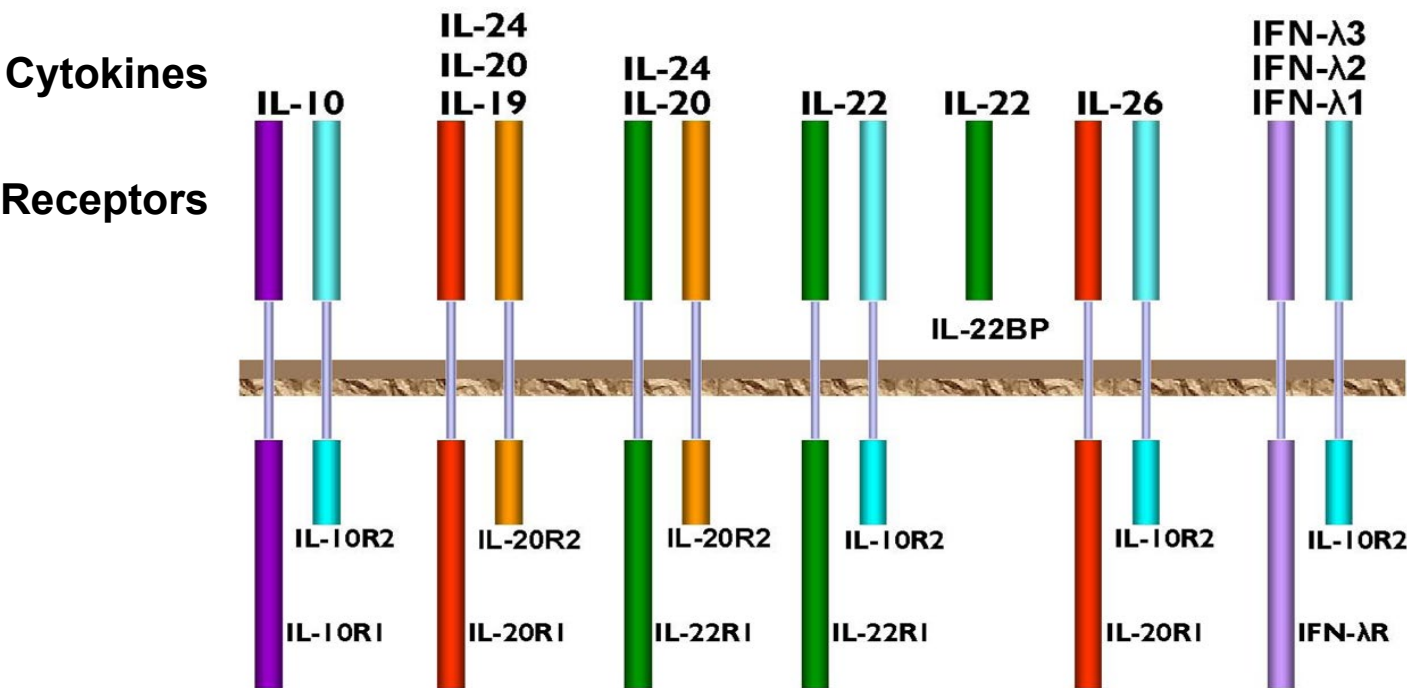
IL-2 Family

One receptor,
multiple partners



IL-10 Family

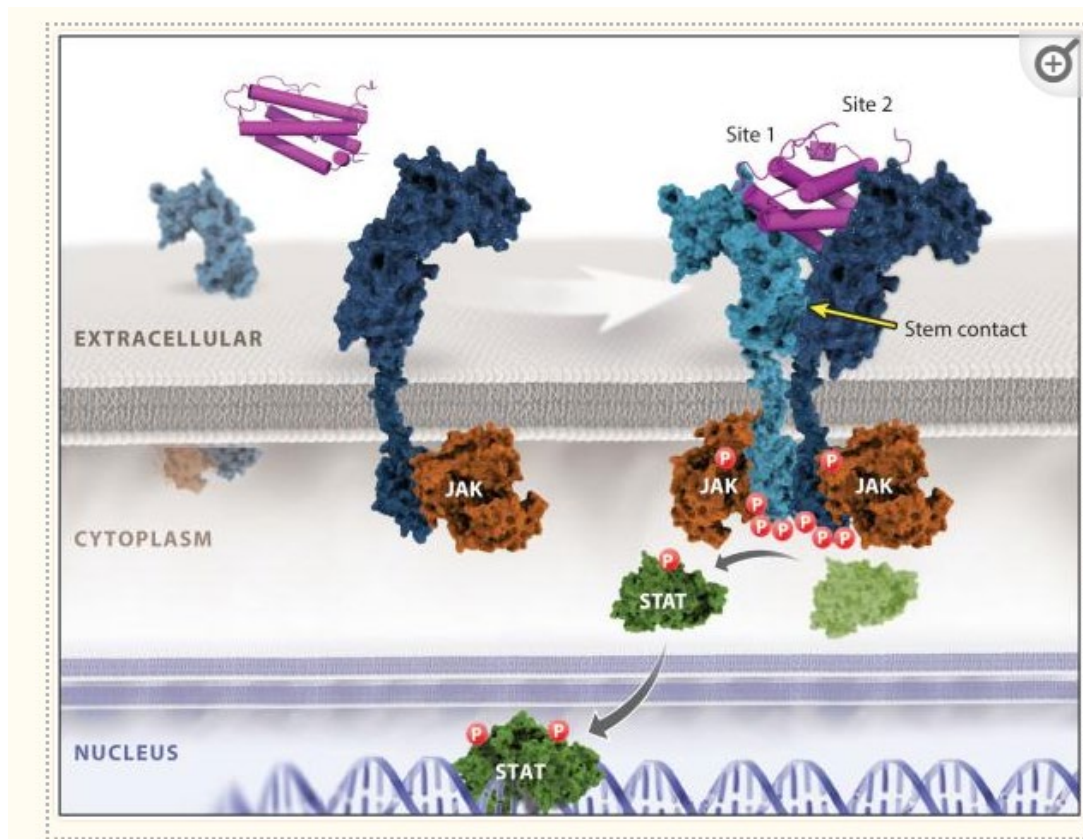
Multiple receptors
heterodimerize with
multiple partners



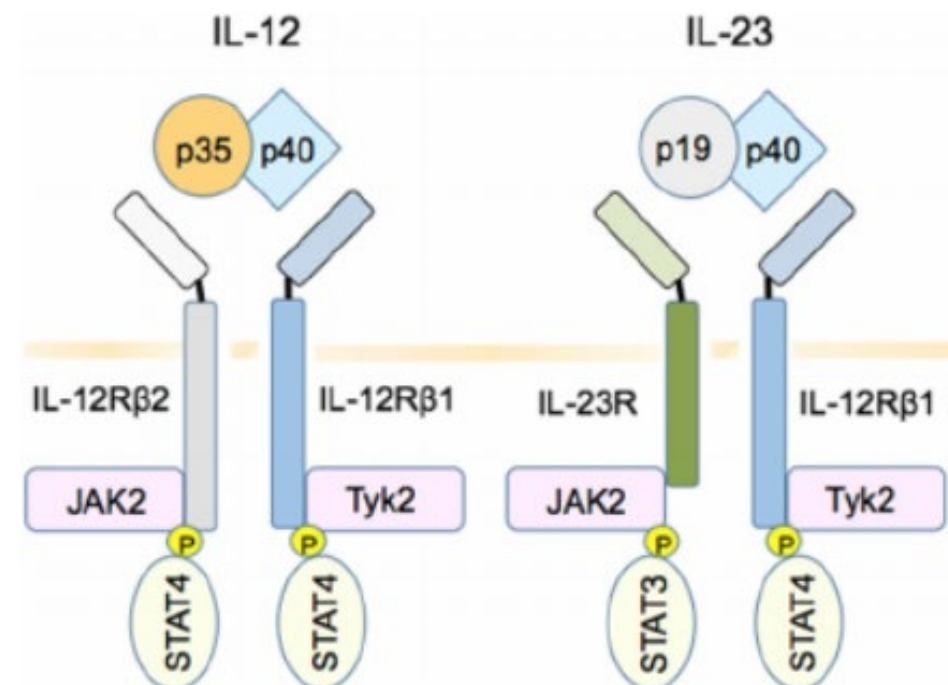
Sources: Top image - Liao *et al.* (2011) *Curr Opin Immun* 23(5): 598-604. Lower image - Zdanov A. (2010) *Cytokine Growth Factor Rev* 21(5): 325-30.

Cytokines are comprised of homomeric or heterodimeric subunits that bind to different receptor subunit combinations

Homomeric Ligands (e.g. IL-2, IL-4)



Heterodimeric Ligands (e.g. IL-12, IL-23)

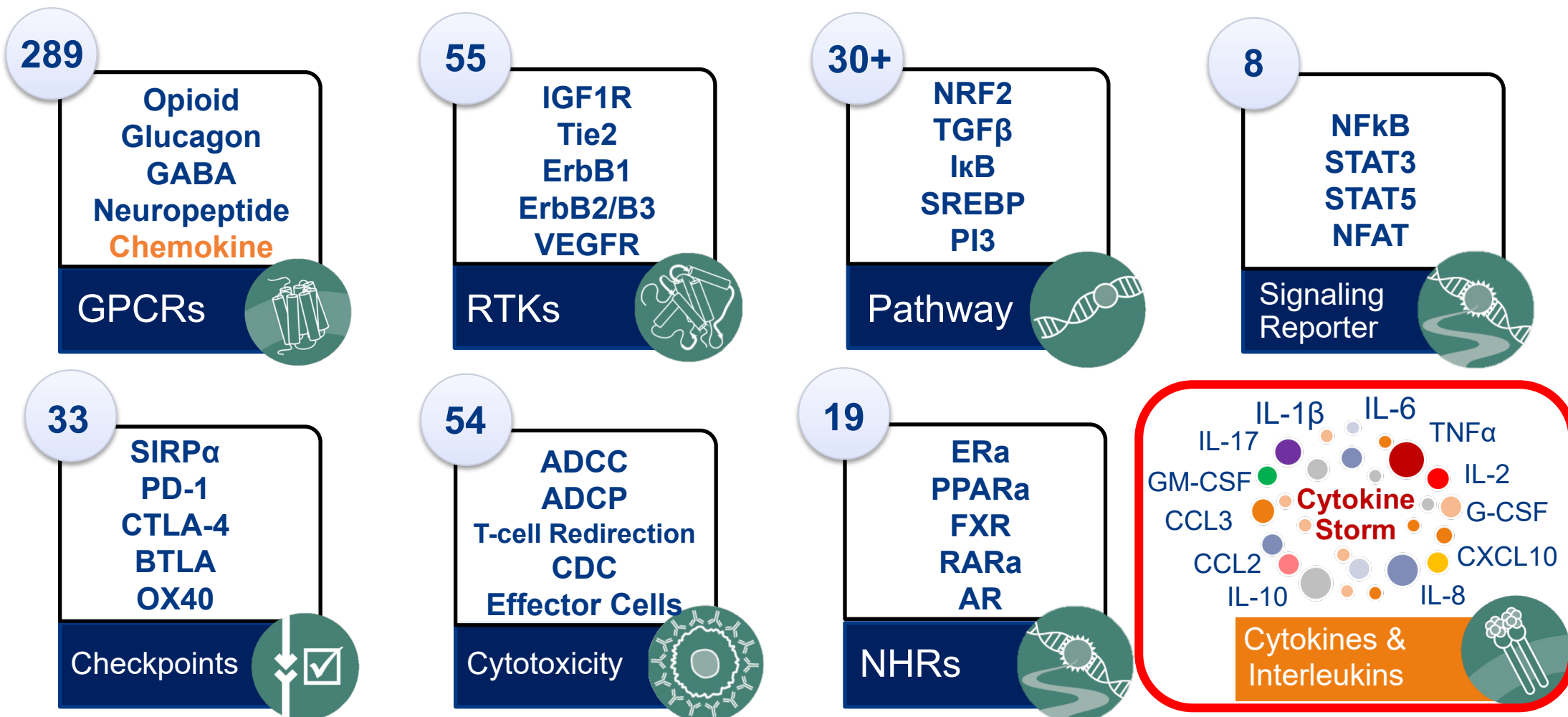


Source (left image): Spangler *et al.* 2015. *Ann Rev Immunol* 33: 139-167.

Source (right image): Choi *et al.* 2015. *Clin Rev Allergy Immunol* 49(3):327-32

Eurofins DiscoverX® Large Menu Cell-Based Assays for Discovery, Potency, & NAb Assay Development

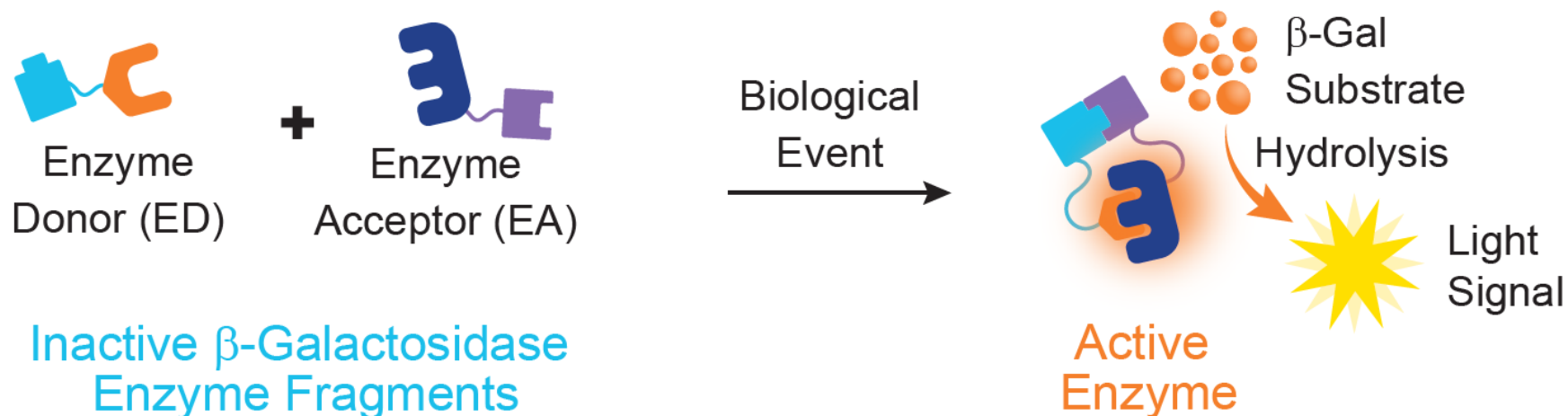
>1,500+ Cell lines and membrane preps to support bioassay development for major drug target classes



NAb = neutralizing antibody

Eurofins DiscoverX[®] Enzyme Fragment Complementation (EFC) Technology

Enabling technologies with a flexible platform based on a split β -galactosidase enzyme



Homogenous

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

Robust

- Large signal-to-background
- Gain of signal assay design
- High assay accuracy, precision and reproducibility

Qualified and Validated

- ICH-based Qualification
- Validated at GMP CRO
- Screened billions of data-points

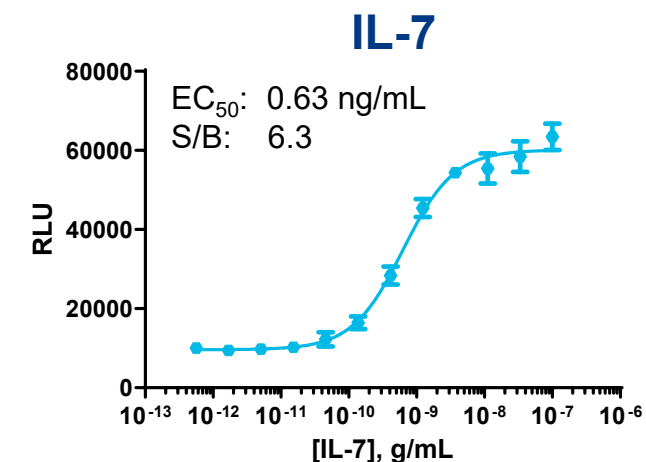
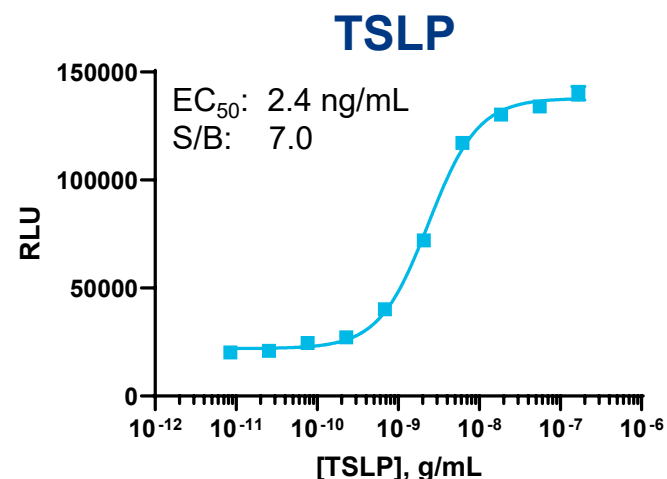
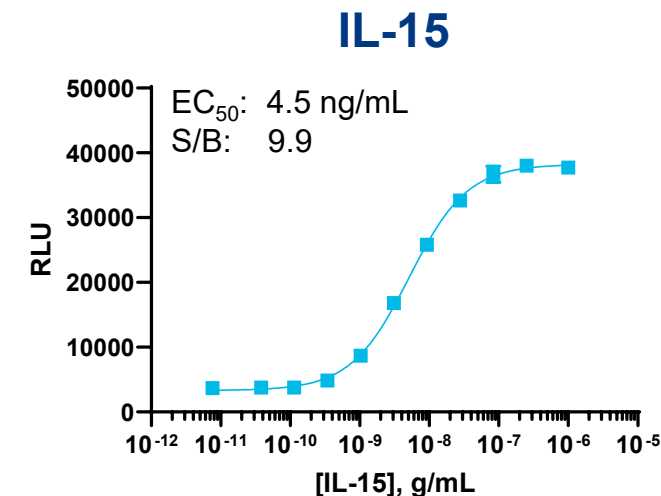
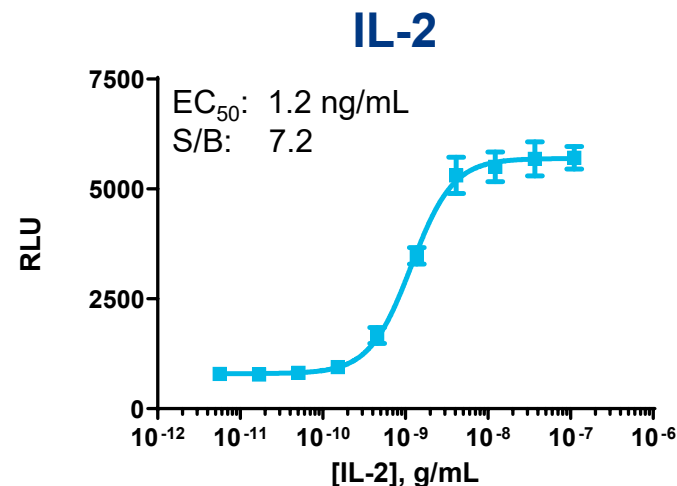
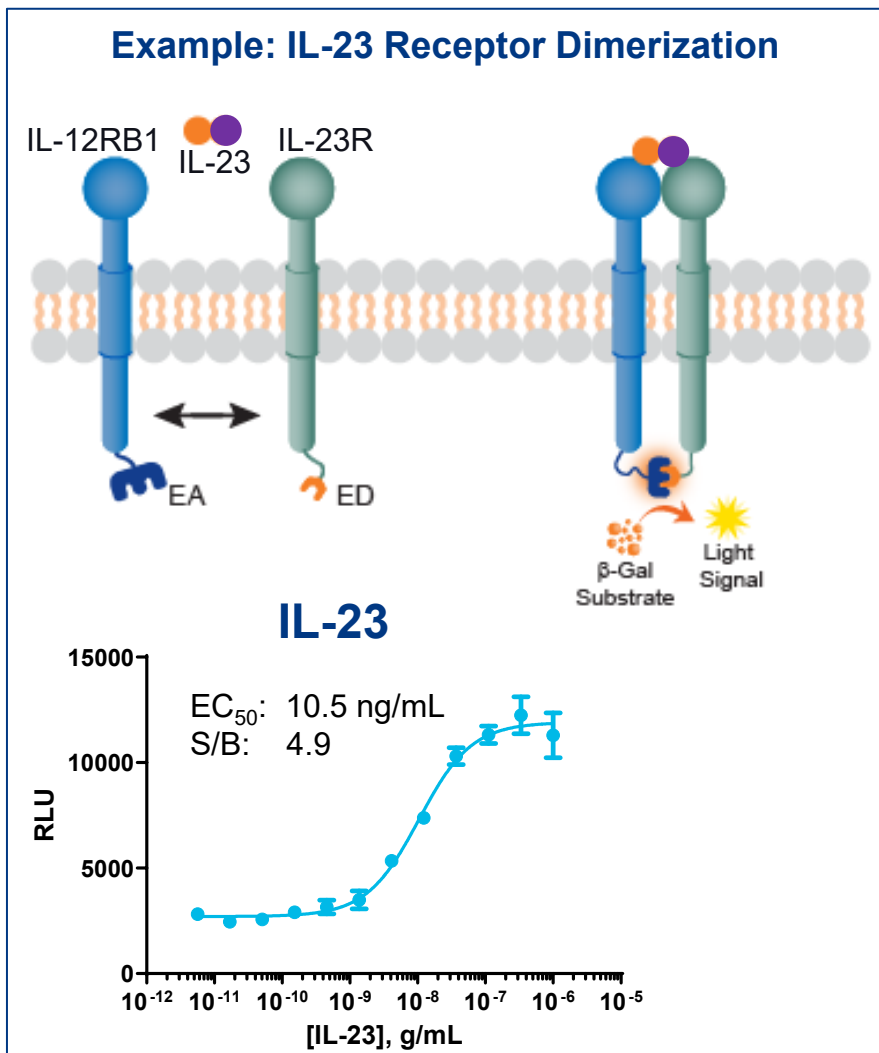
Easy to Transfer

- Assay protocol is similar across platform
- Detailed user manuals
- Method transfer support

ICH = International Council of Harmonization; GMP = Good Manufacturing Practices; CRO = Contract Research Organization

Receptor Dimerization Assay Design

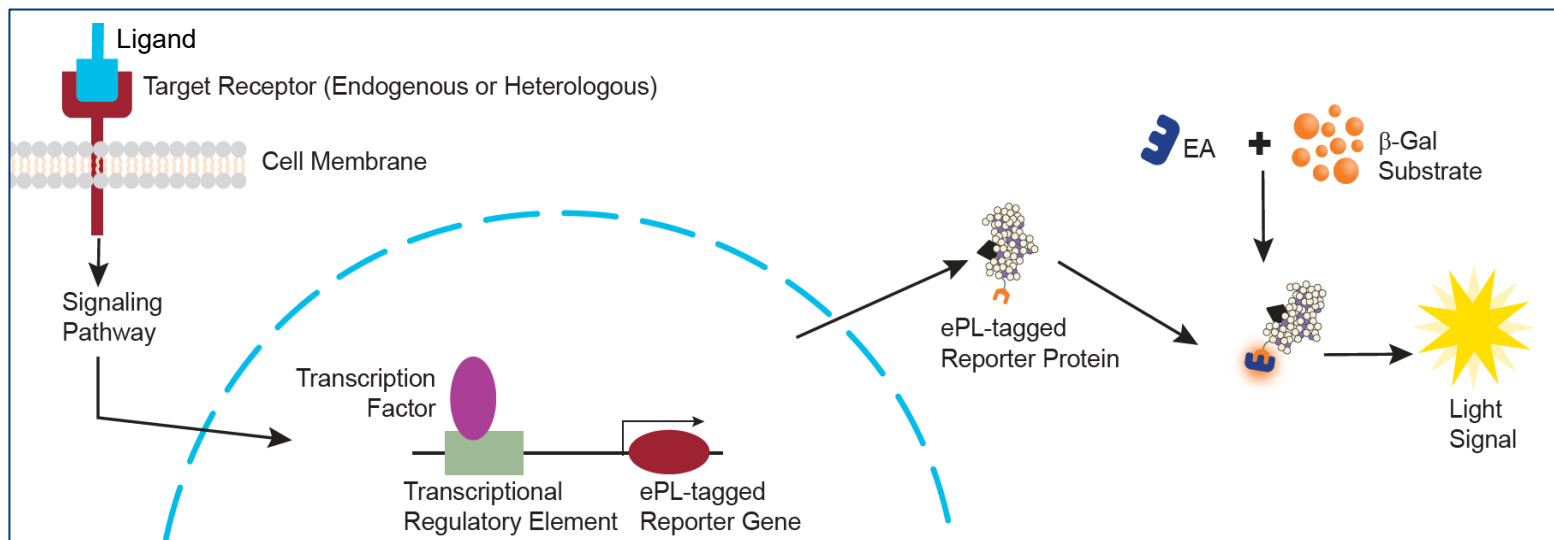
PathHunter® receptor dimerization assays quantify a discrete early step in receptor signaling (proximal)



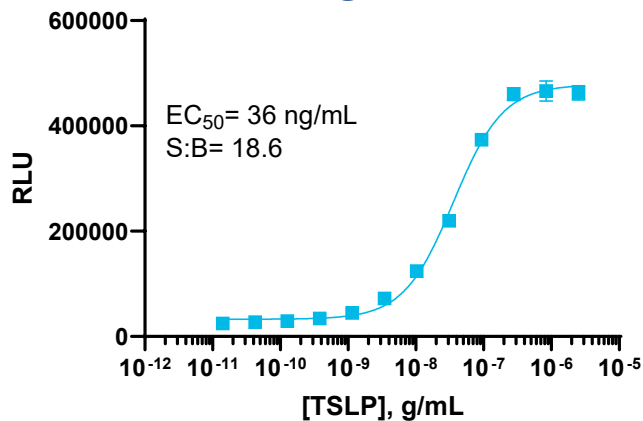
EC₅₀ = Effective Concentration at 50% of ligand S/B = signal-to-background (assay window)

Signaling Reporter Assay Design

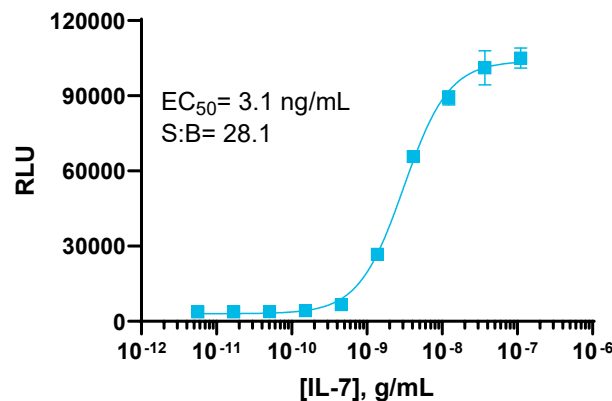
PathHunter® signaling reporter assays quantify a distal receptor signaling event (transcriptional activation)



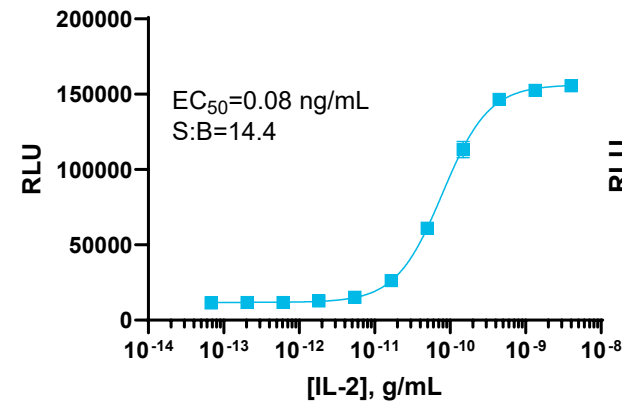
TSLP



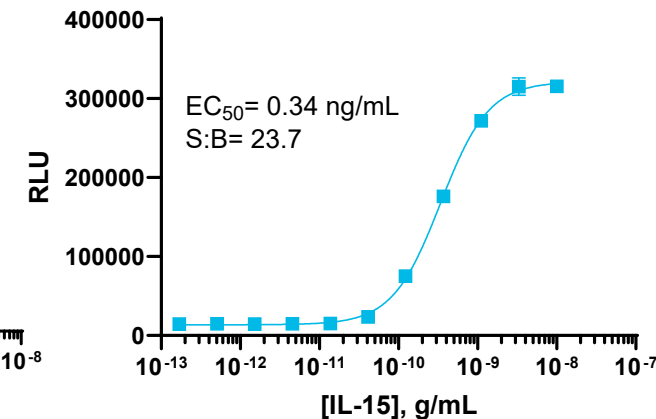
IL-7



IL-2



IL-15

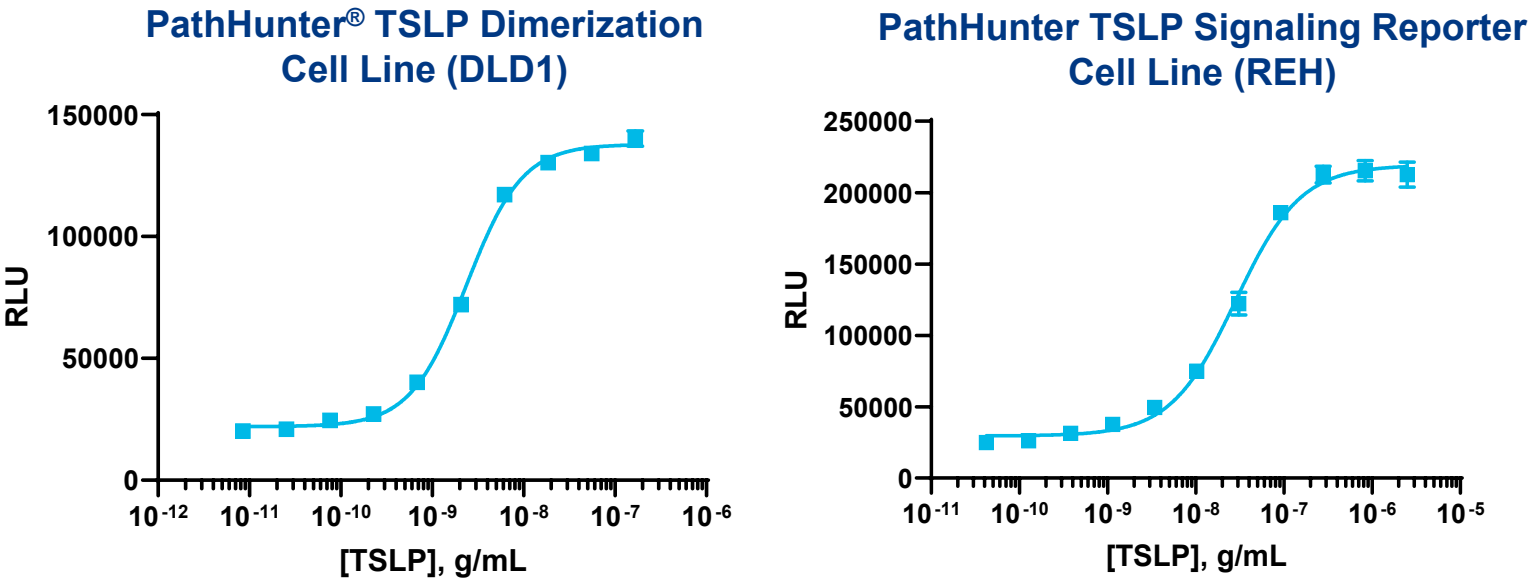


Comparison of Receptor Dimerization vs. Signaling Reporter Assays

Assay window and sensitivity to TSLP (ligand) in different assay formats

The assay window in response to TSLP is slightly larger in the reporter assay.

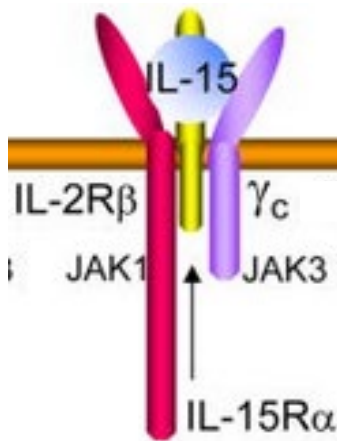
The EC₅₀ is 10-fold lower in the dimerization assay.



Assay Format →	Dimerization Cell Line	Signaling Reporter Cell Line
Cell Background	DLD1	REH
Receptor Subunits: <ul style="list-style-type: none">IL7RACRLF2	<ul style="list-style-type: none">Engineered (Truncated)Engineered (Truncated)	<ul style="list-style-type: none">EndogenousEngineered (Full-length)
Assay Window (S/B)	7.0	8.5
EC ₅₀ (ng/mL)	2.4	28.4
Incubation Time (hours)	16	16

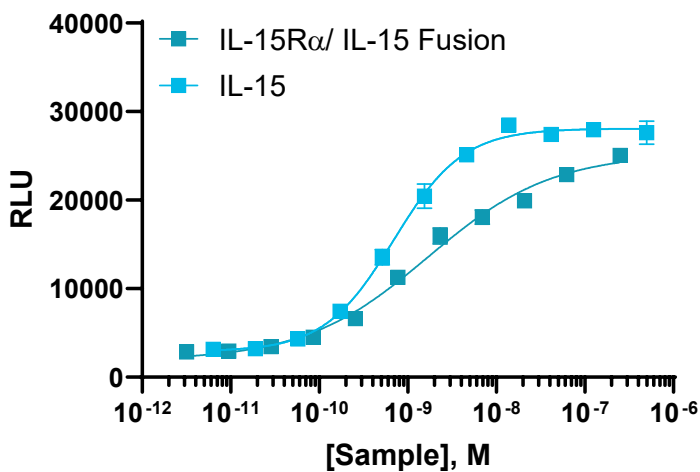
Comparison of Receptor Dimerization vs. Signaling Reporter Assays

Assay window and sensitivity to IL-15 (ligand) in different assay formats

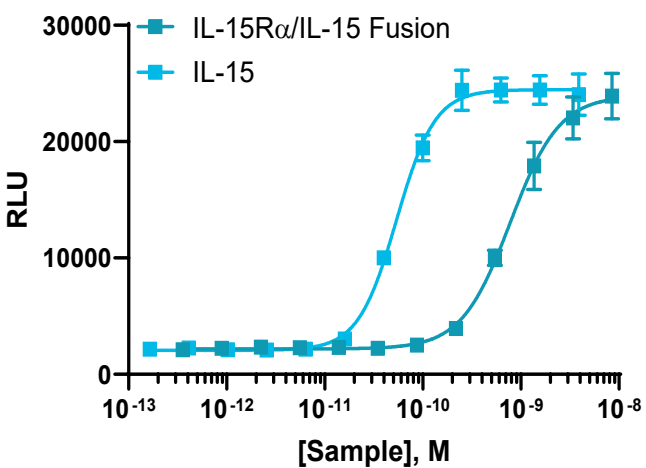


The assay window is greater in the reporter assay and exhibits greater potency (lower EC₅₀).

PathHunter® IL-15 Bioassay (Dimerization)



PathHunter IL-15 Bioassay (Reporter)

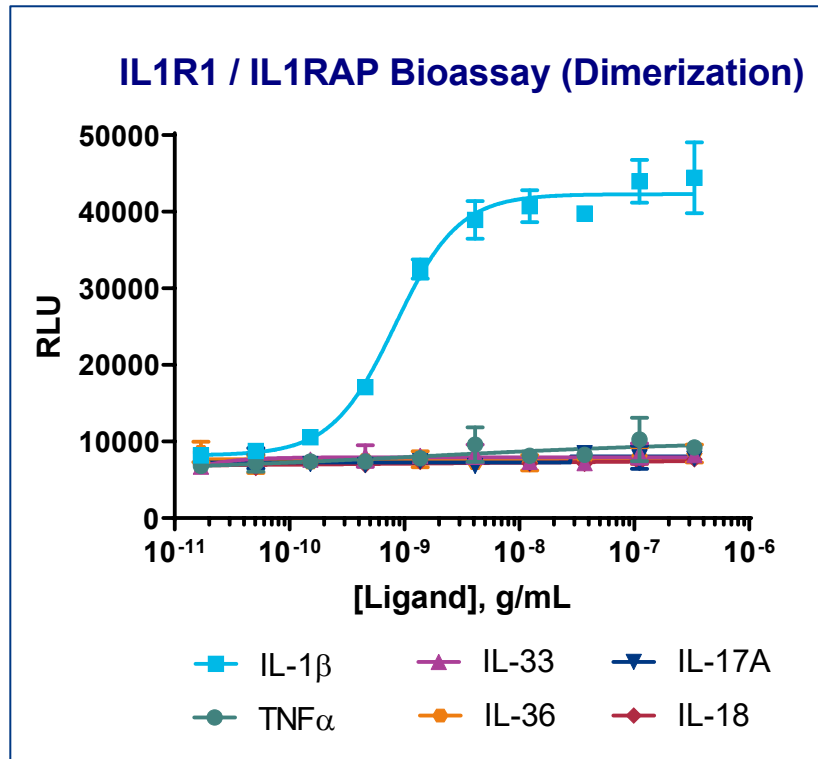


Assay Format →	Dimerization Bioassay	Reporter Bioassay
Cell Background	U2OS	REH
Receptor Subunits: <ul style="list-style-type: none">IL2RBIL2RG	<ul style="list-style-type: none">Engineered (Truncated)Engineered (Truncated)	<ul style="list-style-type: none">Engineered (Full-length)Engineered (Full-length)
Assay Window (S/B): IL-15/IL-15Ra-IL-15 fusion protein	8.8 / 8.7	14.7 / 14.7
EC ₅₀ (nM): IL-15/IL-15Ra-IL-15 fusion protein	0.69 / 1.7	0.054 / 0.78
Incubation Time (hours)	18-20	16-18

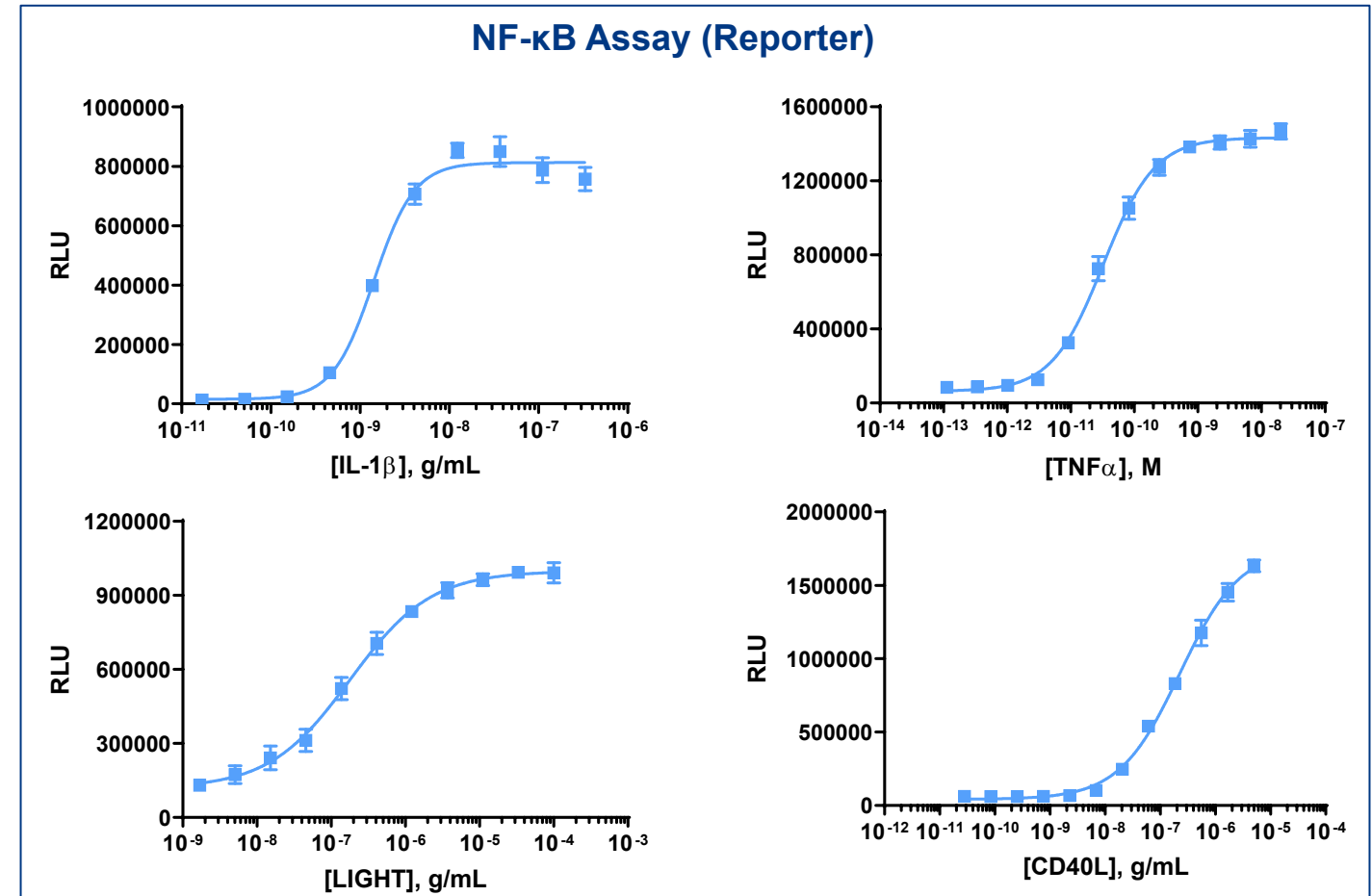
Receptor Dimerization vs. Signaling Reporter Assays

Evaluating Ligand Specificity

Dimerization assay responses are specific, while reporter assays can respond to multiple cytokine signaling pathways



Because of the receptor combination used, the assay response is cytokine-specific

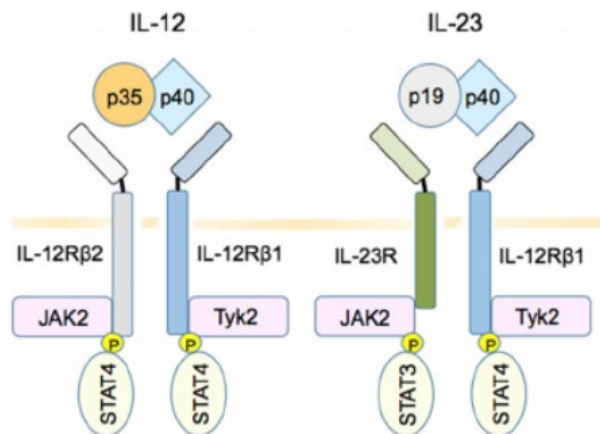


The same reporter gene can be activated by many different cytokines/ligands

Validated Therapeutic Targets for Crohn's Disease, Plaque Psoriasis, & Psoriatic Arthritis

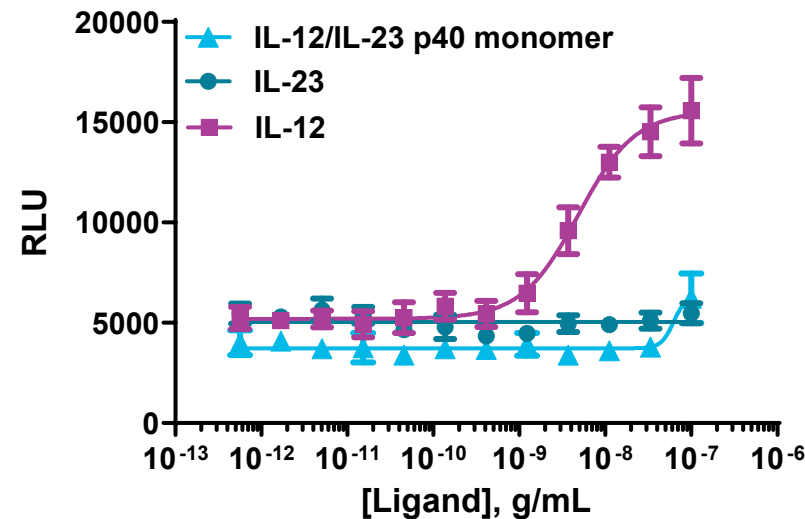
IL-12 and IL-23 dimerization assay specificity

High specificity observed for cognate ligand in each assay compared to little or no specificity observed with p40 monomer (a subunit shared by both ligands).



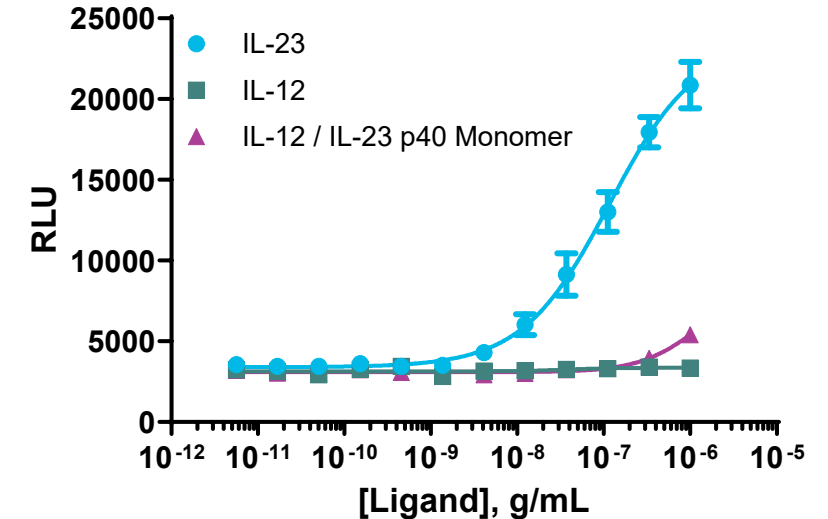
IL-12 Receptor

PathHunter® IL-12RB1/IL-12RB2 Cell Line



IL-23 Receptor

PathHunter IL-23/IL-12RB1 Cell Line

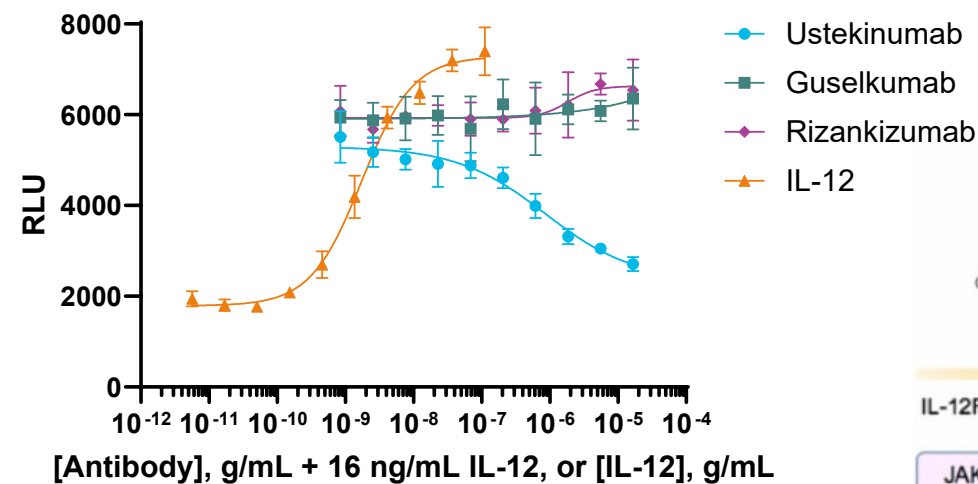


Therapeutic Targeting p40 Subunit of IL-23 & IL-12 Is Active in Both Assays

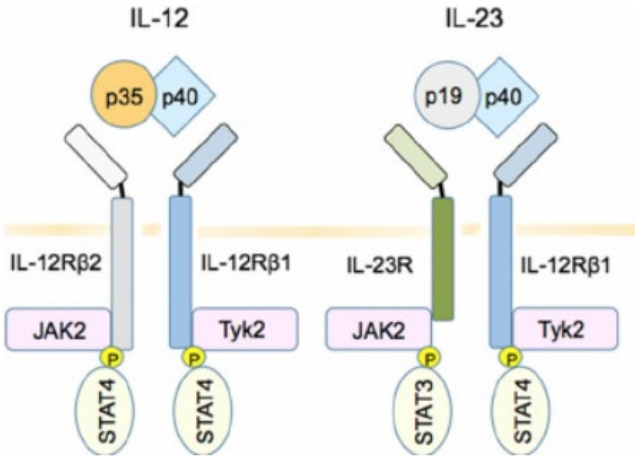
Similar potency for key therapeutic antibodies targeting p19 subunit of IL-23 in the IL-23 receptor assay, but only Ustekinumab is active in IL-12 receptor assay

IL-12 Receptor

PathHunter® IL-12RB1/IL-12RB2 Cell Line



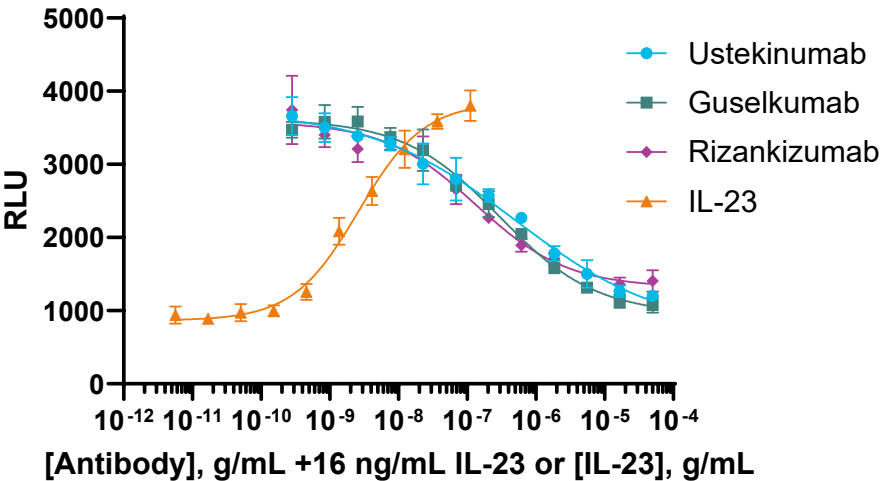
Antibody/Cytokine	Assay Window (S/B)	EC ₅₀ (ng/mL)
Ustekinumab	2.0	857
Guselkumab	-	-
Rizankizumab	-	-
IL-12	3.8	1.7



Ustekinumab (Stelara®) = anti-p40
Guselkumab (Tremfya®) = anti-p19
Rizankizumab (Skyrizi®) = anti-p19

IL-23 Receptor

PathHunter IL-23/IL-12RB1 Cell Line



Antibody/Cytokine	Assay Window (S/B)	EC ₅₀ (ng/mL)
Ustekinumab	3.0	511
Guselkumab	3.3	298
Rizankizumab	2.8	133
IL-23	4.0	2.6

All registered trademarks are the property of their respective owners.

Signaling Reporter Assay Measure IL-4 Activation of Both Type I & Type II IL-4 Receptor Complexes

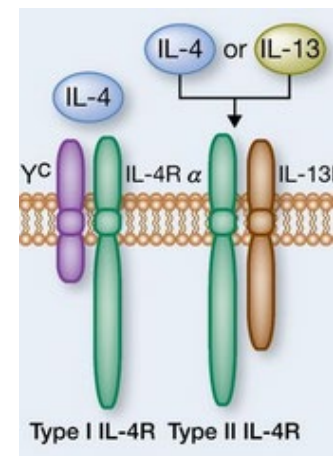
IL-4 receptor (IL-4R) complexes signaling via STAT6

- Type I IL-4R
 - IL-4R + IL-2RG (γ C)
 - Expressed in immune cells and binds only IL-4
- Type II IL-4R
 - IL-4R + IL-13R α
 - Expressed in epithelial cells and binds both IL-4 and IL-13

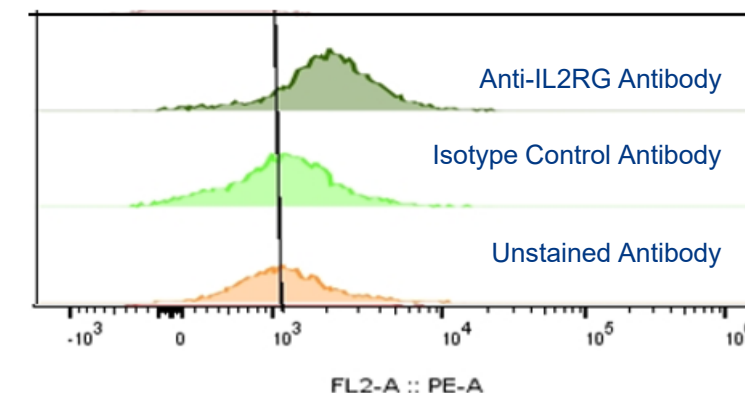
The PathHunter® IL-4/IL-13 signaling reporter assay was used to quantify IL-4 and IL-13 signaling

- Flow cytometry data indicates IL2RG (γ C) receptor is also expressed in the cell line

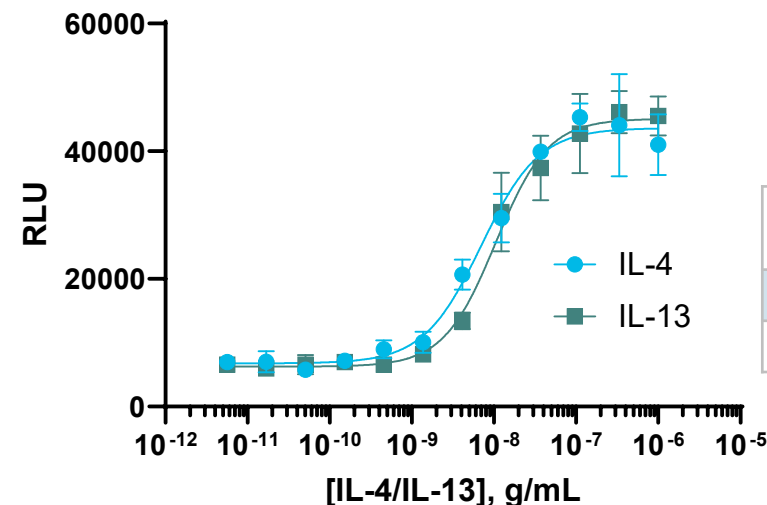
IL-4 results measure combined signaling from both Type I and Type II receptor complexes



Flow Cytometry Analysis



PathHunter IL-4/IL-13 Signaling Reporter Cell Line

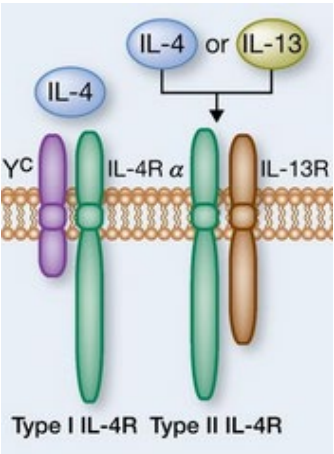


Cytokine	Assay Window (S/B)	EC ₅₀ (ng/mL)
IL-4	6.3	10.0
IL-13	7.7	61.4

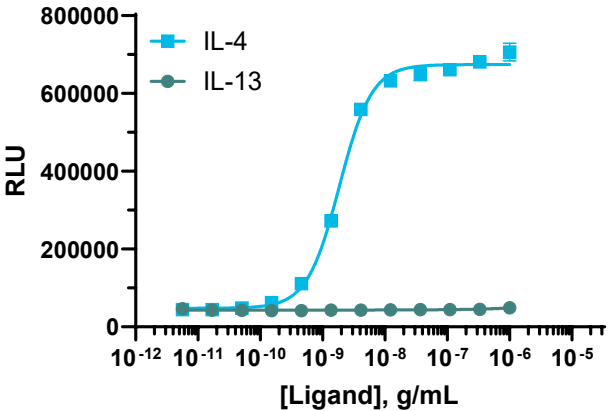
IL-4R Dimerization Assays Are Specific for Either Type I or Type II Receptor Complexes

IL-4 Receptor Dimerization Cell Lines can be used to measure signaling specific for each IL-4 receptor complex

- IL-4 variants that induce exclusively Type I– or Type II–dependent responses could preserve the benefits of IL-4 immunotherapy, but with reduced side effects.

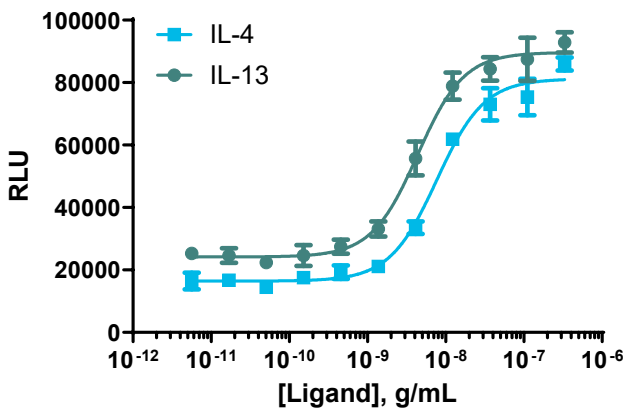


Type I IL-4R: Inflammation
PathHunter® IL-4R/IL-2RG Dimerization Cell Line



Cytokine	S/B	EC ₅₀ (ng/mL)
IL-4	16	1.9
IL-13	-	-

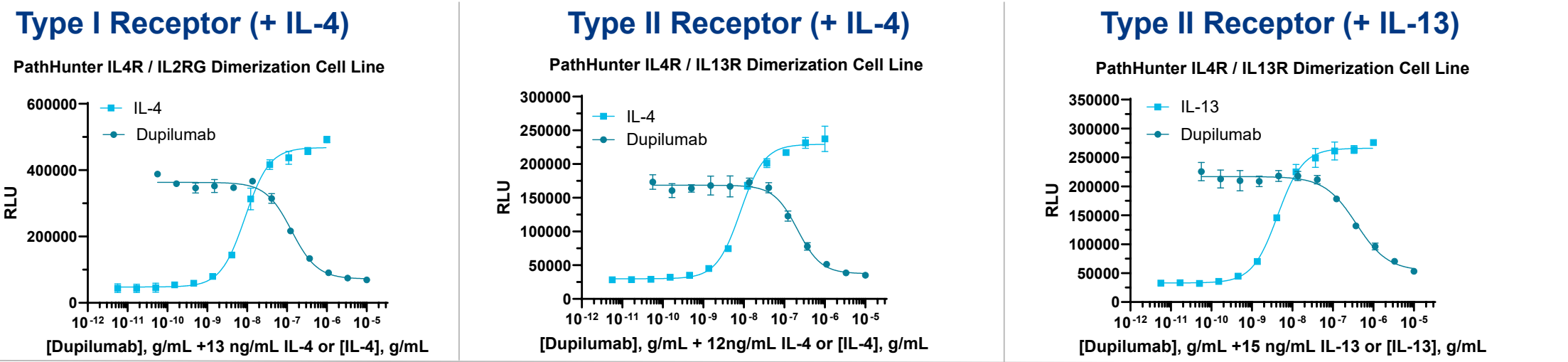
Type II IL-4R: Oncology & Inflammation
PathHunter IL-4R/IL-13R Dimerization Cell Line



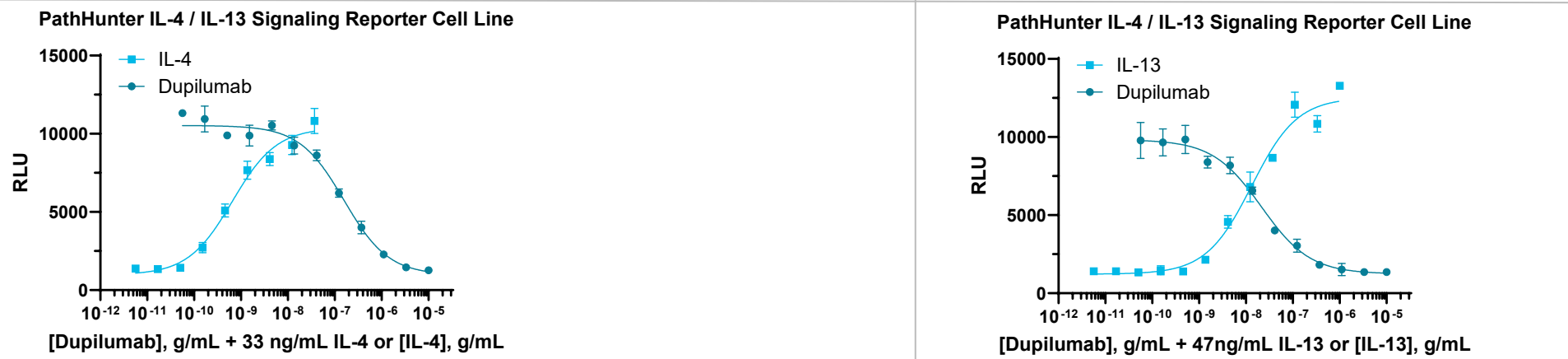
Cytokine	S/B	EC ₅₀ (ng/mL)
IL-4	5.2	7.7
IL-13	3.8	4.4

Dupilumab, an Approved Therapeutic for Eczema & Asthma, Blocks IL-4 & IL-13 Signaling through Both Type I & II IL-4 Receptors

Receptor Dimer



Reporter



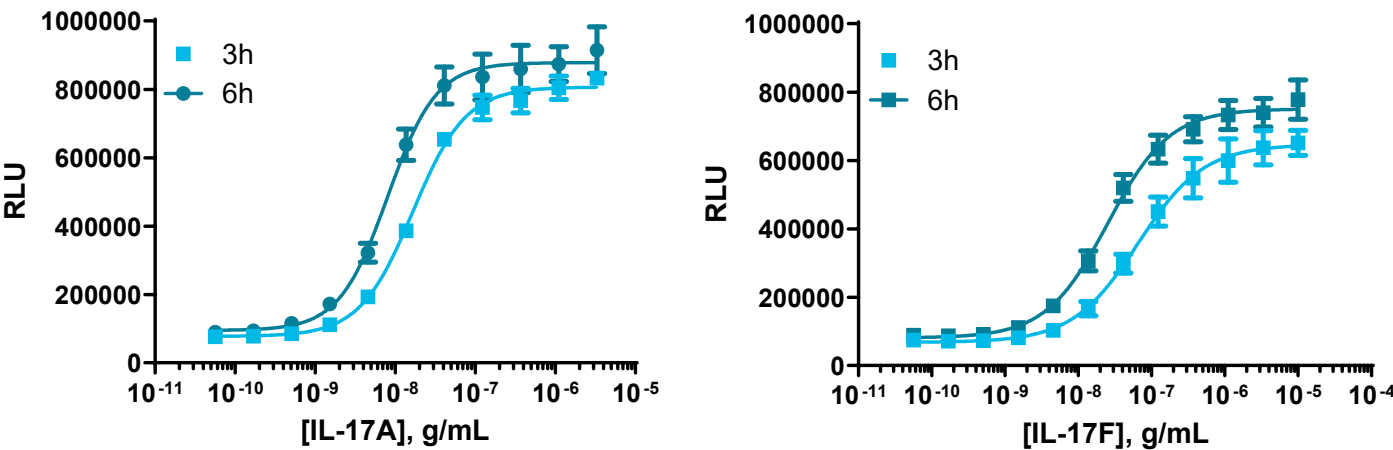
	IL-4		Dupilumab+IL-4		IL-13		Dupilumab+IL-13	
Cell Line	S/B	EC ₅₀ (ng/mL)	S/B	IC ₅₀ (ng/mL)	S/B	EC ₅₀ (ng/mL)	S/B	IC ₅₀ (ng/mL)
IL4R-IL13R Dimerization	8.3	8.4	4.9	210	8.4	4.8	4.2	378
IL4R-IL2RG Dimerization	11.2	8.9	5.6	130	NA	NA	NA	NA
IL-4/IL-13 Signaling Reporter	11.4	0.65	8.9	157.5	9.5	13.2	7.7	21.4

Secukinumab Inhibits IL-17A-mediated Dimerization

Cosentyx® (Secukinumab), which binds to IL-17A, is the first anti-IL-17A therapeutic approved for treatment of plaque psoriasis and psoriatic arthritis

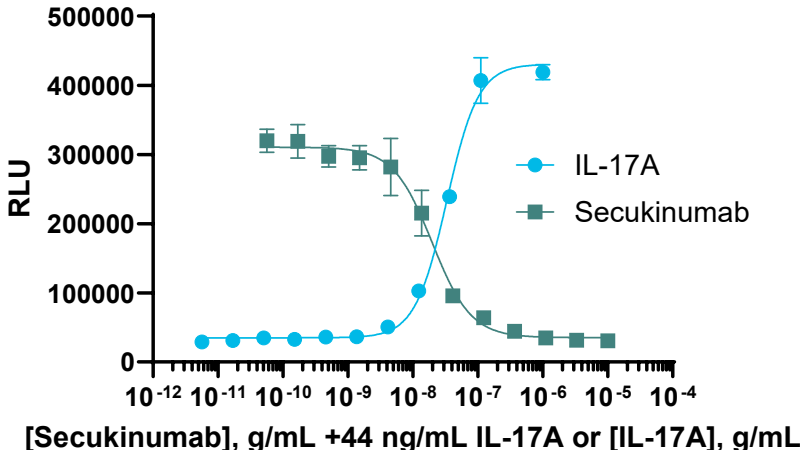
PathHunter® IL-17RA/IL-17RC Dimerization Cell Line Time Course & Inhibition Experiments

Time Course with Ligand (IL-17A and IL-17F)



Cytokine (time point)	Assay Window (S/B)	EC ₅₀ (ng/mL)
IL-17A, 3h	11.6	16.6
IL-17A, 6h	11.6	8.1
IL-17F, 3h	8.8	65.4
IL-17F, 6h	8.8	25.2

Inhibition with Secukinumab



Antibody/Cytokine	Assay Window (S/B)	EC ₅₀ (ng/mL)
IL-17A	14.5	33.3
Secukinumab	10.4	19.8

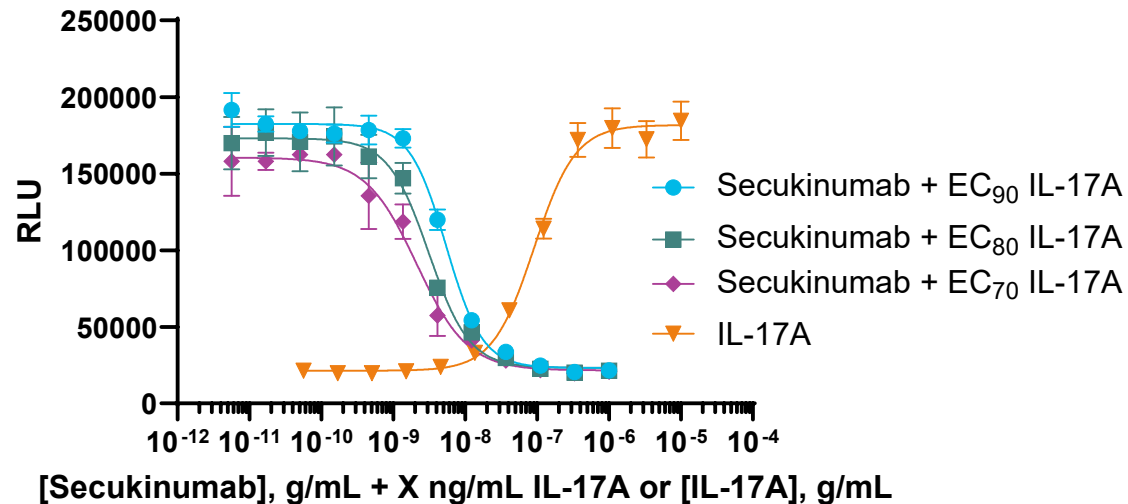
All registered trademarks are the property of their respective owners.

Receptor Dimerization Assays Maintain Robust Activity in the Presence of Normal Human Serum

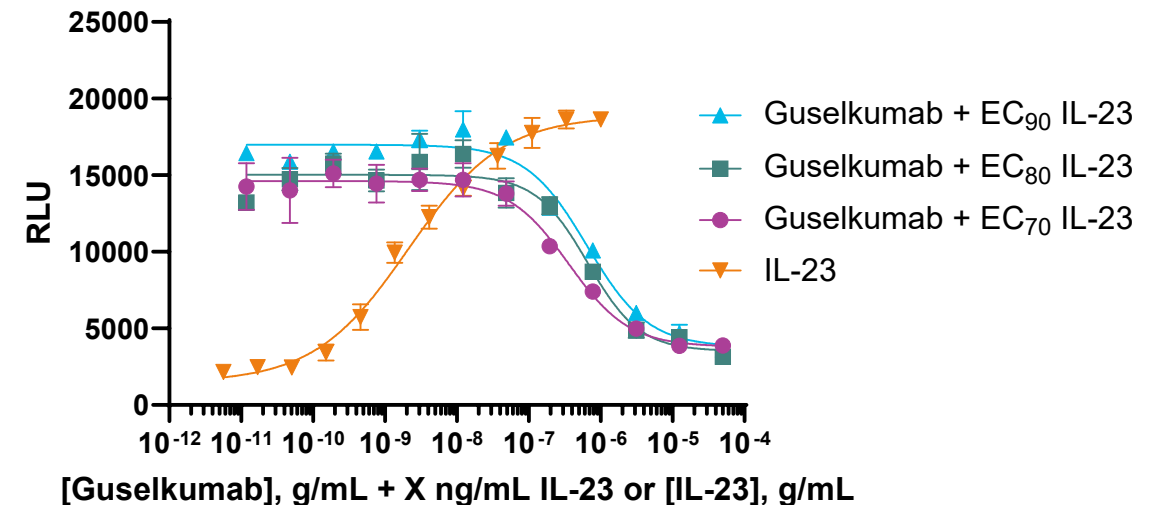
Dimerization assays produce robust agonist and antagonist responses in the presence of 10% normal human serum (NHS), which is required for quantifying the presence of anti-drug neutralizing antibodies (NAb) in human serum samples.

Dimerization Assay Experiments with Samples Prepared in 10% Pooled NHS

PathHunter® IL-17RA/IL-17RC Dimerization Cell Line

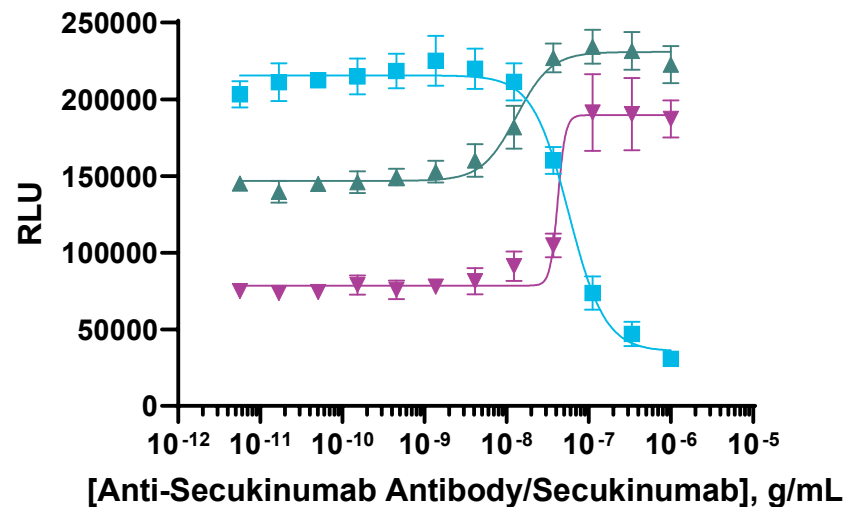


PathHunter IL-23/IL-12RB1 Dimerization Cell Line



Dimerization Assay Experiment with Samples Prepared in 10% Pooled NHS

PathHunter® IL-17RA/IL-17RC Dimerization Cell Line



- Secukinumab+40n g/mL IL-17A
- ▲ Anti-Secukinumab+50ng/mL Secukinumab +40 ng/mL IL-17A
- ▼ Anti-Secukinumab+100ng/mL Secukinumab +40 ng/mL IL-17

Antibodies/Cytokine	Assay Window (S/B)	EC / IC / NC ₅₀ (ng/mL)
IL-17A	9.5	8.4
Secukinumab +40ng/mL IL-17A	6.6	57
Anti-Secukinumab+50ng/mL Secukinumab+40ng/mL IL-17A	1.5	13
Anti-Secukinumab+100ng/mL Secukinumab+40ng/mL IL-17A	2.5	42

FDA guidelines specify a minimum sensitivity of 1 µg/mL for detection of NAb in human serum

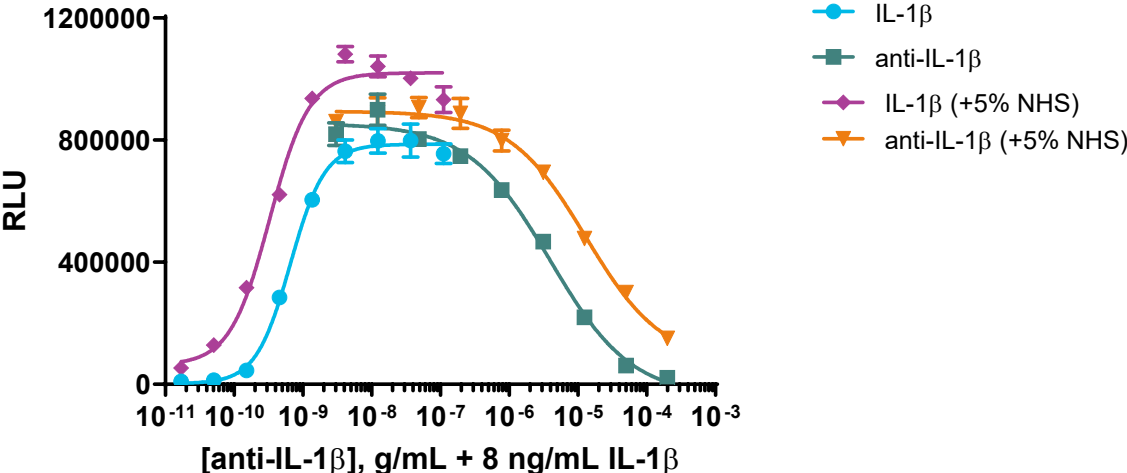
An anti-Secukinumab antibody exhibits potent neutralizing activity of Secukinamb IL-17A inhibition and demonstrates application of the assay for Nab antibody screening of patient serum.

The IL-1 β Reporter Assay Is More Sensitive to the Presence of NHS Format than the Dimerization Assay

Attenuated assay windows and increased IC₅₀ are observed with the IL-1 β reporter assay format in the presence of NHS while dimerization assay results are similar in the presence or absence of NHS

IL-1 β Reporter Assay

U2OS NF- κ B-Reporter Cell Line
CP5, 5K/well; 6hr ligand @ 37°C

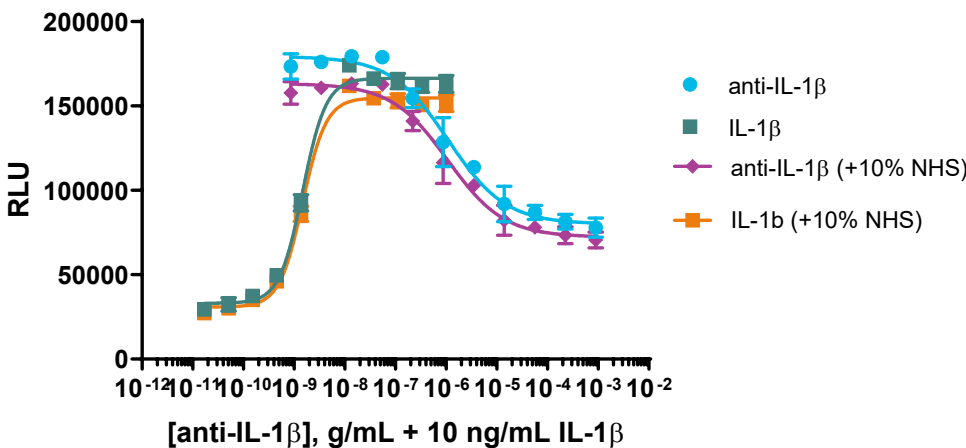


Sample	PBS		+ 5% NHS	
	S/B	EC/ IC ₅₀ , ng/mL	S/B	EC/ IC ₅₀ , ng/mL
IL-1 β (agonist mode)	50.1	0.66	19.5	0.33
Anti-IL-1 β + 8 ng/mL IL-1 β	40.3	4,445	5.7	13,020

Significant loss of assay window and increased IC₅₀ with anti-IL-1 β drug in 5% NHS

IL-1 β Receptor Dimerization Assay

U2OS ILR1A / IL1RAP Bioassay
Part #: 93-1032Y3; Lot #: 17C2411
CP5; O/N recovery; 30' premix @ R/T; 16h ligand @ 37°C



Sample	PBS		+ 10% NHS	
	S/B	EC/ IC ₅₀ , ng/mL	S/B	EC/ IC ₅₀ , ng/mL
IL-1 β (agonist mode)	6.0	1.49	5.5	1.49
Anti-IL-1 β + 10 ng/mL IL-1 β	2.2	1,104	2.2	1,082

No change in assay window or IC₅₀ with anti-IL-1 β drug in 10% NHS

Development of Cytokine Assays for Relative Potency, Characterization & Lot Release Applications

Reflective of MOA of the drug/molecule

Stability-indicating

Relative potency range (50%-150%)

Repeatable, robust, precise, and accurate

Reproducible

- Intra-lot (plate-to-plate)
- Inter-lot (lot-to-lot)
- Analyst-to-analyst

Long-term support – Dedicated well characterized cell banks

Supply agreements

Traceability and quality (audit-ready)

ICH-based Assay Qualification



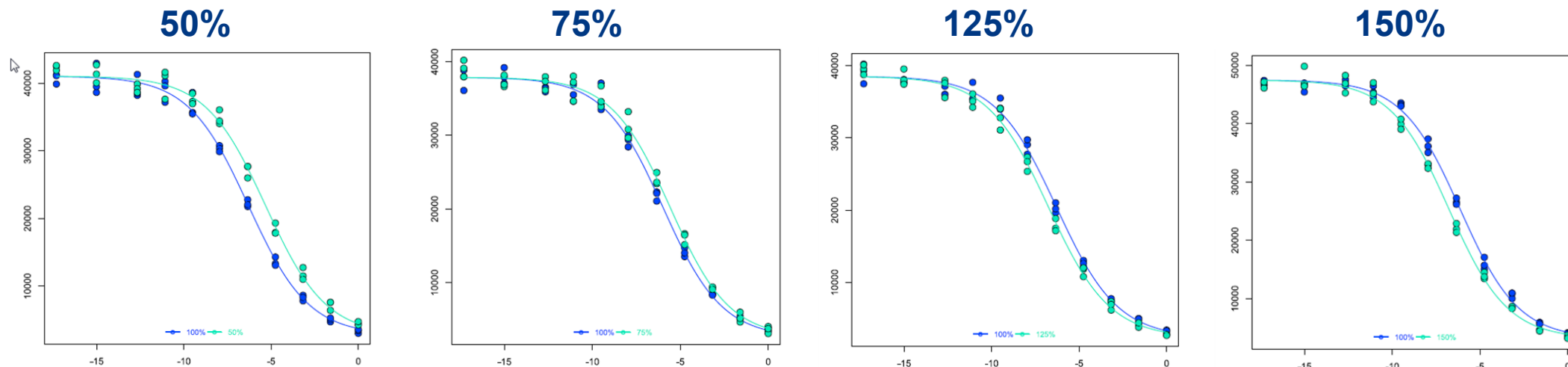
MOA = mechanism of action; ICH = International Council for Harmonization

Relative Potency Analysis for the Tocilizumab (IL-6) Bioassay

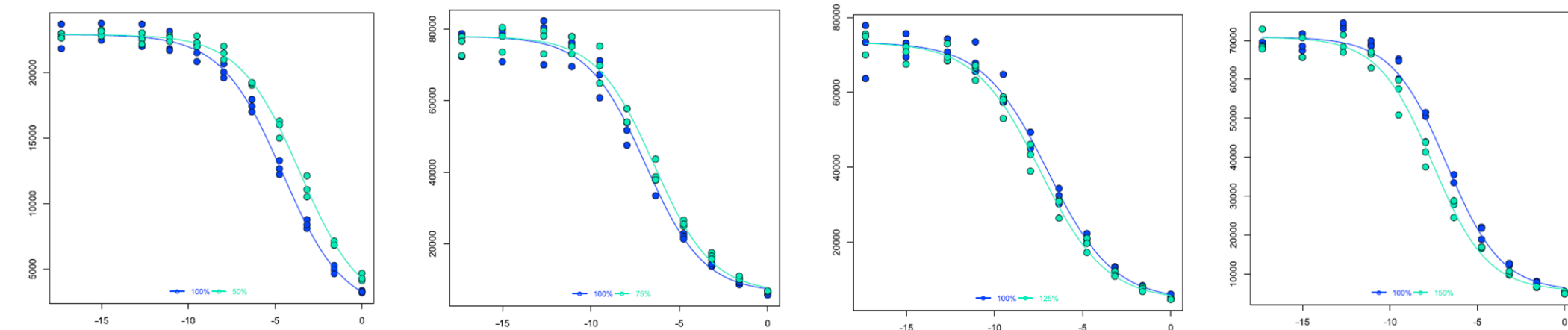
Relative potency analysis for 4 nominal concentrations of tocilizumab

ACTEMRA®
tocilizumab

Analyst 1



Analyst 2



● Reference
● Sample

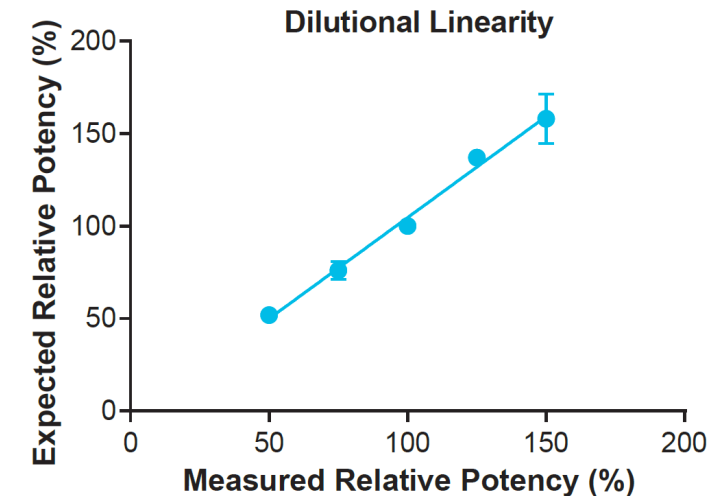
RP values calculated in PLA 3.0. All registered trademarks are the property of their respective owners.

Tocilizumab (IL-6) Bioassay Demonstrates Good Accuracy, Intermediate Precision, & Dilutional Linearity



Qualification of Actemra® using the Tocilizumab Bioassay

Experiment #	Analyst #	Expected RP (%)	Observed RP (%)	Average RP (%)	% RSD	% Recovery
1	1	150	155.5	158.1	8.44	105.4
2	1		163.2			
3	1		141.1			
4	2		172.7			
1	1	125	138.6	137.1	1.66	109.7
2	1		137.4			
3	1		138.6			
4	2		133.8			
1	1	75	74.1	76.1	6.34	101.5
2	1		83			
3	1		75.4			
4	2		71.9			
1	1	50	54.8	52	4.96	103.9
2	1		50.6			
3	1		53.3			
4	2		49.1			



Accuracy: 105.1%

Intermediate Precision: ≤ 8.4%

Dilutional Linearity: $R^2 = 0.9927$

All registered trademarks are the property of their respective owners.

Functional, cell-based assays available for most targets in cell line or assay ready formats (eXpress™ & bioassay kits)

IL-1-like

IL-1
IL-18
IL-33

IL-28-like

IL-28
IL-29

IL-17-like

IL-17
IL-25

IL-4-like

IL-3
IL-4
IL-5
IL-13
CSF2

Non-classified

IL-8
IL-34
CSF1

IL-6/12-like

IL-6
IL-11
IL-12
IL-23
IL-27
IL-31
CSF3

γ-chain utilizing

IL-2R $\alpha\beta\gamma$
IL-2R $\beta\gamma$
IL-4
IL-7
IL-9
IL-15
IL-21
TSLP

IL-10-like

IL-10
IL-19
IL-20
IL-22
IL-24
IL-26

Chemokines

CCR1	CCR10
CCR5	CCR6
CCR4	CCR9
CXCR3	CXCR4
CXCR2	CXCR7
CCR2	CCR3
CCR7	CCR8
CCRL1	CCRL2
CXCR5	CXCR6
CX3CR1	CXCR1

Other Cytokines

TNF α , IFN α , IFN β , IFN γ , CXCRs,
LIF, OSM, G-CSF, GM-CSF, RANK

PathHunter® Cytokine Receptor Assays

- MOA-based dimerization, signaling reporter, and other assay types covering over 80% of targets
- Available in stable cell line or ready-to-use eXpress™ and bioassay kit formats
- Used from discovery through development and into commercial release and stability QC lot testing
- Bioassays are target-based or qualified with a reference ligand or therapeutic

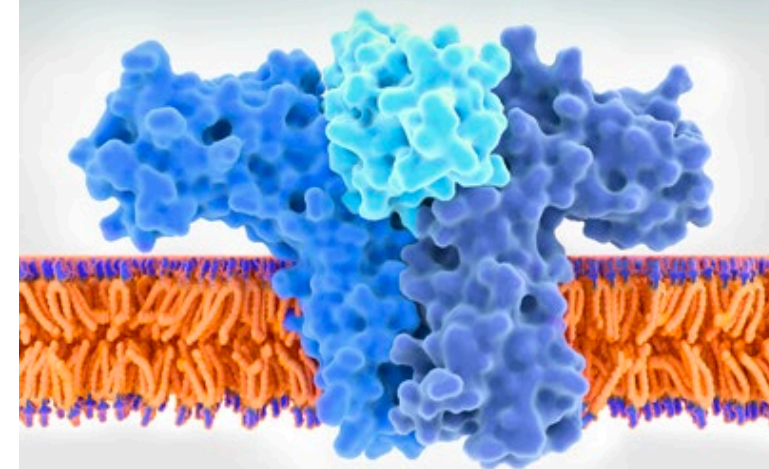
PathHunter Dimerization Assays

- Robust, sensitive, and highly specific assays
- Minimally impacted by the presence of human serum
- Suitable for both relative potency and NAb detection

PathHunter Reporter Assays

- Robust and sensitive but lack receptor specificity
- Provide large assay windows and good sensitivity to ligands
- Suitable for screening and relative potency assays

Cytokine Receptor Product Solutions



Learn more at discoverx.com/target-class/cytokine-receptor/