



Accelerating Development of Therapeutic Interventions for COVID-19

Gaurav Agrawal, Ph.D.

Scientific Development Manager, Eurofins DiscoverX

OUR EXPERTISE
IN YOUR HANDS.
DISCOVER
CONFIDENTLY.

Eurofins DiscoverX is a Global Leader in Supporting Early to Late Stage Drug Discovery

From Discovery to Development to Clinic to Post-Market

20+ years of enabling drug discovery and development programs

10+

Druggable
target classes

1500+

Stable cell line and
membrane preps

20+

core patents

Validated

> 30 Billion data points screened in
assay services with same assays

2000+

Publications across
multiple applications

55+

Qualified and MOA
Based Bioassays

**ICH Based Bioassay
Qualification**

Facilitate downstream validation studies

**Certified CRO
Partners**

Scientific training to enable global
CROs

**Dedicated Scientific
Support**

Experienced team providing scientific support


**20+ Successful
Assay Transfers**

At clients/affiliated CRO sites

50+ Global Programs

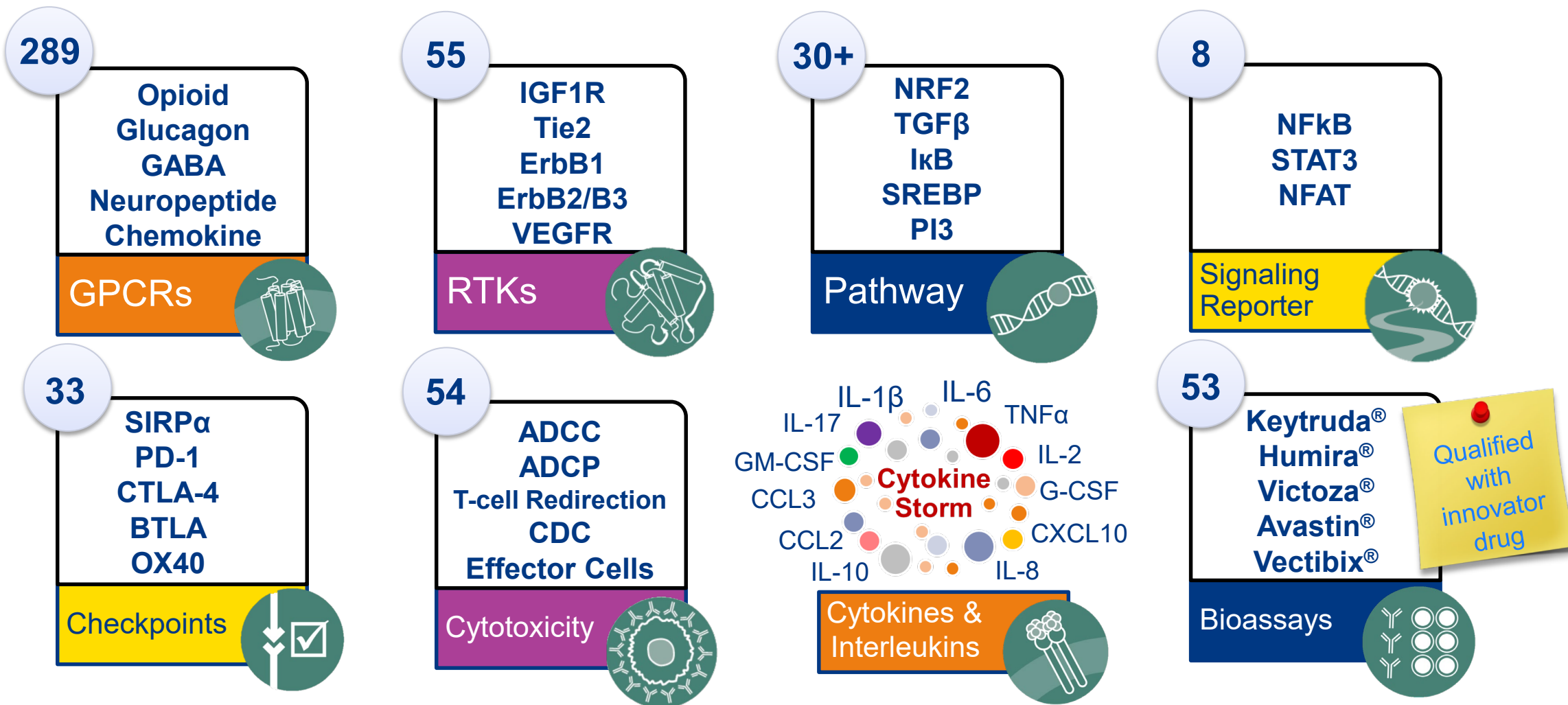
For potency, stability and Nab testing

Global programs using DiscoverX Assays for potency, stability and NAb testing for Drug Release

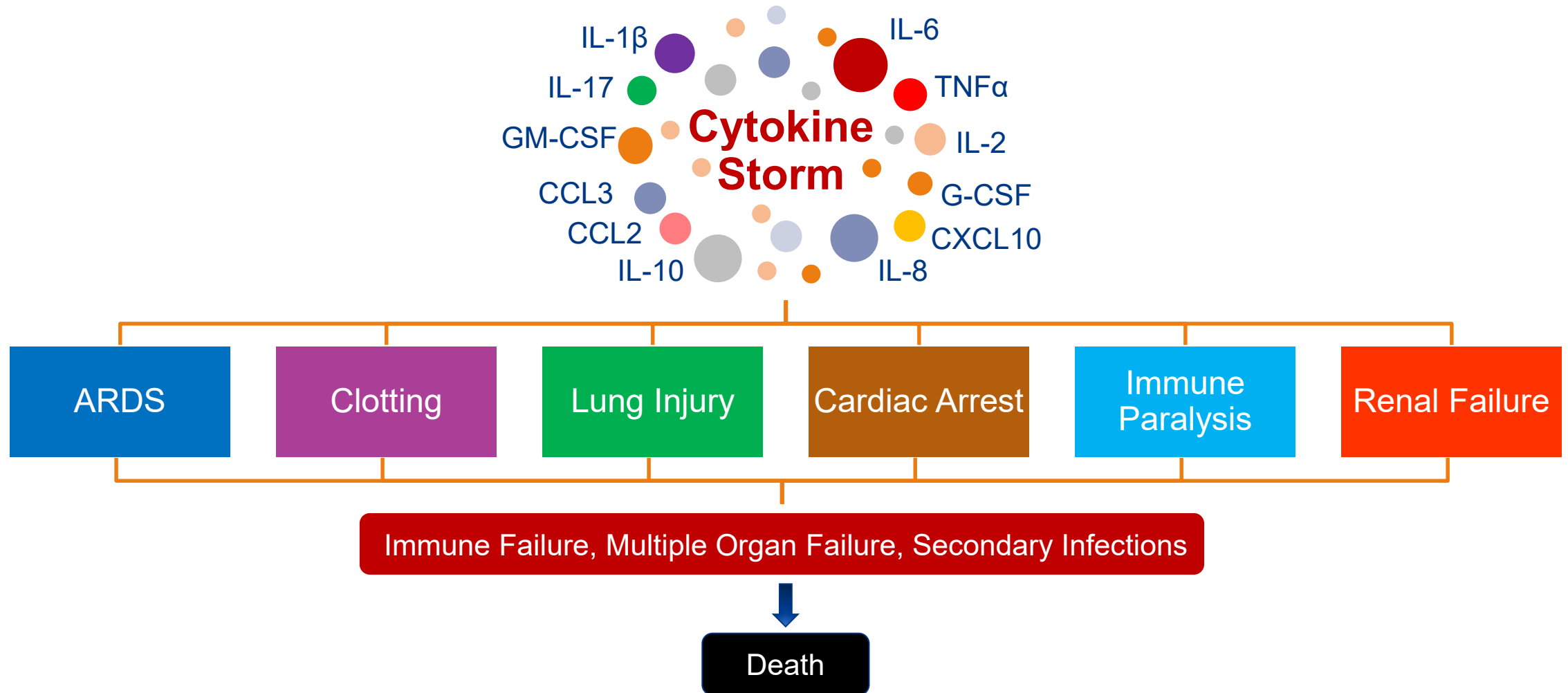
SIRPα Bioassay Multiple Programs in North America for originator biologics Nab assessment programs at CROs	PD-1 Bioassay Multiple Programs in North America and APAC Clinical Development of Biosimilars	VEGF Bioassay Multiple Programs in North America, EU and APAC Clinical Development of Biosimilars	Insulin Bioassay Multiple Programs in and APAC for clinical Development of Biosimilars	CNR1 Bioassay US based Biopharma for originator biologic	GLP1R Bioassay Multiple Programs in North America, EU and APAC Clinical Development of Biosimilars in NA	IGF1R Bioassay Multiple Programs in North America and APAC for originator biologic
Bradykinin Bioassay US based CRO for undisclosed client	PTH1R Bioassay Multiple Programs in North America, EU and APAC Clinical Development of Biosimilars	FGFR Bioassay Nab assessment program for US based Biopharma			ErbB2/ErbB3 Bioassay Multiple Programs in North America, EU and APAC Clinical Development of Biosimilars	Melanocortin Bioassay European Biopharma for an originator biologic
IL-1 Bioassay Multiple Programs in NA for originator biologic	IL-2 Bioassay Multiple Programs in North America and EU for originator biologic	IL-7 Bioassay European Biopharma for originator biologic	IL-10 Bioassay Multiple Programs in North America for originator biologic	IL-31 Bioassay US based pharma for Nab assessment and potency testing programs for originator biologic	RANK Assay Multiple Programs in and APAC for clinical Development of Biosimilars	GM-CSF Bioassay Multiple Programs in EU and NA for originator biologic

Largest Menu of Cell-Based Assays for Discovery, Potency, & NAb Assay Development

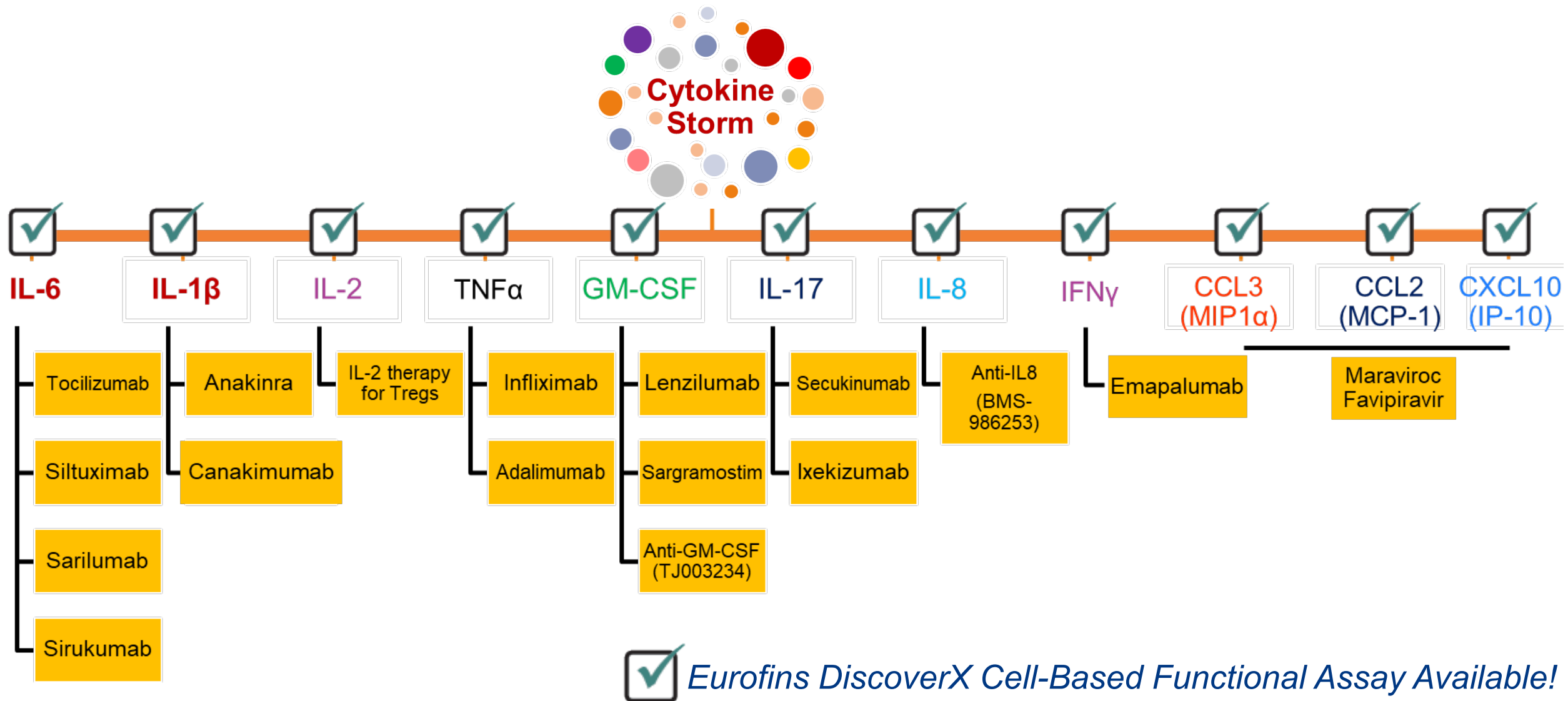
>800 cell lines to support bioassay development for major drug target classes



Cytokine Storm in COVID-19 Pathology

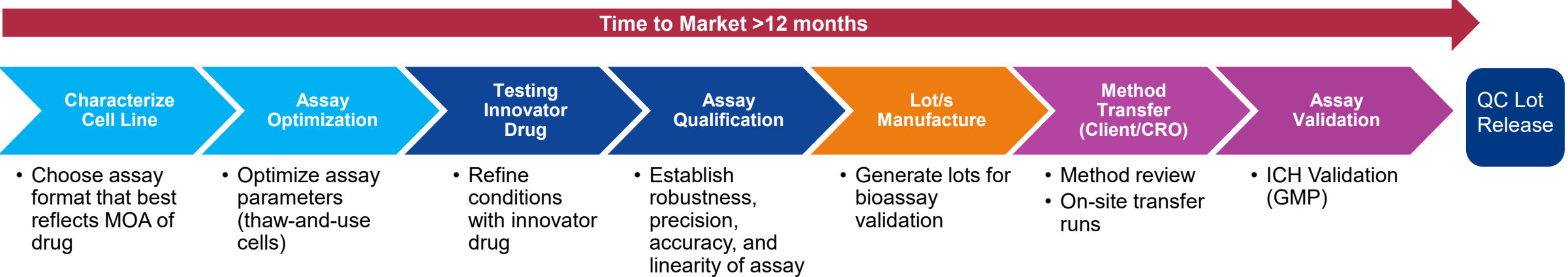


Withstanding the Cytokine Storm in COVID-19: Therapeutic Approach

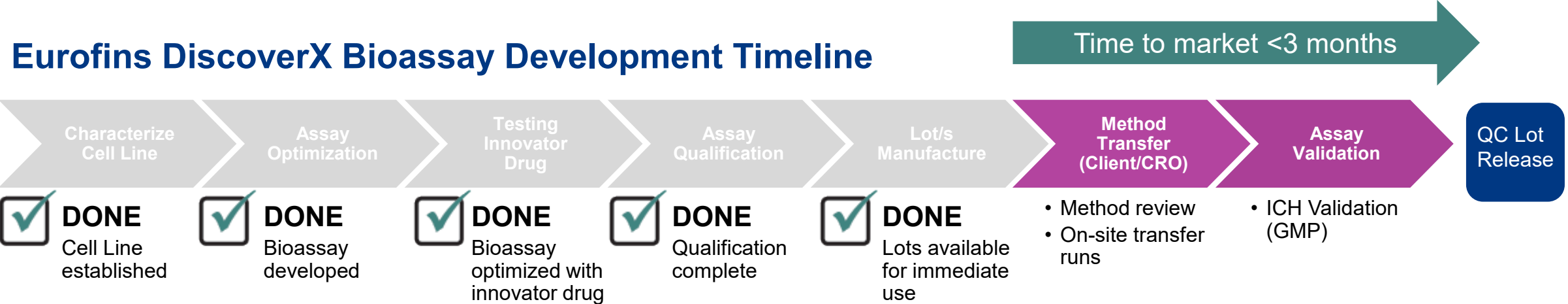


Accelerating Implementation Phase for QC Lot Release with Qualified Bioassays

Typical Bioassay Development Timeline



Eurofins DiscoverX Bioassay Development Timeline



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

Analytical Assays for Therapeutics Targeting Key Proinflammatory Cytokines

1. Qualified Bioassays:

- Case Study with PathHunter® Tocilizumab Bioassay - Qualified with  **ACTEMRA®** *tocilizumab* for therapeutics targeting IL-6 pathway
- IL-1 β - PathHunter® Anakinra Bioassay - Qualified with  **Kineret®** (anakinra)
- GM-CSF - PathHunter® Sargramostim Bioassay – Qualified with **Leukine®** *sargramostim*
A Recombinant GM-CSF–Yeast-Expressed
- TNF α - PathHunter® Adalimumab Bioassay - Qualified with **HUMIRA** *adalimumab*

2. Signaling Reporter Assays

- IL-6
- TNF α
- RANK (TNFSF11)

3. Functional Cell-Based Assays for Small Molecule Inhibitors

- JAK1, JAK2 and JAK3

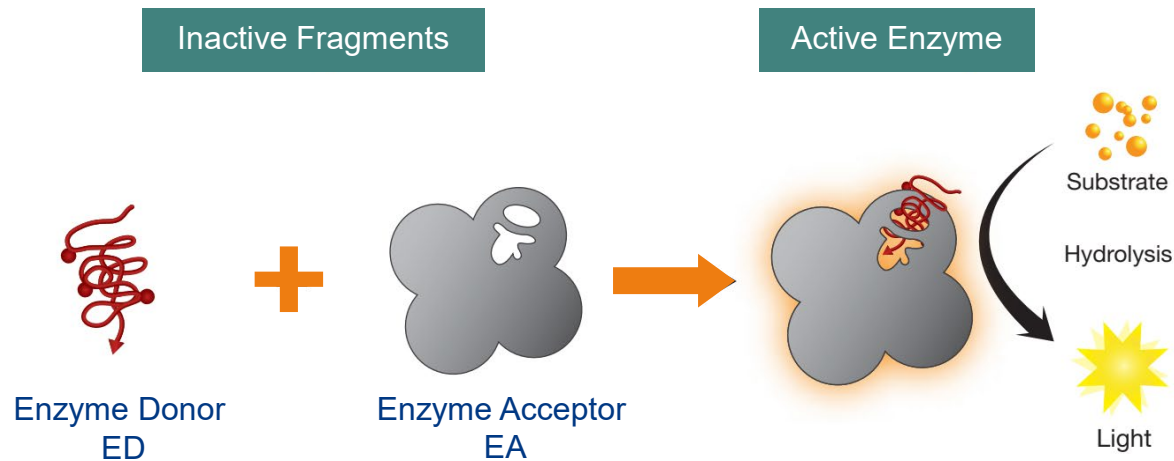
Technology Overview

Enzyme Fragment Complementation



Enzyme Fragment Complementation (EFC) Platform

Enabling Technologies with a Flexible Platform based on a Split β -Galactosidase Enzyme



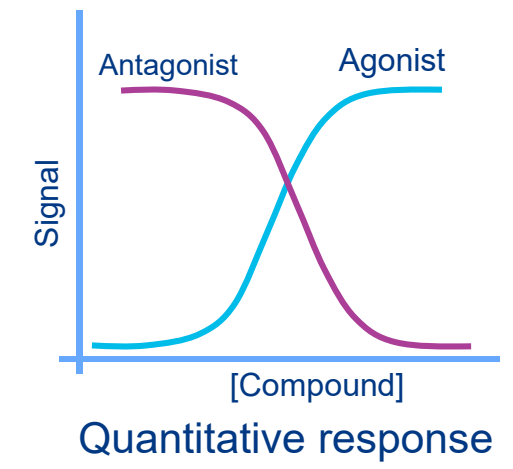
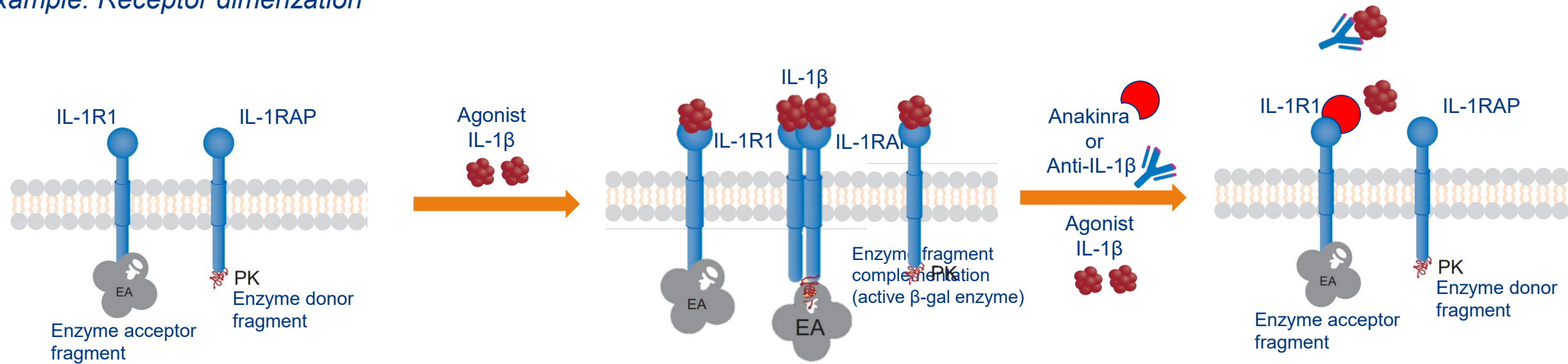
EFC Assay Principle:

- Complementation between two inactive enzyme fragments called the Enzyme Donor (ED) and Enzyme acceptor (EA) results in formation of an active β -gal enzyme.
- The active functional β -gal enzyme hydrolyzes its substrate to produce chemiluminescent signal.

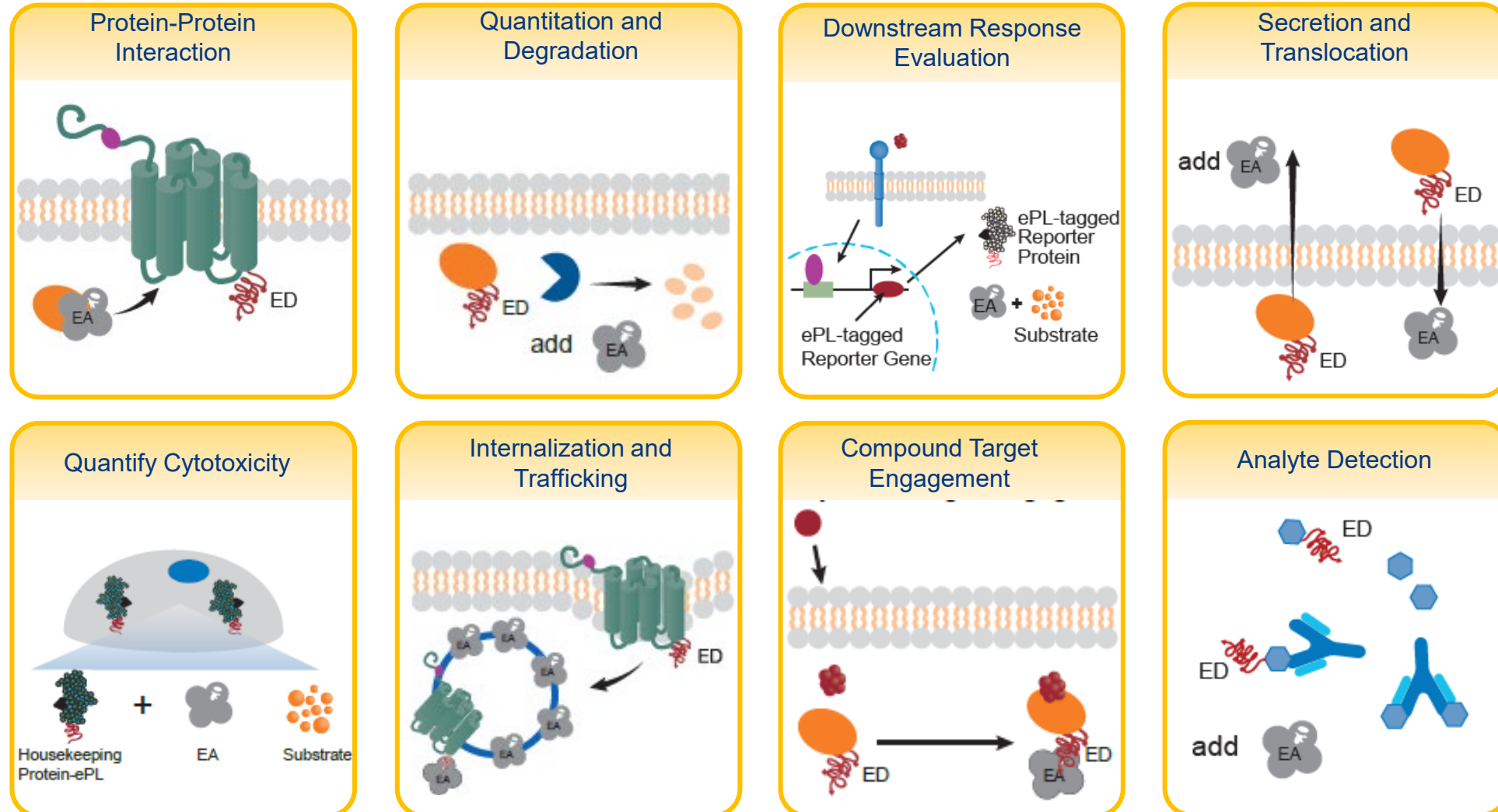
- **Homogenous Format** — Mix-and-read assay format that does not require washing, centrifugation, or filtration
- **Robust** — Enzymatically-amplified assays with a large signal-to-background ratio and high precision with Z' factors >0.7 and lot-to-lot reproducibility
- **Qualified and Validated** — Extensively optimized for hundreds of targets used for screening in billions of data points, and thousands of peer-reviewed publications
- **Scalable** — Easily scalable and HTS-friendly from 96- to 1536-well microplate format

MOA-Based Assays Design using Enzyme Fragment Complementation Technology

Example: Receptor dimerization

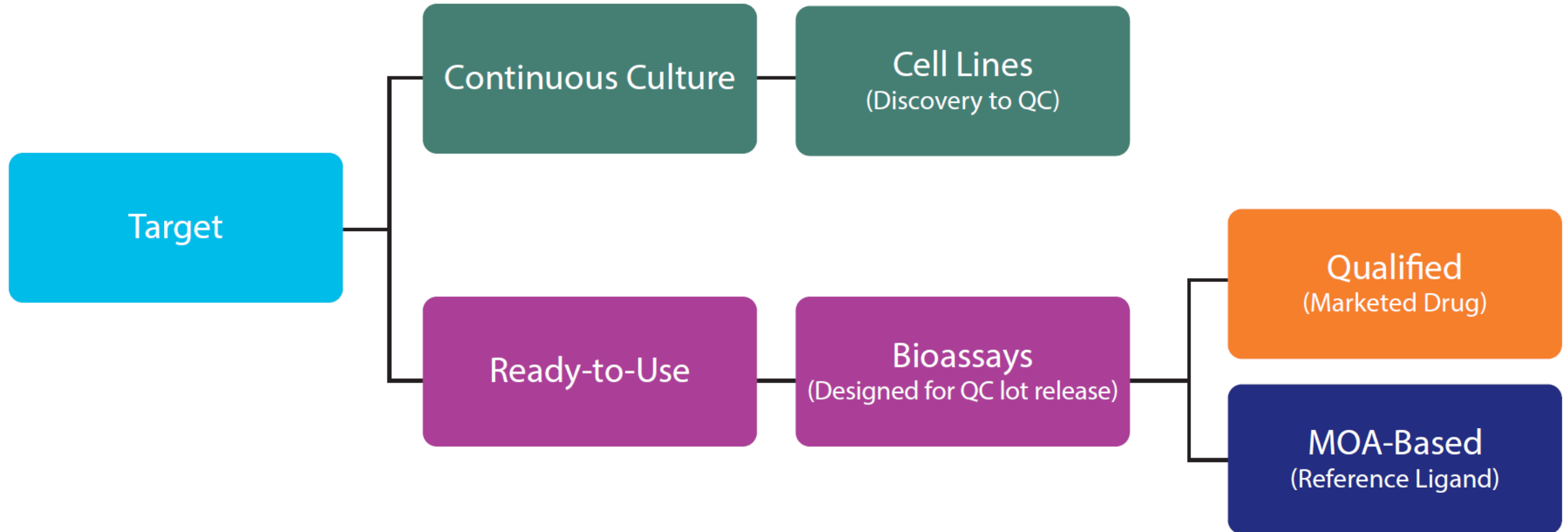


Enzyme Fragment Complementation: a Versatile and Robust Platform for Cell-Based Assays



Flexibility of Formats - Phase-Appropriate Solutions for Every Development Program

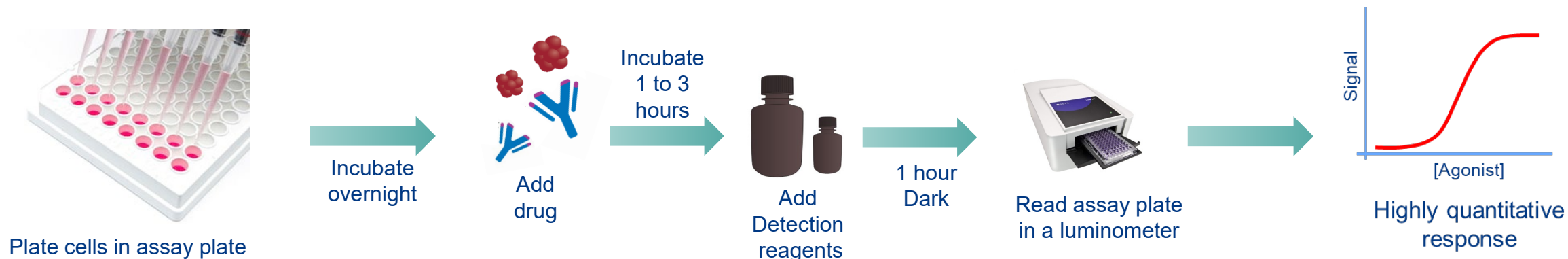
From Continuous Culture to Qualified Bioassays



Eurofins DiscoverX Qualified Bioassays are designed for QC Lot Release Programs

EFC Cell-Based Assays Protocol

Typical assay protocol for stable cell lines and assay ready kits



Simplified, universal assay format

- Easy-to-follow, detailed user manuals
- Rapid, single addition protocol
- Homogeneous, no wash format

Sensitive detection with large S:B ratio

Easily quantified luminescence read-out

- Dose-response curve
- Utilizes any standard plate luminometer

Everything is included!



Bioassay Kit includes all reagents needed for the assay. No need to worry about external sourcing of media reagents that may introduce variability

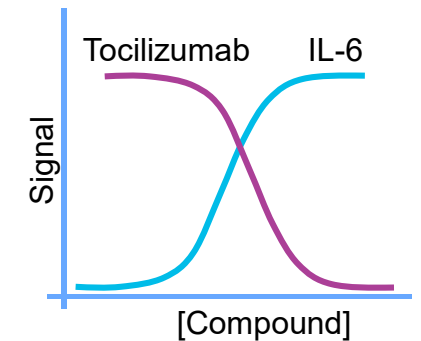
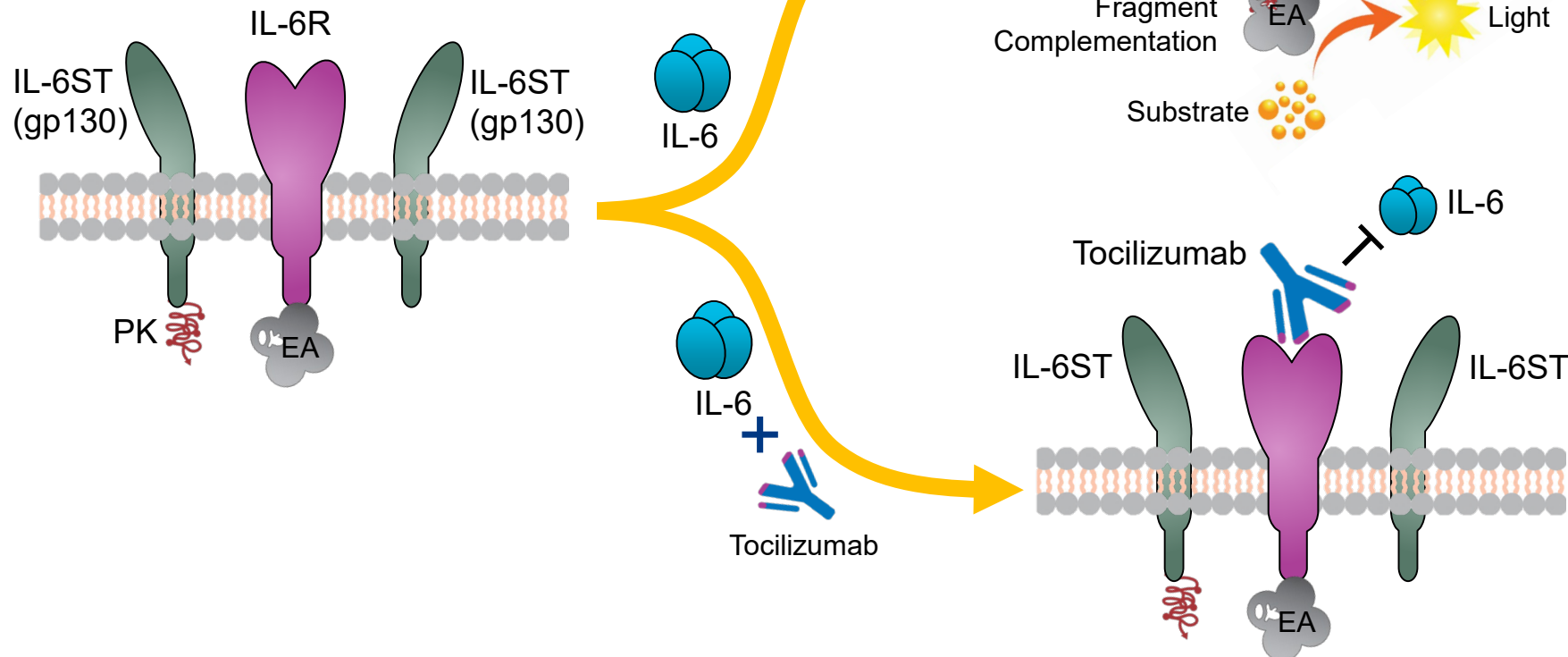
Case Study:

PathHunter[®] Tocilizumab Bioassay - Qualified with Actemra[®] for Therapeutics Targeting IL-6 Activity

PathHunter® Tocilizumab Bioassay - Qualified with Actemra®

MOA-reflective Assay Design: Receptor Dimerization

The assay is designed to detect IL-6 – induced interaction of the IL-6 receptor (IL-6R) with the IL-6 Signal Transducer protein (IL-6ST)



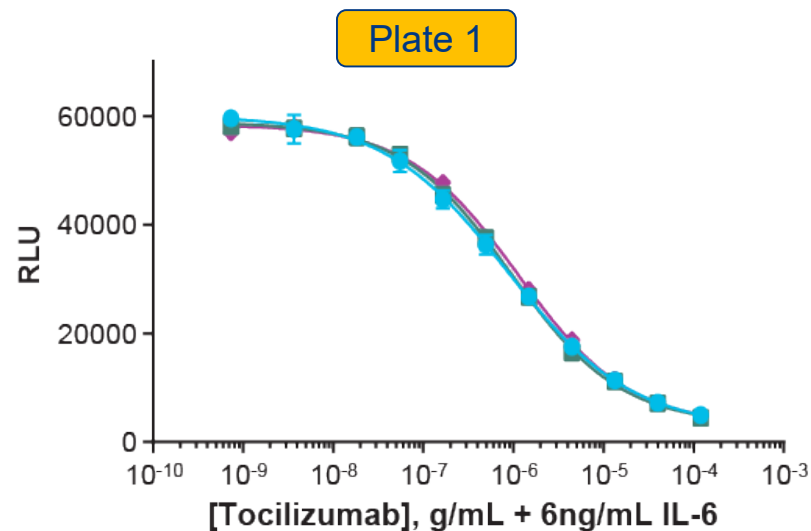
Actemra® is a registered trademark of Chugai Seiyaku Kabushiki Kaisha Corp., a member of the Roche Group

- Assay consistency (%CV) between eight 11-pt DRCs
- Plate uniformity: EC_{80} and IC_{80} (of IL-6 and Tocilizumab) across entire plate
- Plate-to-Plate variability: 3 plates with 11-pt DRCs run on 3 days
- Slope consistency
- Relative potency analysis: Accuracy, precision and linearity of the assay over a range of 50-150% performed by two operators
- Parallel line analysis using PLA3.0 or SoftMax Pro 7.1

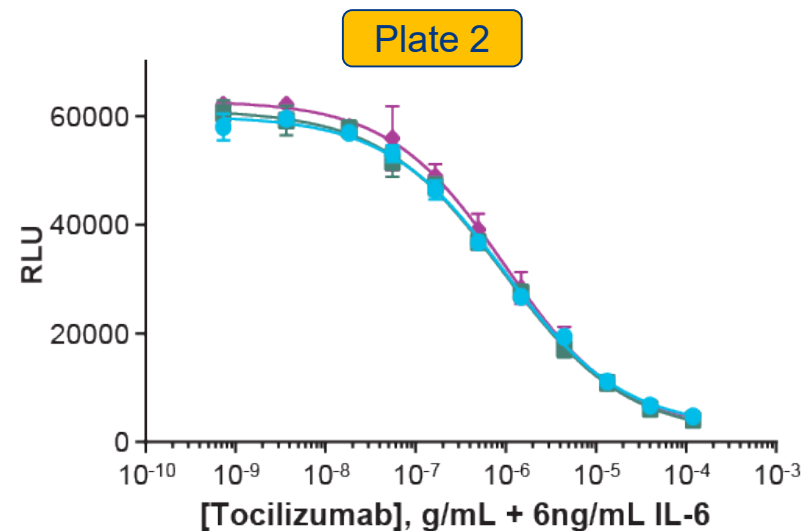


Assessing Intra-Day Plate-to-Plate Variability

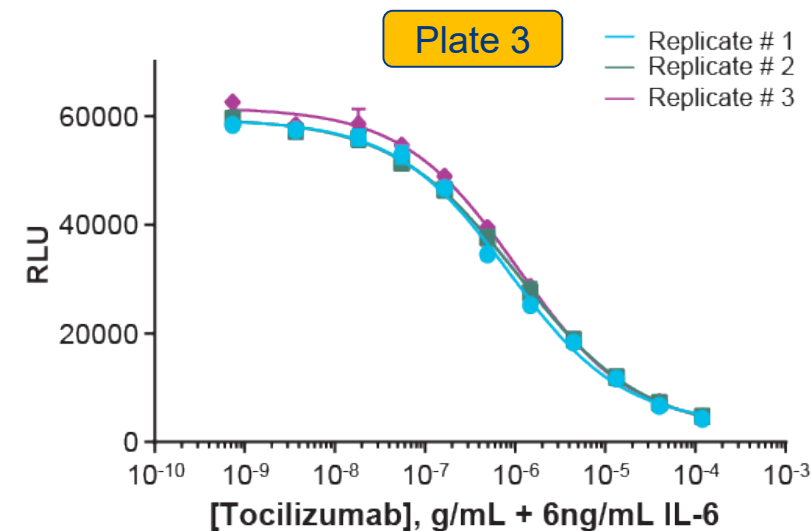
Experiments depicting 3 assay plates with full-plate DRC - Same Day – One operator



Parameter	R1	R2	R3
S/B	12.1	13.7	12.6
IC ₅₀ (ng/mL)	868	935	1,131



Parameter	R1	R2	R3
S/B	13.2	15.2	14.1
IC ₅₀ (ng/mL)	966	943	972



Parameter	R1	R2	R3
S/B	14.5	12.8	13.8
IC ₅₀ (ng/mL)	850	1077	1063

Repeatability and Intermediate Precision (Inter-Plate)

Plate	Sample	S/B	% RSD, S/B	IC ₅₀ , ng/mL	Mean IC ₅₀ , ng/mL	% RSD, IC ₅₀
1	R1	12.1	6.4	868	978	13.9
	R2	13.7		935		
	R3	12.6		1130		
2	R1	13.2	7.1	966	960	1.59
	R2	15.2		943		
	R3	14.1		972		
3	R1	14.5	6.2	850	997	12.8
	R2	12.8		1080		
	R3	13.8		1060		

Intermediate Precision (Inter-Day)

Day	IC ₅₀ , ng/mL	Mean IC ₅₀ , ng/mL	%RSD, IC ₅₀
1	1250	1330	24.6
2	1690		
3	1050		

Intermediate Precision (Inter-Plate): 9.7%

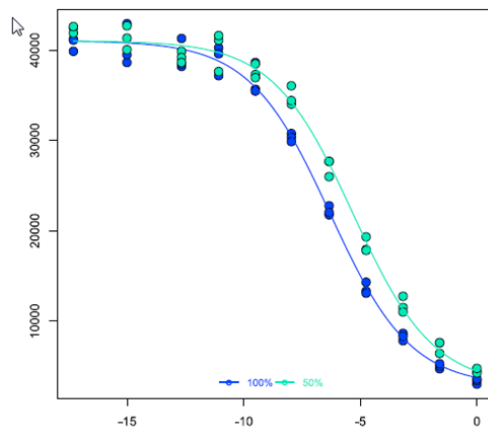
Intermediate Precision (Inter-Day): 24.6%

Parallelism and Relative Potency Estimation

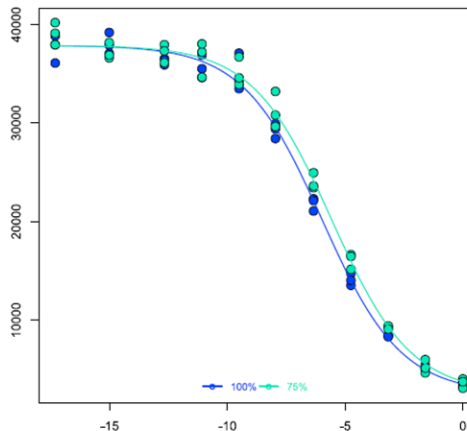
Relative Potency: Parallelism and Potency Estimation (PLA)

Analyst 1

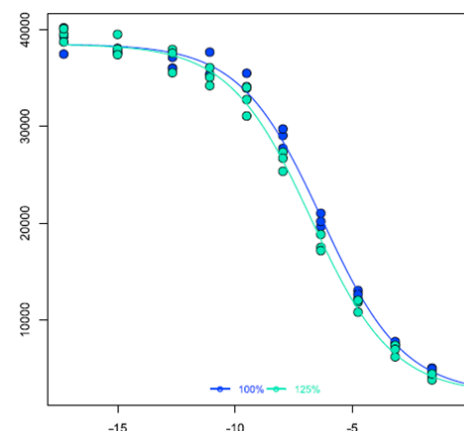
50%



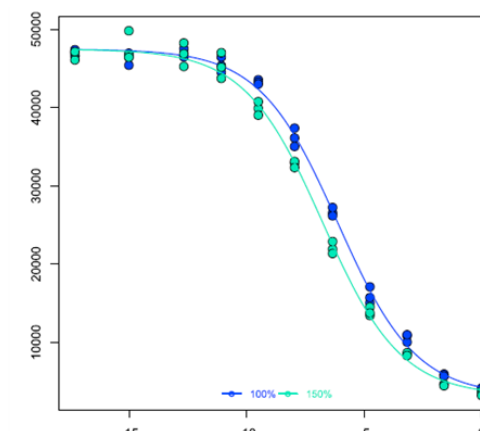
75%



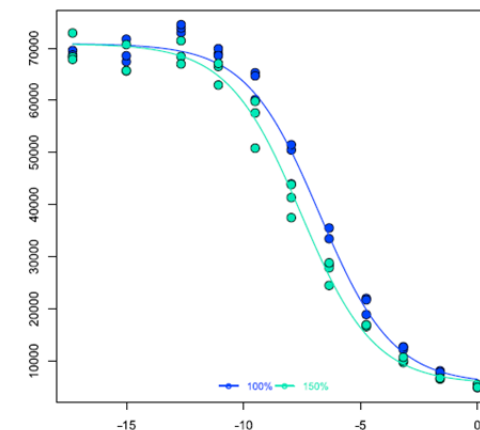
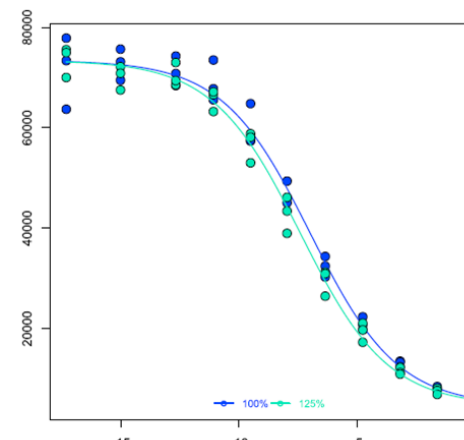
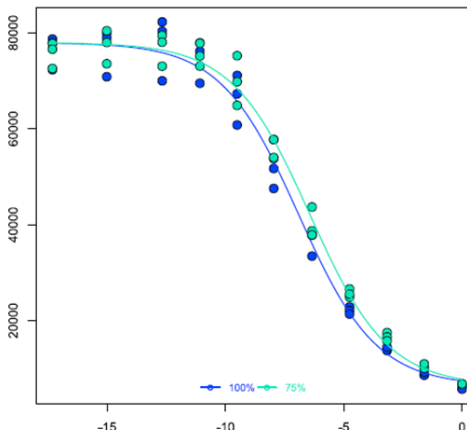
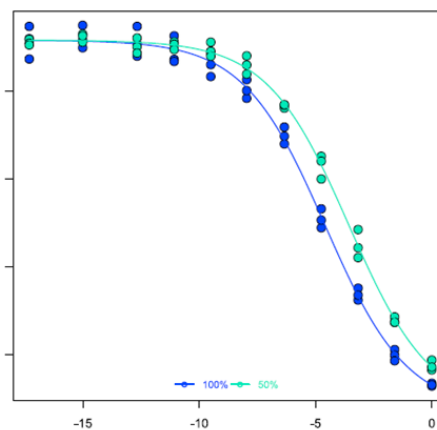
125%



150%



Analyst 2



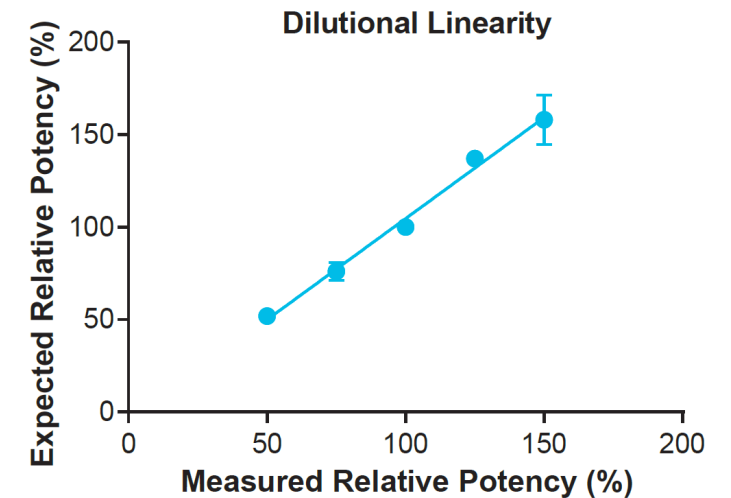
● Reference
● Sample

Note: RP values calculated in PLA 3.0

Tocilizumab Bioassay: Qualification with Actemra®

Summary: Accuracy, Precision and Dilutional Linearity (Single Analyst)

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			



Dilutional Linearity: 99.27%

Accuracy: 105.1%

Precision: 5.4%

Tocilizumab Bioassay is highly accurate, precise and linear!

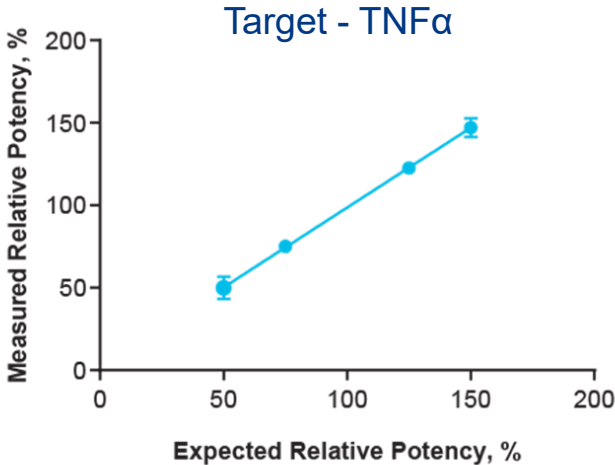
Qualified Bioassays for Precise Potency Determination for Cytokines and Interleukins

ICH Based Assay Qualification

- ✓ Accuracy
- ✓ Precision
- ✓ Range
- ✓ Dilutional Linearity
- ✓ Parallelism
- ✓ Specificity

HUMIRA
adalimumab

Adalimumab Bioassay Qualification



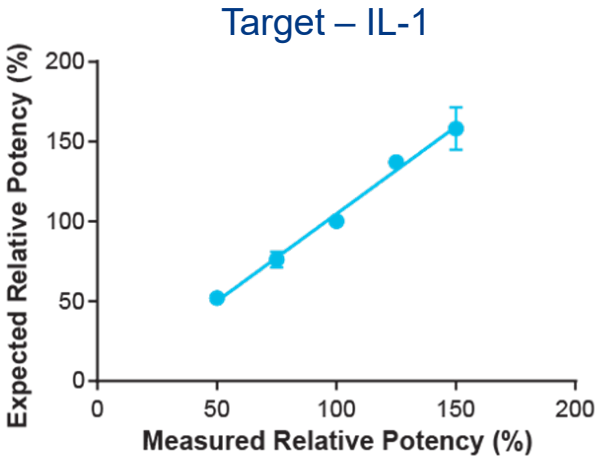
Dilutional Linearity: 99.99%

Accuracy: 105.1%

Precision: 5.65%

Kineret
(anakinra)

Anakinra Bioassay Qualification



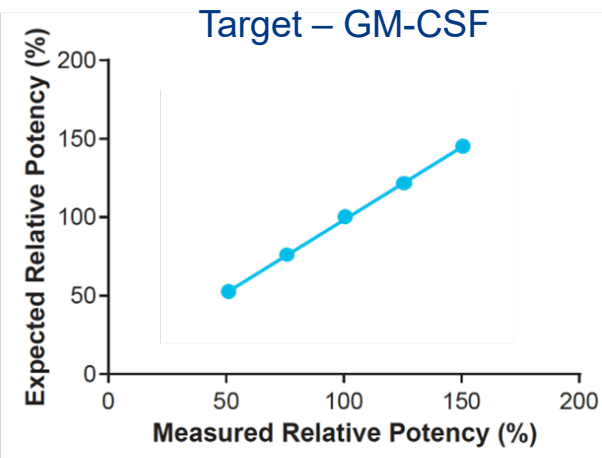
Dilutional Linearity: 96.7%

Accuracy: 99.4%

Precision: 11.4%

Leukine
sargramostim
A Recombinant GM-CSF–Yeast-Expressed

Sargramostim Bioassay Qualification



Dilutional Linearity: 99.56%

Accuracy: 99.1%

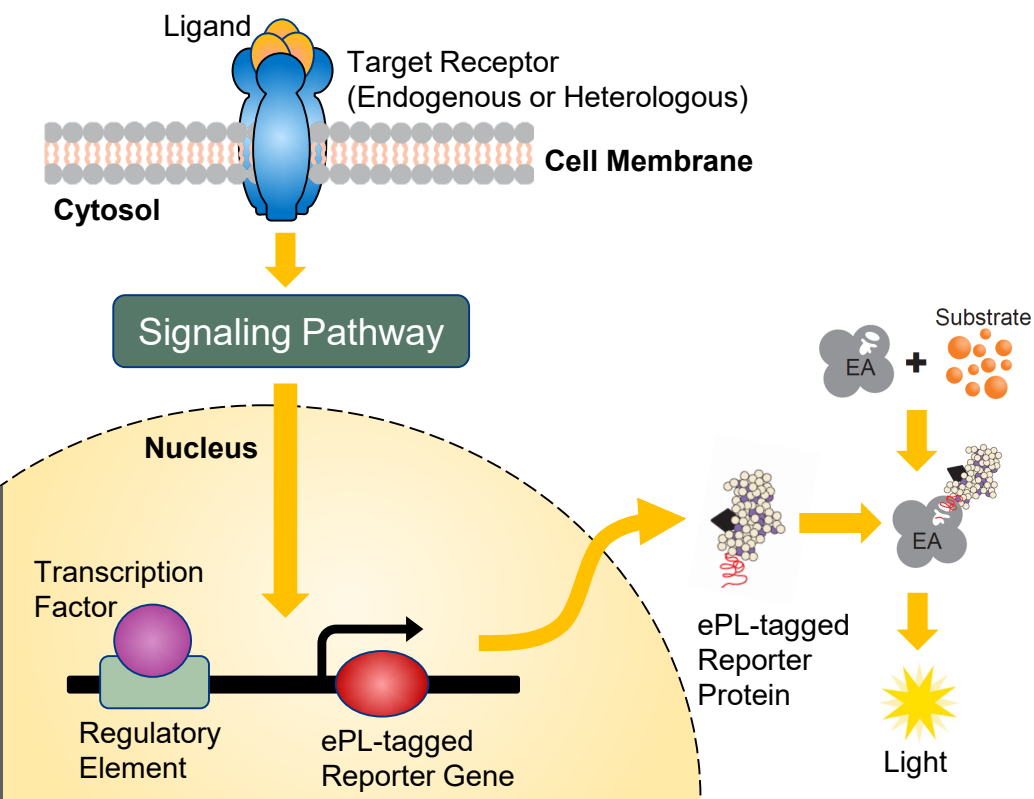
Precision: 3%

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

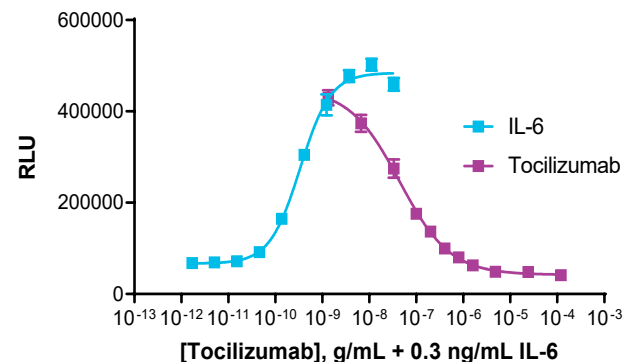
Signaling Pathway Reporter Assays for Cytokines and Interleukins

Assay Design : Signaling Reporter

The assay is designed for quantifying the activation or inhibition of signaling pathways to provide a downstream read-out that is complementary to upstream receptor-based assays to gain a comprehensive understanding of drug molecule's MOA

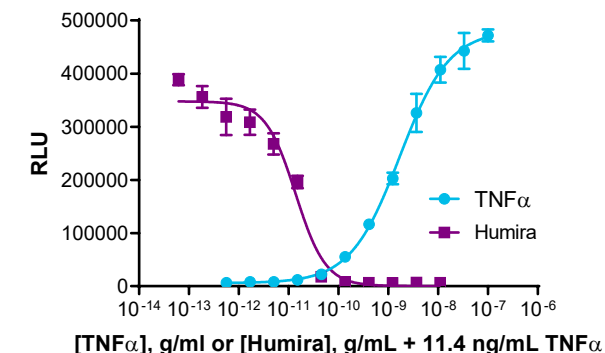


PathHunter® HepG2 STAT3 Pathway Reporter Assay



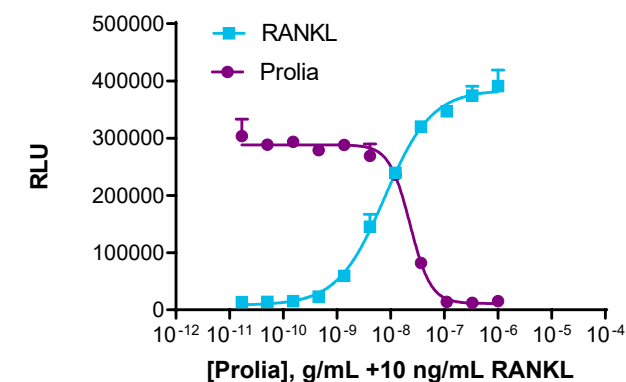
	EC ₅₀ /IC ₅₀	S/B
IL-6	0.33 ng/mL	6.2
Tocilizumab	44 ng/mL	11.6

PathHunter U2OS NF-κB Pathway Reporter Assay



	EC ₅₀ /IC ₅₀	S/B
TNFα	1.7 ng/mL	72.3
Humira	13 pg/mL	64

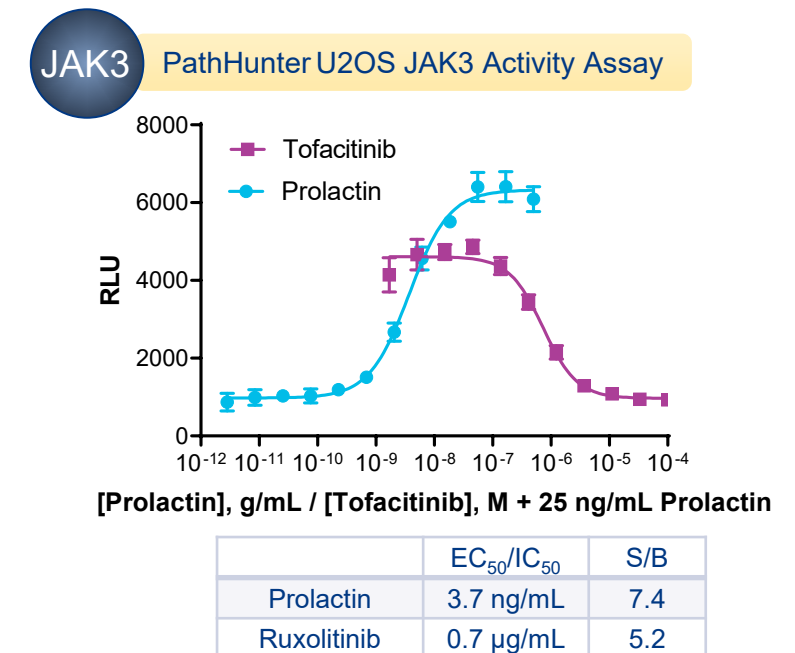
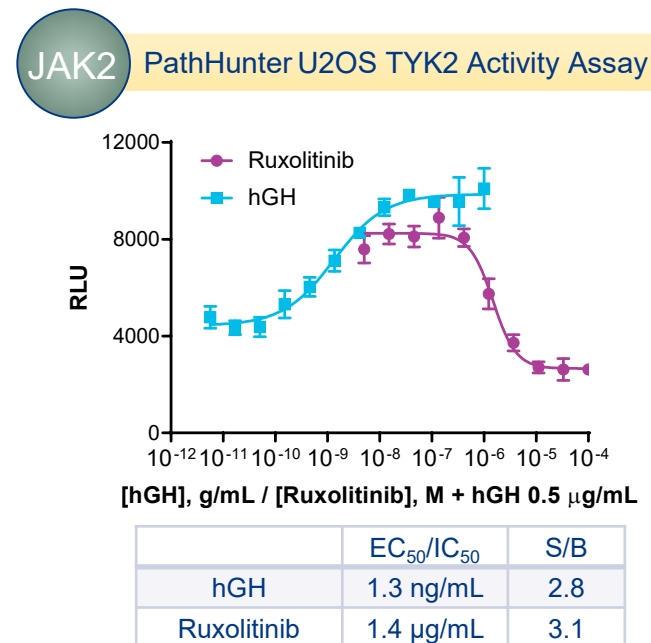
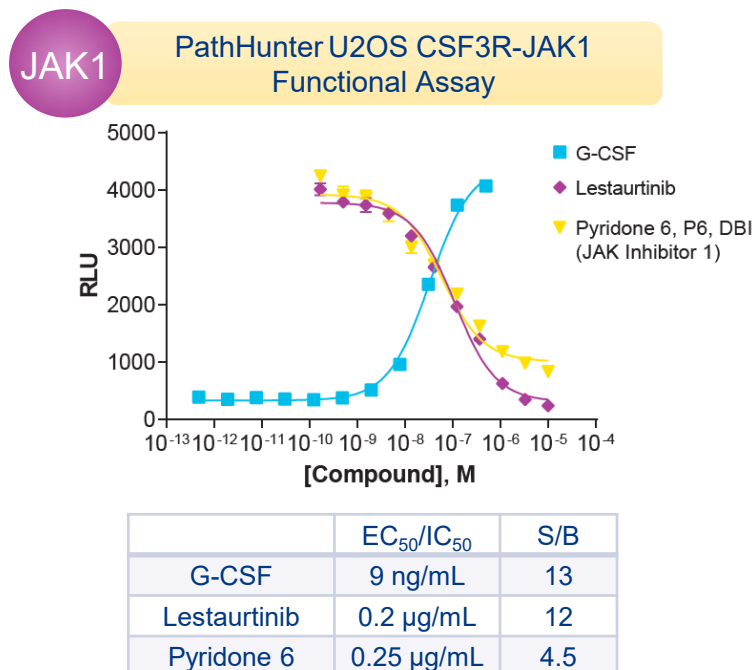
PathHunter U2OS RANK NF-κB Pathway Reporter Assay



	EC ₅₀ /IC ₅₀	S/B
RANKL	7.7 ng/mL	27
Prolia	23 ng/mL	21

PathHunter® Cell-Based Functional Assays for Small Molecule JAK Inhibitors

- There are several active, randomized, controlled trials evaluating the therapeutic potential of JAK inhibitors for treatment of COVID-19 (e.g. Ruxolitinib, Baricitinib)
- JAK inhibitors can have a distinct advantage over other immunomodulatory strategies in COVID-19, as they can exert a dual effects: anti-inflammatory (blockade of multiple, pro-inflammatory cytokines at the same time) and anti-viral effects (inhibiting cellular viral endocytosis)

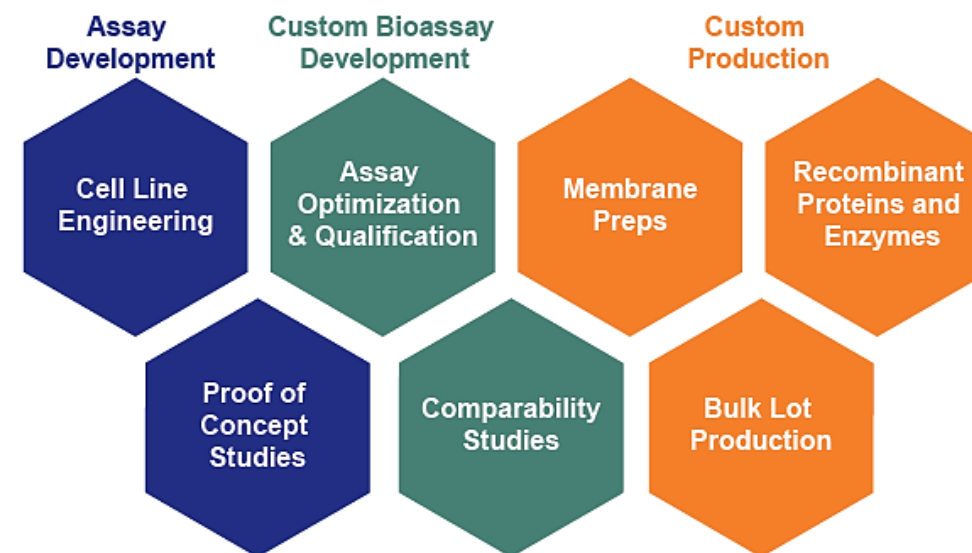




Your Target Biology, Our Expertise — Building a Better Assay Together

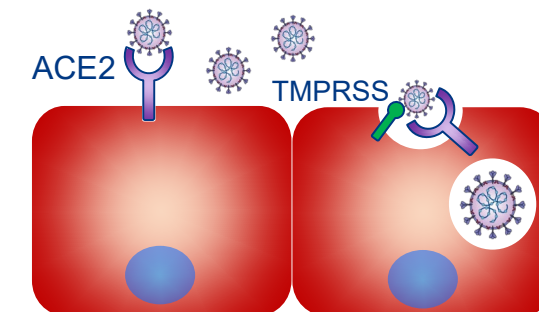
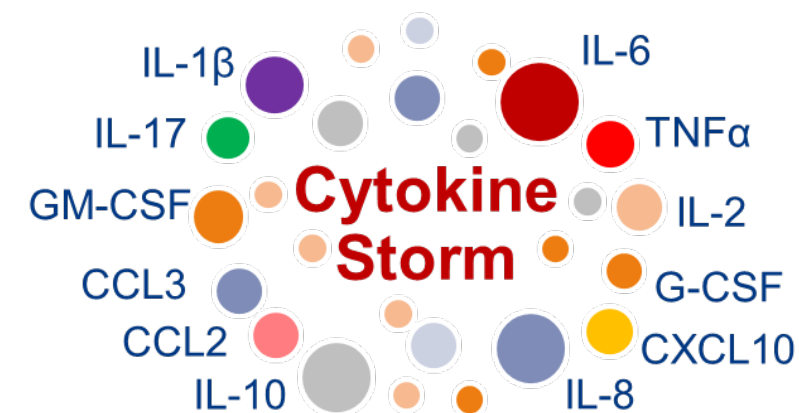
- **Development Expertise** — Decades of cell-based assay development, cell line engineering, and recombinant enzyme development expertise
- **Cell Line Engineering Capability** — Exogenous expression approaches (constitutive vs inducible) or gene editing (e.g. KO/KI with CRISPR/Cas9)
- **Collaborative** — Consultative assay development with regular updates through a dedicated project manager
- **Complete Solution** — Customized assay development with screening and profiling services within the same team

CAD Services Capabilities



Accelerate your drug discovery program with qualified, ready-to-use cell-based interleukin and cytokine assays for Covid-19 research

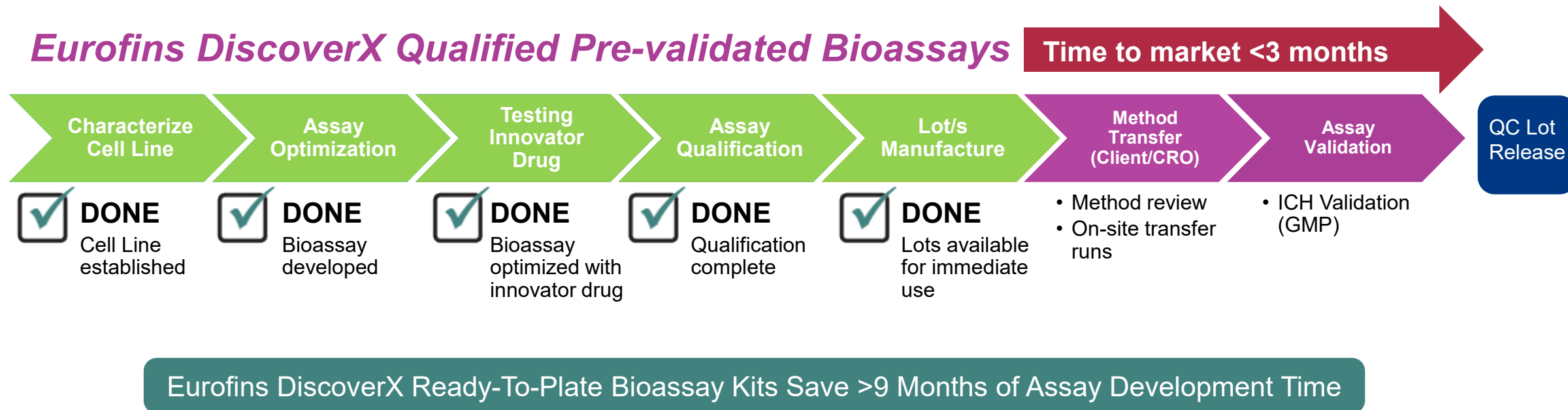
- **Target-Specific** – MOA-reflective, functional assays for drugs targeting pro-inflammatory cytokines, such as IL-1, IL-6, IL-7, GM-CSF and TNF α , all of which are implicated in COVID-19 pathology
- **Qualified Assays** – Robust, reproducible off-the-shelf assays qualified with innovator drugs
- **Complete Solutions** – Custom Assay Development enabling overexpression of other COVID-19-relevant receptors such as ACE2, TMPRSS2, etc.



Visit discoverx.com/covid-19 to learn more

Accelerating Implementation Phase for QC Lot Release with Qualified Bioassays

Typical Bioassay Development Timeline



Eurofins DiscoverX is a Global Leader in Supporting Early to Late Stage Drug Discovery

From Discovery to Development to Clinic to Post-Market

20+ years of enabling drug discovery and development programs

10+

Druggable
target classes

1500+

Stable cell line and
membrane preps

20+

core patents

Validated

> 30 Billion data points screened in
assay services with same assays

2000+

Publications across
multiple applications

55+

Qualified and MOA
Based Bioassays

**ICH Based Bioassay
Qualification**

Facilitate downstream validation studies

**Certified CRO
Partners**

Scientific training to enable global
CROs

**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