Construction Structure Construction Construc

Implementing MOA-Reflective Cytotoxicity Assays Using Readyto-Use KILR[®] Target & Effector Cells from Screening to Lot Release

Andrew Green Sr. Business Development Manager Eurofins DiscoverX

OUR EXPERTISE IN YOUR HANDS. DISCOVER CONFIDENTLY.

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



The Eurofins Discovery PRODUCTS COMPANY

From Discovery to Development to Clinic to Post-Market

20+ Years of Enabling Drug Discovery and Development Programs

Image: constraint of the second sec		10+ Druggable target classes	1500 Stable cell line a membrane prep	+ 20+ and os		ents	2000+ Publications across multiple applications	
			55+ Qualified & MOA-based bioassays	Validated >30 Billion Data Points screened in assay services with same assays			S 1	3 Certified CRO Partners Scientific training to enable global CROs
ICH-Based Bioassay Qualification Facilitate downstream validation studies	Dedicated Scientific Experienced to scientific supp	d Support eam providing ort	20+ Successful Assay Transfers At clients/affiliated CRO sites		ful fers RO sites	70+ (For poter testing	Glo ncy,	bal Programs stability and NAb



FDA approves 100th monoclonal

NEWS · 05 MAY 2021

DiscoverX

The Eurofins Discovery PRODUCTS COMPANY

Therapeutic mAbs are the fastest growing class of biological therapeutics, and over 100 mAbs are already approved by the FDA

Many of these therapeutic antibodies work by activating antibody-dependent cell-mediated cytotoxicity (ADCC) or other effector-mediated functions (ADCP, CDC, T-cell Redirection)

Starting from Phase I and even earlier, functional assays to assess effector function (e.g. ADCC) should be performed to evaluate therapeutic antibody potency and in other cases for safety



antibody product Thirty-five years on from the FDA's approval of a first monoclonal antibody, these biologics account for nearly a fifth of the agency's new drug approvals each year.

Therapeutic antibodies to induce ADCC approved in the US						
Antibody	Туре	Target				
Rituximab/other similar anti-CD20	Chimeric IgG1	CD20				
Trastuzumab	Humanized IgG1	HER2				
Alemtuzumab	Humanized IgG1	CD52				
Cetuximab	Chimeric IgG1	EGFR				
Ofatumumab	Human IgG1	CD20				
Pertuzumab	Humanized IgG1	HER2				
Dinutuximab	Chimeric IgG1	GD2				

Challenges with Existing Cytotoxicity Assays

🛟 eurofins

DiscoverX

The Eurofins Discovery PRODUCTS COMPANY



Enzyme Fragment Complementation (EFC) Technology

eurofins DiscoverX

The Eurofins Discovery PRODUCTS COMPANY



KILR[®] Cytotoxicity Assay Overview ADCC Example

🛟 eurofins

The Eurofins Discovery PRODUCTS COMPANY



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

🛟 eurofins

The Eurofins Discovery PRODUCTS COMPANY



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

🛟 eurofins

The Eurofins Discovery PRODUCTS COMPANY



Ready-to Use Format Easy-to-Use KILR[®] Raji ADCC Bioassay Kit Protocol



The Eurofins Discovery PRODUCTS COMPANY



*Effector cells sold separately

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



The Eurofins Discovery PRODUCTS COMPANY



Condition	S/B	EC ₅₀ , ng/mL
Continuous Culture	19	2.6
Assay-ready	6.7	2.6

Excellent Concordance between EC₅₀

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

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

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

Content of the Eurofins Discovery PRODUCTS COMPANY



	Sample	Description	S/B	HillSlope	E _{Max}	EC ₅₀ , pg/mL
	hCD20-mab13	Non-fucosylated anti-CD20 (Rituximab)	5.4	1.152	88%	92.8
	hCD20-ga-mab13	Non-fucosylated anti-CD20 (Obinituzimab)	5.2	1.307	84%	46.5
-	hCD20-mab12	Fc Null variant of Rituximab				
	hCD20-mab1	Wild-type Rituximab	4.5	1.270	69%	288
٠	lsotype (IgG1)	Human IgG1 isotype control				

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

Repeatability (Over 3 Days); Single Analyst KILR[®] Raji ADCC Bioassay Kit

eurofins

The Eurofins Discovery PRODUCTS COMPANY



Parameter	Day 1	Day 2	Day 3	% RSD
EC ₅₀ , pg/mL	6.2	7.03	7.4	8.9%
S/B	9.8	10.6	8.3	12.2%
E _{Max} , %	135%	114%	75.8%	27.7%

Excellent inter-day repeatabilty

Consistent Performance Between Analysts (Relative Potency) & eurofins KILR[®] Raji ADCC Bioassay Kit

The Eurofins Discovery PRODUCTS COMPANY

DiscoverX



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

		D	isc	ove	erX
The Eurofins	Discovery	PRODU	ICTS	сом	PANY

eurofins

Nominal RP, %	Analyst	Observed RP, %	Average RP, %	% RSD	Average % Recovery	
	1	156				
	1	179.4				
150	1	177.4	166.6	7.0	111.1%	
	2	166				
	2	154.3				
	1	130.6				
105	1	131.2	107.2	26	101.8%	
120	1	126	127.3	3.0		
	2	121.4				
	1	82		13.0	106.5%	
	1	116				
100	1	112.7	106 5			
100	1	98.4	100.5			
	1	117.8				
	2	112.2				
	1	70.4		4.9		
75	1	76.6	73.9		98.5%	
10	1	77.3	10.0			
	2	71.1				
50	1	47.8				
	1	45.8			96.9%	
	1	40.9	48.5	13.2		
	2	58.4				
	2	49.4				



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

Multiple MOAs Supported by KILR® Cytotoxicity Assays



The Eurofins Discovery PRODUCTS COMPANY



KILR ADCP Assay Concept and Workflow

Content of the Eurofins Discovery PRODUCTS COMPANY







Sample	S/B	E _{Max}	EC ₅₀ , ng/mL
Rituximab	6.2	83%	2.44

ADCP Observed for Diverse Antigens in Solid and Heme Cancer Models

60-

HER2+ Breast Cancer

Trastuzumab

100·

-



100-

eurofins

Cetuximab Human IgG1 80-80 Human IgG1 40-60-% ADCP % ADCP 60. %ADCP 40-20-**40** 20-20 0-0 -20+ 10-11 10-10 10-9 10-8 10-7 10-6 10-5 10-4 10-11 10-10 10-9 10-8 10-7 10-6 10-5 **10**-11 **10**-10 10-9 10-8 10-7 -20-[Antibody], g/mL [Antibody], g/mL [Antibody], g/mL CD38+ B Cell Model HER2+ Gastric Cancer HER2- Breast Cancer 80 - Antibody X 100· Trastuzumab 100--Human IgG1 Human IgG1 60 80. 75-% ADCP 60-**40** % ADCP % ADCP 50-40-**20**· 25-20-0 **10**-11 **10**⁻¹⁰ 10⁻⁹ **10**-8 **10**-7 10⁻¹⁰ 10⁻⁹ 10⁻⁵ 10⁻⁷ 10⁻⁶ 10⁻⁵ 10⁻⁴ 10-11 10-10 10 10-8 10-7 10-6 10-5 -20--20--25-[Daratumumab], g/mL [Antibody], g/mL [Antibody], g/mL

HER2- Breast Cancer

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

18

10-6

10-6

Higher Percentage of Phagocytosis Seen with Plate-Based, KILR[®] Assay Compared to Flow-Based Assay Format

eurofins DiscoverX

The Eurofins Discovery PRODUCTS COMPANY



Evaluate ADCP Activity of Therapeutics with IgG1 and IgG4(S228P) Fc Regions



The Eurofins Discovery PRODUCTS COMPANY



Results indicate developers using an IgG4(S228P) Fc should evaluate ADCP activity of their therapeutics, preferably in multiple donors

mediated cytotoxicity applications such as:

🔅 eurofins

DiscoverX

The Eurofins Discovery **PRODUCTS COMPANY**





Visit discoverx.com/kilr-bioassays to learn more

ADCC, ADCP, CDC, T-cell Redirection and CAR-T KILR ADCC Bioassays allows discrimination of differences in Fab and Fc regions to enable rank ordering of antibodies

KILR[®] cytotoxicity platform offers MOA-based assays for variety of cell-

KILR ADCC Bioassays, when used in combination with KILR CD16 Effector Cells, produce highly accurate and precise data suitable for use in relative potency assays

KILR ADCP Bioassays provide a robust, MOA-reflective readout for ADCP in a simple, plate-based assay format that minimizes macrophage handling

KILR ADCP Bioassays have utility for evaluating ADCP activity of therapeutics with IgG1 and IgG4(S228P) Fc regions