

***New cell based assays and tools for epigenetics,  
biologics, metabolic markers and more.....***

Craig Malcolm: Product Manager Cell Analysis & Proteomics

Cell-Based Assay Tour - March 2014

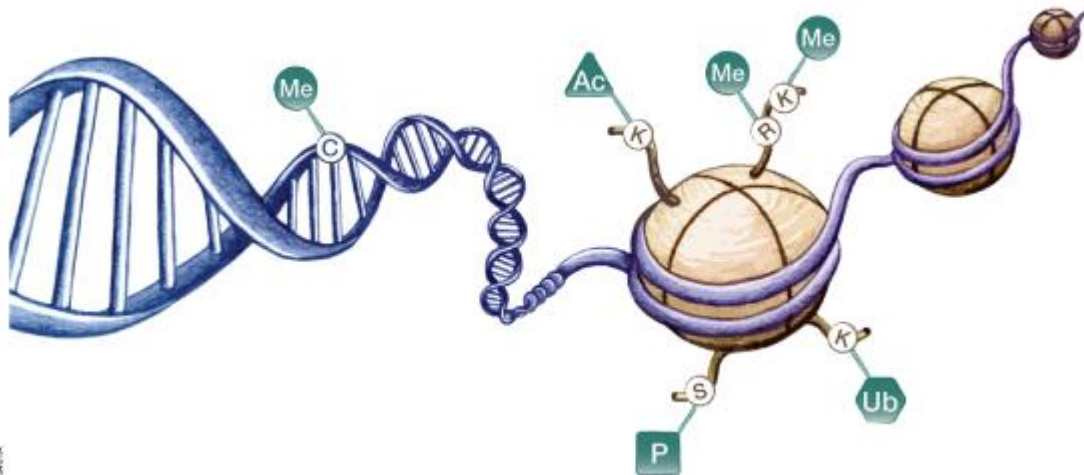
## *Outline*

- Brief outline of Promega's epigenetics portfolio
  - new isozyme-selective cell-based histone deacetylase assays
- Cell based assays for Biologics characterisation
  - Reporter gene bioassays for antibody-dependent cellular cytotoxicity (ADCC)
  - Additional reporter gene assays for other biologic therapeutic mAb's
- Recently launched suite of luminescence read-out metabolite assays
- New assays for oxidative stress (ROS-Glo)

## ***Epigenetics - definition***

- heritable changes in gene expression that arise from changes in chromosomes without alteration of DNA sequence
- changes occur throughout all stages of development or in response to environmental factors such as exposure to toxins or chronic stress and are implicated in diseases such as cancer
- Epigenetic mechanisms of gene regulation, which collectively make up the epigenome, include modifications to DNA and histone components of nucleosomes as well as expression of noncoding RNAs (ncRNAs)
- modifications can affect gene accessibility to DNA-binding and regulatory proteins such as methyl-CpG-binding proteins, transcription factors, RNA polymerase II and other components of the transcriptional machinery, ultimately altering transcription patterns, often in tissue- and cell-specific ways

## Typical epigenetic modifications



Epigenetic mechanisms involved in regulation of gene expression

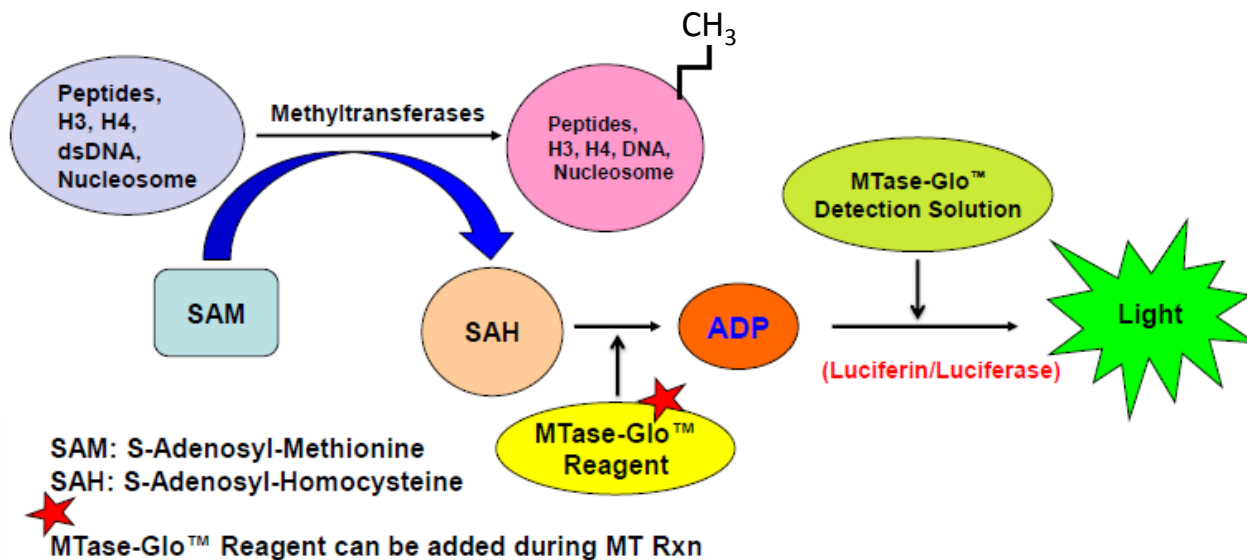
Cytosine residues within DNA can be methylated, and lysine and arginine residues of histone proteins can be modified. Me = methylation, Ac = acetylation, P = phosphorylation, Ub = ubiquitination.

## ***Promega's epigenetics portfolio - I***

- **Non-cell based assays**
- **MethylEdge™ Bisulfite Conversion System**
  - efficient DNA conversion and recovery with reduced template fragmentation using a protocol that can be completed in less than two hours
  - kit includes includes desulphonation and cleanup. Component
  - no additional cleanup kit required
- **Methylation-Specific Restriction Enzymes** *eg.* HpaII, MboI, MspI, and Sau3AI
  - valuable tools for studying DNA methylation patterns
  - coupled with robust DNA purification and PCR reagents - allow for detection of methylation-specific events

## Promega's epigenetics portfolio - II

- Methyltransferase (MTase-Glo) – *in development*
- Universal assay
- Based on production of s-adenosylhomocysteine (SAH)
- Designed for characterising the effects of compounds on DNA, protein and small molecule methyltransferases



## ***Promega's epigenetics portfolio - III***

- **Cell-based & biochemical assays** - simple luminescence-based assays for
- **histone deacetylase (HDAC) & sirtuin (SIRT) activities**
  - HDAC-Glo™ Assays are used to determine histone deacetylase activity in cell-based or biochemical formats
  - flexible, sensitive bioluminescent assays provide a highly predictive, easy-to-use method for accurate and efficient inhibitor profiling.
- **DUB-Glo™ Protease Assay**
  - homogeneous, bioluminescent assay that measures the activity of numerous deconjugating enzymes including deubiquitinating (DUB), deSUMOylating (SEN) and deneddylating (NEDP) proteases
  - these proteases reverse the protein modification by ubiquitin and ubiquitin-like proteins (Ubl proteins) and thus are integral components in the complex mechanisms of post-translational protein regulation in eukaryotes.

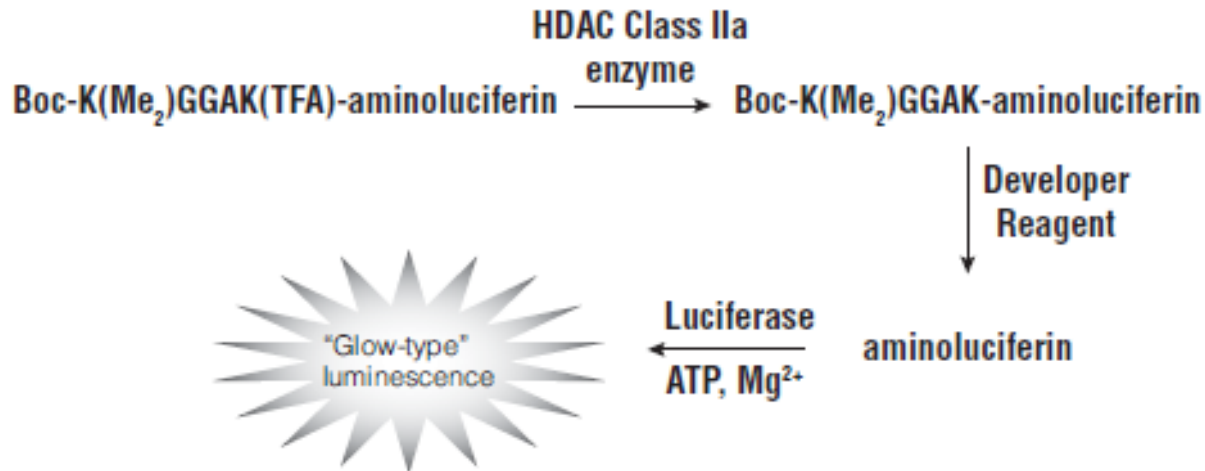


## ***New HDAC-Glo™ Class IIa and 2 selective assays***

- HDAC-Glo™ Class IIa and 2 selective assays launched 31/01/14
- single-reagent-addition, homogeneous, luminescent assays that measure the relative activity of histone deacetylase (HDAC) Class IIa and Class I enzyme 2, respectively, from cells, extracts or purified enzyme sources
- assays use an isozyme-selective acetylated, live-cell-permeant, luminogenic peptide substrate that can be deacetylated by HDAC activities
- deacetylation of the peptide aminoluciferin substrate is measured using a coupled enzymatic system in which a protease in the Developer Reagent cleaves the deacetylated peptide from the aminoluciferin substrate, releasing aminoluciferin
- aminoluciferin is quantified in a reaction using Ultra-Glo™ Recombinant Luciferase (firefly)
- signal from the assay reaction can be measured within 15–45 minutes after reagent addition with no sample manipulation
- HDAC-mediated luminescent signal is persistent, with a half-life of greater than 2 hours, allowing batch processing of multiwell plates.



## HDAC-Glo™ Class IIa - mechanism



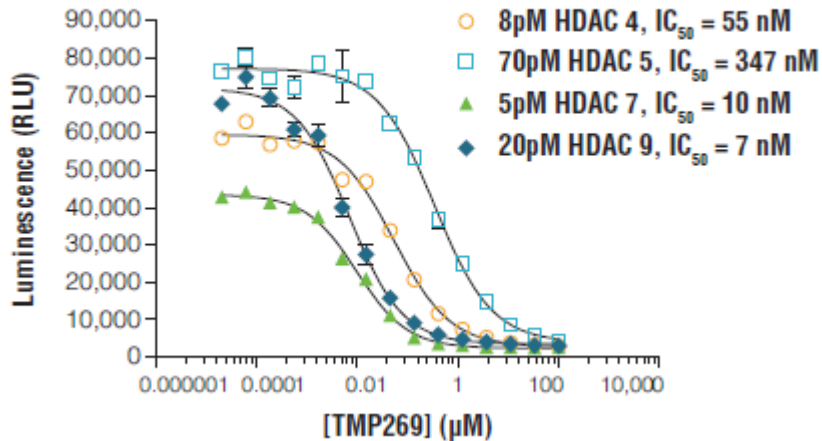
**Note:** The three enzymatic events occur in a coupled, nearly simultaneous reaction that is proportional to deacetylase activity.

### Assay Advantages

- **Simple Measurement of Deacetylating Activities:** Single-reagent-addition, homogeneous "add-mix-measure" protocol.
- **Sensitive:** The assay provides 100-fold or better sensitivity than comparable fluorescence methods.
- **Utility:** The assay may be used with recombinant enzyme sources or in a cell-based format.
- **Fast Data Acquisition:** Collect maximal signal in as few as 20 minutes with persistent, "glow-type" steady state signal half-life.

## HDAC-Glo Class IIa data

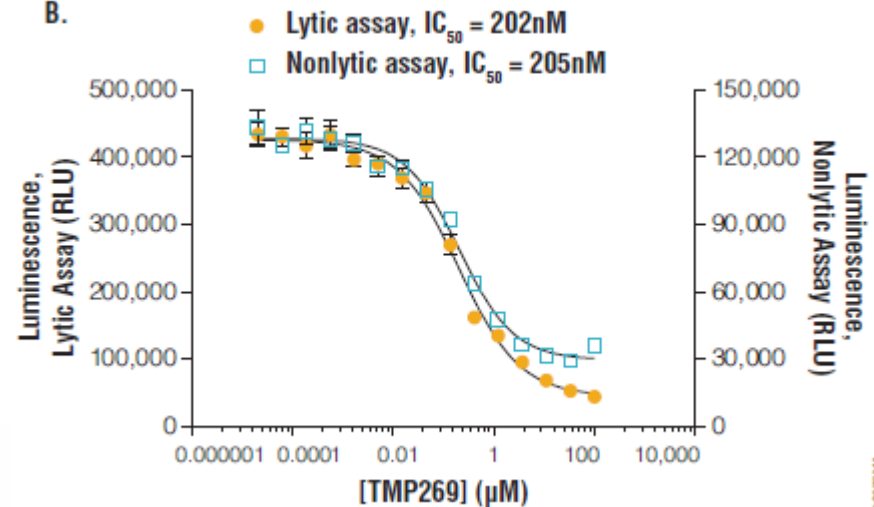
A.



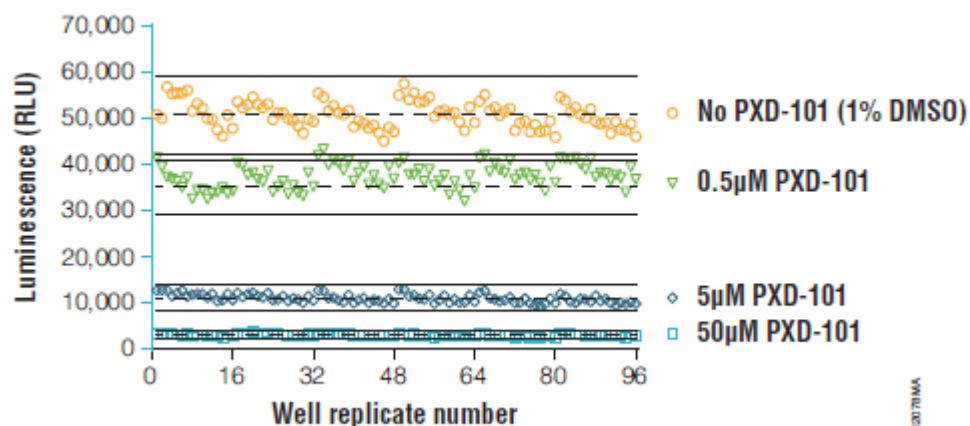
- cell-based assay format
- compare lytic/non-lytic assay
- using K562 cells as HDAC IIa source
- 2,500 cell/well (lytic)
- 10,000 cells/well (non-lytic)
- 384-well format
- assay volume 40μl

- biochemical assay format
- compare HDAC 4, 5, 7 and 9
- 384-well format
- assay volume 20μl

B.



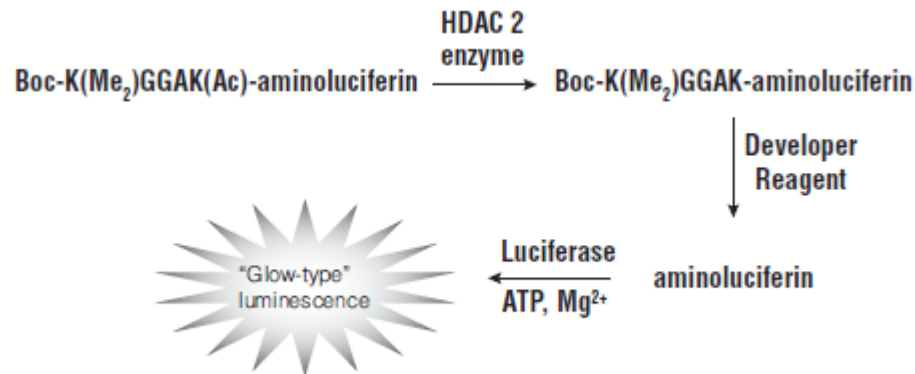
# *Z' analysis in 384-well format (HDAC-Glo™ IIa assay)*



Biochemical assay format using purified HDAC 7 & PXD-101 inhibitor

Compound Test Condition (n = 96)	Signal:Background	% CV	Z'
no PXD-101 (1% DMSO)	N.A.	5.6	N.A.
0.5µM PXD-101	1.4	5.5	0.10
5µM PXD-101	4.6	8.3	0.72
50µM PXD-101	16.7	8.7	0.81

## HDAC-Glo™ Class 2 - mechanism

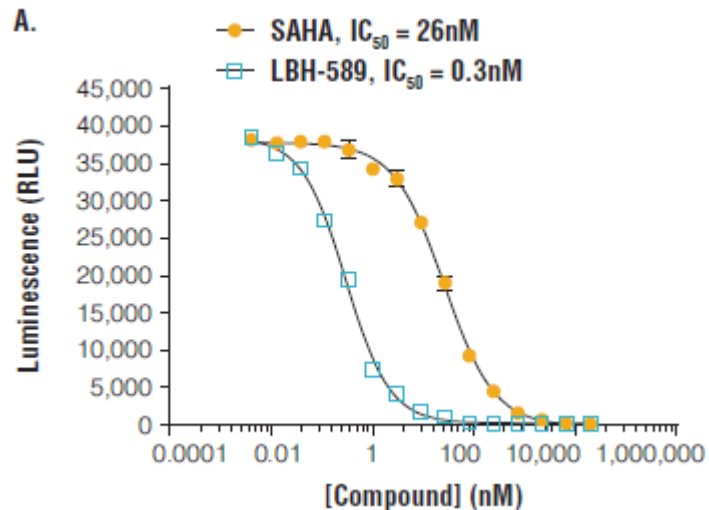


**Note:** The three enzymatic events occur in a coupled, nearly simultaneous reaction that is proportional to deacetylase activity.

### Assay Advantages

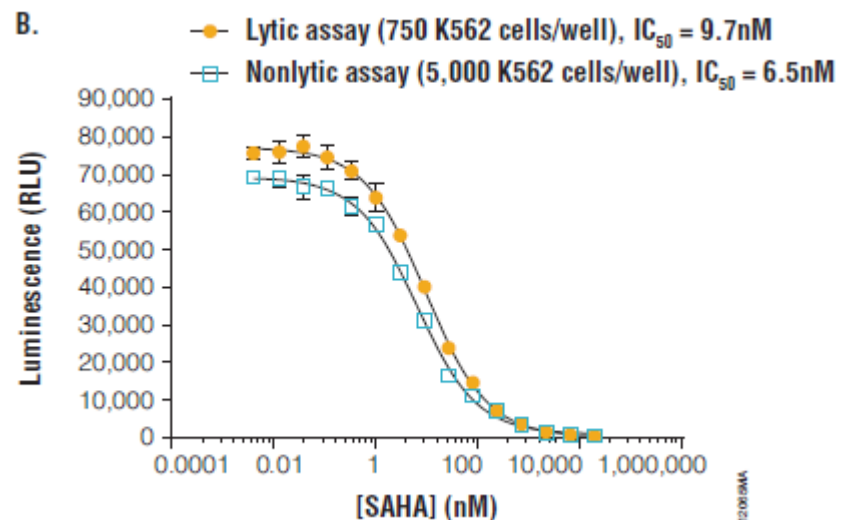
- **Simple Measurement of Deacetylating Activities:** Single-reagent-addition, homogeneous "add-mix-measure" protocol.
- **Sensitive:** The assay provides 100-fold or better sensitivity than comparable fluorescence methods.
- **Utility:** The assay may be used with recombinant enzyme sources or in a cell-based format.
- **Fast Data Acquisition:** Collect maximal signal in as few as 20 minutes with persistent, "glow-type" steady state signal half-life.

## HDAC-Glo Class 2 data

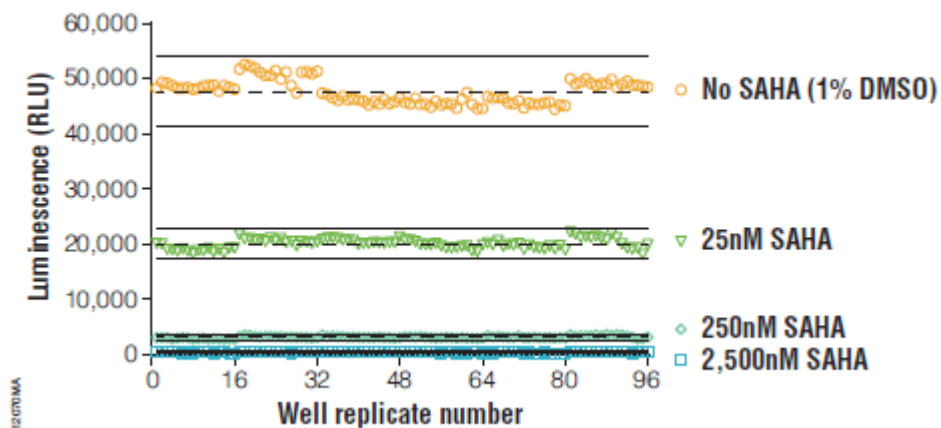


- cell-based assay format
- compare lytic/non-lytic assay
- using K562 cells as HDAC 2 source
- 750 cells/well (lytic)
- 5,000 cells/well (non-lytic)
- 384-well format
- assay volume 40 $\mu$ l

- biochemical assay format
- Recombinant HDAC 2
- Inhibitors SAHA & LBH-589
- 384-well format
- assay volume 20 $\mu$ l



## ***Z' analysis in 384-well format (HDAC-Glo™ 2 assay)***



Biochemical assay format using purified HDAC 2 & SAHA inhibitor

Compound Test Condition (n = 96)	Signal:Background	% CV	Z'
no SAHA (1% DMSO)	N.A.	4.5	NA
25nM SAHA	2.4	4.4	0.67
250nM SAHA	15.1	6.1	0.84
2,500nM SAHA	125.0	11.1	0.86



***Tools for biologics – ADCC reporter assays and other cell-based assays for mAb characterisation***



## 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

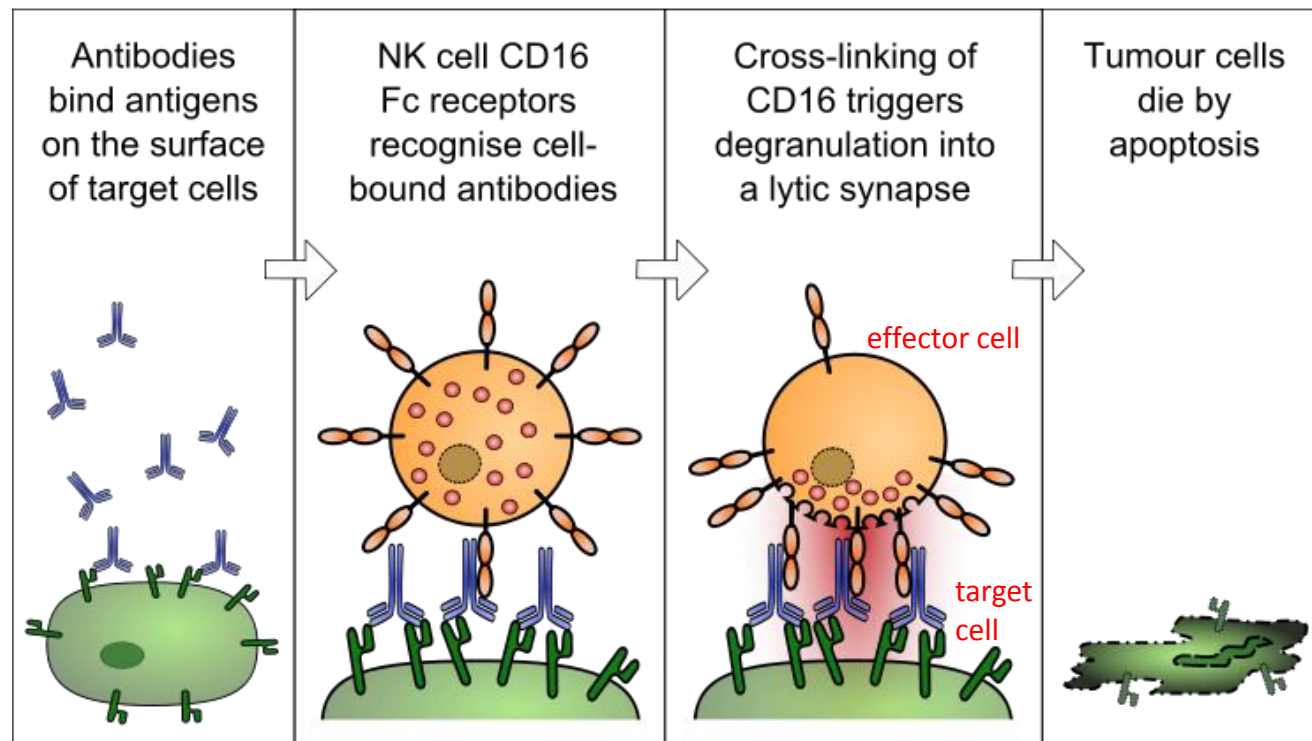


Image source: Wikipedia

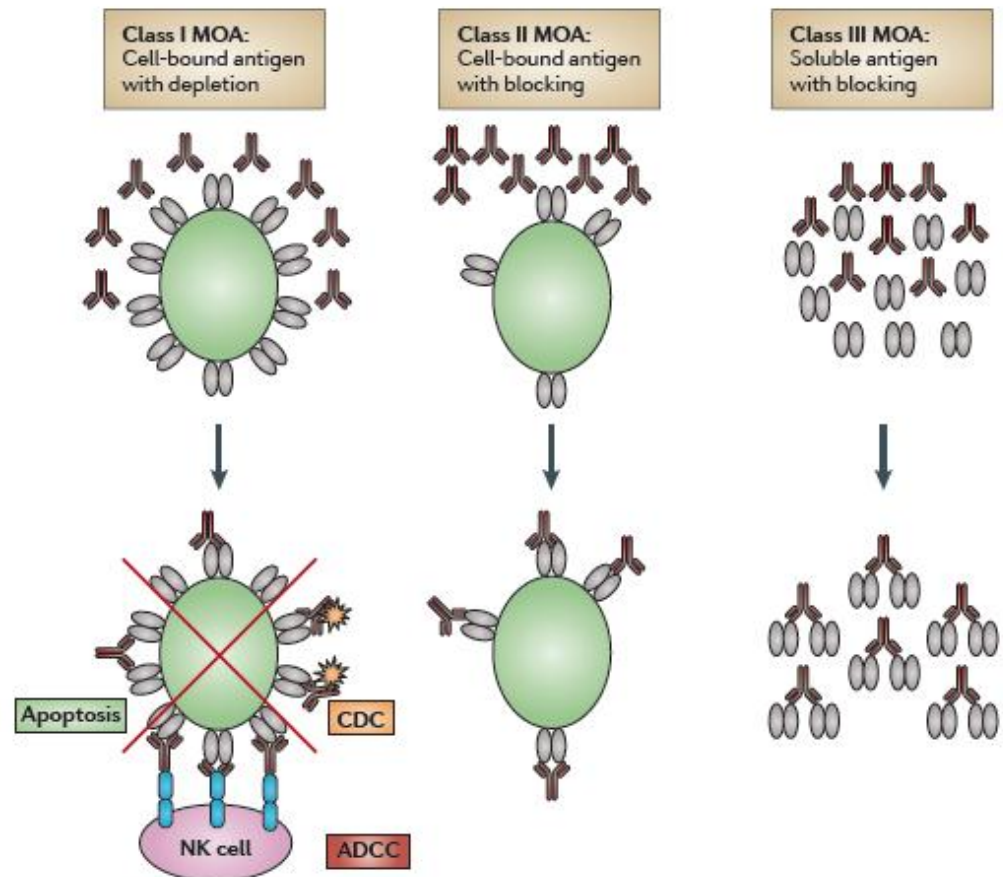
# Mechanism of action (MOA) for mAb - definitions

## Classification of therapeutic antibodies based on their putative mechanisms of action

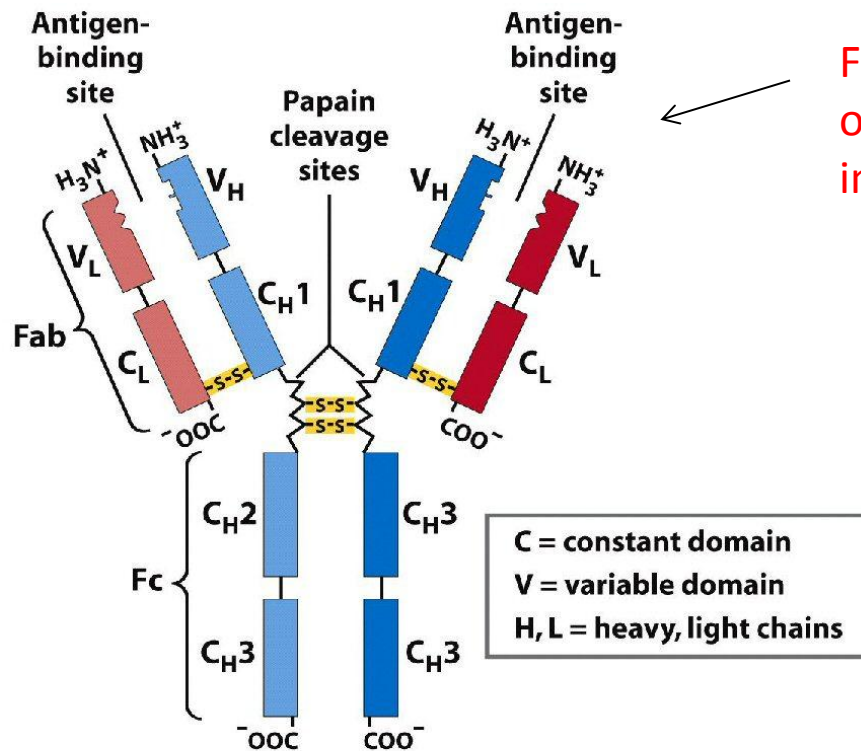
**Class I:** mAbs bind cell surface antigens and the Fc effector function is part of the MOA

**Class II:** mAbs bind cell surface antigens but MOA does not involve Fc effector function

**Class III:** mAbs bind and neutralize soluble antigens, blocking ligand from binding to its cognate receptor (e.g., Avastin)



# Antibody structure



Fab domain - binds to antigen expressed on target cells *eg.* cancer cells, virus infected cells

Fc domain – binds to Fc receptor (*eg.* FcγRIIR, CD16) on immune system effector cells

Figure 5-21a  
Lehninger Principles of Biochemistry, Fifth Edition  
© 2008 W. H. Freeman and Company

# Traditional ADCC Assays

## Effector cells

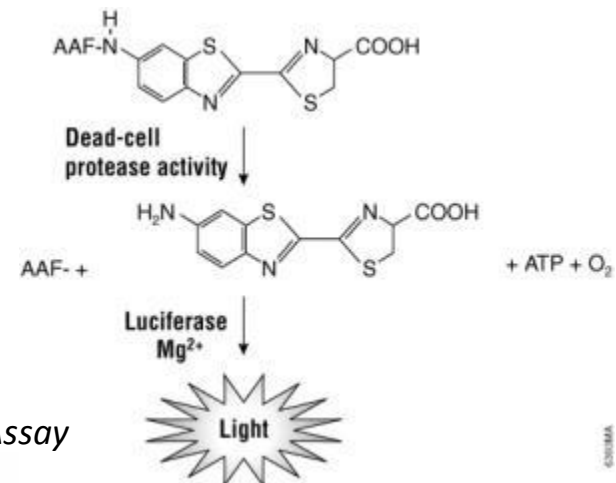
- PBMCs (peripheral blood mononuclear cells) ← difficult to isolate & purify, variable yields & efficacy
- Natural killers cells (NK) from PBMCs
- NK cell lines (eg. NK92) ← stability of Fc receptor expression an issue

## Target cells

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

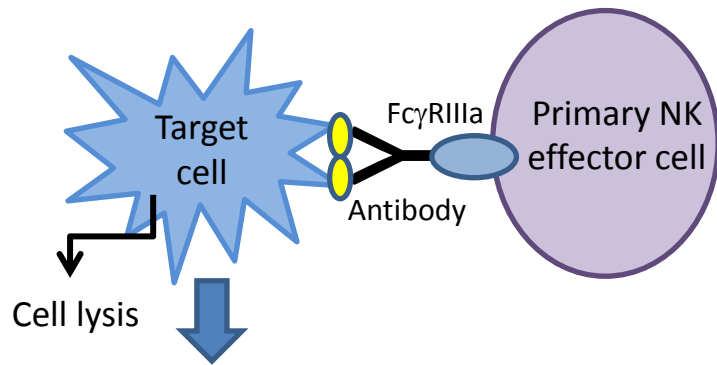
additional assay step

CytoTox-Glo™ Cytotoxicity Assay



# Classic ADCC assay vs ADCC reporter bioassay

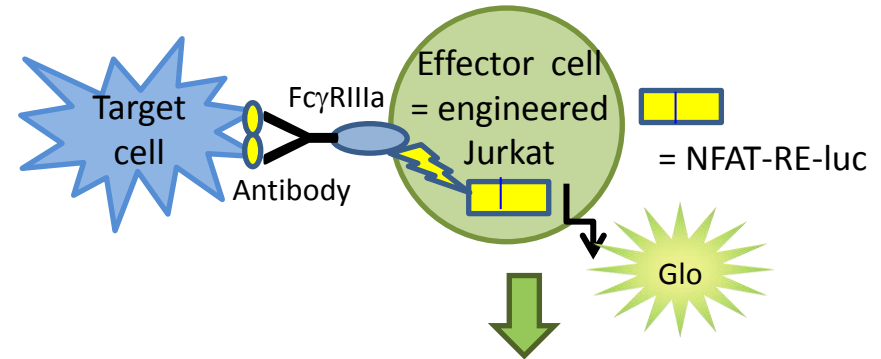
## Classic ADCC assay



**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



## Scientific basis of ADCC reporter bioassay

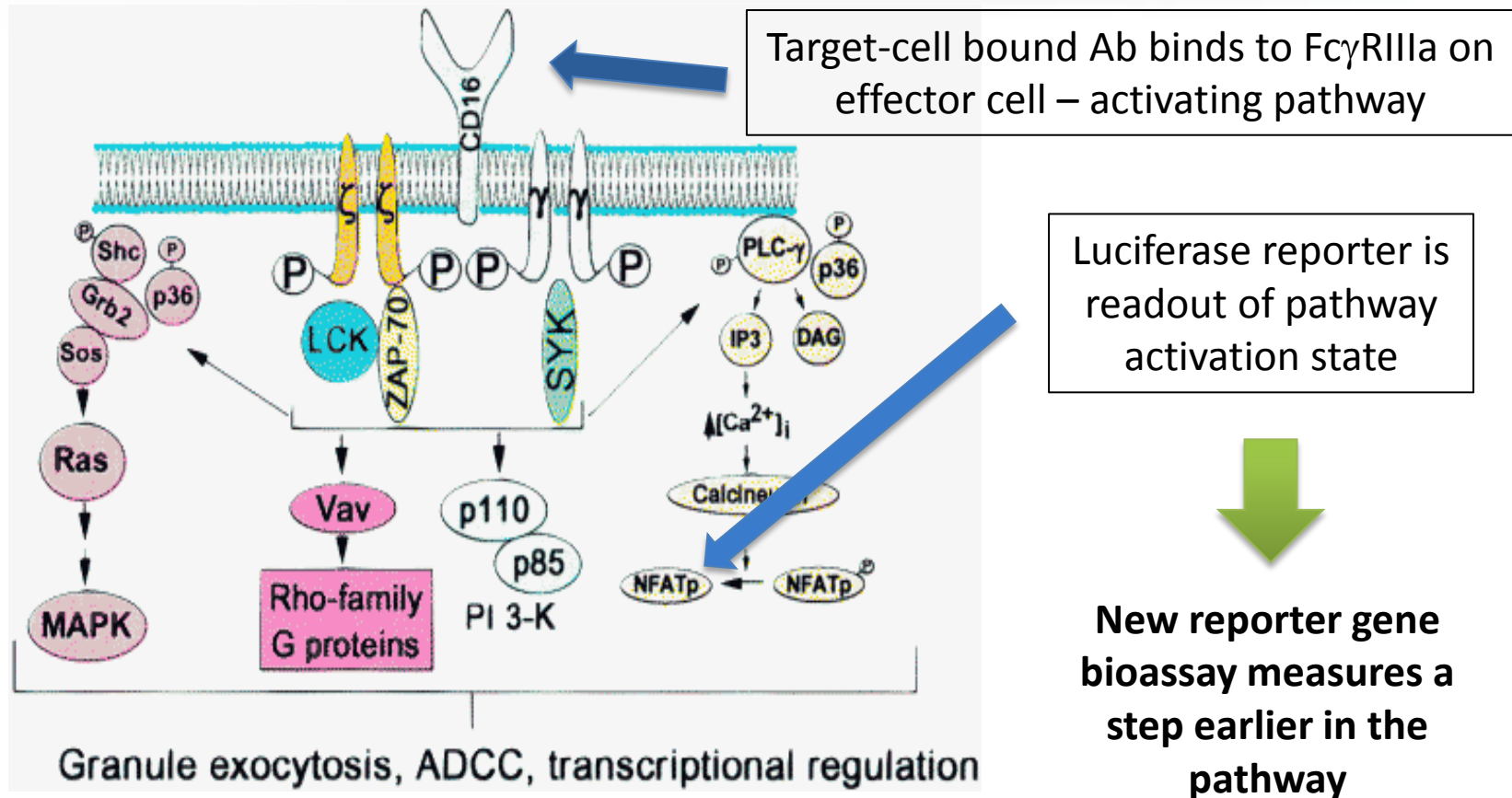
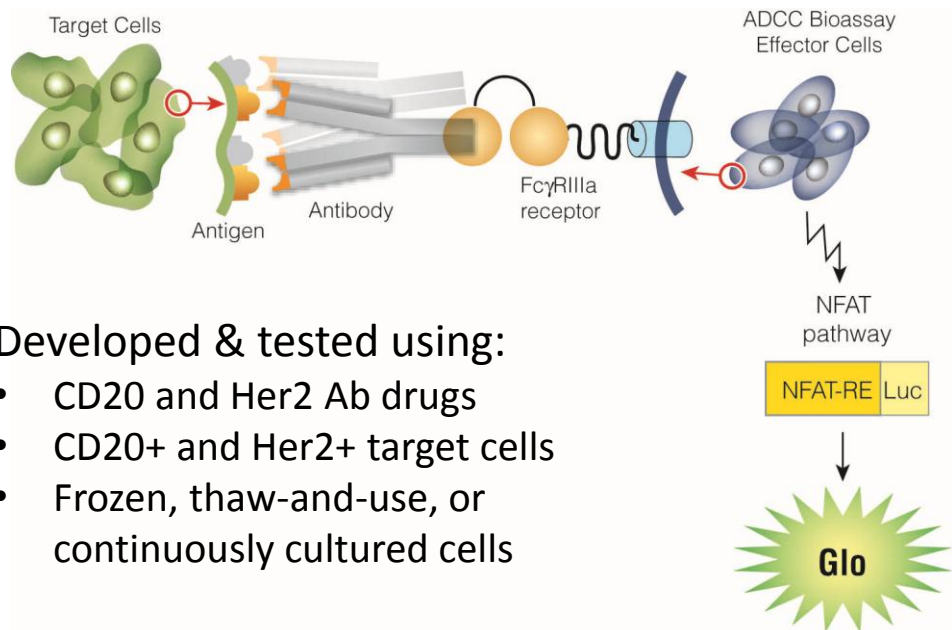


Image source: Leibson-PJ, *Immunity* 1997

# 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

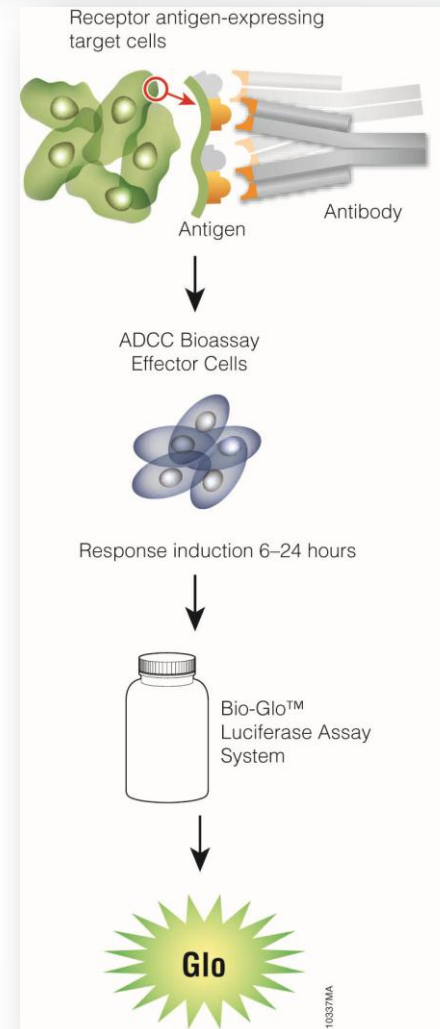
1. Effector cells are engineered to express FcγRIIIa (V158) and NFAT-RE luc2 luciferase
2. 'Cells as reagents' (thaw-and-use)
3. Homogeneous assay format – simple 'add-mix-read' bioluminescent assay
4. Optimized and robust assay reagents and protocol
5. Performance characteristics that meet needs of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



# ADCC reporter bioassay protocol

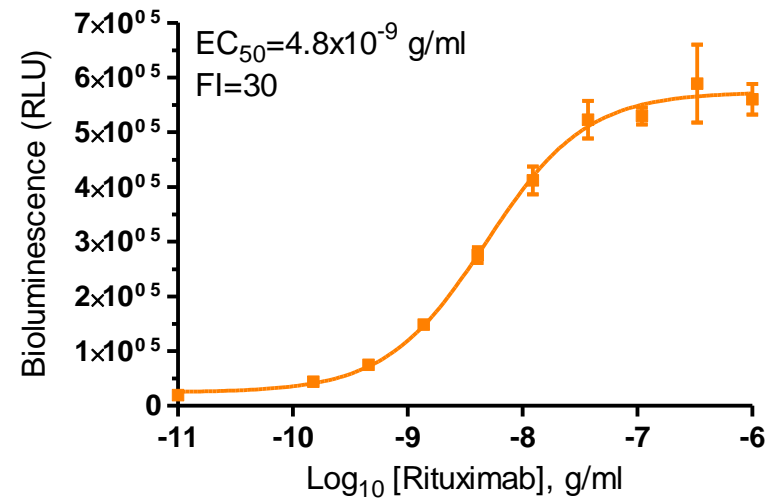
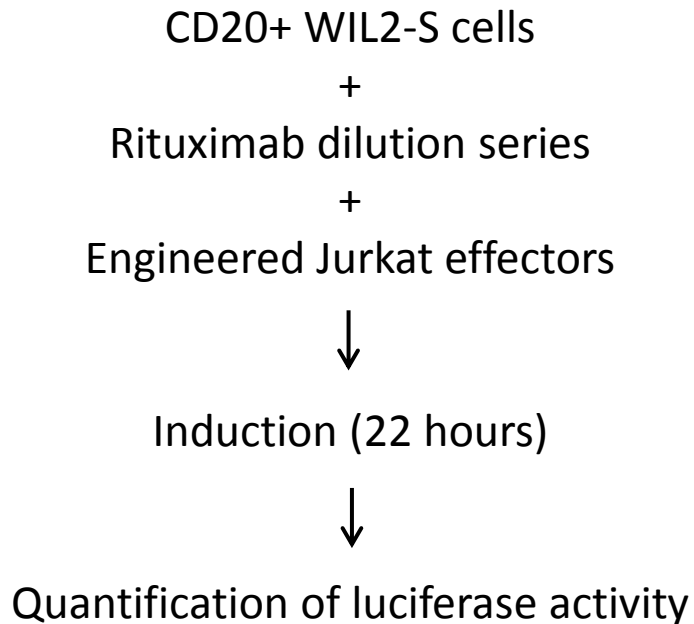
## *Single day bioassay – no cell preparation!*

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 – typical results

## Assay protocol:



## Specifics

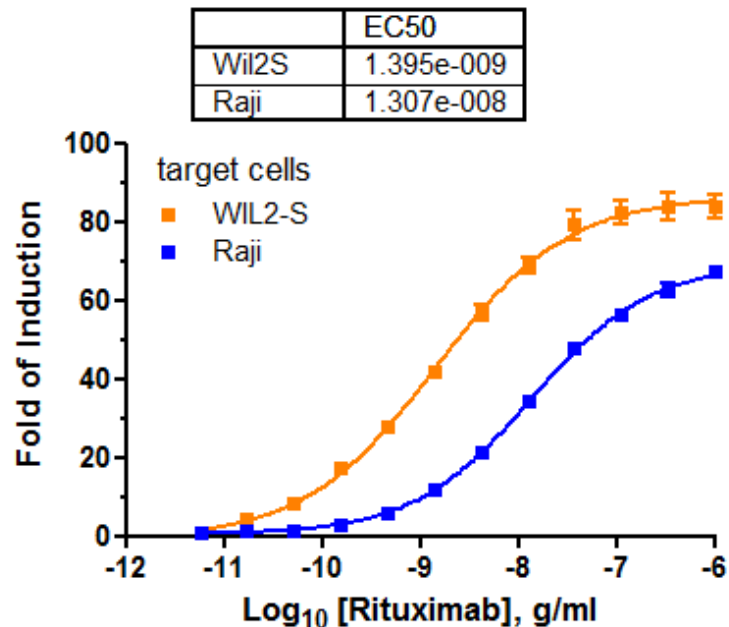
- **E:T ratio = 6:1**  
(150k effector cells:25k WIL2-S target cells, per well)

## Use of different target cells

*Suspension or adherent target cells can be used*

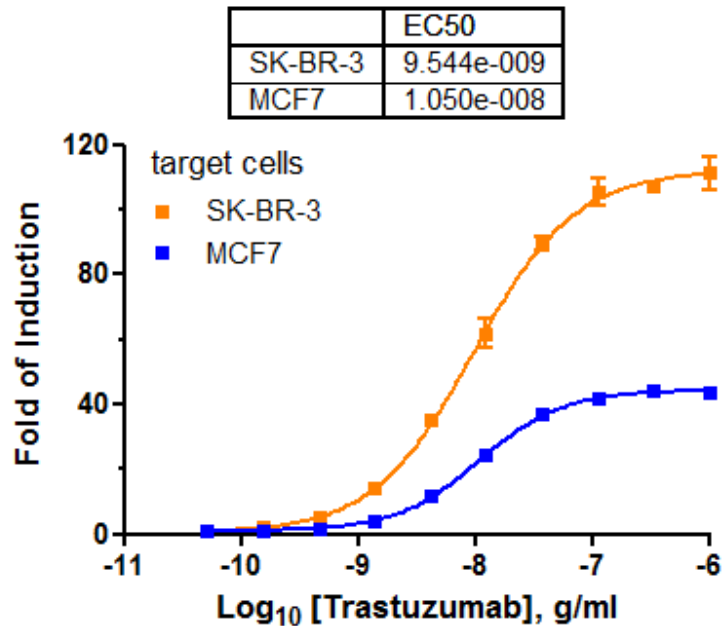
### Rituximab (anti-CD20)

**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**



## *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

# ***ADCC reporter bioassay is specific and reflective of the mechanism of action of the biologic material***

Target cells, effector cells and specific antibody

■ Wil2-S, Jurkat/NFAT-luc+FcγRIIIa, Rituximab

No Target cells

● NO Wil2-S, Jurkat/NFAT-luc+FcγRIIIa, Rituximab

No Effector cells or no FcγRIIIa

▲ Wil2-S, Jurkat/NFAT-luc (NO FcγRIIIa), Rituximab

▲ Wil2-S, NO Jurkat/NFAT-luc+FcγRIIIa, Rituximab

No antibody or non-specific antibody

▼ Wil2-S, Jurkat/NFAT-luc+FcγRIIIa, NO Rituximab

▼ Wil2-S, Jurkat/NFAT-luc+FcγRIIIa, Trastuzumab

**Assay signal is dependent on:**

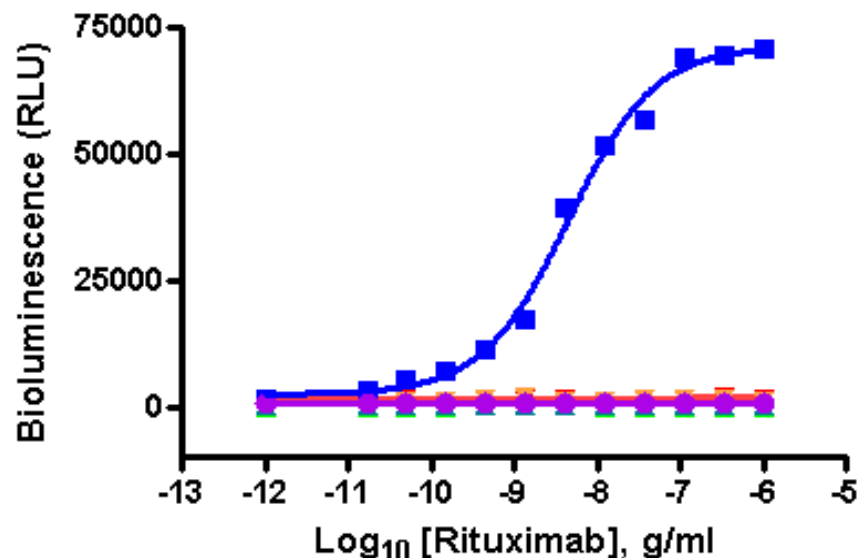
Presence of Target cells

+

Presence of FcγRIIIa receptor

+

Appropriate specific antibody



**Reflects correct MOA**

# Bioassay characteristics - ICH guideline Q2 [R1]

## Validation of Analytical Procedures:

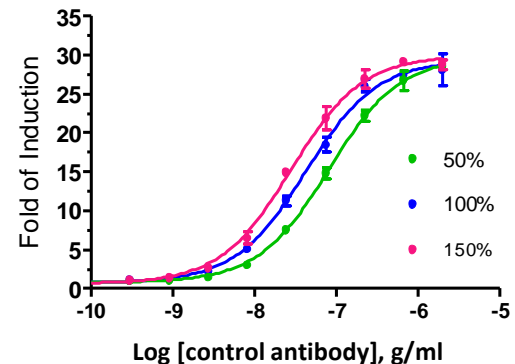
- Accuracy
- Precision:
  - ✓ Repeatability (intra-assay precision)
  - ✓ Intermediate precision (day to day, analyst-to analyst)
  - ✓ Reproducibility (lab to lab)
- Specificity
- Linearity
- Range
- Robustness

**Assay is well controlled  
(precise, accurate, robust,  
reproducible)**

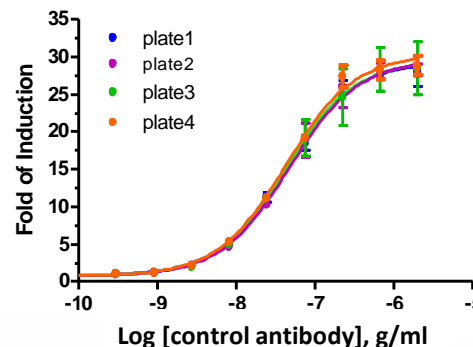
### Design:

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

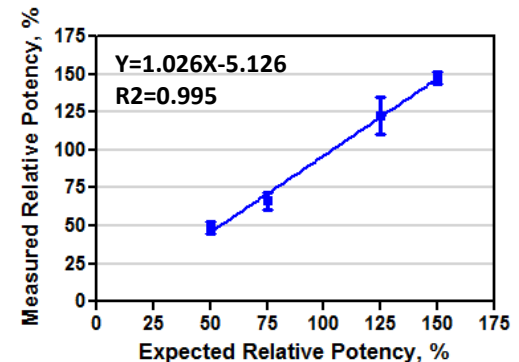
## Relative potency



## Repeatability

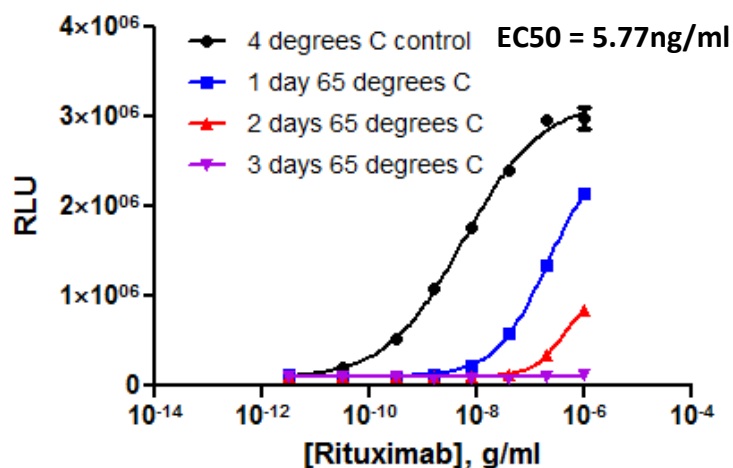


## Linearity

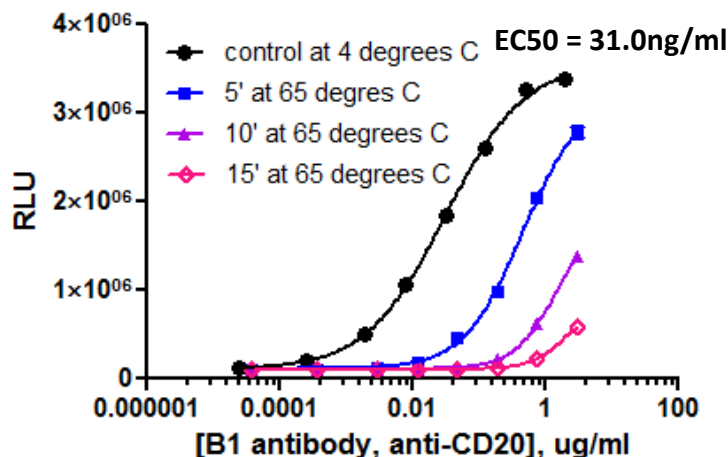


# Stability indicating for Fc effector function – activity of heat-treated antibody drugs

## Rituximab

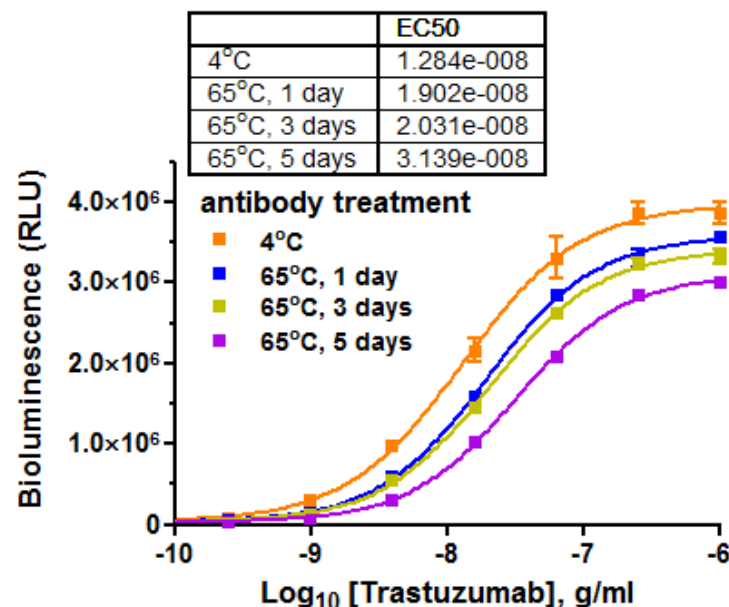


## Tositumomab



## Activity of heat-treated antibody drugs

## Trastuzumab



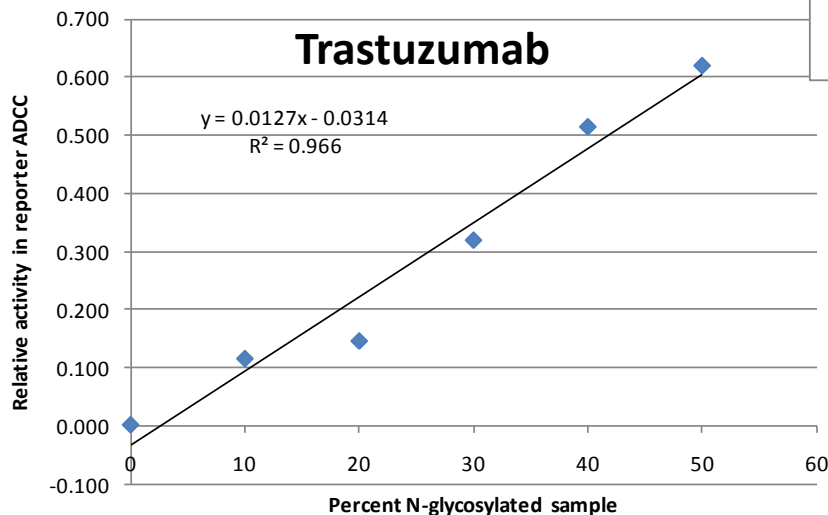
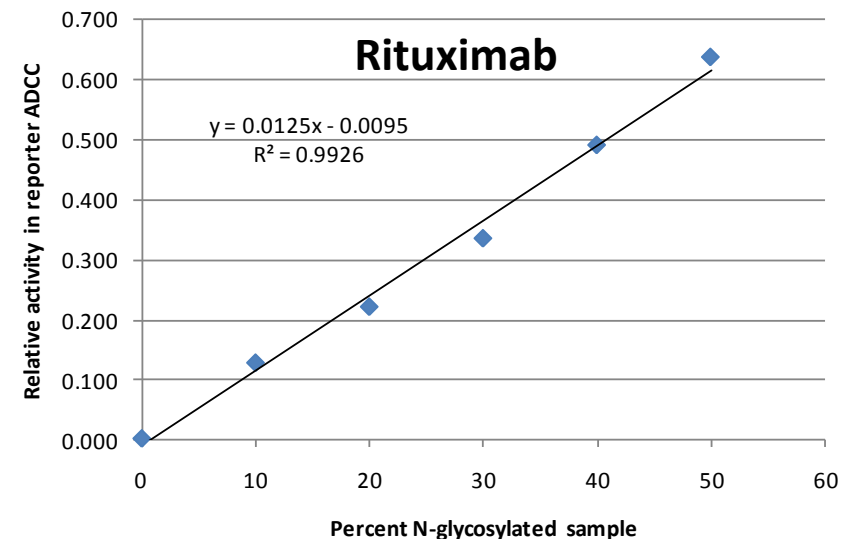
Assay is stability-indicating & usable as a QC lot-release assay



# 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

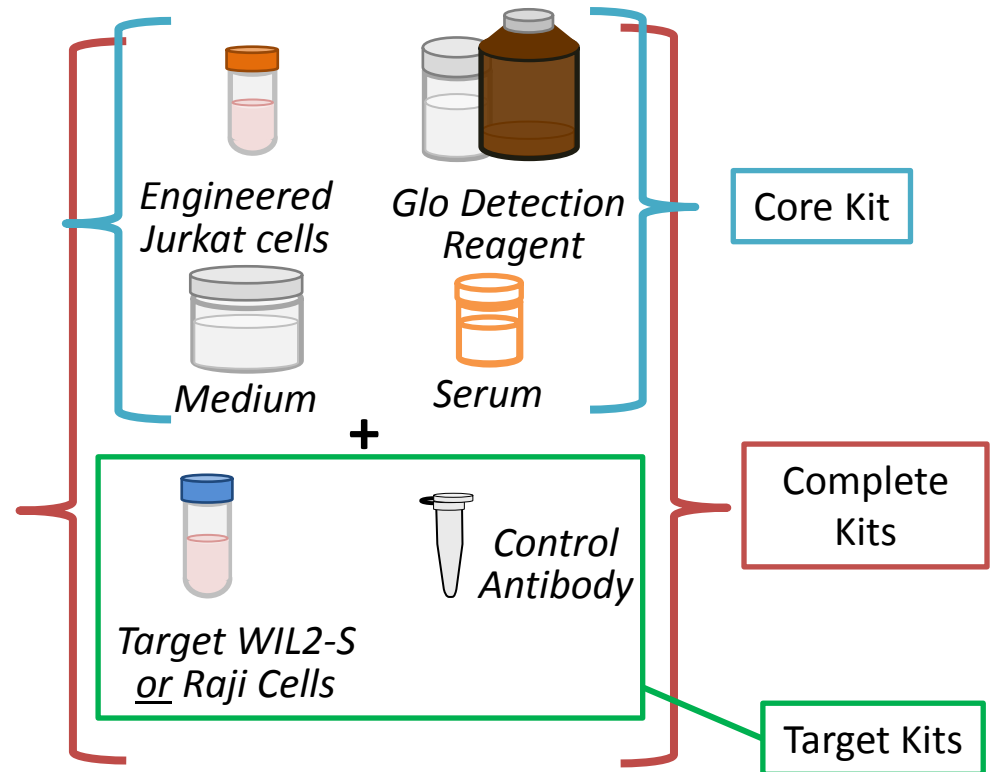


*Small differences in Fc effector activity in ADCC pathway activation are easily distinguished in the ADCC reporter bioassay*

# ADCC reporter 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:  
Includes control antibody and target cells plus effector/target cells and reagents
3. Target Kits: Control antibody and target cells



**Note:** the ADCC Bioassay Effector Cells alone are also available for propagation and banking under a license agreement

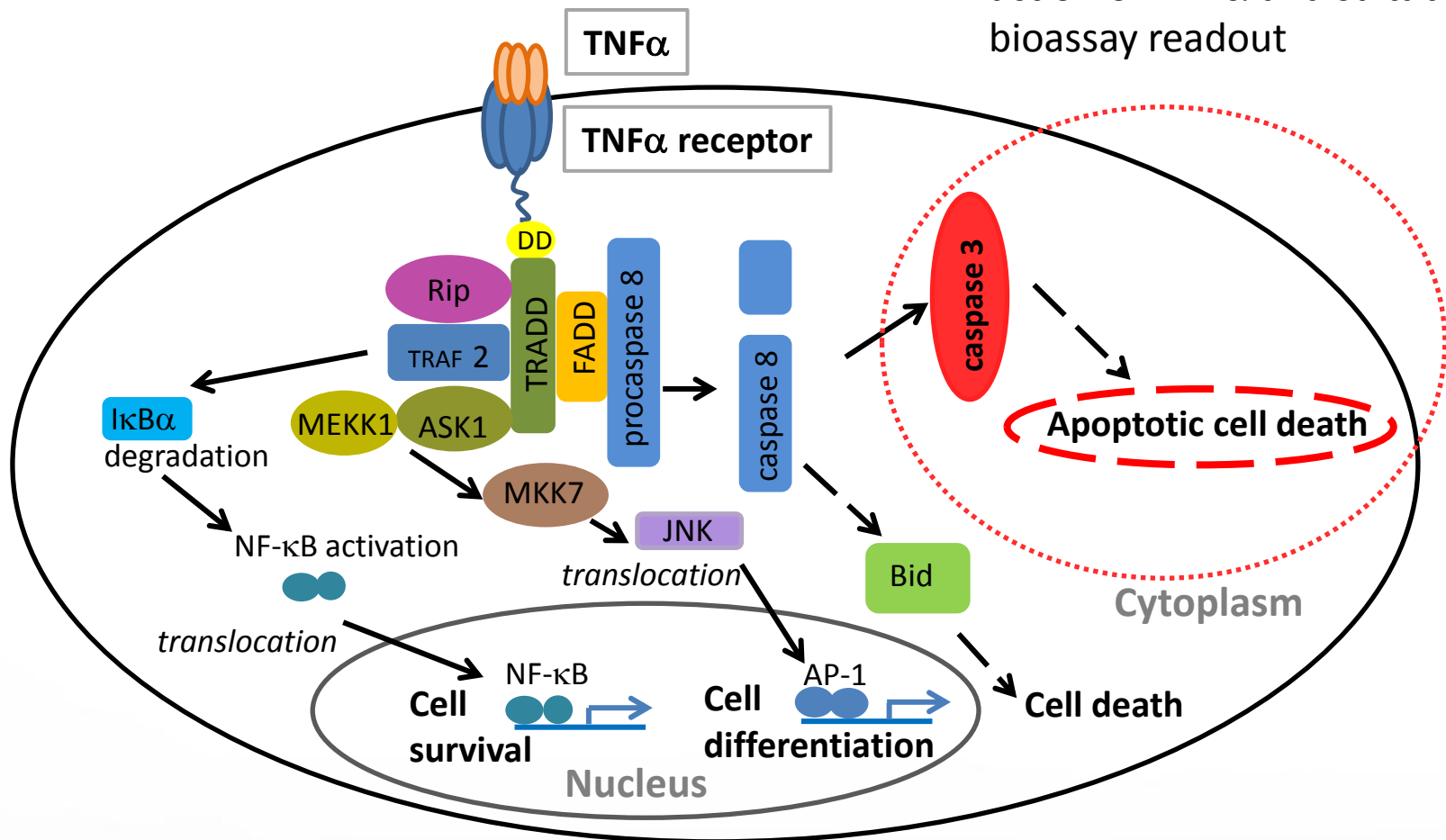
## ***Other cell-based assays for therapeutic mAb characterisation***

<b>Assay</b>	<b>Cell type</b>	<b>Antibodies</b>	<b>Readout</b>
TNF $\alpha$ blocker assay	U937 cells	Adalimumab, Infliximab	Caspase 3/7 activation
VEGF blocker assay	HEK 293	Avastin	NFAT/ <i>luc2P</i> reporter assay
CTLA-4 receptor blocker assay	Jurkat	Yervoy	IL2 promoter/ <i>luc2P</i> reporter assay

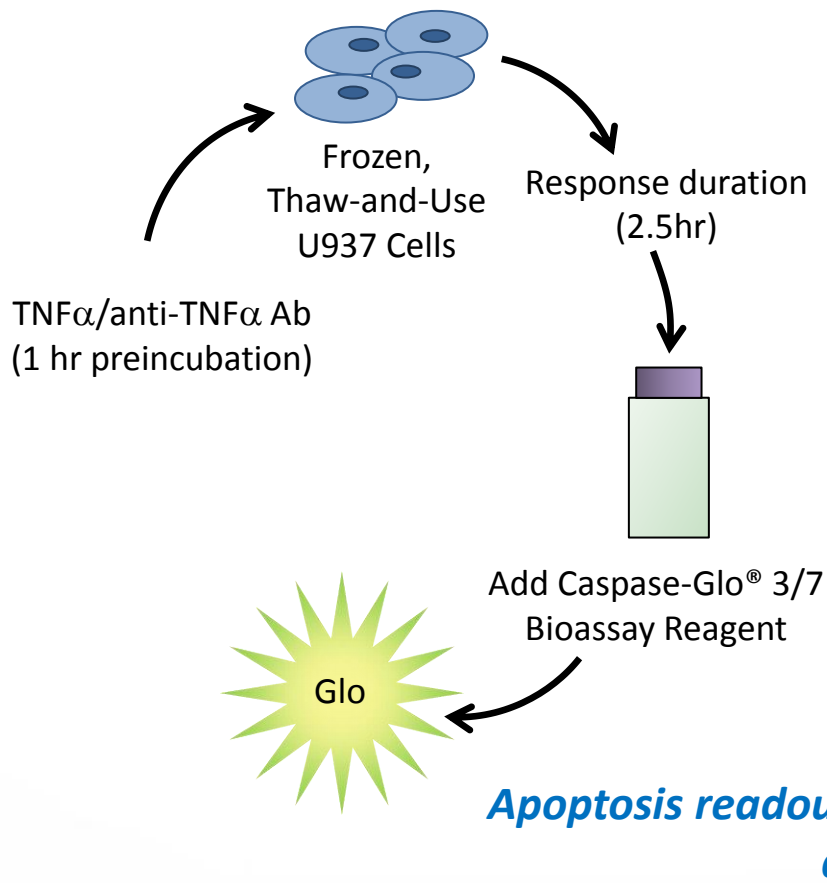
These are just examples – more assays for biologics are being added all the time....

# *TNF $\alpha$ blocker assay*

Apoptosis is a mechanism of action of TNF $\alpha$  and suitable as a bioassay readout



## ***Bioluminescent TNF $\alpha$ blocker apoptosis bioassay***

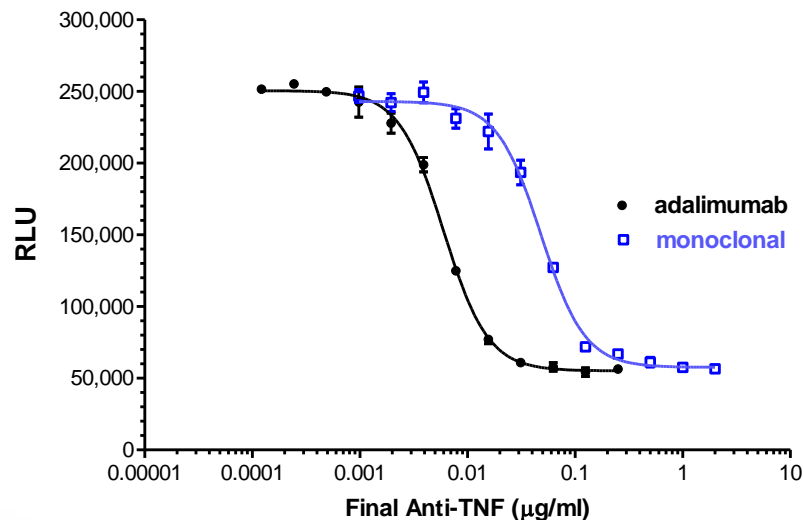


### **Protocol:**

1. TNF $\alpha$  and anti-TNF $\alpha$  Ab are co-incubated; the TNF $\alpha$  and anti-TNF $\alpha$  mix is added to thaw-and use U937 cells in assay medium.
2. Response is induced in as little as 2.5hr.
3. Caspase-Glo<sup>®</sup> 3/7 Assay is added and luminescence is measured immediately.

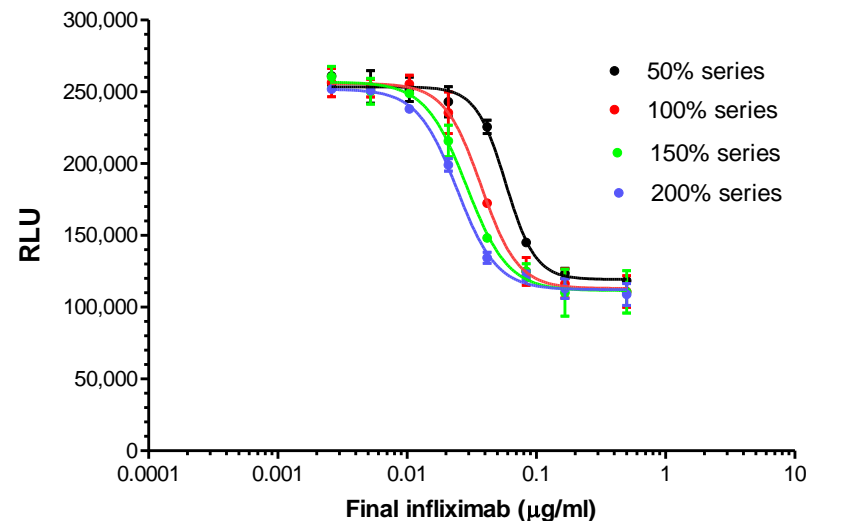
# ***TNF $\alpha$ blocker apoptosis bioassay can be used to quantify activity of on-market TNF $\alpha$ blocker drugs***

**A. Fully human **adalimumab**** demonstrates significant improved potency in TNF $\alpha$  blocking biological activity compared with a mouse mAb.



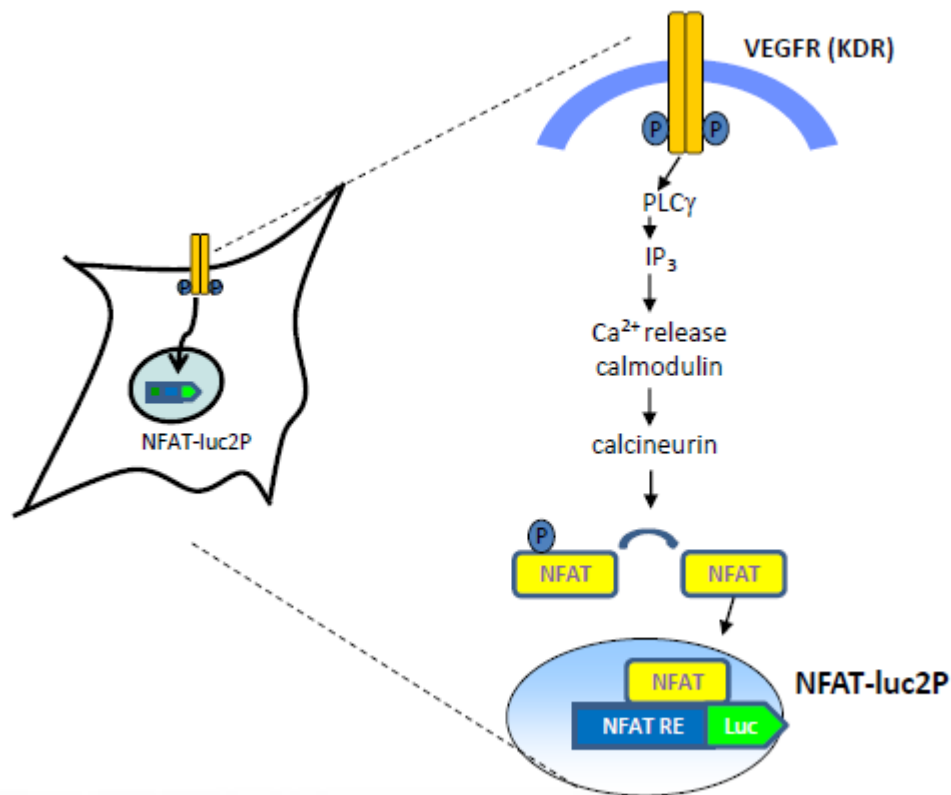
	adalimumab	monoclonal
EC50	0.005997	0.04813

**B. Good potency discrimination is obtained using a series of **infliximab** preparations of different theoretical potencies.**



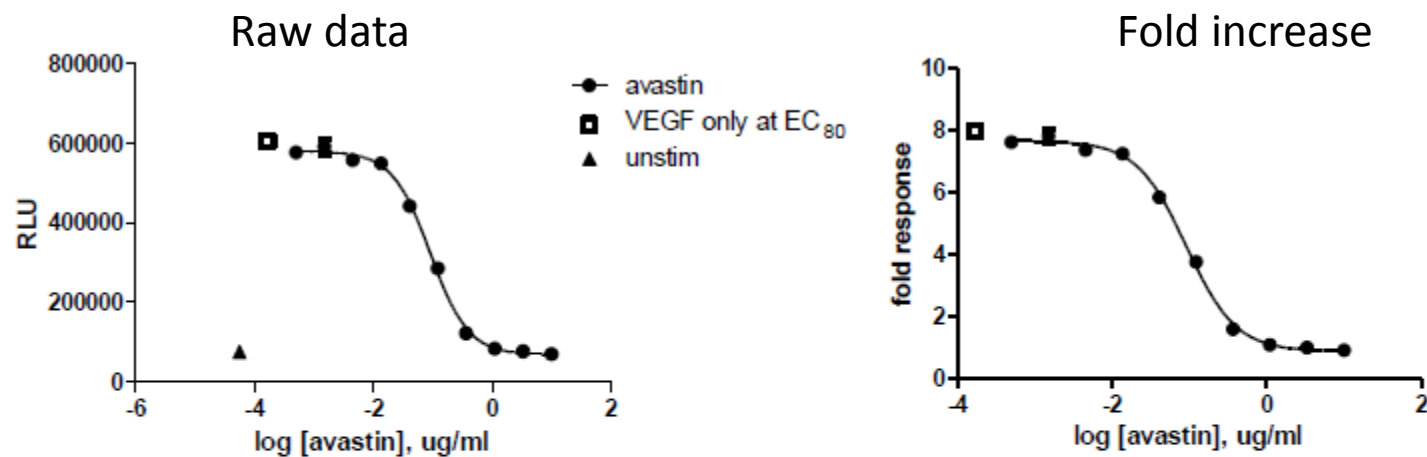
	50% series	100% series	150% series	200% series
EC50	0.05798	0.03743	0.02890	0.02437

## *VEGF blocker assay principle*





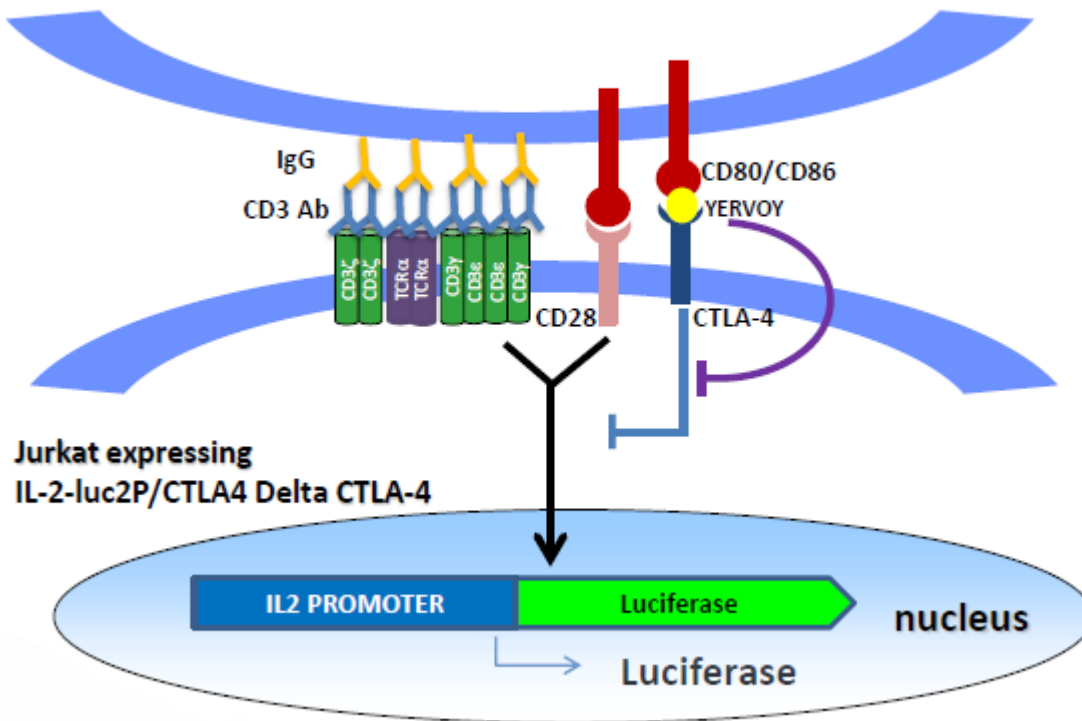
# Avastin titration



Incubate cells with Avastin titration for 30 minutes prior to adding VEGF at EC<sub>80</sub> concentration (14 ng/ml)  
 IC<sub>50</sub> of Avastin = 92 ng/ml

## Yervoy/CTLA-4 assay principle

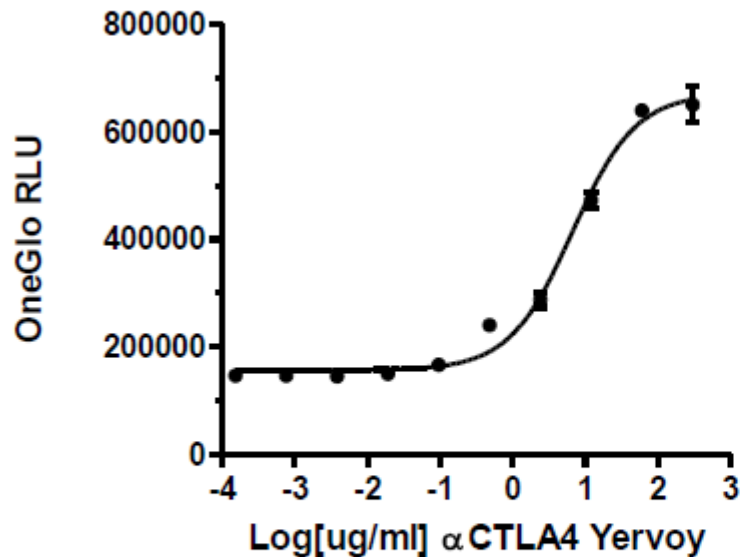
### Antigen Presenting Cell Raji



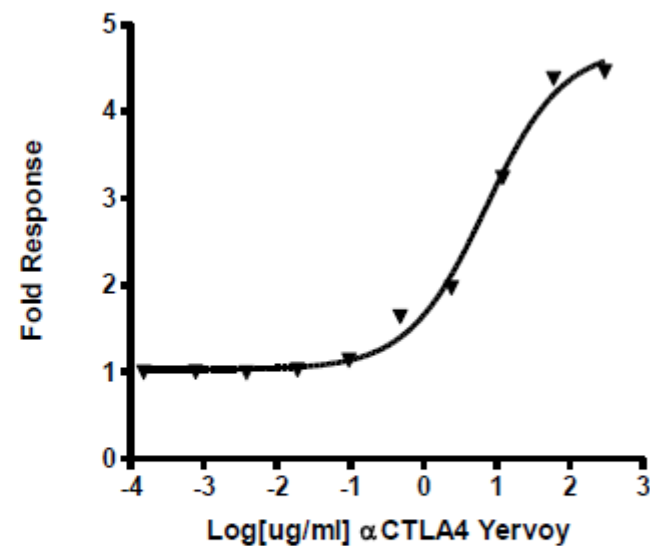
- Activation of T cells and cytokine release such as IL2 production require CD3 antibody induced TCR activation and CD80/CD86 co-stimulatory signals to CD28.
- CTLA-4 receptor expression sequester ligands away from CD28, thus preventing co-stimulatory signals from being received.
- Yervoy Antibody blockade of CTLA-4/CD80/CD86 interactions promote T cell activation

# *Jurkat/IL2-luc2P/CTLA-4 Delta stable cell line performance*

Raw data



Fold increase

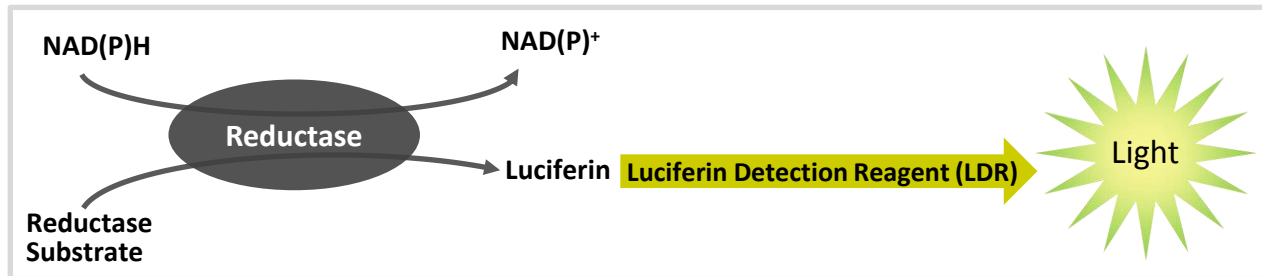


Ec50 = 6.64ug/mL

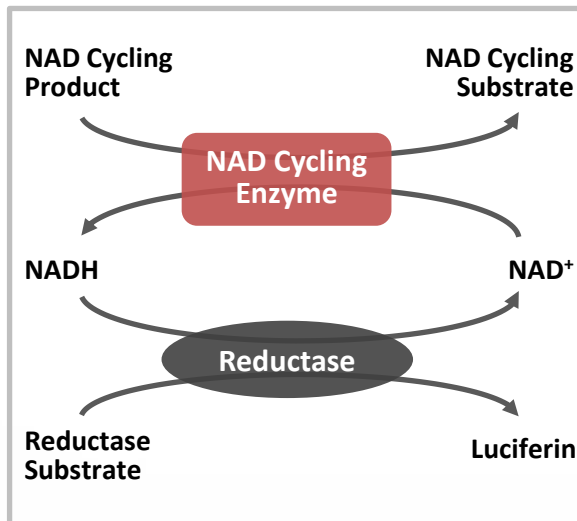
***New assays for metabolic markers***

# *A novel Proluciferin Substrate was Combined with Specific Cycling Enzymes to Develop Three Assays*

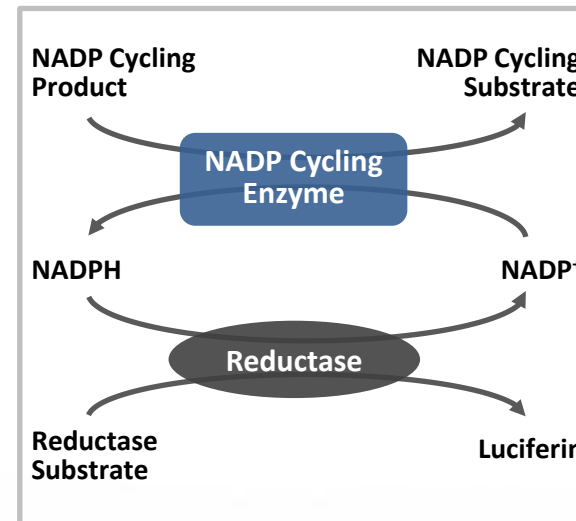
**NAD(P)H-Glo™** Detection System- detects reduced forms **NADH** and **NADPH**



**NAD/NADH-Glo™ Assay** – detects non-phosphorylated forms



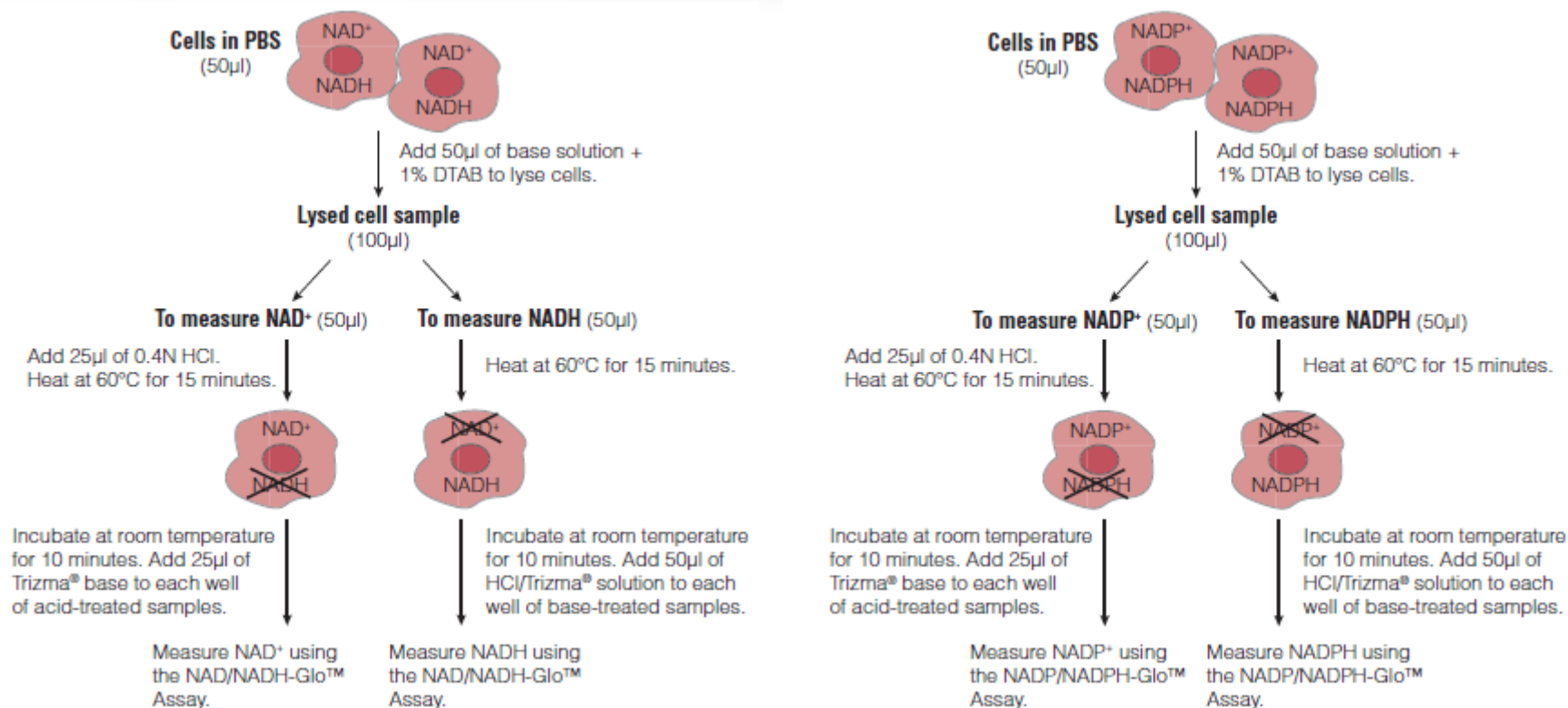
**NADP/NADPH-Glo™ Assay** – detects phosphorylated forms



## ***Features of the assays: sensitivity , specificity and wide assay window***

	NAD(P)H-Glo Detection system	NAD/NADH-Glo Detection system	NADP/NADPH-Glo Detection system
<b>Limit of Detection (LOD)</b>	25nM (625fmol/25μl)	1nM (25fmol/25μl)	1nM (25fmol/25μl)
<b>Linearity</b>	25nM – 50μM	1 – 500nM	1 – 500nM
<b>Signal-to-background (S/B max)</b>	~400	~250	~250
<b>Cells/well for Total dinucleotides</b>	NA	500 -25,000	500 – 12,000
<b>Cells/well for Individual dinucleotides</b>	NA	2,000 -100,000	2,000 – 100,000

# *Sample preparation for measuring NAD<sup>+</sup> /NADH or NADP/NADPH individually*



Both assays employ an sample acidification or heat-treatment (60°C) step to selectively destroy a specific metabolite

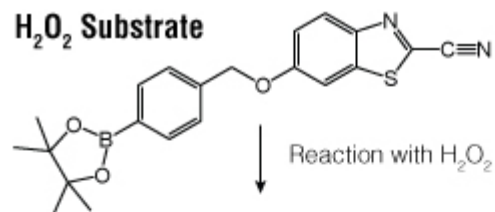


## ***ROS-Glo™ $H_2O_2$ Assay***

## ***ROS-Glo™ H<sub>2</sub>O<sub>2</sub> assay***

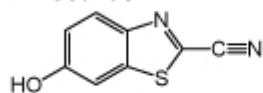
- ROS-Glo™ H<sub>2</sub>O<sub>2</sub> Assay is a homogeneous, fast and sensitive bioluminescent assay that measures the level of hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>), a reactive oxygen species (ROS), directly in cell culture or in defined enzyme reactions
- A derivatised luciferin substrate is incubated with sample and reacts directly with H<sub>2</sub>O<sub>2</sub> to generate a luciferin precursor
- Addition of ROS-Glo™ Detection Solution converts the precursor to luciferin and provides Ultra-Glo™ Recombinant Luciferase to produce light signal that is proportional to the level of H<sub>2</sub>O<sub>2</sub> present in the sample.

# **ROS-Glo™ mechanism & protocol**



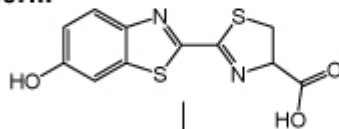
Reaction with H<sub>2</sub>O<sub>2</sub>

**Luciferin Precursor**



ROS-Glo™ Detection Solution

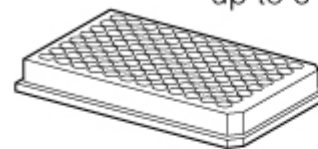
**Luciferin**



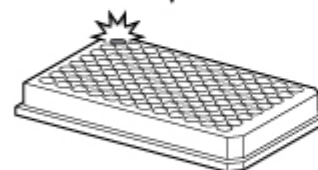
**Light**

11359MA

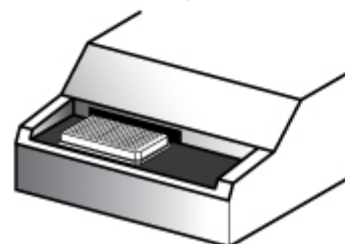
Treat samples.  
Add H<sub>2</sub>O<sub>2</sub> Substrate  
Solution. Incubate for  
up to 6 hours.



Add ROS-Glo™ Detection  
Solution. Incubate for  
20 minutes.

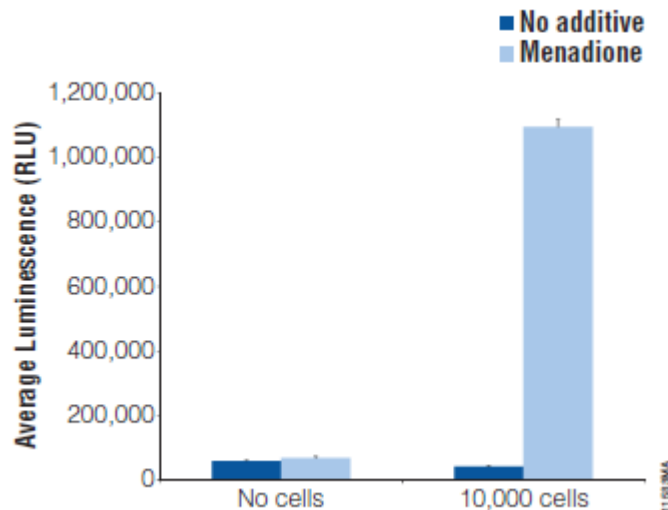


Read luminescence.



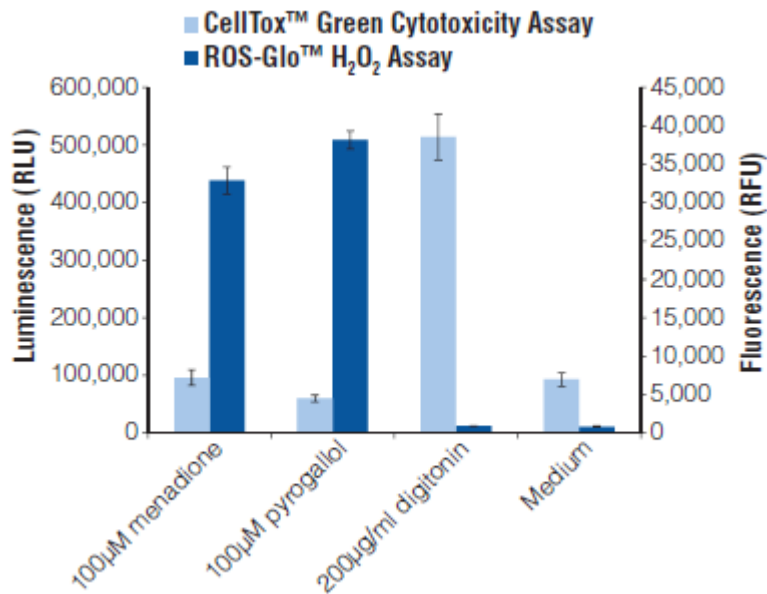
11360MA

## ***ROS-Glo™ assay data***



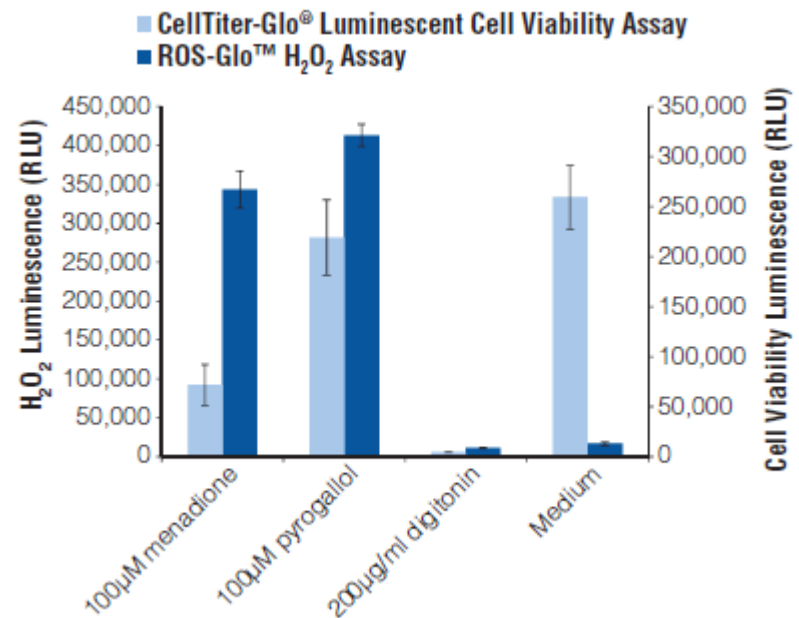
- Hep G2 cells
- 96-well plate format
- 20 $\mu$ M menadione
- 10,000 cells/well
- 2 hour incubation with menadione
- Add ROS-Glo™ Detection Solution
- Luminescence read after 20 minutes

## ROS-Glo™ - multiplexing



Multiplexing a fluorescent cytotoxicity readout with luminescence ROS-Glo™ to obtain both readings from the same population of cells

(Hep G2 plated at 2,000 cells/well in 384-well plate)



After 2 hours compound treatment, ROS-Glo™ assays are conducted on media aliquoted into a separate plate followed by addition of CellTiter-Glo viability reagent to original plate.

(Hep G2 plated at 2,000 cells/well in 384-well plate)

***Questions??***