

Accelerating Potency Bioassay Implementation from Development to QC Lot-Release & Ensuring Long-term Assay Reproducibility

Alpana Prasad, Ph.D.

Director, Innovations and Strategic Partnerships

Well Characterized Biologics & Biological Assays November 1, 2022

OUR EXPERTISE IN YOUR HANDS, DISCOVER CONFIDENTLY.

Eurofins DiscoverX[®] Is a Global Leader in Cell-Based Assays



For Screening, Profiling, Characterization, Potency, & NAb Applications

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 pate	ents Pub	000+ lications across multiple ications
			55+ Qualified & MOA-based bioassays	Validated >30 Billion Data Points Screened in assay services with same assays		4 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		70+ Global Programs For potency, stability and NAb testing		

Bioassay Implementation Workflow

A Holistic Approach Supporting Programs from Development to Post-Market

🔅 eurofins

DiscoverX

The Eurofins Discovery PRODUCTS COMPANY



Enzyme Fragment Complementation (EFC) Robust Platform for Interrogating Different MOAs

Simple, add-and-read assay format Homogenous · Easily quantified luminescence read-out Inactive Active β-Gal **Enzyme Fragments** Enzyme Sensitive · Enzymatically-amplified assay Substrate Detection High precision & accuracy Hydrolysis · Large dynamic range Light Robust Enzyme Enzyme · High assay reproducibility Signal Donor (ED) Acceptor (EA) Easy to Similar workflow across platform Transfer · Compatible with most readers **Protein/Protein Interactions Quantify Cytotoxicity** Internalization & Trafficking **Compound Target Engagement Receptor Dimerization Downstream Response Evaluation Quantification & Degradation** ED-tagged Reporter ePL-tagged Housekeeping Reporter Gene Protein-ED 🔨 ED

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

Constructions DiscoverX

Examples of Target Biology Translated into Assay Design

eurofins DiscoverX

The Eurofins Discovery PRODUCTS COMPANY



Ready-to-Use Qualified Bioassay Method From Assay Development to Transfer

The Eurofins Discovery PRODUCTS COMPANY **Qualified Bioassays MOA-based Bioassays**

Assay Selection & Testing Innovator Drug Assay Optimization Assay Qualification Cell Line Characterization Choose assay format Optimize assay Refine conditions with Establish robustness, that best reflects MOA of parameters (thaw-&-use originator drug or clinical precision, accuracy & linearity of assay cells only) molecule drug Optimize with marketed drug Select relevant MOA Culture & freeze conditions Robustness: 11-point DRC or reference standard • Plate uniformity: EC₈₀/IC₈₀ Establish stable cell line Cell density Dilution series across entire plate Retrovirally engineered Assay reagents & buffers Establish suitability of assay Reproducibility and · Ligand incubation time Cell backgrounds to determine relative potency intermediate precision: 3 Cell plating time Functional response plates with 11-point DRC Determine range and Assay windows Assay plate type linearity of the assay over over 3 days Passage stability Incubation time with 50%-150% with single Second analyst: Determine detection reagents Sterility analyst (at least 3 range and linearity over Assess sensitivity to matrix independent runs) 50% -150% of assay (normal human serum) Confirm manufacturability Establish QC release **Bioassay kit &** Transfer Multiple lots criteria for production lots detailed protocol method Scale-up (~500/lot)

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

🔅 eurofins

DiscoverX



Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

All registered trademarks are the property of their respective owners

Assays Implemented in Global Biologics Programs

(Characterization, Stability Studies, NAb Detection & Potency Testing in Lot Release)

The Eurofins Discovery PRODUCTS COMPANY

🛟 eurofins

Anti-SIRPα	Anti-PD1	BTLA	Anti-VEGF	IGF1R	GLP1R	CNR1	BRDKBR2	INSRb
Multiple programs in NA & EU for originator biologics	Multiple programs in NA & APAC for originator & biosimilars	US-based program for originator biologic	Multiple programs in NA, EU & APAC for biosimilars	Multiple programs in NA and APAC for originator biologic	Multiple programs in NA, EU & APAC for originator & biosimilars	US-based program for originator biologic	US-based CRO	Multiple programs in NA, EU & APAC for biosimilars
Potency & NAb	Potency	Not disclosed	Potency	Potency	Potency	Potency	Potency	Potency
IL-17R	IL-2R	IL-31R	11 C	*	5. 	MCxR	FSHR	PTHR1
EU biopharma for biosimilar	Multiple programs in NA and EU for originator biologic	US-based pharma for originator biologic		And and a		Multiple programs in NA & EU for originator biologic	APCA-based program for biosimilar	Multiple programs in NA, EU & APAC for biosimilars
Not disclosed	Not disclosed	Potency & NAb		15		Not disclosed	Potency	Potency
IL-10R	IL-23R	IL-7R				CALCRL/RAMP3	CXCR4	CXCR2
Multiple programs in NA for originator biologic	US-based biopharma for originator biologic	EU-based program for originator biologic		- North America: El L = Euron		NA-based pharma for originator biologic	EU-based pharma for originator biologic	US-based program for originator biologic
Potency	Potency	Potency		– North America, EO – Europ	e, AFAC - Asia Facilic	Potency	NAb	Not disclosed
TSLPR	CSF2R	CSF1R	RANK	FGFR	ErbB2/ErbB3	GM-CSF	AXL	CD16 Effectors
EU-based pharma for originator biologic	Multiple programs in NA & EU	Multiple programs in NA & EU	Multiple programs in NA & APAC for biosimilars	US-based biopharma for originator biologic	Multiple programs in NA, EU & APAC for biosimilars	Multiple programs in NA & EU for originator biologic	US-based program for originator biologic	Multiple programs in NA & EU for ADCC
Potency	Potency	Potency	Potency	NAb	Potency	Potency	Not disclosed	Potency

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

Qualified, off-the-shelf bioassays

DiscoverX



Case Study: KILR® Raji Bioassay for Therapeutics Driving ADCC Mechanism

Challenges with Existing Cytotoxicity Assays Addressed by KILR[®] Platform

🛟 eurofins |

DiscoverX

The Eurofins Discovery PRODUCTS COMPANY

What assay development scientists face	Radioactivity or fluorescent dye-based assays are inherently leaky, less robust & produce high background	KILR assays are non-radioactive and not leaky	Little to no background	High assay background
	Low assay throughput restricts screening and characterization efficiency	KILR assays are high- throughput compatible (384-well)	Ease of implementation from screening to lot release	[Antibody]
JLJL	Surrogate assays fail to reflect the MOA. These are <i>"Predictive" and not "Reflective"</i>	KILR assays are end- point assays and specifically measure target cell killing	Physiologically- reflective of <u>TRUE MOA</u>	KILL Effector Cell ADCC
	High donor-to-donor variability associated with effector cells	KILR CD16 Effector Cells are single donor- derived	<u>Eliminate</u> donor variability	(S 100 + Uesp () Appropriate S 50 - 2 - 1 - 0 - 1 - 2 - 3 Log (Anti-CD20 Ab) (ng/mL)

KILR[®] Cytotoxicity Assay for ADCC Overview

🛟 eurofins 🛛

The Eurofins Discovery PRODUCTS COMPANY



KILR[®] Assay ADCC Demonstrated Using a Variety of Antibodies, Antigens, and Target Cell Types



Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

eurofins DiscoverX

The Eurofins Discovery PRODUCTS COMPANY

Eliminating Donor Variability in ADCC Assays Using KILR[®] CD16 Effector Cells

🔅 eurofins

The Eurofins Discovery PRODUCTS COMPANY



Continuous Culture vs Ready-to-Assay Cell Formats KILR[®] Raji ADCC Bioassay Kit Performance

Assay Ready

Continuous Culture

Condition S/B EC₅₀, ng/mL **Continuous Culture** 2.6 19

6.7

2.6

14

Assay Ready

formats

Raw Luminescence (RLU) 150000-150000-100000 100000 RLU RLU 50000-50000-10-12 10-11 10-10 10-9 10-8 10-7 10-6 10-5 10-12 10-11 10-10 10-9 10-8 10-7 10-6 10-5 [Anti-hCD38 antibody], g/mL [Anti-hCD38 antibody], g/mL %Killing (%ADCC activity) 60-60-40 40-% ADCC % ADCC 20 20-10⁻¹¹ 10⁻¹⁰ 10⁻⁹ 10⁻⁸ 10⁻⁷ 10⁻⁶ 10⁻⁵ 10-11 10-10 10-9 10-8 10-7 10-6 10-5 [Anti-hCD38 antibody], g/mL [Anti-hCD38 antibody], g/mL -20--20-

Condition	E _{Max}	EC _{₅0} , ng/mL
Continuous Culture	45%	2.6
Assay Ready	51%	2.6

Excellent Concordance between EC₅₀ of the two

Comparable % Killing (% ADCC) with KILR Raji Bioassay Cells Relative to Continuous Culture

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

eurofins



Suitable for Screening and Rank Ordering of Antibodies KILR[®] Raji ADCC Bioassay Kit



Isotype (IgG1) Human IgG1 isotype control -- -- -- --

KILR Raji Bioassay allows discrimination of differences in Fab and effector regions of related antibodies

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands – Discover Confidently

Effectors used: KILR CD16 Effector Cells

eurofins

DiscoverX

The Eurofins Discovery PRODUCTS COMPANY

15

Repeatability (over 3 Days); Single Analyst

Raw Luminescence (RLU) 120000-200000-200000 150000 150000-80000 RLU RLU RLU 100000 100000 40000 50000 50000 0-0-0+ 10-12 10-10 10-14 10⁻⁸ **10**-16 10⁻¹⁶ 10-14 10-12 10-10 1(10-16 10-14 10-12 10-10 10-8 [Rituximab], g/mL {Rituximab], g/mL [Rituximab], g/mL % Killing (% ADCC activity) 150 150 100-100-⁻ % ADCC 75-2 100 9 ADCC 8 50 % ADCC 50-50 50· 25 0-0-0-10-12 10-12 10-10 10-12 10-10 10-16 10-14 10⁻¹⁰ 10-8 10-16 10-14 10-8 10-14 10-16 10 {Rituximab] g/mL {Rituximab], g/mL {Rituximab], g/mL

Day 2

Day 3

Parameter Day 1 Day 2 Day 3 % RSD EC_{50} (pg/mL) 7.03 8.9% 6.2 7.4 S/B 9.8 8.3 12.2% 10.6 E_{Max} (%) 135% 114% 75.8% 27.7%

Excellent inter-day repeatability

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

Effectors used: KILR CD16 Effector Cells

KILR® Raji ADCC Bioassay Kit

Day 1

🔅 eurofins DiscoverX

The Eurofins Discovery PRODUCTS COMPANY

Consistent Performance Between Analysts (Relative Potency) & eurofins KILR[®] Raji ADCC Bioassay Kit The Eurofins Discovery PRODUCTS COMPANY



Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

Effectors used: KILR CD16 Effector Cells

17

DiscoverX

Phase-Appropriate Method Qualification Suitability of KILR[®] Raji Bioassay Cells for Relative Potency Assays

Nominal RP, %	Analyst	Observed RP, %	Average RP, %	% RSD	Average % Recovery
	1	156			
	1	179.4	166.6	7.0	111.1%
150	1	177.4			
	2	166			
	2	154.3			
	1	130.6			
105	1	131.2	407.0	0.0	101.8%
125	1	126	127.3	3.6	
	2	121.4			
	1	82			
	1	116	106.5	13.0	106.5%
100	1	112.7			
100	1	98.4			
	1	117.8			
	2	112.2			
	1	70.4			
75	1	76.6	73.9	4.9	98.5%
	1	77.3			
	2	71.1			
	1	47.8			
	1	45.8		13.2	96.9%
50	1	40.9	48.5		
	2	58.4			
	2	49.4			



🔅 eurofins

Parameter	Value	Specification
Accuracy (Average % Recovery)	103%	100% ± 20%
Repeatability	14.2%	≤20%
Intermediate Precision	≤13.2%	≤20%
Linearity (R ²)	0.9926	≥0.95

Assay demonstrates very good accuracy, repeatability, intermediate precision, and dilutional linearity

Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands - Discover Confidently

Effectors used: KILR CD16 Effector Cells

The Eurofins Discovery PRODUCTS COMPANY

DiscoverX

18

Ready-to-Use Format 🛟 eurofins Simple and Fast KILR® Raji ADCC Bioassay Kit Protocol DiscoverX The Eurofins Discovery PRODUCTS COMPANY <8 hours Thaw Wash Plate Target Count Cells Add Detection Add Effector Reagent Antibody Cells* Cryopreserved 1 hour @RT KILR Target Cells **Detect Target Opsonize Target** Cell Death Cells with Antibody 4 hours @37°C Bioassay Kit Includes Required Reagents for the Assay Pati-surface at-content 1 Mar CofA a eurofins **Cell Plating** Certificate Cryopreserved Detection **Assay Plates User Manual** of Analysis Reagent Cells Reagent *Effector cells sold separately

Same KILR® Target Cells Enable Assessment of Multiple MOAs

🛟 eurofins |

The Eurofins Discovery PRODUCTS COMPANY

ody-Dependent			
lar Phagocytosis	Complement-Dependent Cytotoxicity	T-cell Redirection (with T-cell Engagers)	Chimeric antigen receptor T-cells
pendent phagocytosis sosomal degradation	Activated complement system leads to formation of membrane attack complex (MAC)	Bi-specific antibody engages T-cell with cancer cells for killing	Engineered T-cells recognize and kill cancer cells
Macrophage Macrophage	Complement C1q Target Antigen MAC KILR Target Cell	Activated T-Cell CD3 BITE CD19 CD19 CD19 KILR Target Cell	Activated T-Cell
	ar Phagocytosis spendent phagocytosis vosomal degradation	Cytotoxicity Activated complement system leads to formation of membrane attack complex (MAC)	Outy-Dependent lar Phagocytosis Outprement-obpendent Cytotoxicity I-cell Redirection (with T-cell Engagers) pendent phagocytosis rsosomal degradation Activated complement system leads to formation of membrane attack complex (MAC) Bi-specific antibody engages T-cell with cancer cells for killing Macrophage Image: Complement Cruption of membrane attack complex (MAC) Image: Complement Cruption of membrane attack complex (MAC) Image: Complement Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC) Image: Cruption of membrane attack complex (MAC)

Largest Menu of Cell-based Assays for Potency & NAb Assay Development



>1500+ Cell Lines and Membrane Preps to support bioassay development for major drug target classes



Drug Discovery & Development Phase-Appropriate Solutions | Our Expertise in Your Hands – Discover Confidently

All registered trademarks are the property of their respective owners 21

Options to Accelerate Your Bioassay Development Plan

Content of the Eurofins Discovery PRODUCTS COMPANY

Catalog	 30 ready-to-use bioassays 1500+ cell lines
Customize	 Optimize & qualify with your clinical molecule New MOA/background
Create	Parental Cell LinesRetroparticles
Visit <i>discoverx.com</i> /	bioassays to learn more

Come visit us at our booth! We look forward to discussing your program requirements.