

***ADCC Reporter Bioassay:  
A Novel, Bioluminescent Cell-Based Assay for Quantifying  
Fc Effector Function of Antibodies***

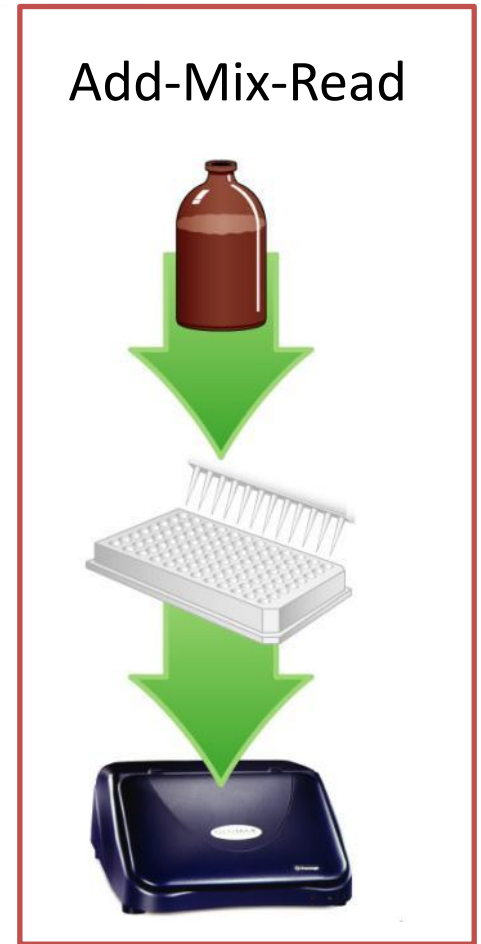
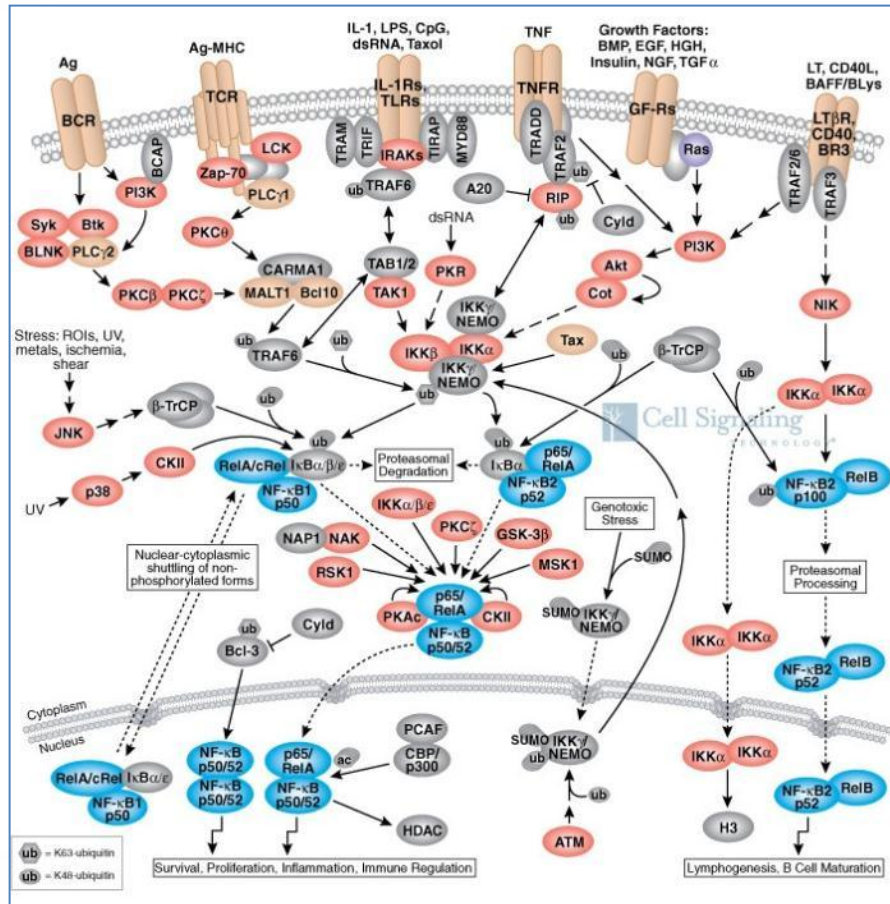
Richard Somberg, Ph.D.

October 2012

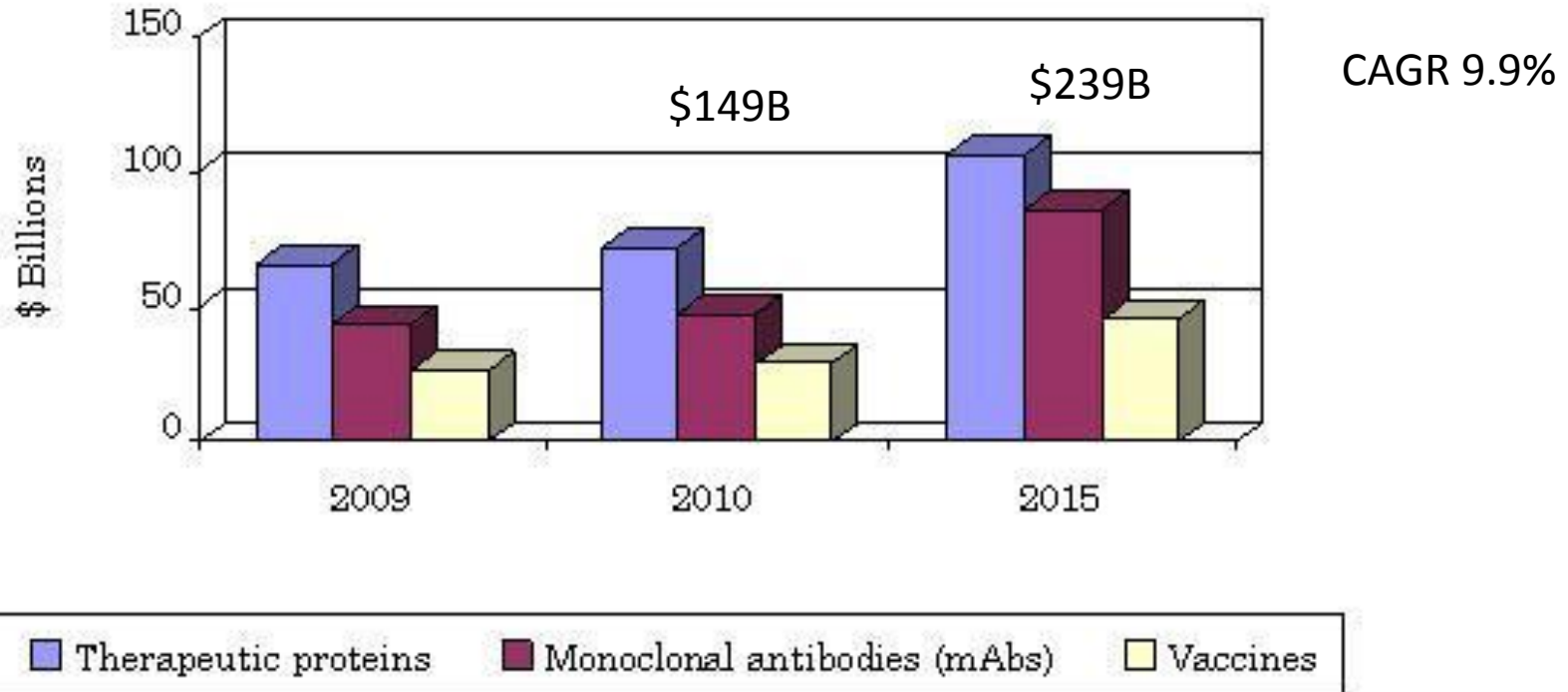
## Outline

- Introduction to ADCC – *Problem with classic ADCC assays*
- Principle of the ADCC Reporter Bioassay
- “Cells as reagents” – *Frozen, thaw-and-use format*
- Performance – *Specific, Linear, Precise, Accurate, Reproducible, as well as Potency & Stability Indicating*
- Testing Ab variants – *Glycosylation & Fucosylation*
- Commercial formats - *Kits & Cell Propagation Model*

# Complicated Biology — Simple Assay



## Global Biologics Market & Forecast

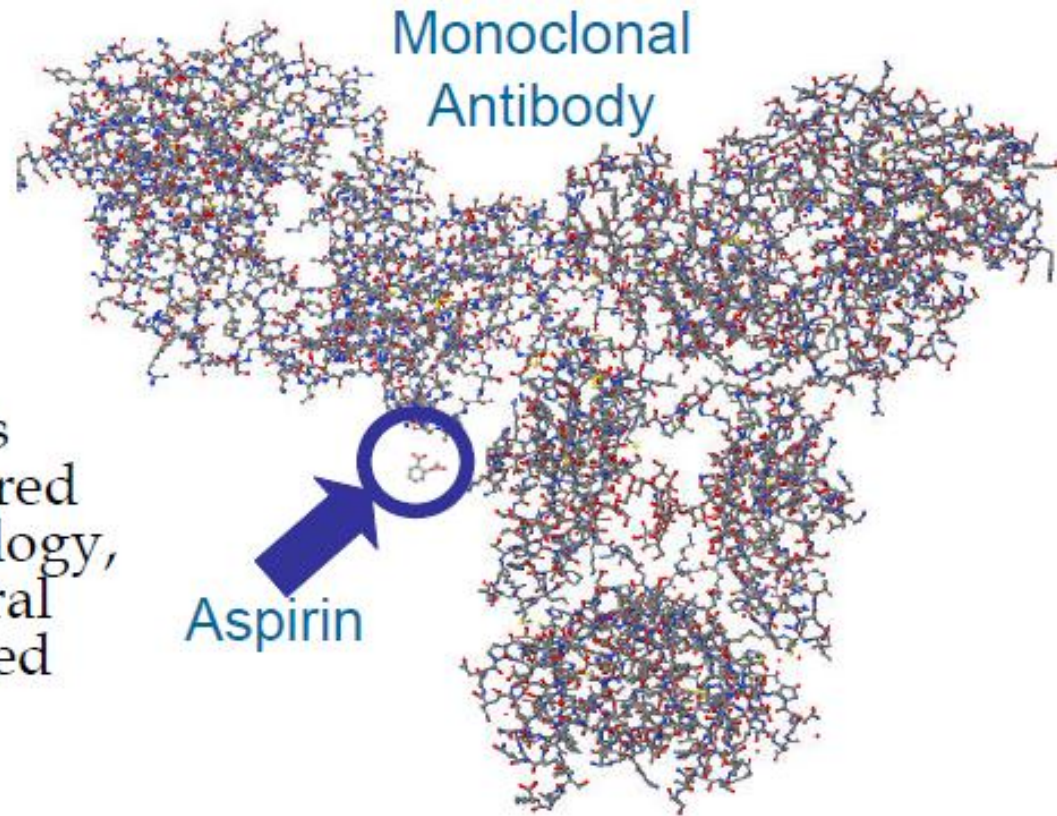


- Monoclonal antibodies (mAb) ~ 1/3<sup>rd</sup> of total biologics market
- mAb = \$48 billion in 2010, expected \$86 billion by 2015 (CAGR) of 12.4%

Source: BCC Research

## ***Biological Product – Monoclonal Antibody (mAb)***

- Biological products are **generally** produced using a living system or organism.
- Biological products may be manufactured through biotechnology, derived from natural sources, or produced synthetically.



Source: FDA webinar, Feb 15, 2012

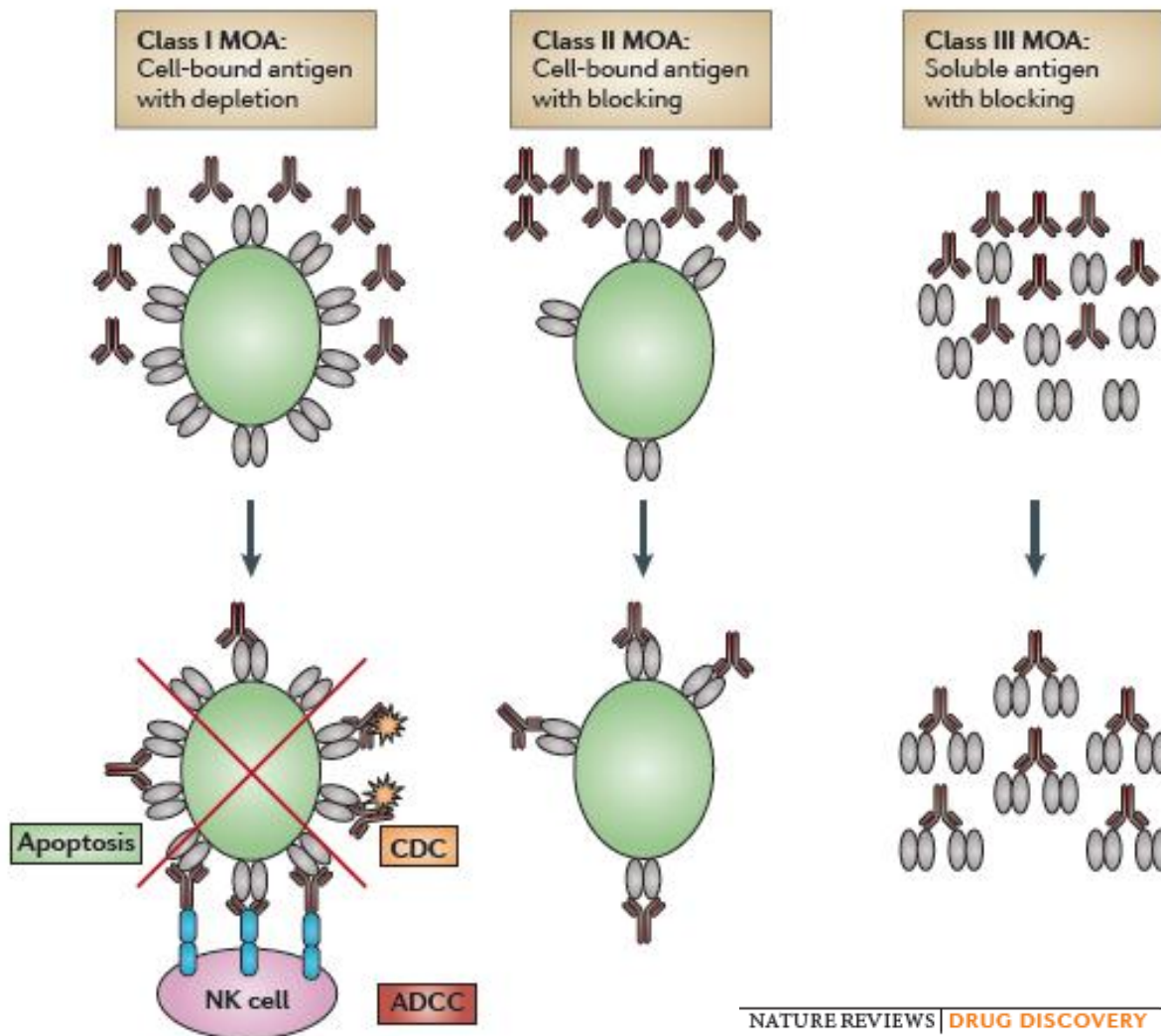
## *An Ideal Bioassay...*

- Reflective of the mechanism of action (MOA) of the biological product
- Well controlled (precise, accurate, robust, reproducible)
- Stability-indicating
- Usable as a QC lot-release assay

Modified from Chana Fuchs (DMA/CDER)

*In this webinar, we will demonstrate how the novel ADCC Reporter Bioassay fulfills each of these elements*

# Mechanism of Action (MOA) for mAb



# ***Introduction to ADCC***



## What is ADCC?

**Antibody-dependent cell-mediated cytotoxicity (ADCC)** is the main MOA of antibodies through which virus-infected or other diseased cells are targeted for destruction by components of the cell-mediated immune system, such as NK cells

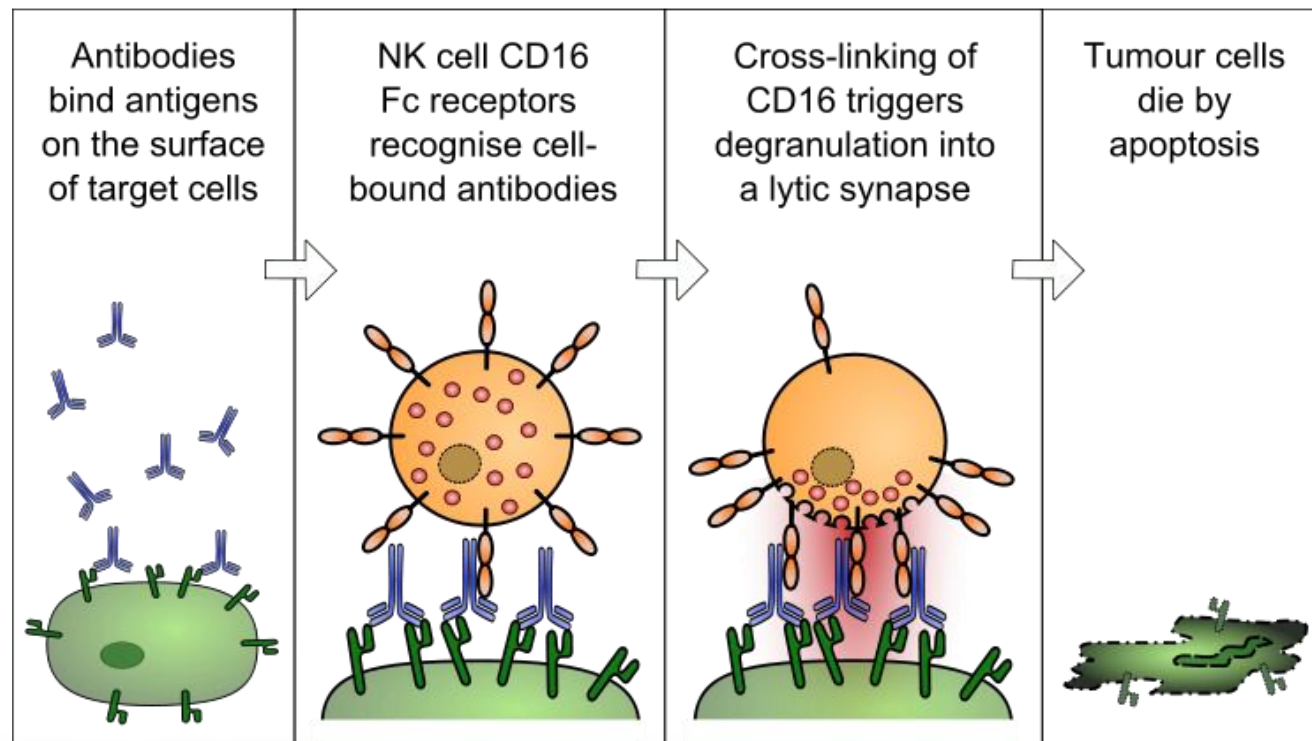


Image source: Wikipedia

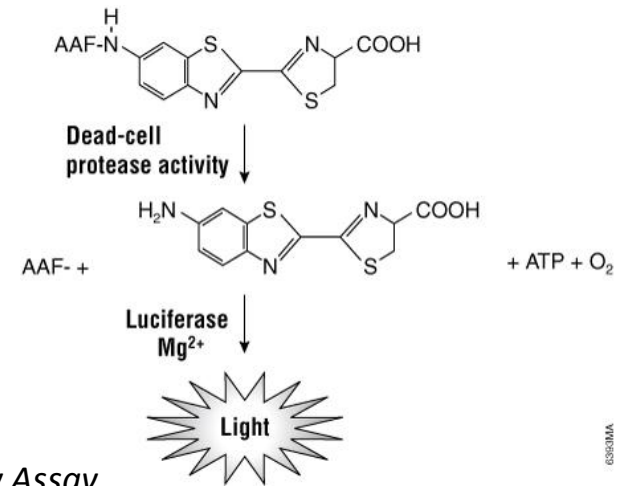
# Classic ADCC Assays

## Effector cells

- PBMCs (peripheral blood mononuclear cells)
- NK from PBMCs
- NK cell lines

## Target cells

- Load with chromium-51 or Eu
- Monitor cell lysis (LDH, Calcein AM, GAPDH, CytoTox-Glo™)



*CytoTox-Glo™ Cytotoxicity Assay*

# ***The Problem with Current ADCC assays***

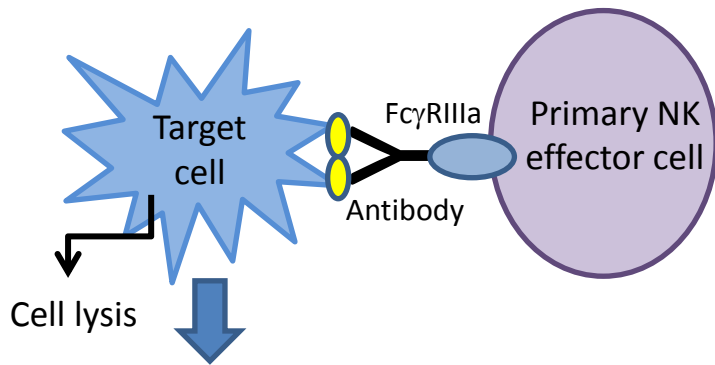
## ***Case Study – ADCC Challenge***

- A company acquired a late-stage mAb drug and needed to switch the manufacturing cell line and process to fit into their standard process
- FDA requested that an assay examining the mAb mechanism of action (ADCC) be used to demonstrate similarity between process change
- Classic ADCC assays had poor reproducibility, high variability, and were not suitable for use
- The company developed a reporter assay with low variability, high reproducibility, and used it to successfully demonstrate similarity and make the manufacturing cell line change

***Solution: A Better ADCC Bioassay***

# Classic ADCC assay vs ADCC Reporter Bioassay

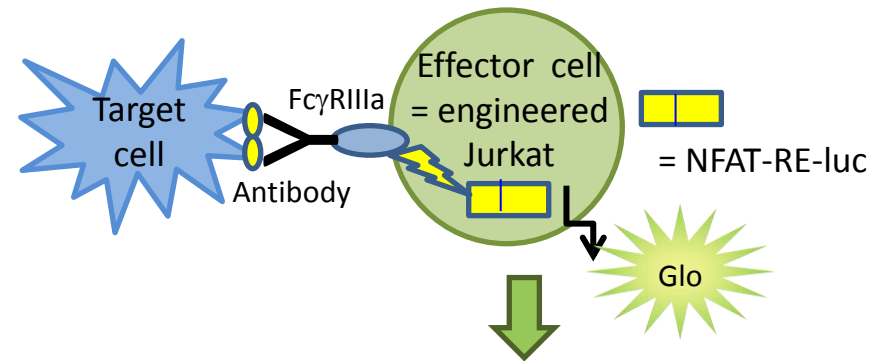
## Classic ADCC assay



**Signal is from target cell**

- High variability of assay - mainly due to primary NK cells
- Spontaneous lysis of target & effector cells results in high background

## Reporter-based ADCC bioassay



**Signal is from effector cell**

- Reduced variability by replacing NK cells with genetically engineered stable cell line
  - FcγRIIIa (V158)
  - NFAT-RE luc2
- Improved bioassay performance with robust reagents and assay design

# ***Principle of the ADCC Reporter Bioassay***

# Scientific Basis of ADCC Reporter Bioassay

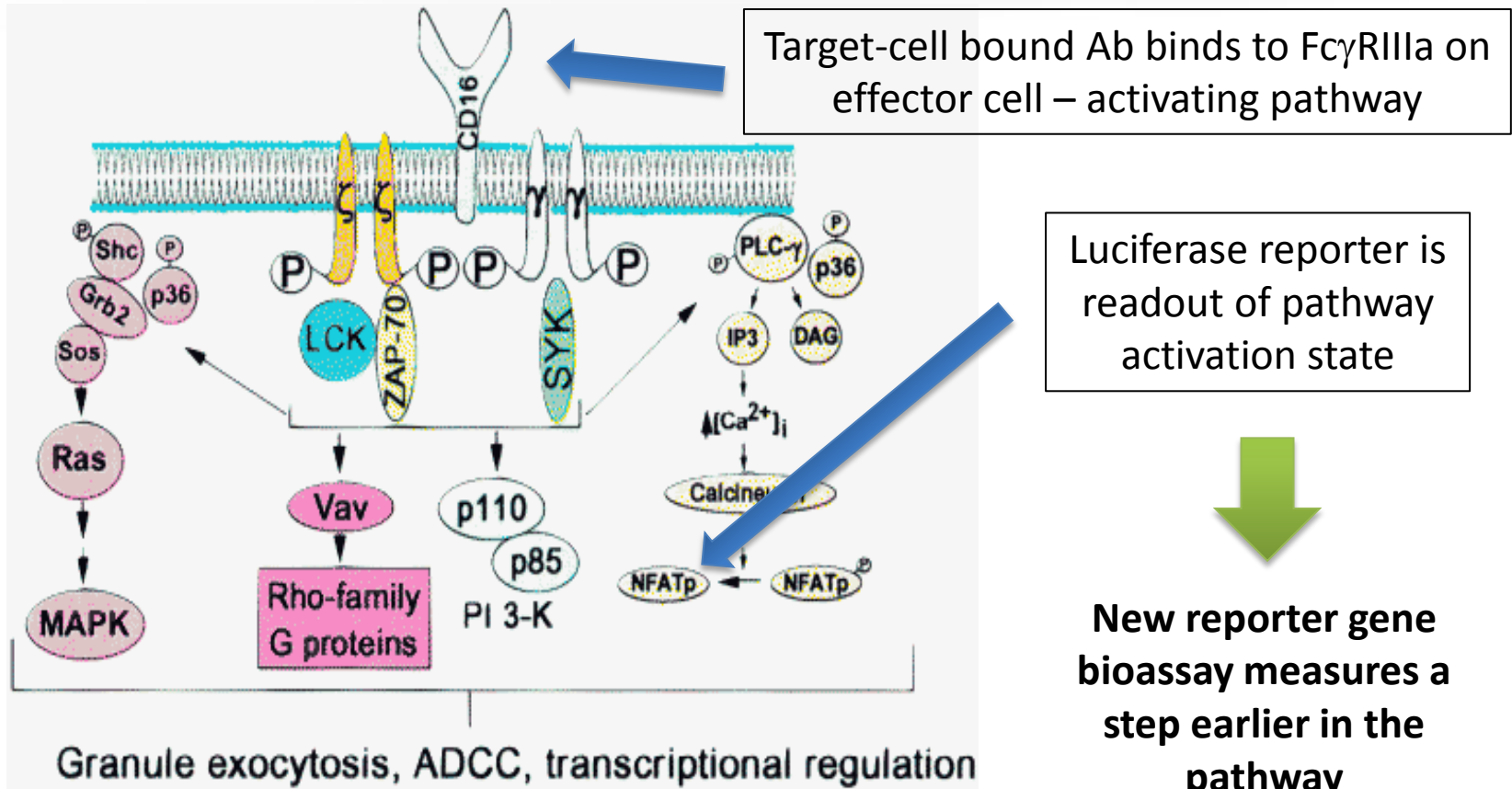
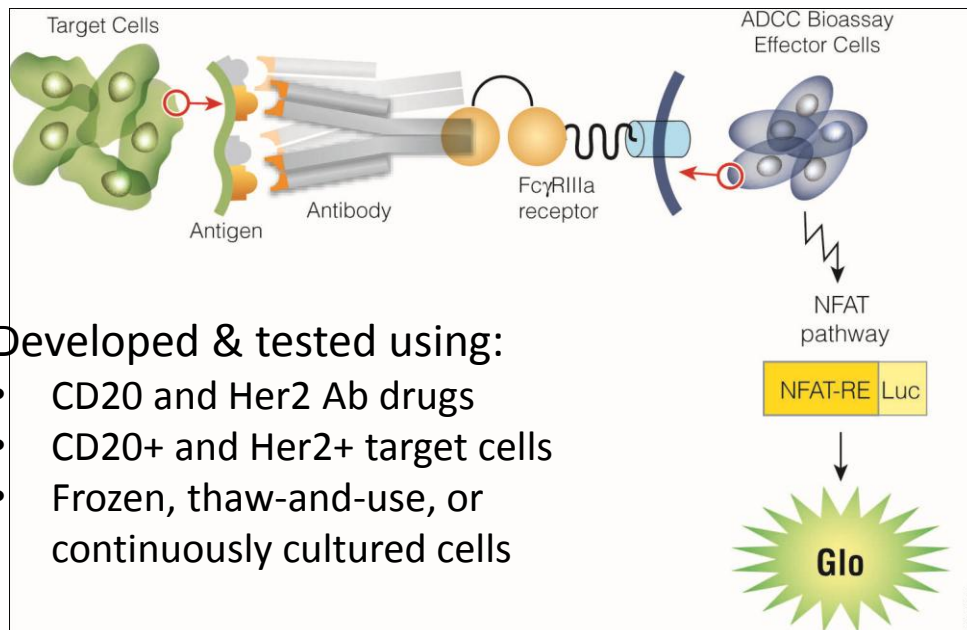


Image source: Leibson-PJ, *Immunity* 1997



# ADCC Reporter Bioassay - Development

## Low Variability NFAT-RE Luciferase bioassay



Developed & tested using:

- CD20 and Her2 Ab drugs
- CD20+ and Her2+ target cells
- Frozen, thaw-and-use, or continuously cultured cells
- Extensive 'alpha' evaluations:
  - tested in multiple global biopharma & biotechs
  - tested in multiple systems

1. Effector cells are engineered to express FcγRIIIa (V158) and NFAT-RE luc2 luciferase
2. 'Cells as reagents' (thaw-and-use)
3. Homogeneous assay format – simple 'add-mix-read' bioluminescent assay
4. Optimized and robust assay reagents and protocol
5. Performance characteristics that meet needs of stability testing, lot release and Ab characterization

# Bioassay Characteristics - ICH Guideline Q2 [R1]

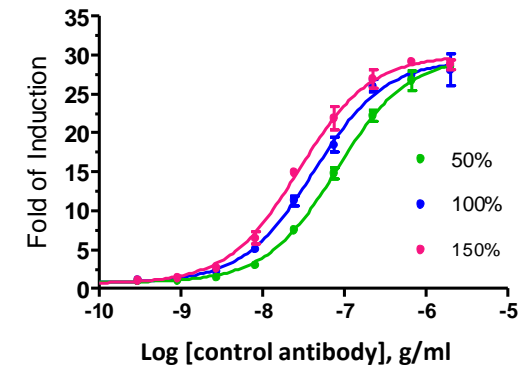
## Validation of Analytical Procedures

- Accuracy
- Precision:
  - ✓ Repeatability (intra-assay precision)
  - ✓ Intermediate precision (day to day, analyst-to analyst)
  - ✓ Reproducibility (lab to lab)
- Specificity
- Linearity
- Range
- Robustness

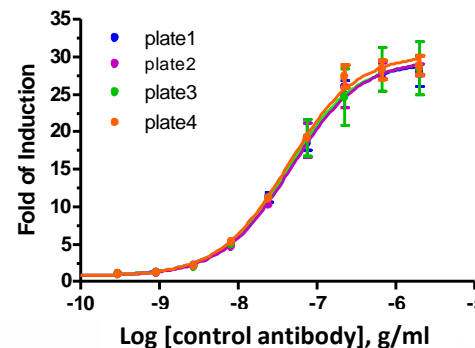
**Design:**

- Two analysts
- Three days
- Four plates per day
- ✓ 100% vs 50%
- ✓ 100% vs 75%
- ✓ 100% vs 125%
- ✓ 100% vs 150%

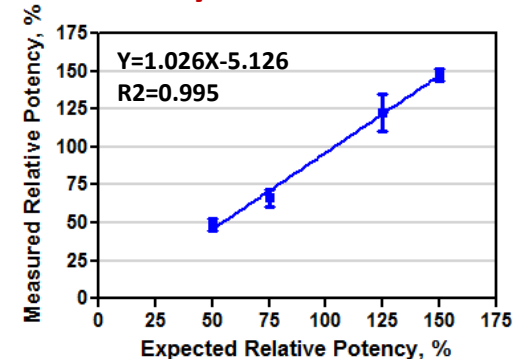
### Relative potency



### Repeatability



### Linearity

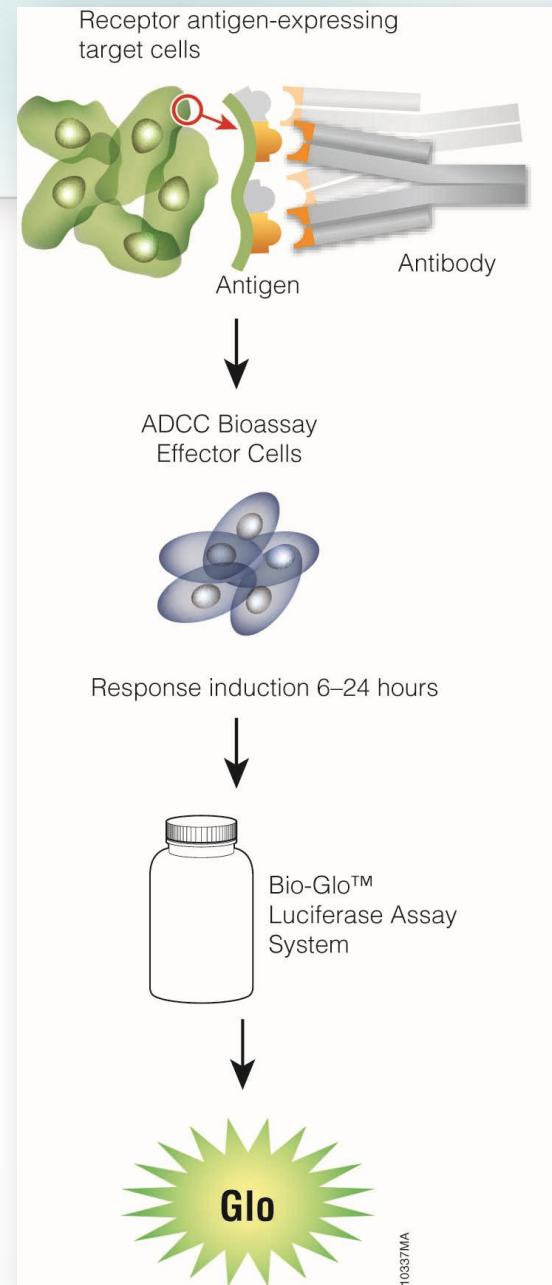


# ***Simple Protocol***

# ADCC Reporter Bioassay Protocol

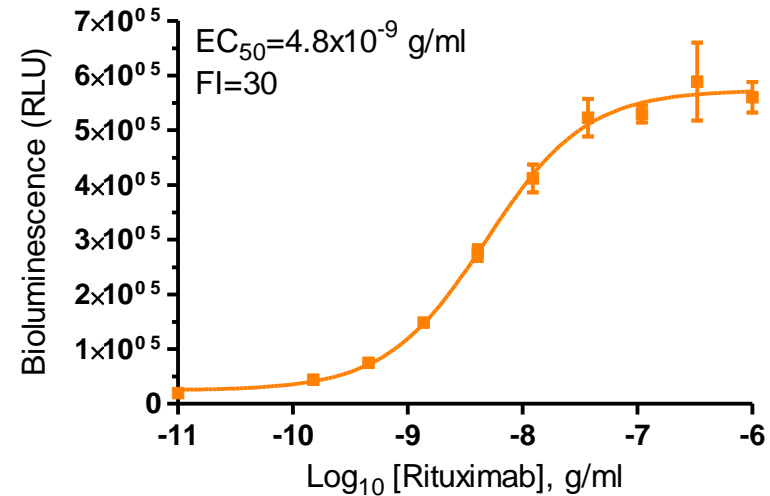
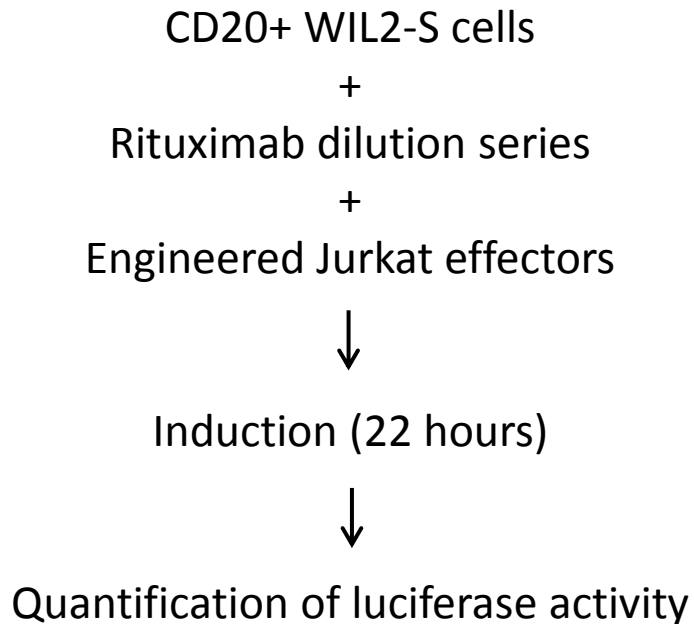
## Single day bioassay

1. Incubate control, reference or test antibody with target cells.
2. Add engineered effector cells containing:
  - FcγRIIIa (V158)
  - NFAT-RE luc2 luciferase
3. Incubate to allow for pathway activation (as short as 6 hours).
4. Add luciferase detection reagent and measure luminescence.



# ADCC Reporter Bioassay – Initial Results

## Assay protocol:



## Specifics

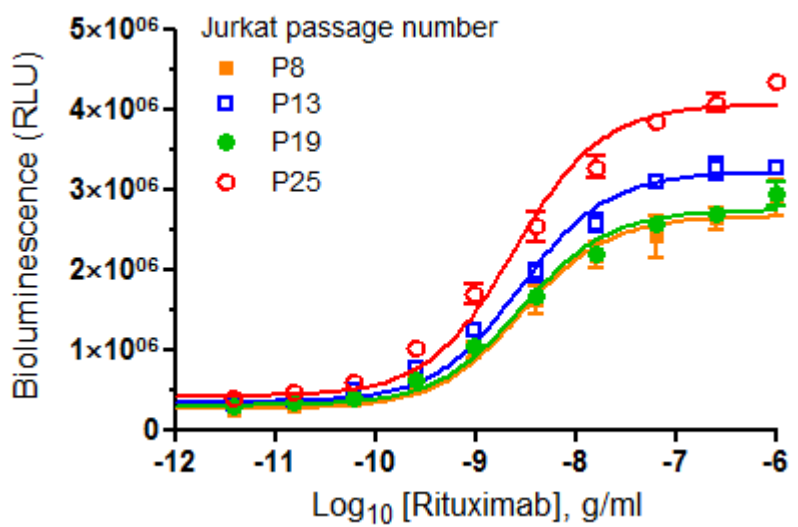
- **E:T ratio = 6:1**  
(150k effector cells:25k WIL2-S target cells, per well)

***Cell Selection and Frozen,  
Thaw-and-Use Format***

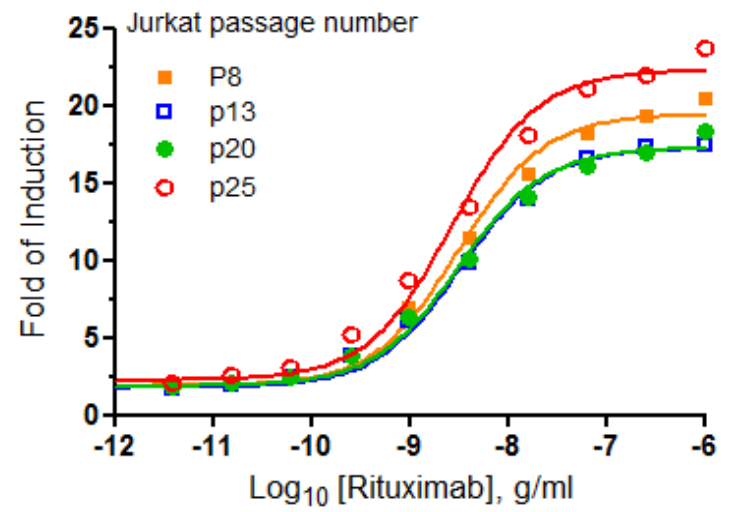
# Engineered Effector Cell Clone Selection

Clone selection based on maximizing RLU, fold induction, and passage stability

### Bioluminescence



### Fold Induction



	P8	p13	p20	p25
EC50	3.094e-009	3.211e-009	3.077e-009	2.665e-009

- E:T ratio = 7.5:1
- 6hr induction
- Bio-Glo™ Luciferase Assay System

# Cells as Reagents

## Frozen, Thaw-and-Use Cells

1. Human cell lines
  - Developed as Thaw-and-Use for immediate use in bioassay
  - Designed to give good recovery and robust response upon thawing
2. Thaw-and-Use format
  - Cell propagation conditions & defined freezing protocol control assay performance for a consistent bioassay response
  - No pre-culturing prior to assay means less variability introduced
  - Indefinite storage
  - Identical cells in bioassay, day-to-day
3. Minimizes pre-assay planning, time & labor
  - Ample cell banks provide long-term supply

**→ No cell culture required with cells in frozen, thaw-and-use format**



## Complete QC on Cells

### Production cell batches are rigorously tested:

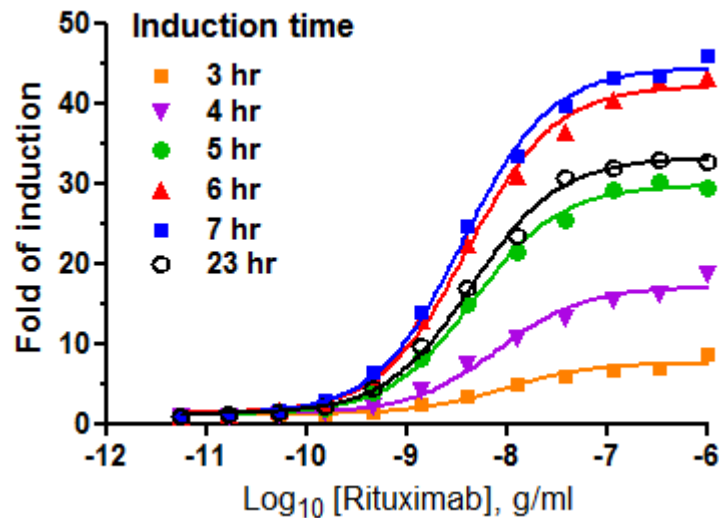
- STR analysis – cell ID profile (human)
- CO1 analysis (cytochrome oxidase) – test for presence of species (human and other potential contaminants)
- Cell doubling time under propagation conditions
- Mycoplasma (Hoechst and direct culture)
- Sterility
- Cell density
- Cell viability after thaw
- Fill volume
- ADCC Reporter Bioassay ( $EC_{50}$  and fold induction)



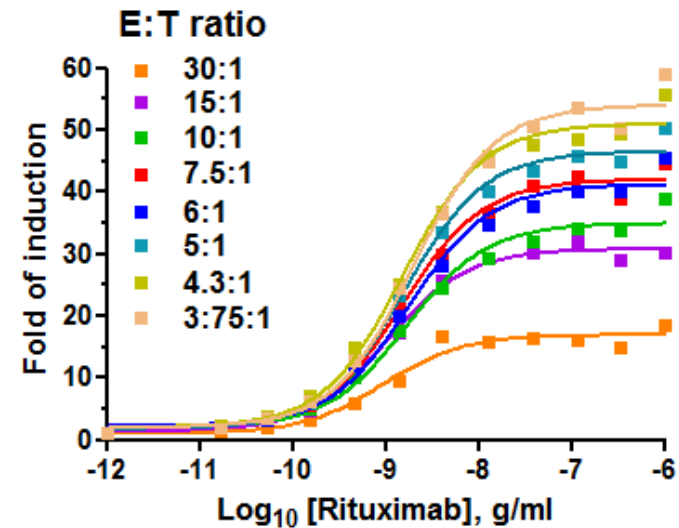
# ***Optimization Studies***

# Critical Assay Parameters

## Induction time



## E:T ratio with constant Effector cell number



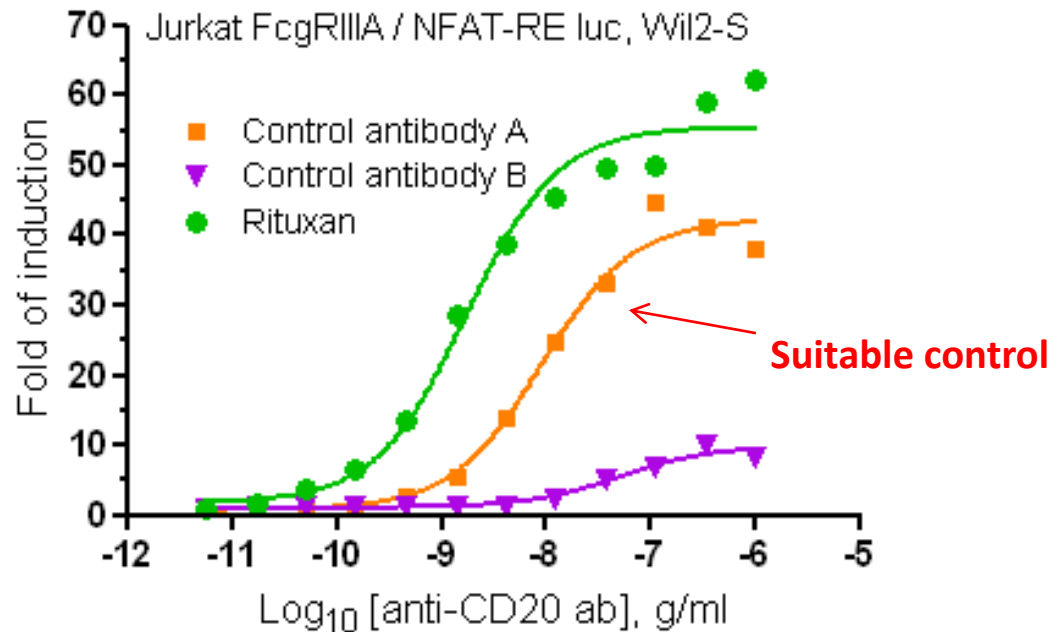
## Other parameters tested:

- Assay buffer: serum concentration, use of low IgG serum
- Cell numbers per well
- Pre-plating and incubation time: target cell plating, antibody/target cells incubation
- Assay plates: White flat, V- or U-bottom plates

# Selection of Control Antibody

## Requirements:

- $EC_{50}$  close to the range of biologic Ab drugs
- Good fold induction
- Good stability



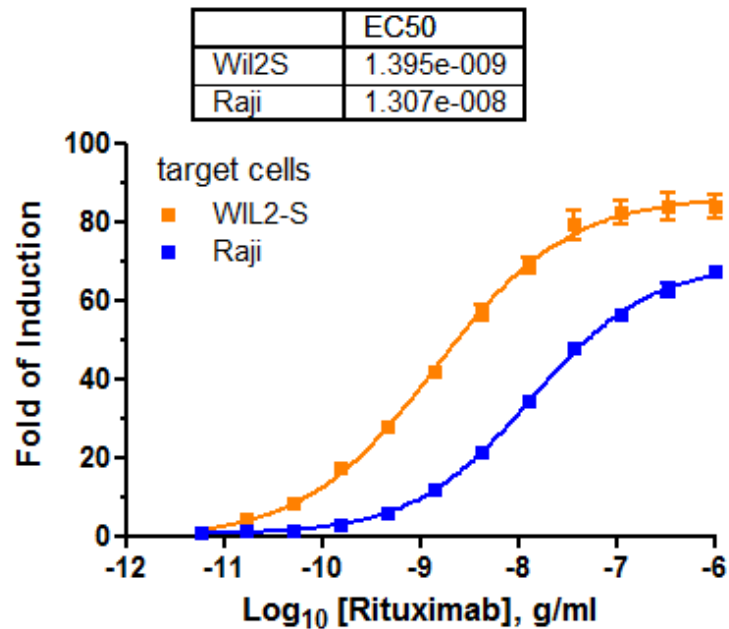
Anti-CD20	Rituximab	Control Ab (A)	Control Ab (B)
$EC_{50}$ (g/ml)	$1.7 \times 10^{-9}$	$8.9 \times 10^{-9}$	$47.3 \times 10^{-9}$
Fold Induction	62	38	8

# Use of Different Target Cells

*Suspension or adherent target cells can be used*

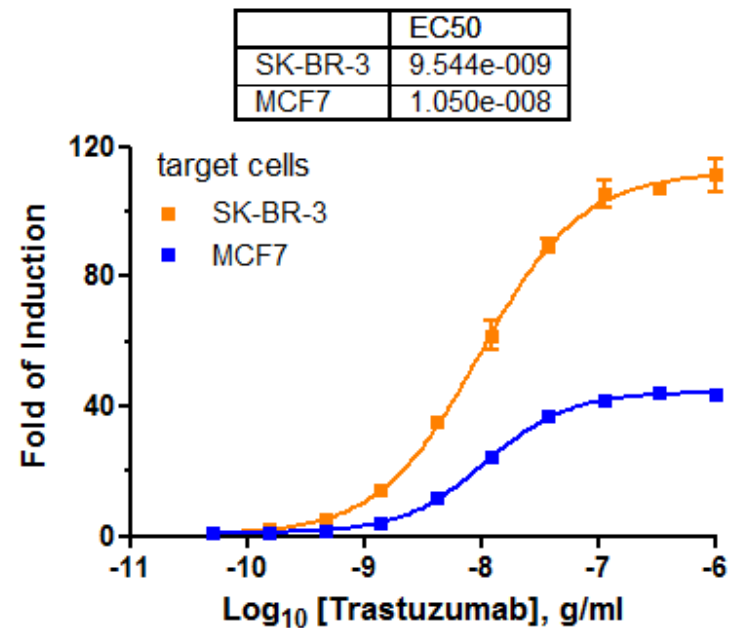
## Rituximab (anti-CD20)

**CD20<sup>+</sup> B cell lines (suspension) as target cells**



## Trastuzumab (anti-Her-2)

**Her2<sup>+</sup> breast cancer cell lines (adherent) as target cells**

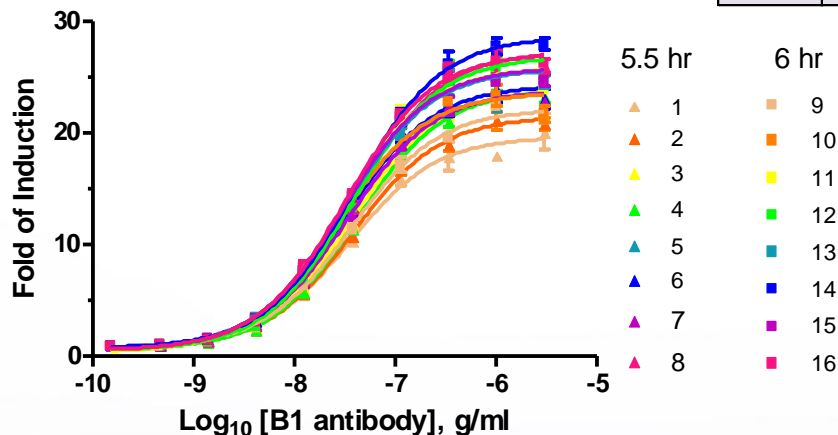


# Bioassay Development: Optimization Using DOE

## Variables:

1. Induction time
2. Target/Ab pre-incubation
3. Effector cell number
4. Target cell number

run	induction time hr	Target cell / Ab	Jurkat cell	Target cell
		incubation time(mins)	plating number (K)	plating number (K)
1	5.5	30	75	10
2	5.5	30	75	12.5
3	5.5	30	90	12
4	5.5	30	90	15
5	5.5	45	75	10
6	5.5	45	75	12.5
7	5.5	45	90	12
8	5.5	45	90	15
9	6	30	75	10
10	6	30	75	12.5
11	6	30	90	12
12	6	30	90	15
13	6	45	75	10
14	6	45	75	12.5
15	6	45	90	12
16	6	45	90	15



## Outputs & Results:

Good response (fold induction) = 19-27

Good (low) L-term values = 0.1-0.2\*

\* a measure of assay precision around the EC50 determination  
(log width of the 95% confidence interval around logEC50)

# *Performance*

## *Qualification Studies*

- **Parallelism** and measurement of **potency** relative to the reference antibody
- **Linearity & accuracy** of observed versus expected potencies across the desired working range of potencies
- **Precision**
  - intra-assay
  - intermediate (inter-assay) precision
- **Specificity** to show response is dependent on specific antibody and the presence of target cells and FcγRIIIa on effector cells, and not other components
- **Stability-indicating** to show the bioassay is capable of detecting loss of structural integrity of an antibody

*These qualification studies are critical to demonstrate a useful and effective ADCC bioassay*



# ADCC Reporter Bioassay is Specific

Target cells, effector cells and specific antibody

■ Wil2-S, Jurkat/NFAT-luc+F $\gamma$ R11a, Rituximab

No Target cells

● NO Wil2-S, Jurkat/NFAT-luc+F $\gamma$ R11a, Rituximab

No Effector cells or no F $\gamma$ R11a

▲ Wil2-S, Jurkat/NFAT-luc (NO F $\gamma$ R11a), Rituximab

▲ Wil2-S, NO Jurkat/NFAT-luc+F $\gamma$ R11a, Rituximab

No antibody or non-specific antibody

▼ Wil2-S, Jurkat/NFAT-luc+F $\gamma$ R11a, NO Rituximab

▼ Wil2-S, Jurkat/NFAT-luc+F $\gamma$ R11a, Trastuzumab

## Assay signal is dependent on:

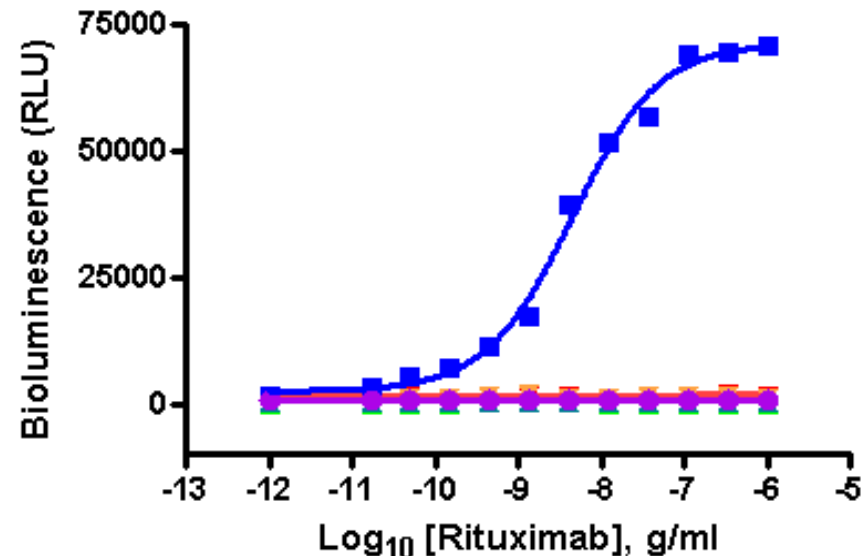
Presence of Target cells

+

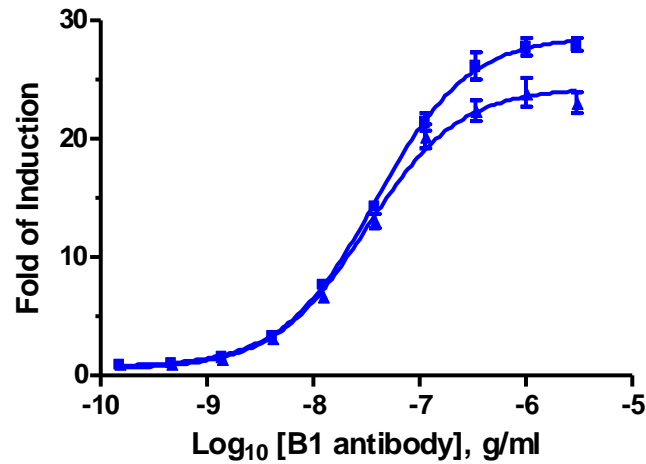
Presence of F $\gamma$ R11a receptor

+

Appropriate specific antibody

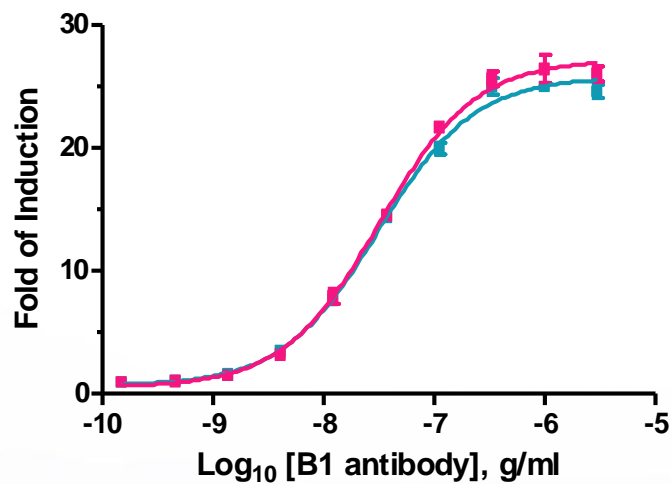


# ADCC Reporter Bioassay is Robust



## Time of induction

Run	Induction time	EC <sub>50</sub>
1	6.0 hr	3.15x10 <sup>-8</sup> g/ml
2	5.5 hr	3.83x10 <sup>-8</sup> g/ml

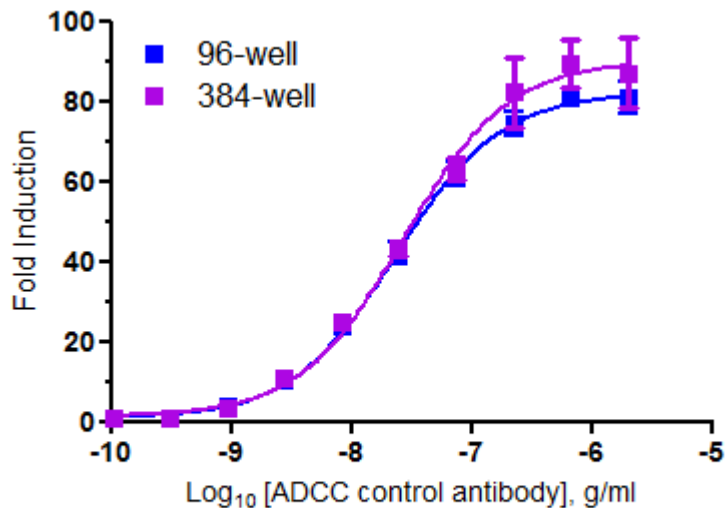


## E:T ratio and cell # per well

Run	E:T ratio	E cell #	T cell #	EC <sub>50</sub>
1	7.5:1	75k	10k	3.09x10 <sup>-8</sup> g/ml
2	6:1	90k	15k	3.83x10 <sup>-8</sup> g/ml

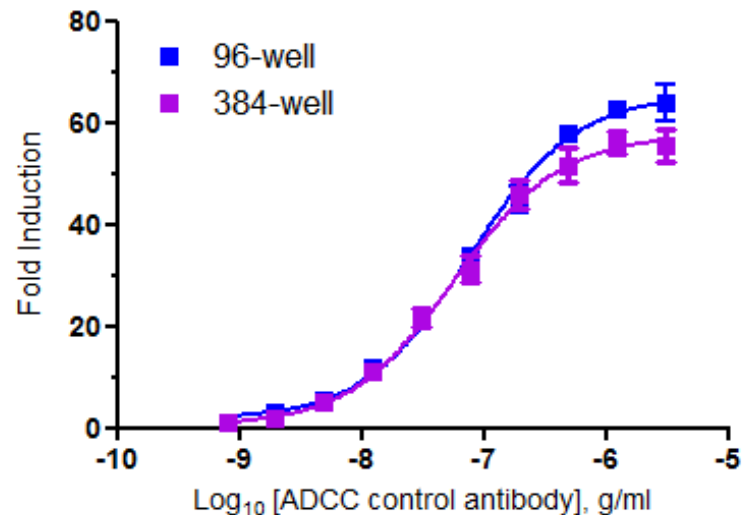
# Miniaturization to 384-well Plates

## WIL2-S target cells



	96-well	384-well
EC50	2.286e-008	2.651e-008

## Raji target cells



	96-well	384-well
EC50	7.434e-008	5.583e-008

Assay volume per well	Target cells	Antibody	Effector cells	Bio-Glo™
96-well plate	25µl	25µl	25µl	75µl
384-well plate	5µl	5µl	5µl	15µl

# Assay Qualification Results

Bioassay uses frozen-thaw-and-use cells for both effector cells and WIL2-S target cells

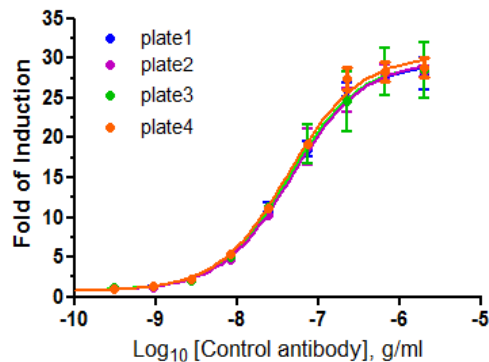
## Design:

- Three days
- Four plates per day
- ✓ 100% vs 50%
- ✓ 100% vs 75%
- ✓ 100% vs 125%
- ✓ 100% vs 150%

## Representative plate layout

	1	2	3	4	5	6	7	8	9	10	11	12	Plate1
A													
B		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		100%
C		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		50%
D		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		100%
E		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		50%
F		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		100%
G		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		50%
H													

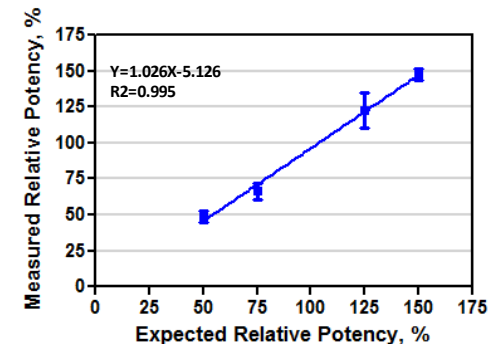
## Repeatability



**Precision** = average of RSD (%) = 7.3%  
**Accuracy** = average of Recovery (%) = 95.8%

	Antibody Sample	Measured Potency (%)	Mean Potency (%)	SD %	Recovery (%)	RSD (%)
day 1	50%	48.5	48.9	3.9	97.7	7.9
		45.2				
		52.9				
day 2	75%	63.1	66.4	5.9	88.5	8.9
		62.9				
		73.2				
day 3	125%	112.1	123.0	12.3	98.4	10.0
		136.3				
		120.5				
day 1	150%	148.8	147.6	3.6	98.4	2.4
		150.4				
		143.6				

## Linearity



*Good repeatability, accuracy, precision and linearity were obtained*

# Assay Qualification Results with Raji Target Cells

## Analyst 1

Day	Antibody Sample	Measured Potency (%)	Mean Potency (%)	SD (%)	Recovery (%)	CV (%)
1	50%	49.9	51	0.7	102	1.4
2		51.3				
3		50.5				
1	75%	78.9	76	5.11	101	6.7
2		78.8				
3		70				
1	125%	118.6	117	1.19	94	1
2		116.9				
3		116.3				
1	150%	143.2	145	3.91	97	2.7
2		142.5				
3		149.6				

**Precision: 2.95%**

**Accuracy: (recovery average): 98.5%**

**Linearity:**  
 $Y=0.922x+5.0$

## Analyst 2

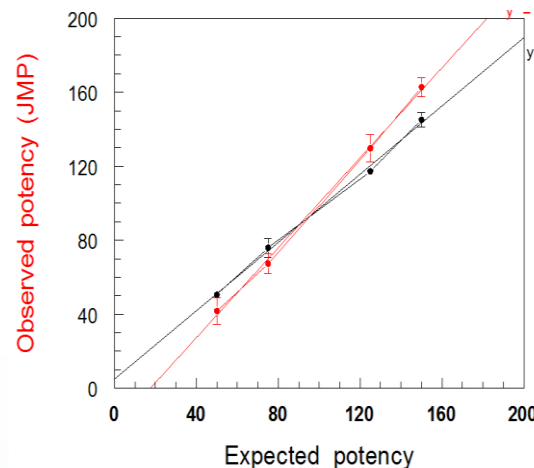
Day	Antibody Sample	Measured Potency (%)	Mean Potency (%)	SD (%)	Recovery (%)	CV (%)
1	50%	38.4	41.8	7.2	83.5	17.2
2		47.2				
3		33.2				
4		48.2				
1	75%	59.6	67.4	5.2	89.9	7.7
2		70.2				
3		69.3				
4		70.5				
1	125%	120	129.7	7.5	103.7	5.8
2		132.3				
3		137.8				
4		128.6				
1	150%	160.2	162.8	5.2	108.5	3.2
2		158.2				
3		162.7				
4		170				

**Precision: 8.47%**

**Accuracy (recovery average): 96.4%**

**Linearity:**  
 $Y=1.22x-21.3$

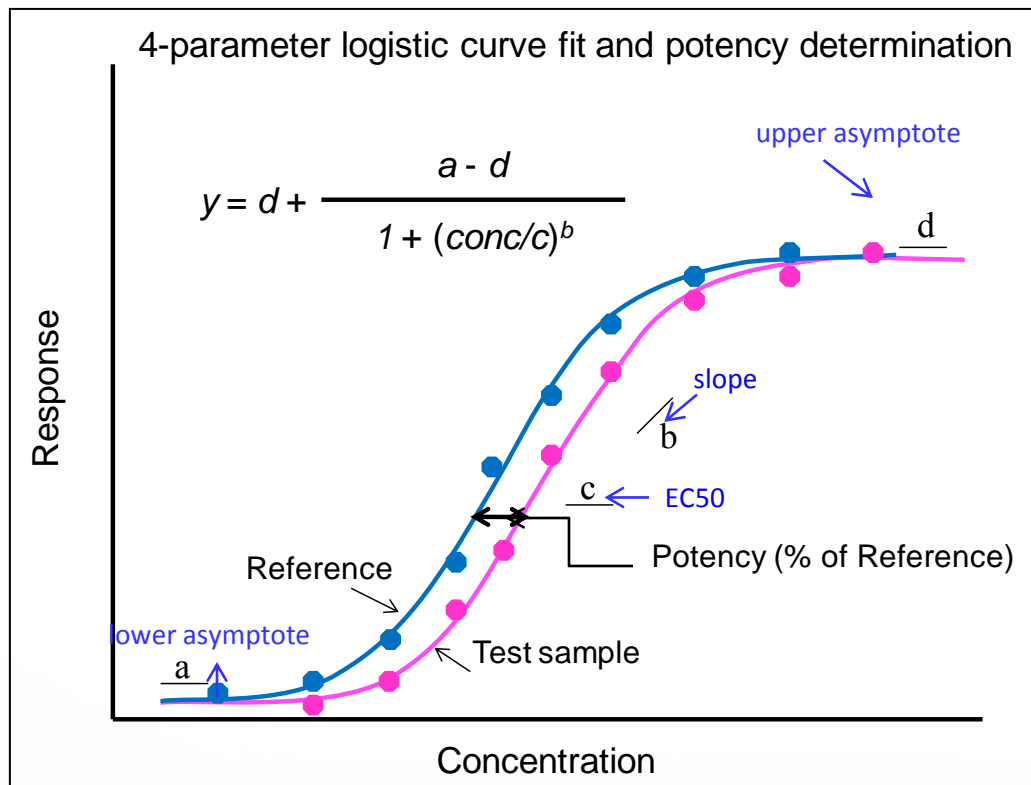
## Linearity



# Potency Determinations Using Quantitative Bioassays

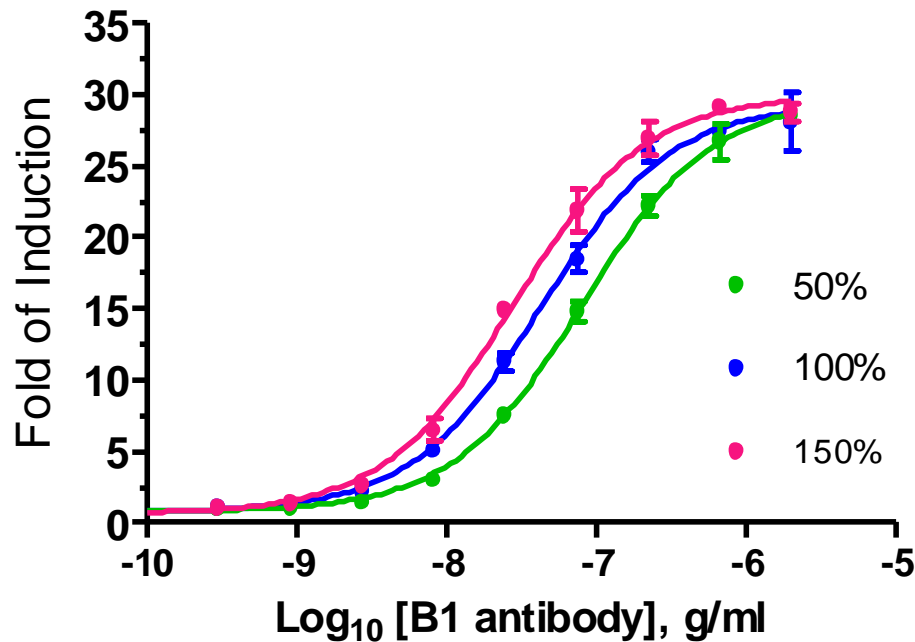
A test sample of unknown biological activity is compared with a reference sample of established biological activity in a dose-response study in the test system.

**The bioassay establishes potency relative to a reference standard**



- Curve fitting and statistical methods determine parallelism
- Parallel curves signify equivalent means of effecting biological activity
- Relative potency is quantified through shift of response along the x-axis
- Slope difference suggests non-equivalent means of effecting response if it falls outside of acceptance criteria; a manufactured lot would fail if this were so

# Measurement of Relative Potency & Parallelism



	100%-1	100%-2	100%-3	100%-4
EC50	4.230e-008	4.693e-008	4.323e-008	4.245e-008

relative potencv    102%    92%    100% control    102%

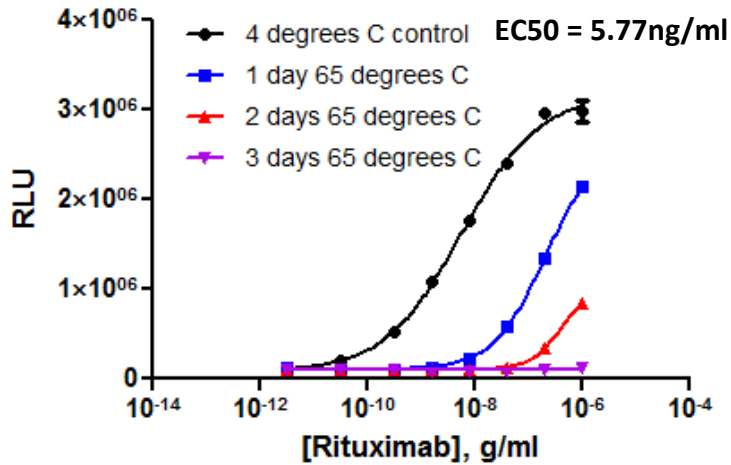
*Parallelism and relative potency determined with JMP Software*

# ***Stability Indicating***



# Stability Indicating for Fc Effector Function

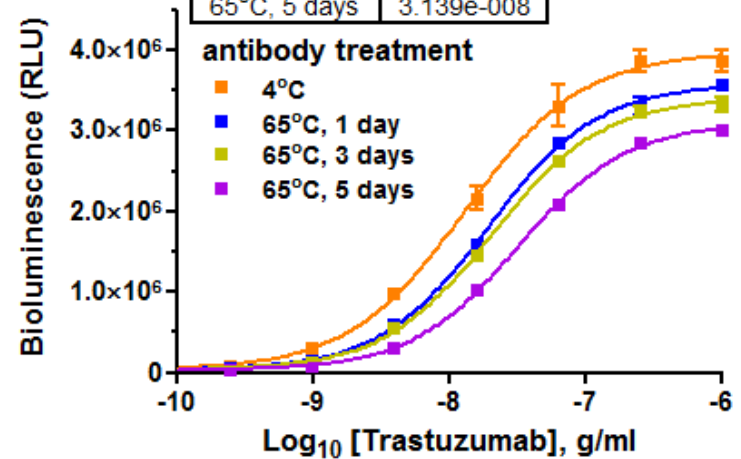
## Rituximab



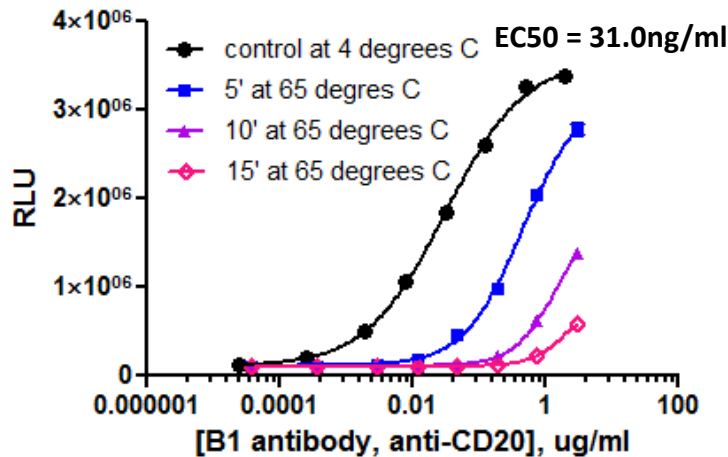
## Activity of heat-treated antibody drugs

## Trastuzumab

	EC50
4°C	1.284e-008
65°C, 1 day	1.902e-008
65°C, 3 days	2.031e-008
65°C, 5 days	3.139e-008

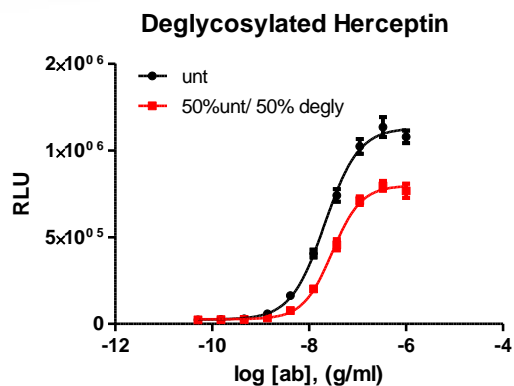


## Tositumomab

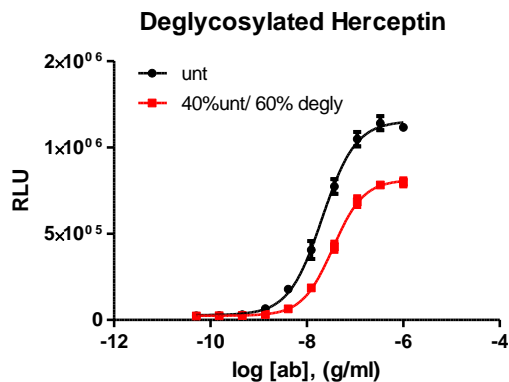


# ***Antibody Variants***

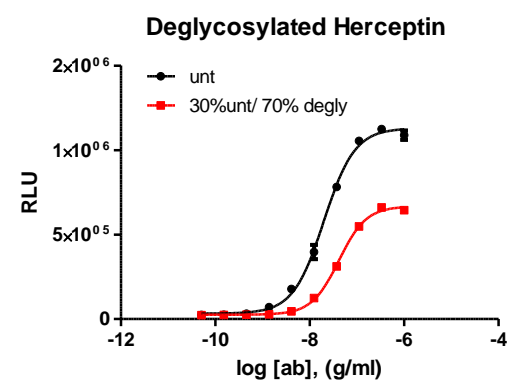
# Analysis of Mixed Glycosylation mAbs



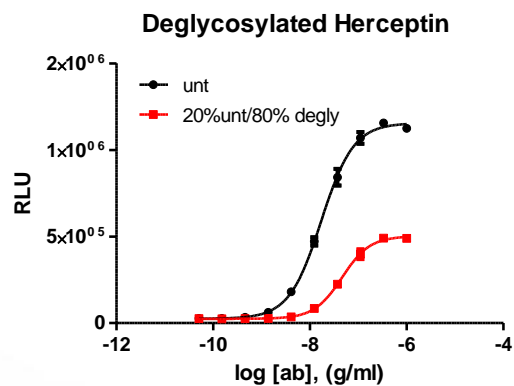
	unt	50%unt/ 50% degly
EC50	2.110e-008	2.990e-008



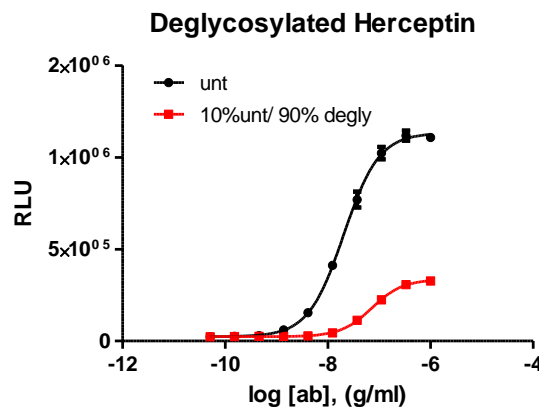
	unt	40%unt/ 60% degly
EC50	2.082e-008	3.486e-008



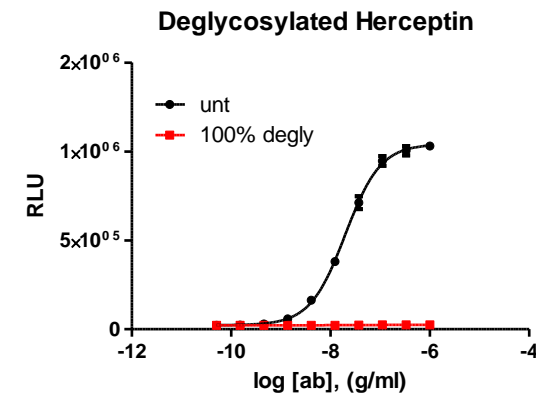
	unt	30%unt/ 70% degly
EC50	2.002e-008	4.153e-008



	unt	20%unt/80% degly
EC50	1.720e-008	4.626e-008



	unt	10%unt/ 90% degly
EC50	2.037e-008	7.174e-008



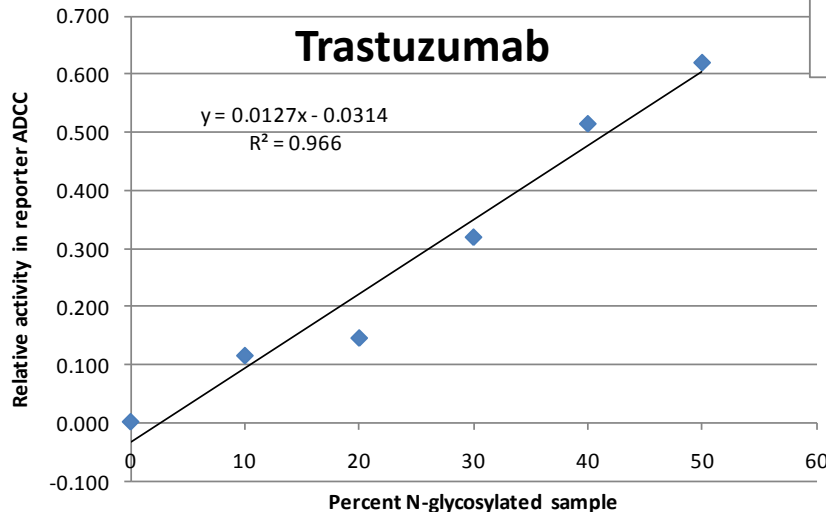
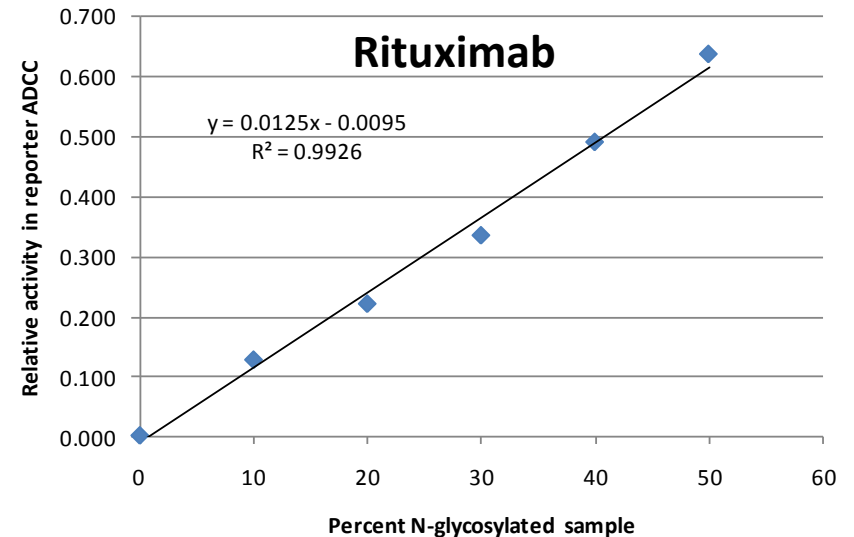
	unt	100% degly
EC50	1.988e-008	3.202e-008

Target cells: SKBR3; Unt = 100% glycosylated

# ADCC Reporter Bioassay Activity Correlates with Amount of Antibody N-glycosylation

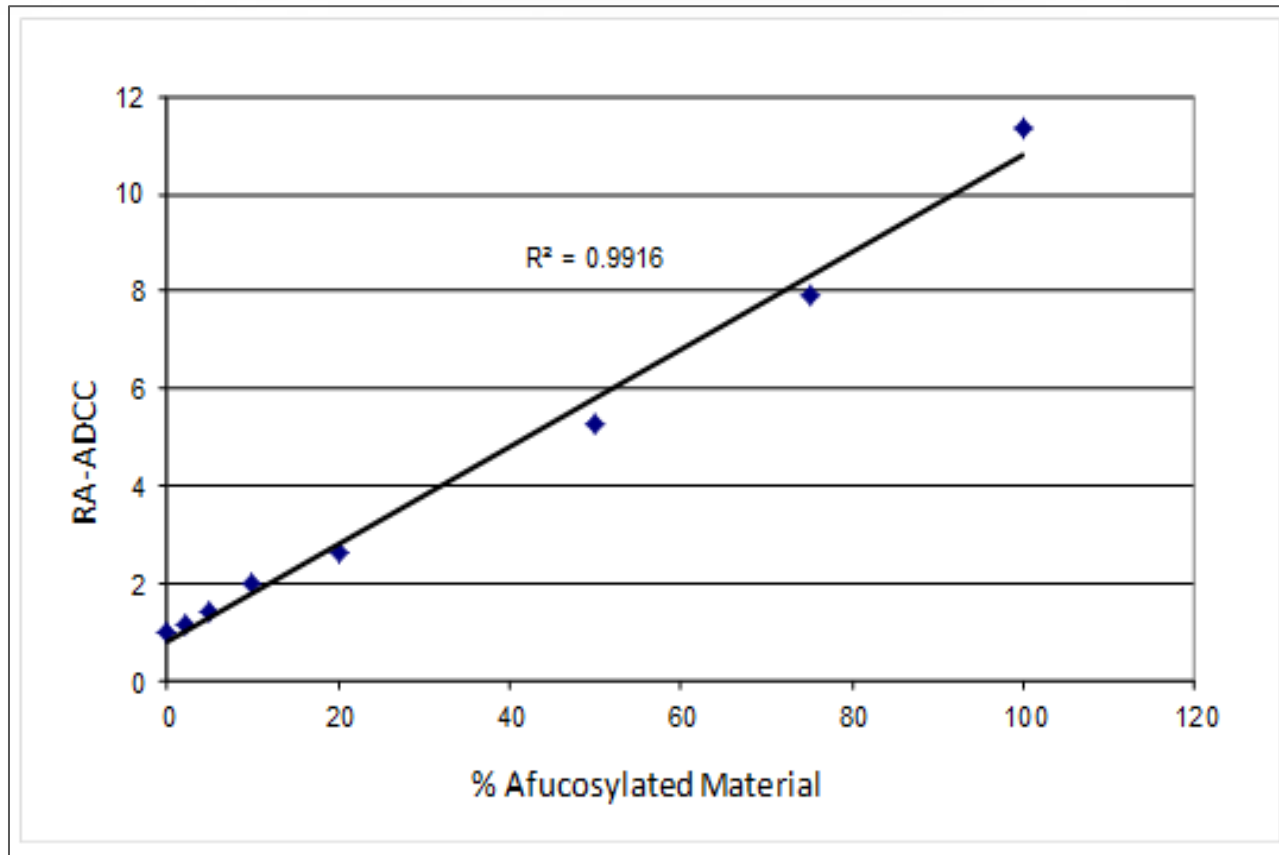
## Rituximab and Trastuzumab:

Linear correlation obtained between percentage of N-glycosylated antibody in blended antibody samples and relative luciferase reporter activity in ADCC reporter bioassay



*Small differences in Fc effector activity in ADCC pathway activation are easily distinguished in the ADCC reporter bioassay*

# *ADCC Reporter Activity Correlates with Amount of Antibody Afucosylation*



Linear correlation shown between percentage of afucosylated antibody in blended antibody samples and relative luciferase reporter activity in ADCC reporter assay

# Bioassay Characteristics - ICH Guideline Q2 [R1]

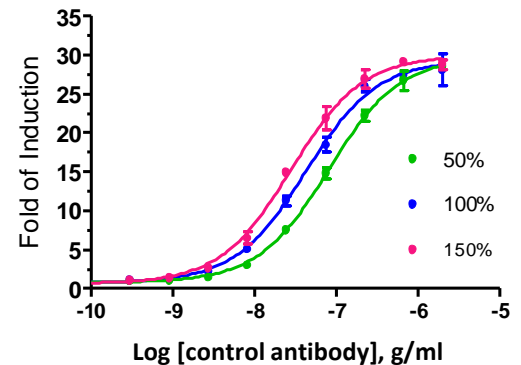
## Validation of Analytical Procedures:

- Accuracy
- Precision:
  - ✓ Repeatability (intra-assay precision)
  - ✓ Intermediate precision (day to day, analyst-to analyst)
  - ✓ Reproducibility (lab to lab)
- Specificity
- Linearity
- Range
- Robustness

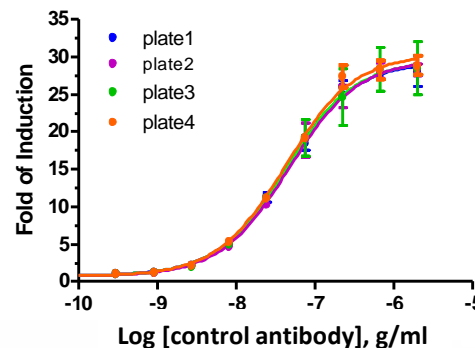
### Design:

- Two analysts
- Three days
- Four plates per day
  - ✓ 100% vs 50%
  - ✓ 100% vs 75%
  - ✓ 100% vs 125%
  - ✓ 100% vs 150%

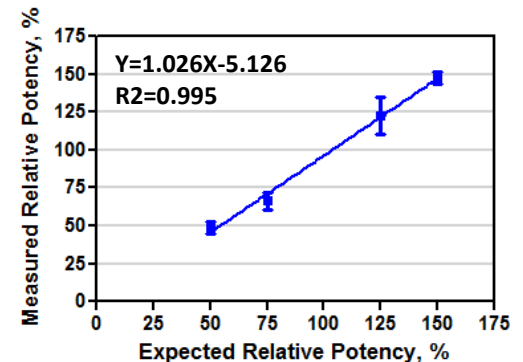
## Relative potency



## Repeatability



## Linearity



# ***External Evaluations***

## *Updates from Clients...*

- Approved manufacturing cell line switch by a pharmaceutical company
- Submitted in an IND filing
- Being developed for lot-release testing
- Charles River Laboratories and Catalent are providing ADCC Reporter Bioassay services
- Adopted by major pharmaceutical companies

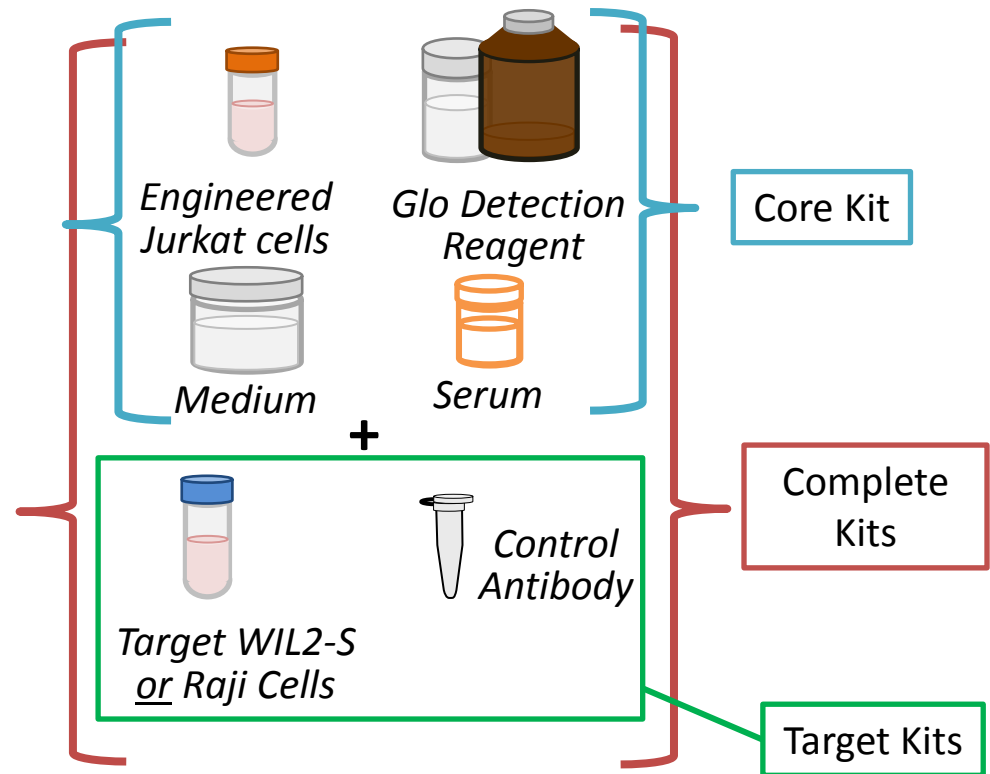


# ***Kit Formats***

# ADCC Reporter Bioassay Kit Configurations

**To be more flexible to research needs, we offer multiple kit formats:**

1. Core Kits:  
1X kit – Cat.# G7017  
5X kit – Cat.# G7018
2. Complete Kits:  
Available as Custom material
3. Target Kits:  
To be available later in year

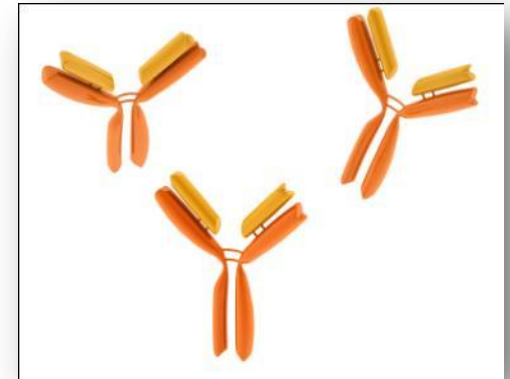


**Note:** the ADCC Bioassay Effector Cells are available for propagation and banking under a unique purchase agreement

# Summary of the ADCC Reporter Bioassay

## Features

- Low variability
- Engineered effector cells to replace primary NK cells (Jurkat FcγRIIIa/NFAT-RE luc2)
- “Cells as reagents”, frozen, thaw-and-use format – consistency & convenience
- Simple & robust protocol & reagents
- Broad applicability in use with multiple target cells – suspension or adherent



## Benefits

- Demonstrates precision, accuracy, linearity, robustness
- Can quantify potency and stability of therapeutic Ab drugs
- Can differentiate biological activity of Fc effector function in ADCC MOA resulting from small changes in Ab glycosylation

## *For more information*



Richard Somberg, PhD  
Strategic Collaborations Manager  
[Richard.somberg@promega.com](mailto:Richard.somberg@promega.com)

Neal Cosby, PhD  
Strategic Marketing Manager  
[Neal.cosby@promega.com](mailto:Neal.cosby@promega.com)

Or

Custom Order Department  
[COD@promega.com](mailto:COD@promega.com)

### ***Acknowledgements:***

Frank Fan, Terry Surowy, Jey Cheng, Rich Moravec, Denise Garvin, Aileen Paguio, Pete Stecha