

# ***Cell Health and Mechanistic Toxicity Assays***

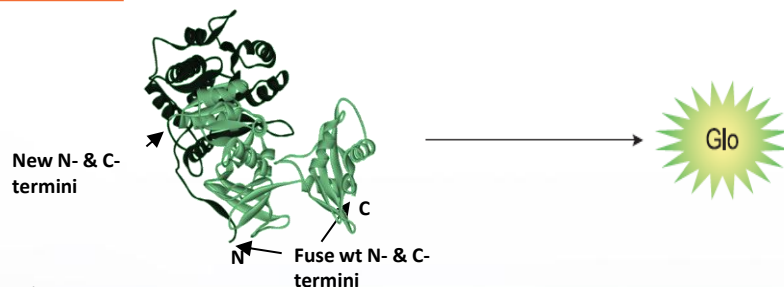
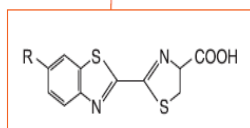
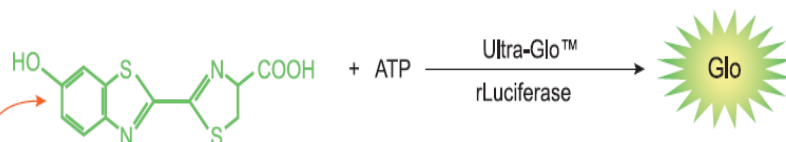
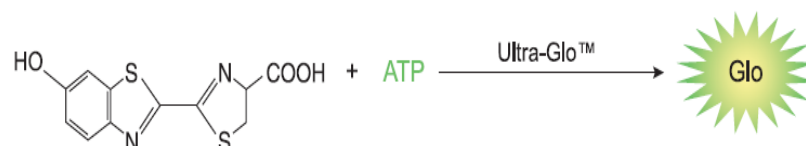
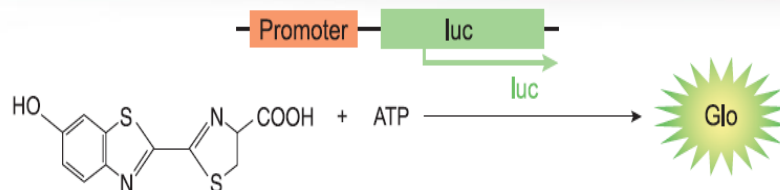
**Frank Fan, Ph.D.  
Director of Research**

**March, 2014**

# *Outline*

- Bioluminescent assays
- Live and Dead Cell Assays
- “Real Time” Assays
- Apoptosis
- Stress Events Leading to Cytotoxicity
- Validation of assays for 3D culture models

# Bioluminescent Assays Developed From Firefly Luciferase Chemistry



- Reporter gene assays
- GPCR & Nuclear Receptor

## Cell Viability

- Kinase assays
- cAMP & PDE assays
- P-glycoprotein assay

## Caspases/proteases

- CYP450 assays
- MAO assays

## GSH assays

- HDAC assays

- GloSensor cAMP
- GloSensor caspase

# ***Cell Health Assays Overview***

## **Viable cells detected using markers of active metabolism**

- **Cellular conversion of indicator dyes (MTT / MTS / Resazurin)**
- **Protease marker**
- **ATP content**
- **Real time viability assay using NanoLuc**

## Dead cells detected using marker of membrane integrity

- LDH release
- Protease release
- Dye uptake / staining

## Apoptosis detected using caspase activities

## Biochemical markers of cell stress leading to cytotoxicity

- Mitochondrial toxicity
- Oxidative stress (ROS and GSH:GSSG ratio)
- NADH

## Luciferase reporters of cell stress pathways leading to cytotoxicity

# Metabolic & Enzymatic Indicators of Cell Viability

## Reagent

### Tetrazolium Reagents

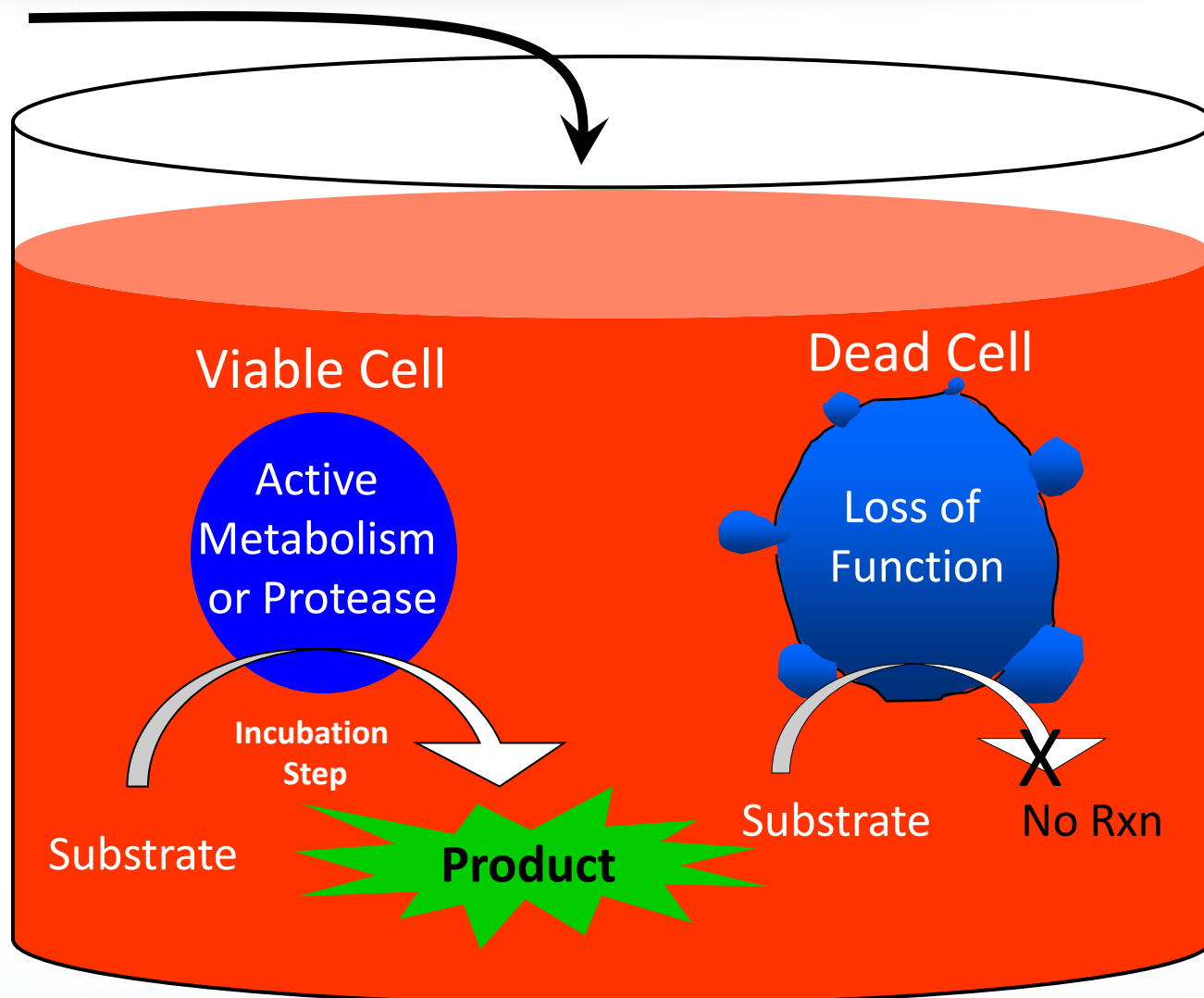
- MTT, MTS, XTT, WST

### Redox Indicators

- Resazurin

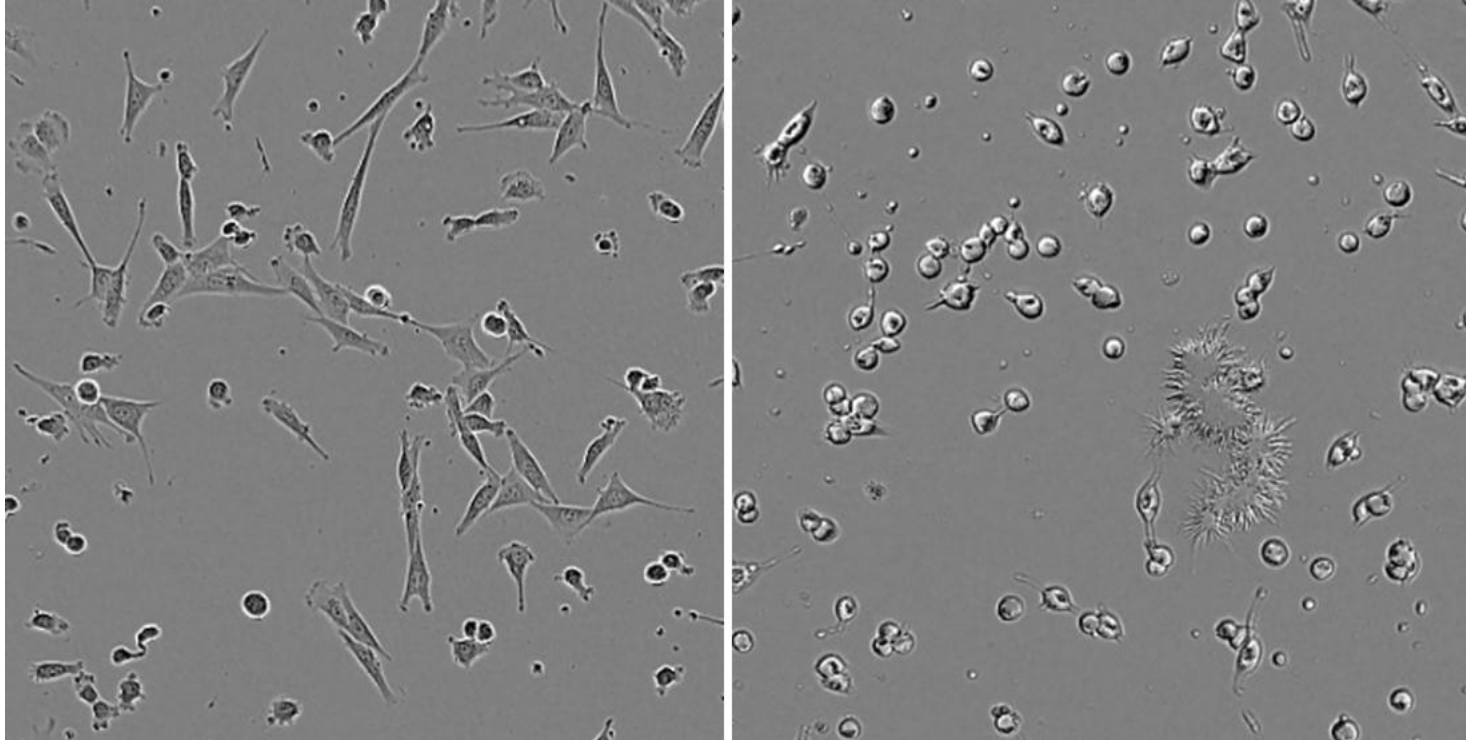
### Enzyme Substrates

- Protease Substrates  
GF-AFC



## ***Balb 3T3 Cells Treated with MTT for 4 Hours***

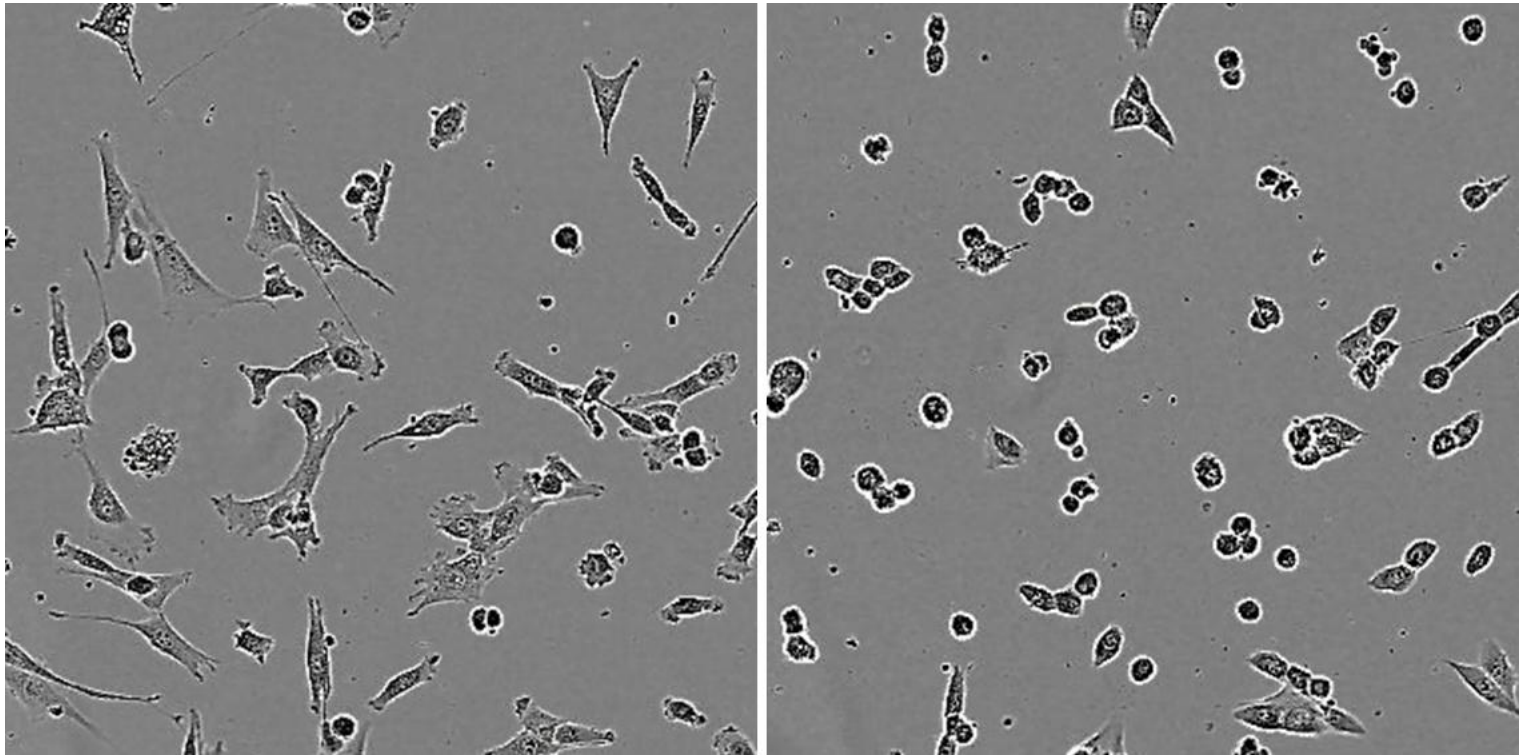
Same field of cells imaged immediately after addition of MTT and after 4 hours incubation.



Images captured by Tracy Worzella using Incucyte instrument from Essen Biosciences.



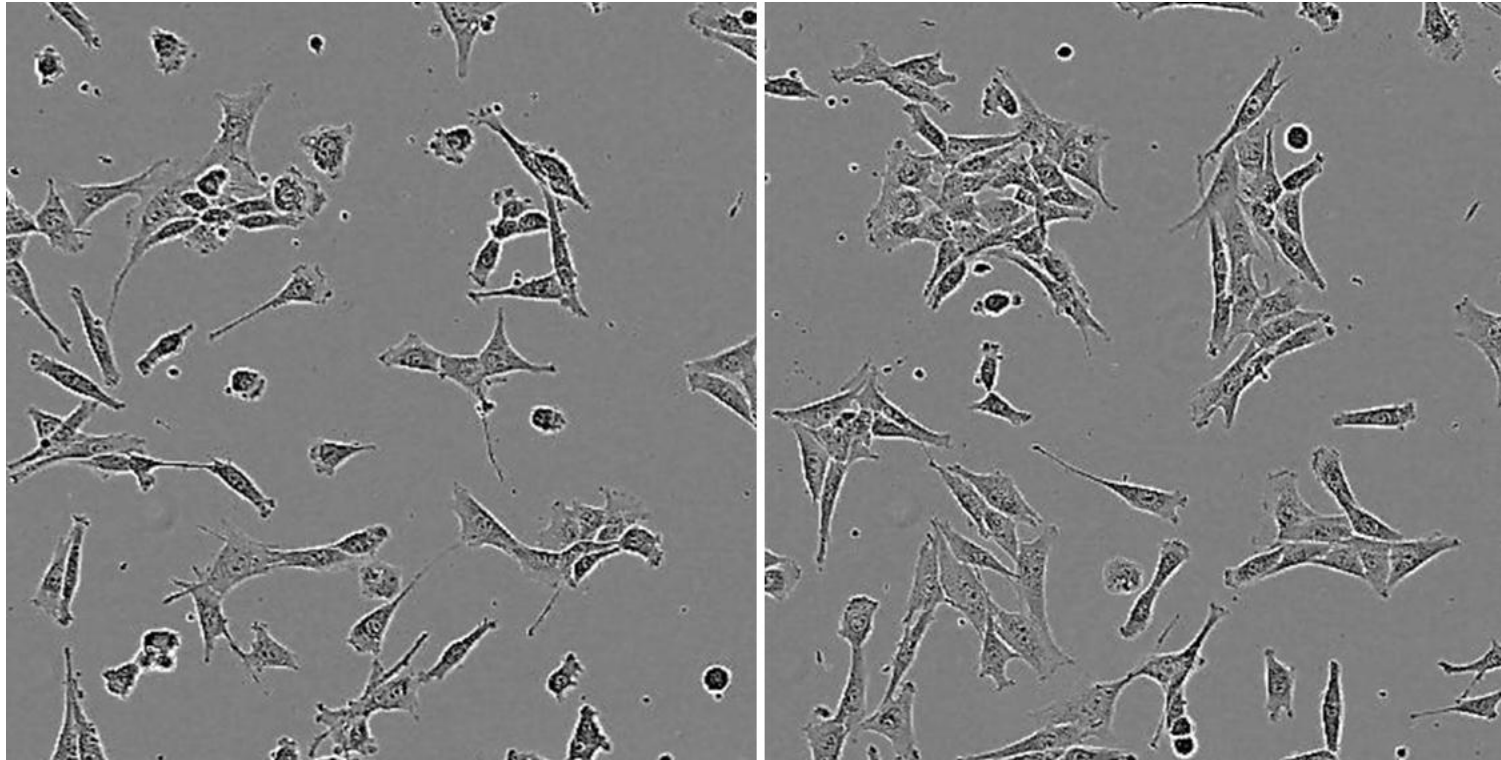
## ***Balb 3T3 Cells Treated with Resazurin for 4 Hours***



Images captured by Tracy Worzella using Incucyte instrument from Essen Biosciences.

## ***Balb 3T3 Cells Treated with GF-AFC for 4 Hours***

4 h exposure used for comparison; but 30 min is usually adequate



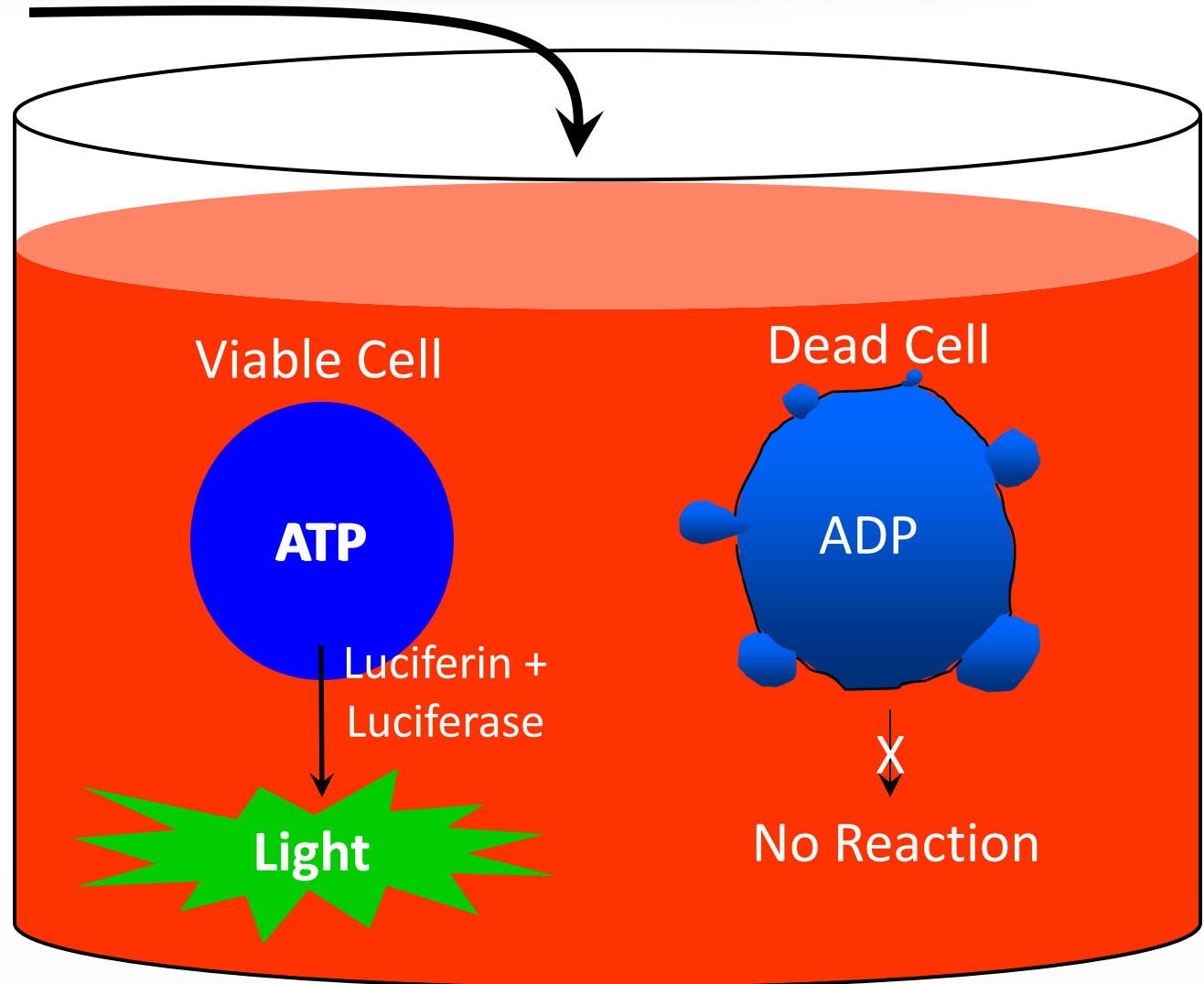
Images captured by Tracy Worzella using Incucyte instrument from Essen Biosciences.



# ATP Assay for Cell Viability (immediate lysis)

CellTiter-Glo<sup>®</sup>  
Reagent

- Lysis Solution
- ATPase Inhibitors
- Luciferin
- UltraGlo Luciferase

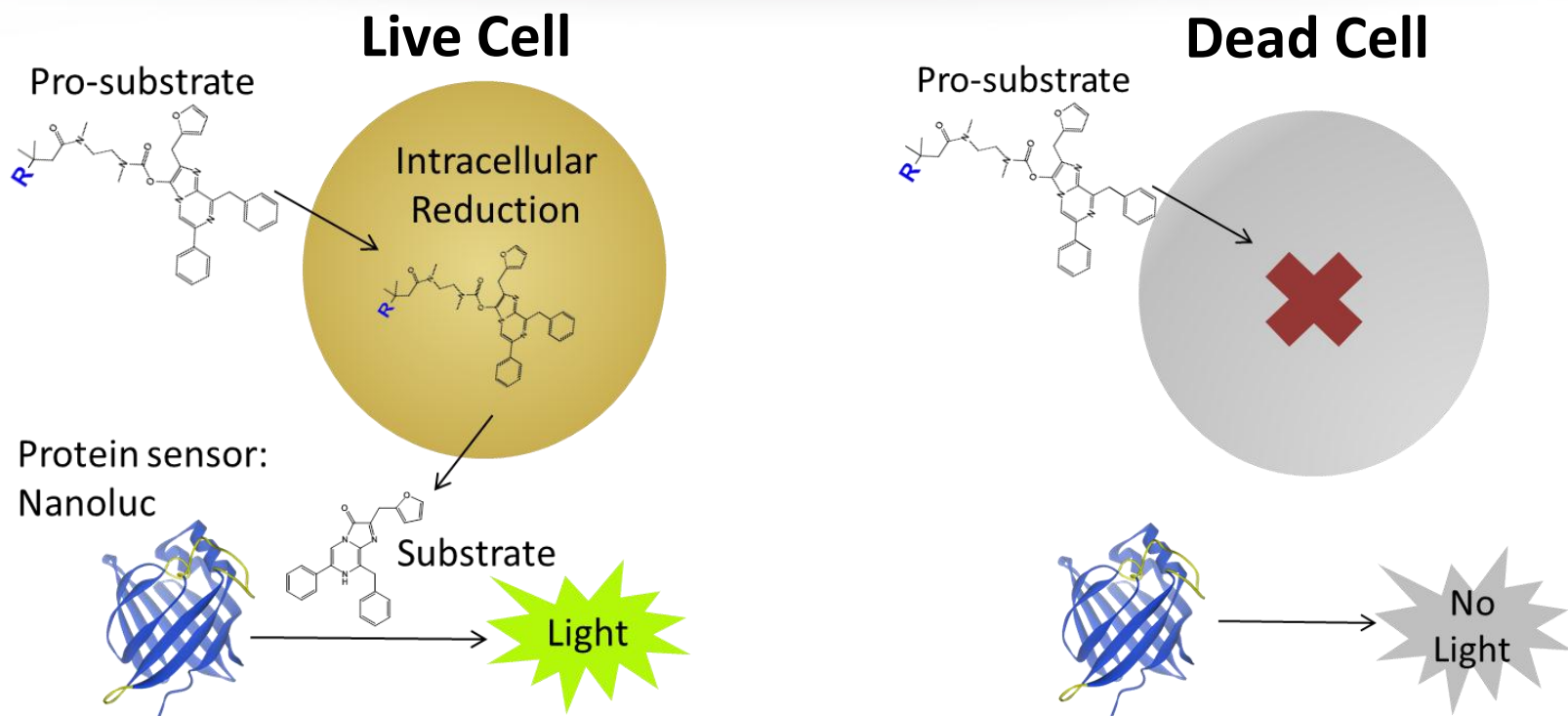


# ***Advantages & Disadvantages of Viability Assays***

<b>Assay</b>	<b>Advantages</b>	<b>Disadvantages</b>
<b>MTT / MTS</b>	Widely used Inexpensive	MTT has 2 step protocol 1-4 hour incubation Interference by reducing compounds Toxic to cells Limited sensitivity
<b>Resazurin</b>	Inexpensive Fluorescent readout Good sensitivity	1-4 hour incubation Interference by reducing compounds Toxic to cells Fluorescence interference
<b>Protease</b>	30 min protocol Cells remain viable Better sensitivity than resazurin Good choice for multiplexing	Fluorescence interference
<b>ATP</b>	10 min protocol Best sensitivity No fluorescence interference Lysis step stops reaction immediately (no incubation with viable cells)	Lytic protocol dictates sequence for multiplexing

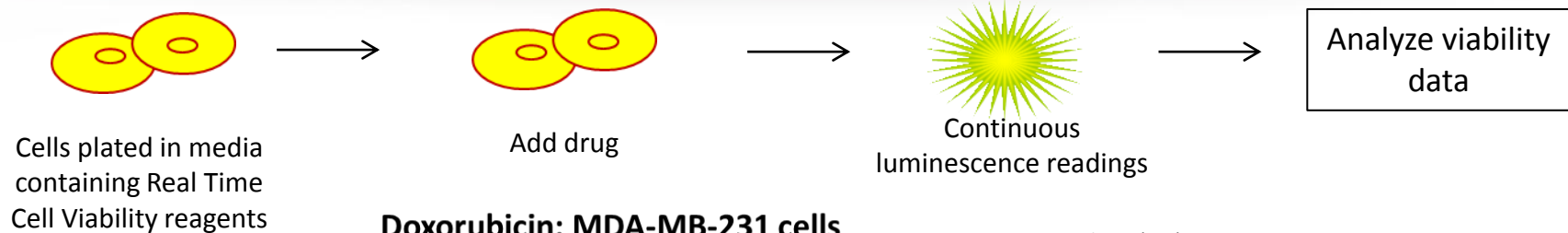
***“Real Time” Cell Viability Assay Based  
on NanoLuc Luciferase  
...in Development***

# Real Time Cell Viability Assay Measures Reducing Potential of the Cell

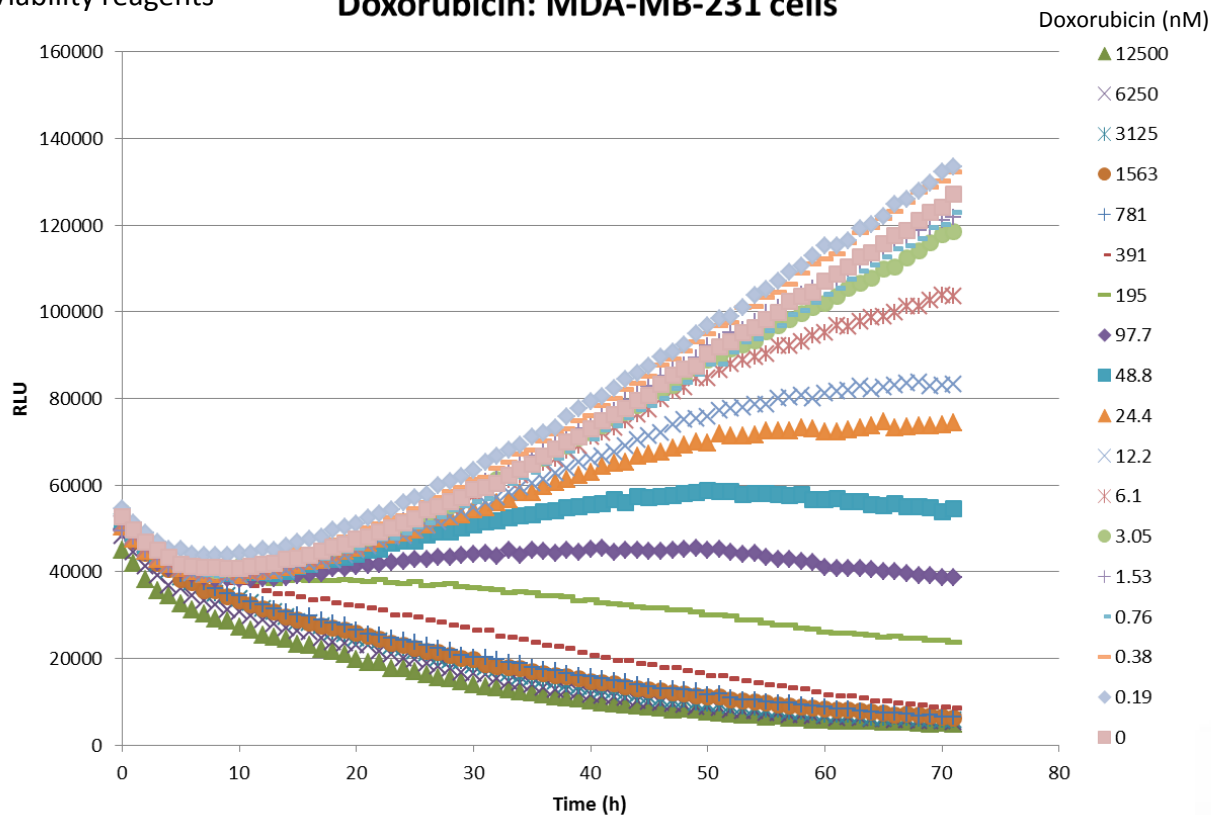


- NanoLuc protein sensor is present outside of the cells
- Pro-furimazine substrate enters the cell and is reduced by the cell to form furimazine
- Furimazine (substrate) diffuses from the cell and is rapidly used by NanoLuc to produce light

# Measure changes in viability over an extended period of time



## Doxorubicin: MDA-MB-231 cells



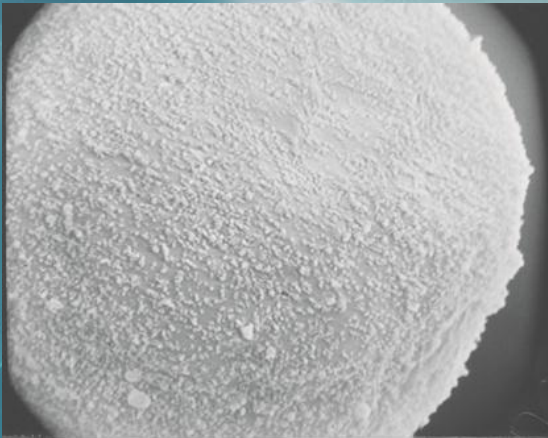
Doxorubicin	
Time (h)	IC50 (nM)
8	ND
16	ND
24	556
32	208
40	116
48	81
56	56
64	38
72	27



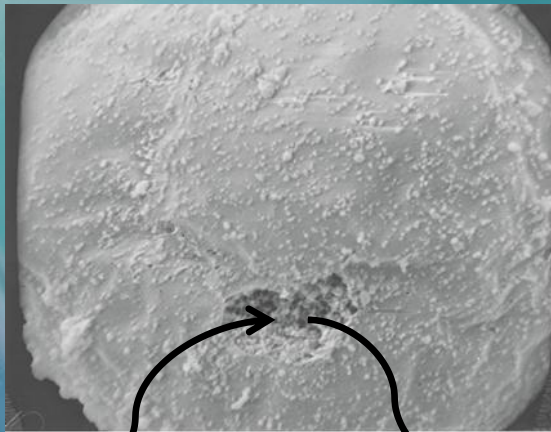
## ***Detecting Dead Cells: Two Basic Approaches***

***The definition of cell viability is based on membrane integrity.***

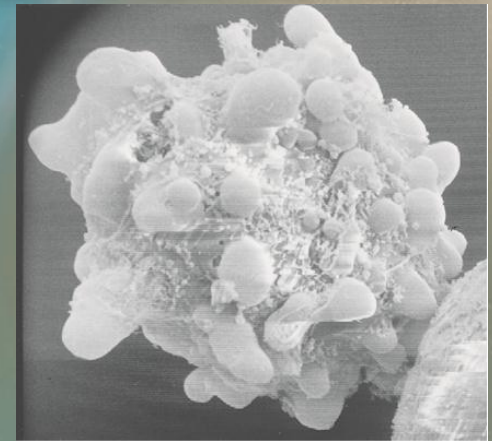
Live



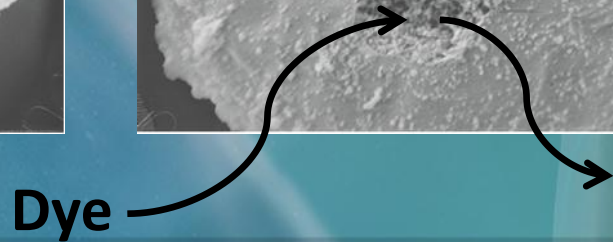
Dead



Apoptotic



Dye



Enzyme  
Marker



# ***Cell Health Assays Overview***

Viable cells detected using markers of active metabolism

- Cellular conversion of indicator dyes (MTT / MTS / Resazurin)
- Protease marker
- ATP content

**Dead cells detected using marker of membrane integrity**

- **LDH release**
- **Protease release**
- **Dye uptake / staining**

Apoptosis detected using caspase activities

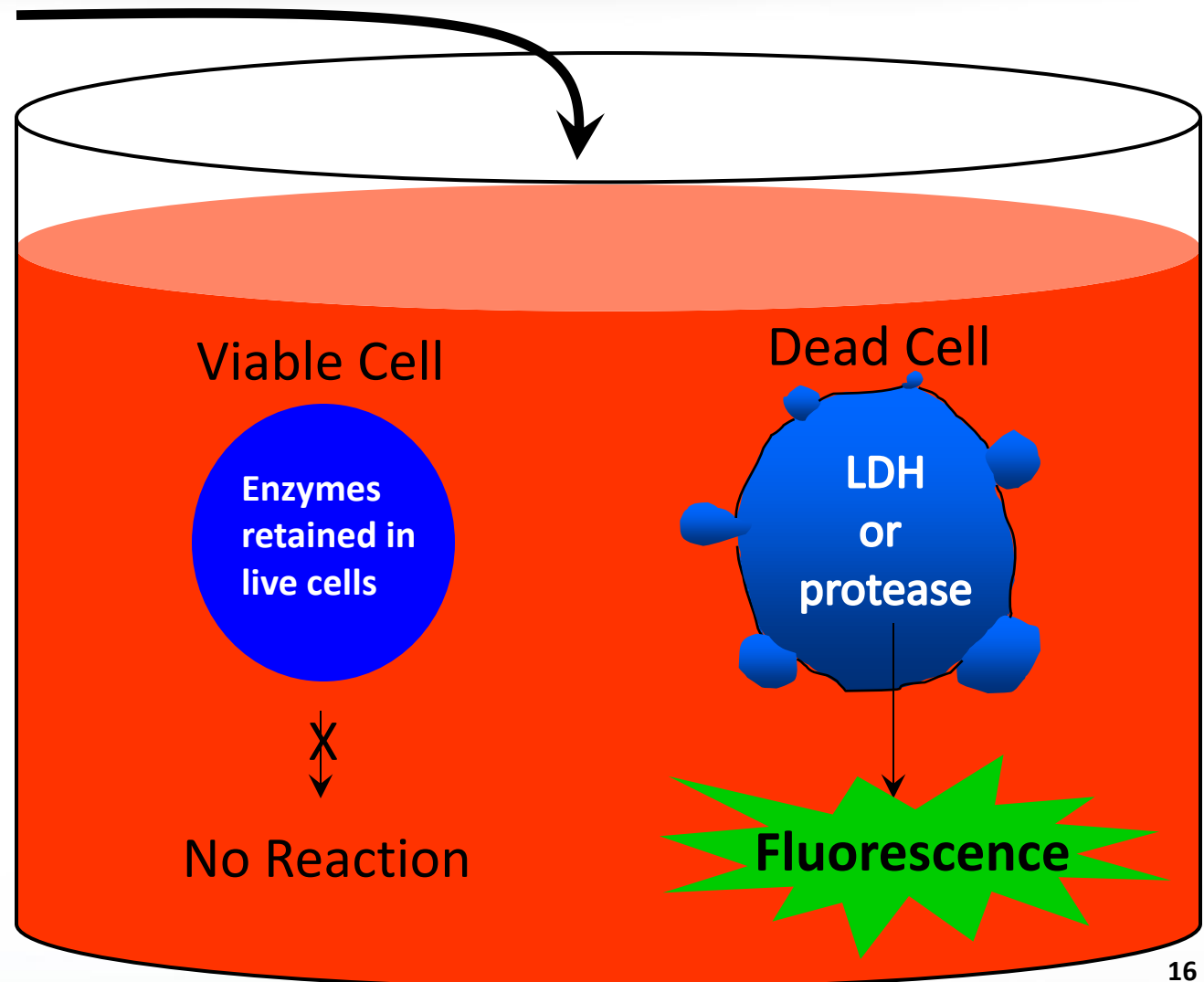
Biochemical markers of cell stress leading to cytotoxicity

- Mitochondrial toxicity
- Oxidative stress (ROS and GSH:GSSG ratio)
- NADH

Luciferase reporters of cell stress pathways leading to cytotoxicity

# Enzyme Marker Release Assay to Detect Dead Cells

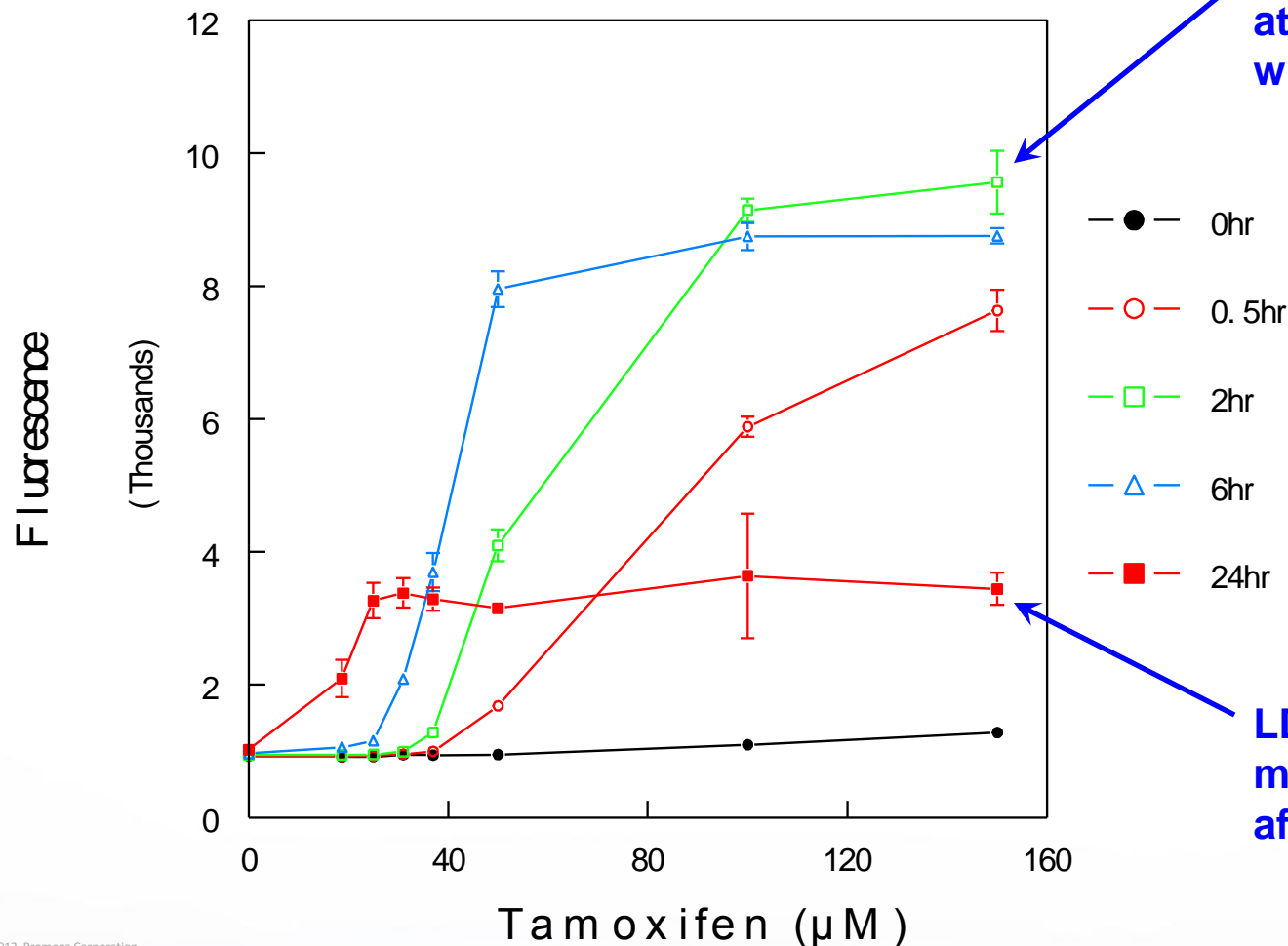
Detection  
Reagent



# *Stability of released enzyme activity in culture medium becomes a limitation*

## **LDH-Release Assay Time Course**

Tamoxifen Treated HepG2 Cells



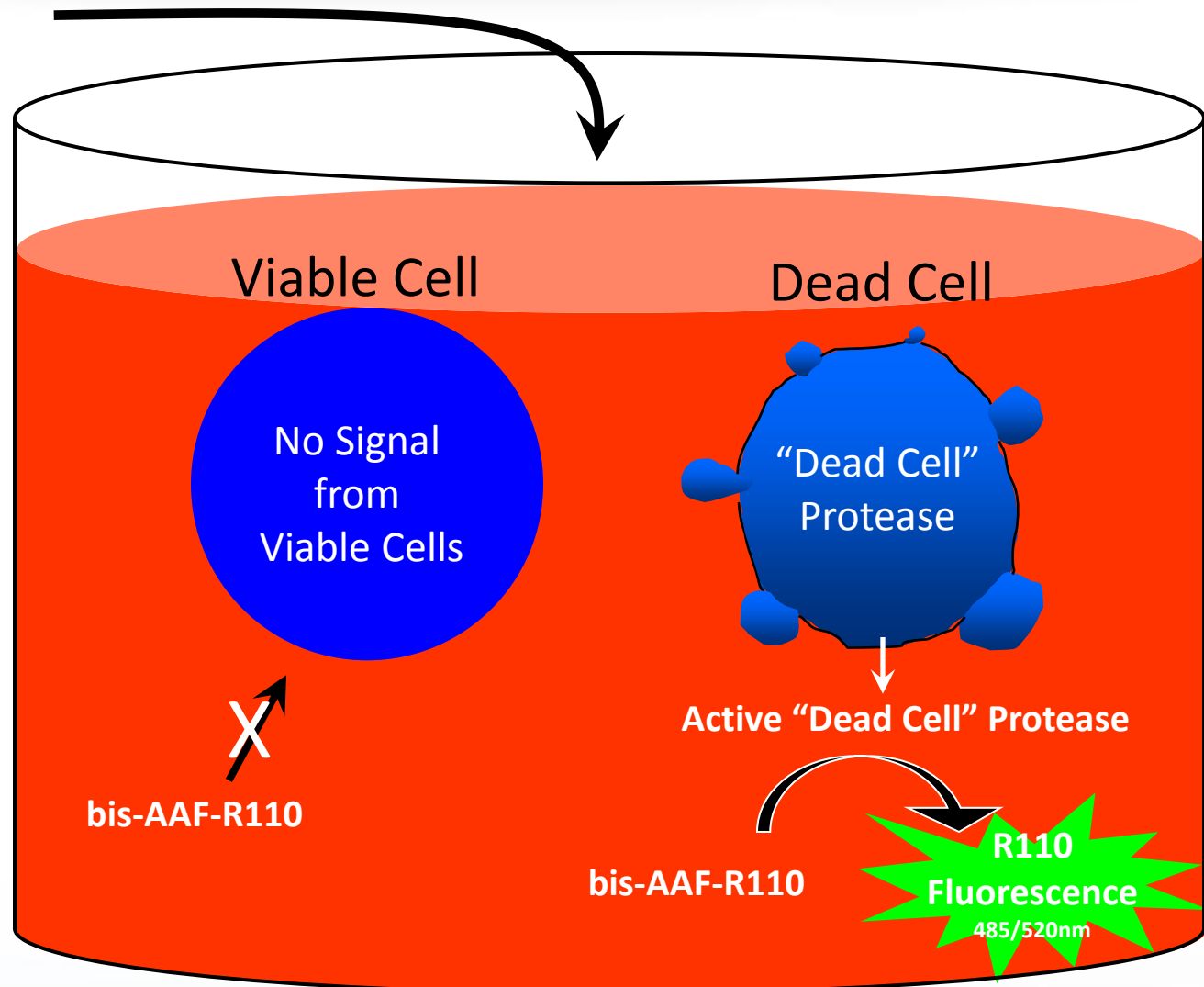
**Cells are all dead at 2 hr treatment with 150µM Tamox**

**LDH activity in medium decreases after 24 hours**

# Dead Cell Protease Assay (Fluorescent)

## Impermeable Protease Substrate (bis-AAF-R110)

- “Dead” Cell protease remains active long after cell death
- Only signal is from “Dead” Cells
- Impermeable substrate can not enter viable cells

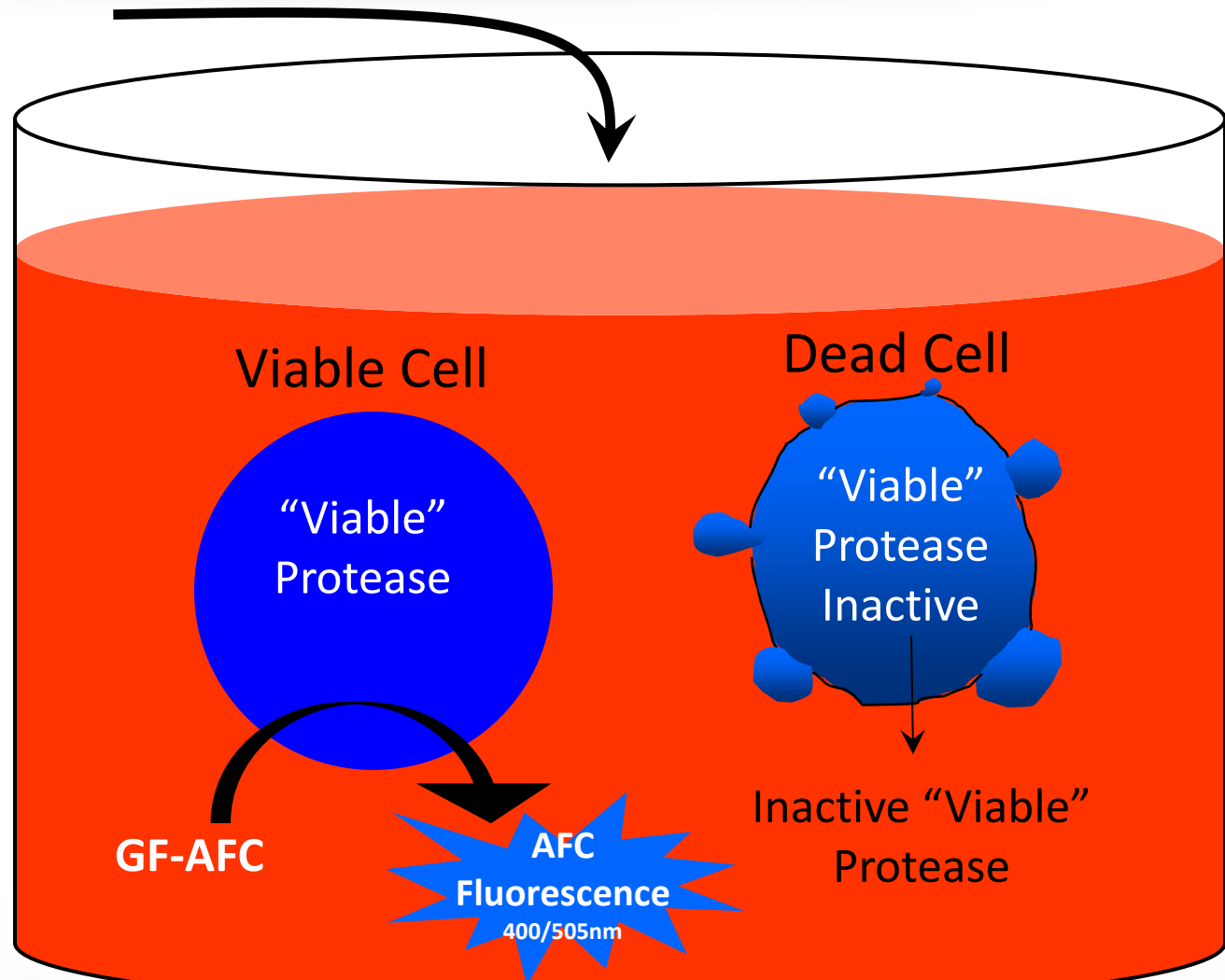




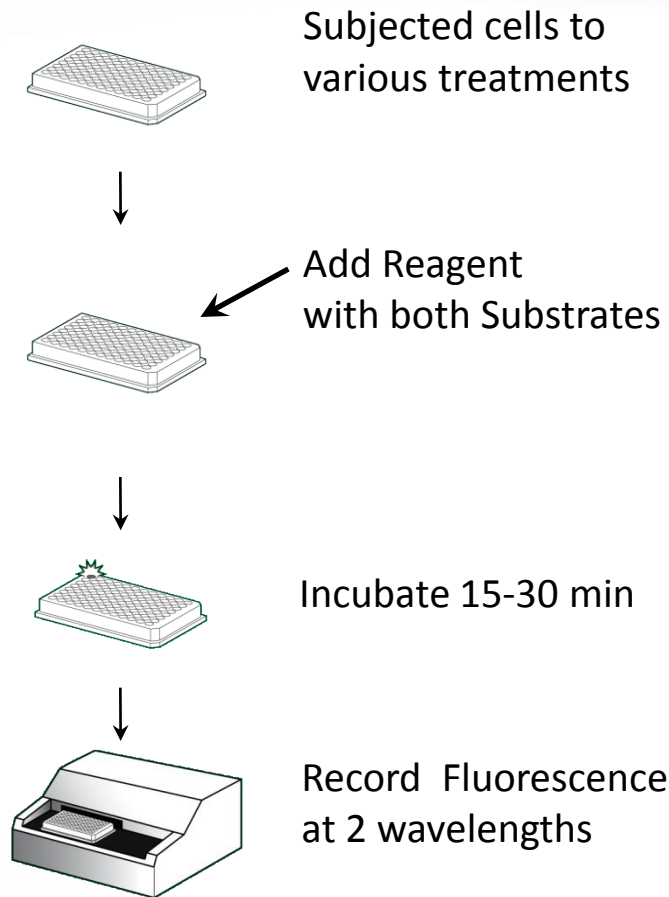
# Viable Cell Protease Assay

## Cell Permeable Protease Substrate (GF-AFC)

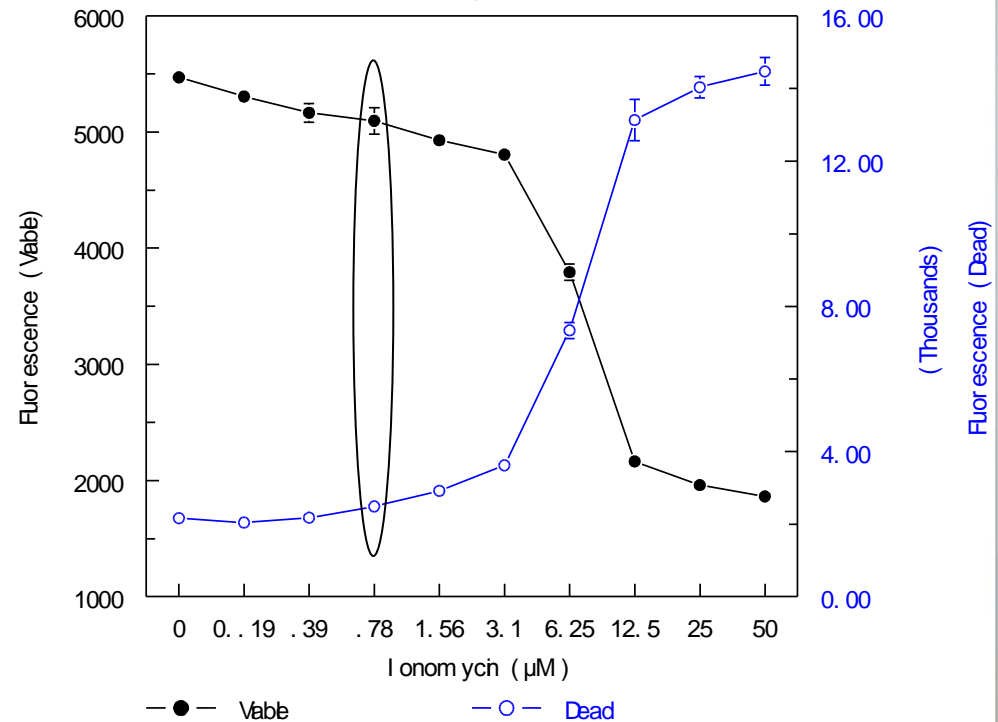
- Viable cells retain protease activity and generate signal
- “Viable” Cell protease becomes inactive upon cell death



# Multiplexing Measurement of Viable Cells & Dead Cells Simultaneously



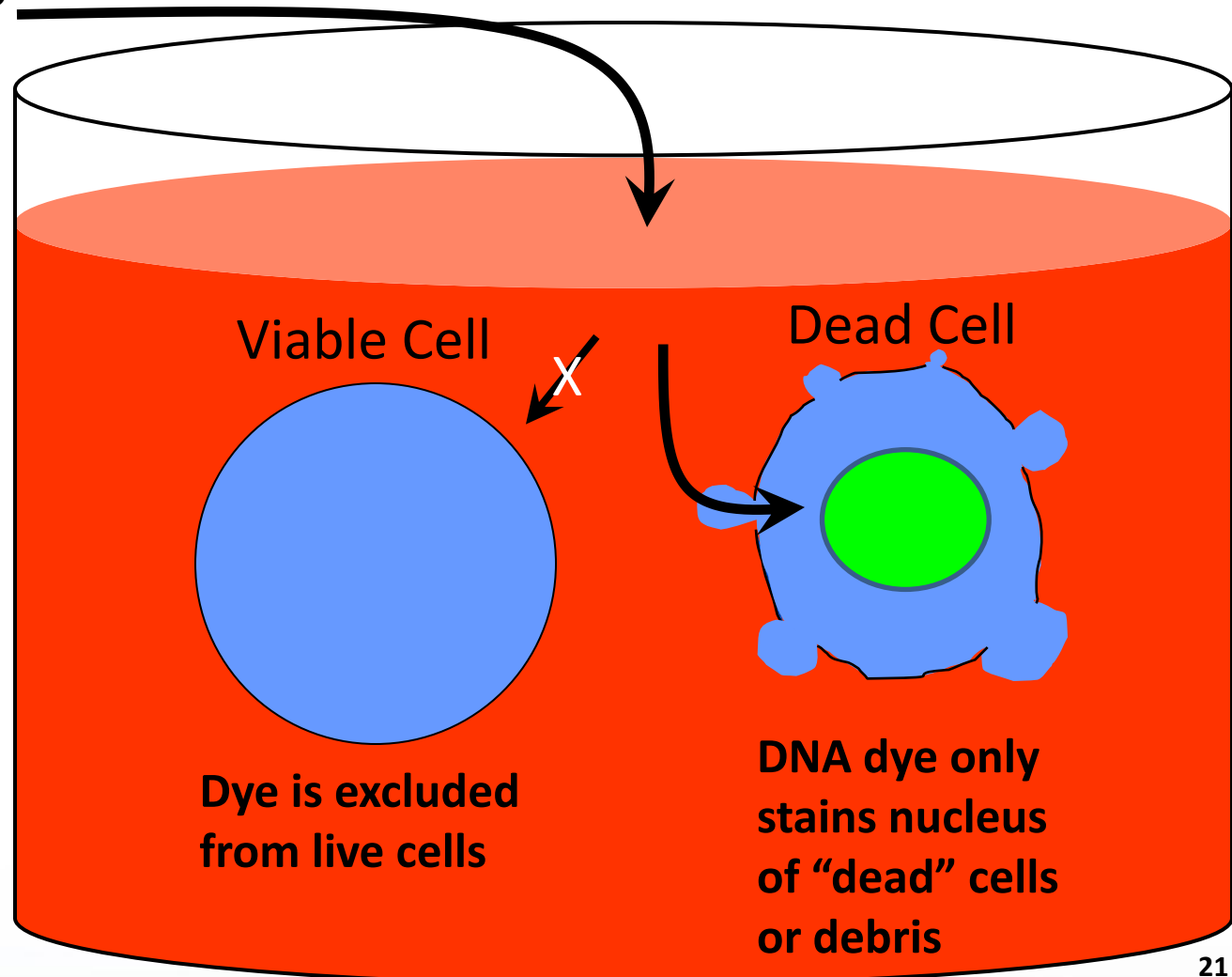
**Multiplex Assay of Viable and Dead Cells by Measuring Protease Activities**



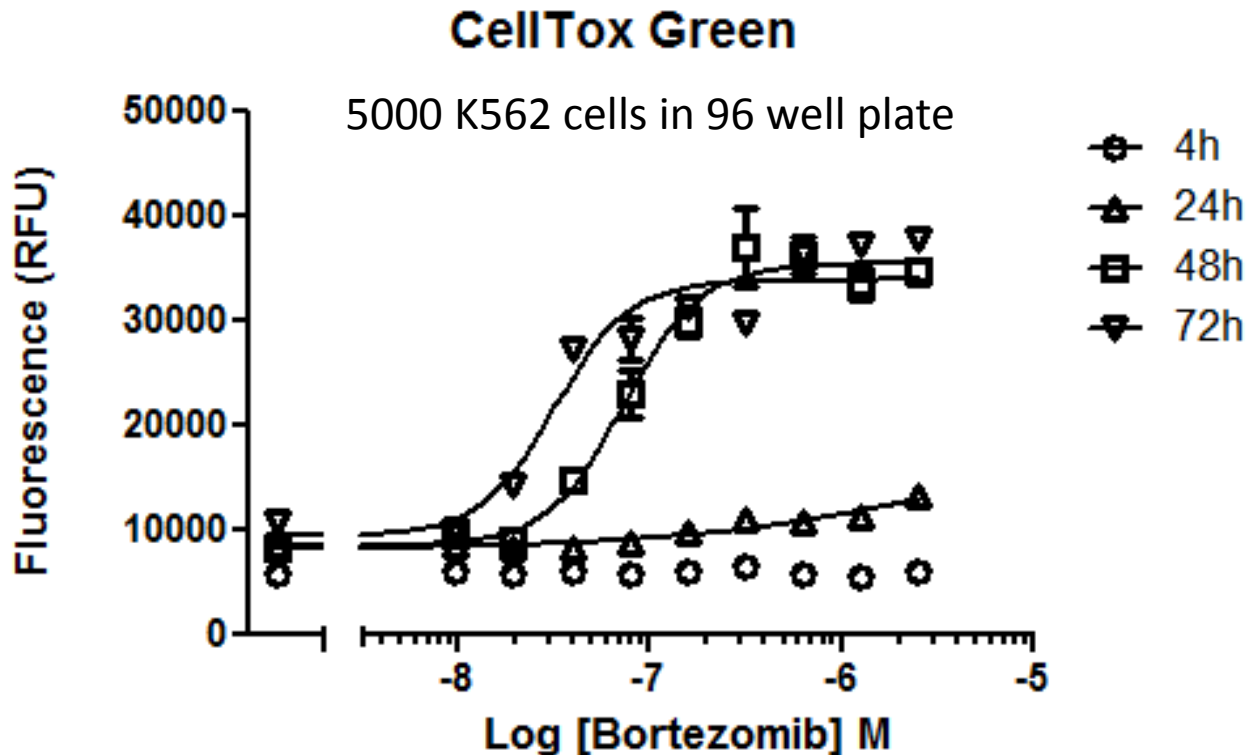
# ***DNA Dye Staining to Detect Dead Cells*** ***(Overcomes some limitations of short half-life markers)***

**Non-permeable  
DNA dye**

**Staining of dead  
cells results in a  
fluorescent signal  
that is stable.**

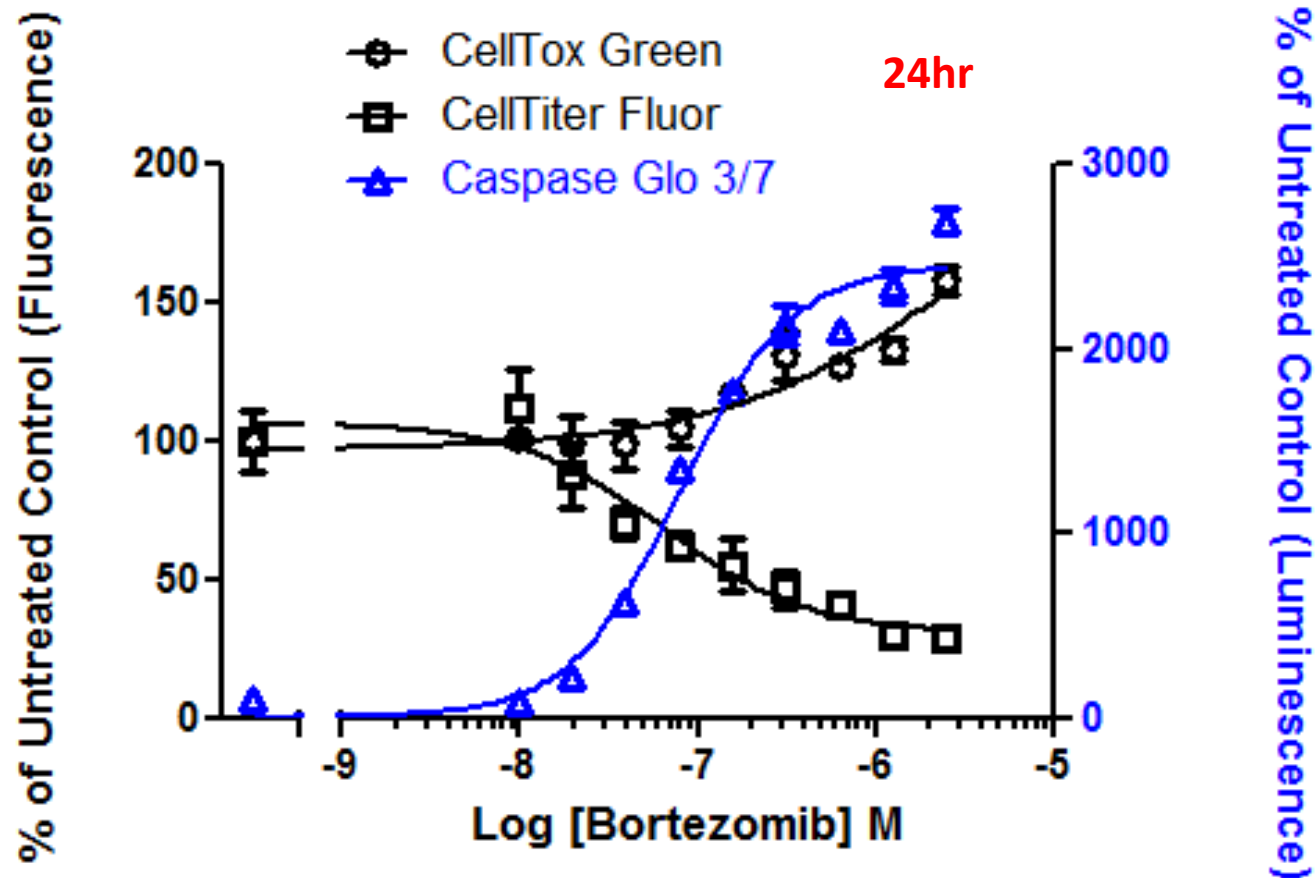


# *Reading the Same Plate Multiple Times to Detect the Onset of Cell Death (on the fly decision)*



First appearance of cell death may trigger further experimentation with the same sample (e.g. How did the cells die? ...apoptosis?)

# *Samples with CellTox Green can be Multiplexed with Cell Viability and Apoptosis Assays*





# ***Advantages & disadvantages of assays to detect dead cells***

Assay	Advantages	Disadvantages
<b>LDH release</b>	<p>Widely used and accepted</p> <p>Absorbance or fluorescent options</p>	<p>Limited sensitivity</p> <p>Limited half-life of LDH in medium</p>
<b>Protease release</b>	<p>Designed for multiplexing</p> <p>More sensitive than LDH</p> <p>Fluorescent reagent is simpler than formulation for LDH assay</p> <p>Fluorescent or luminescent options</p>	<p>Limited half-life of protease marker</p> <p>Fluorescence interference (fluorescent format only)</p>
<b>DNA Staining</b>	<p>Non-toxic / real time assay</p> <p>Staining persists for 72 hours</p> <p>Good choice for multiplexing</p>	<p>Fluorescence interference</p> <p>Less sensitive than amplified protease release assay</p>

# ***Assays to Determine Cell Stress Events Leading to Toxicity***

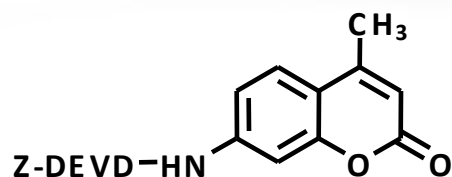
## ***Determining Mechanisms Leading to Cytotoxicity***

Going beyond the standard assays available to detect live or dead cells.

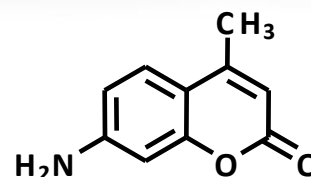
Assay chemistries and approaches to detect...

- Apoptosis
- Oxidative stress (ROS and GSH:GSSG ratio)
- Mitochondrial toxicity
- Genetic reporters to detect stress response pathways

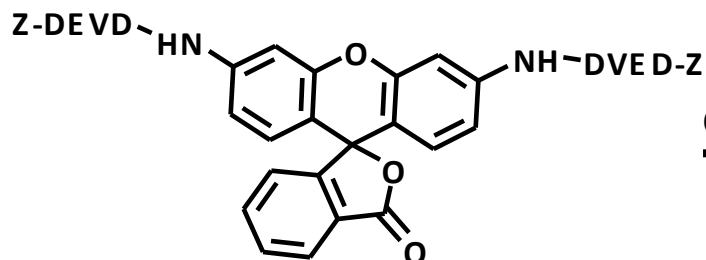
# AMC, R110 and aminoluciferin substrates for measuring caspase activity



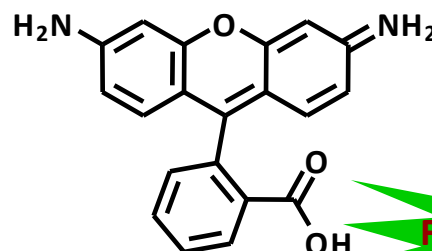
Caspase 3



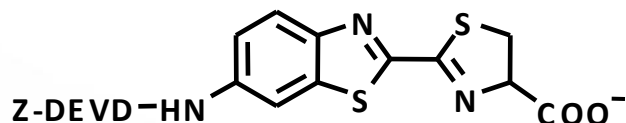
Fluorescence



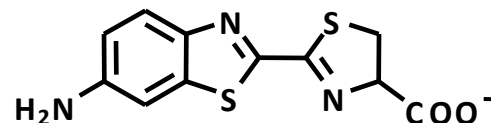
Caspase 3



Fluorescence

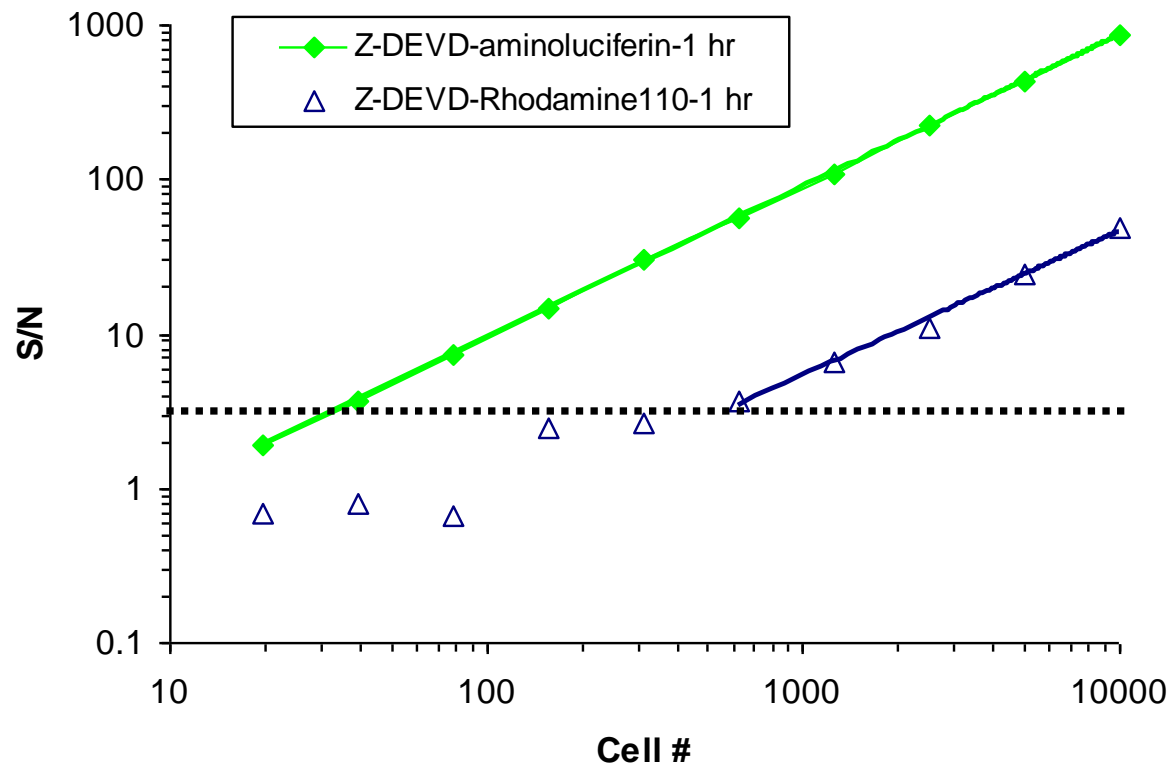


Caspase 3



Luminescence  
Luciferase + ATP

# ***Luminescent assay has better sensitivity and dynamic range***

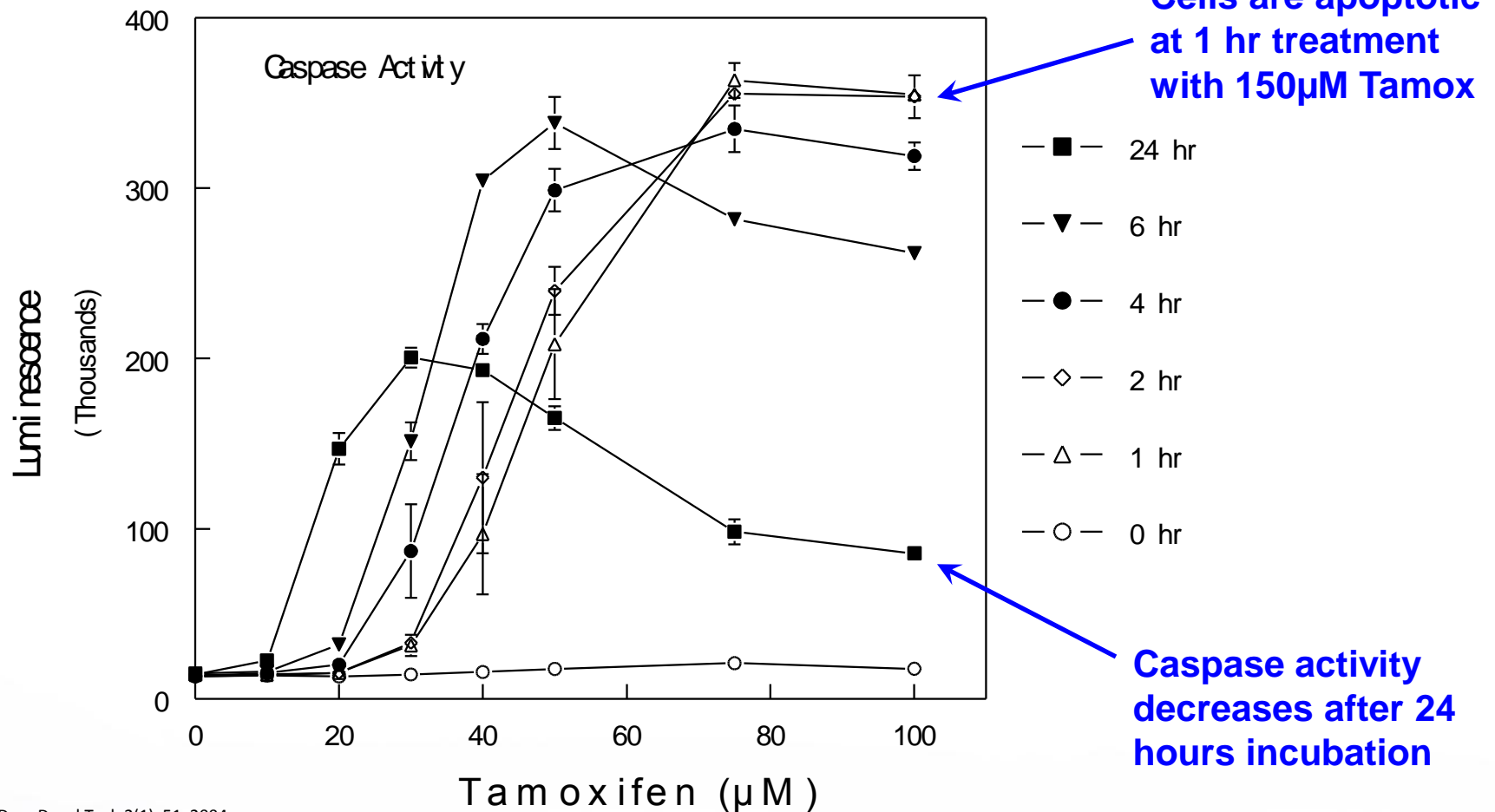


O'Brien, M et al. (2005) Homogeneous bioluminescent protease assays: Caspase-3 as a Model. *J Biomol Screen* **10**: 137.



# *Caspase-Glo<sup>®</sup> 3/7 Time Course Indicates Caspase Activity is Transient*

## Tamoxifen Treatment of HepG2 Cells



# ***Oxidative Stress Assays***

## ***Oxidative Stress Assays***

**Oxidative stress:** an imbalance between the production of reactive oxygen species (ROS) and the cell's capacity to detoxify the ROS or to repair the oxidative damage.

Markers of oxidative stress:

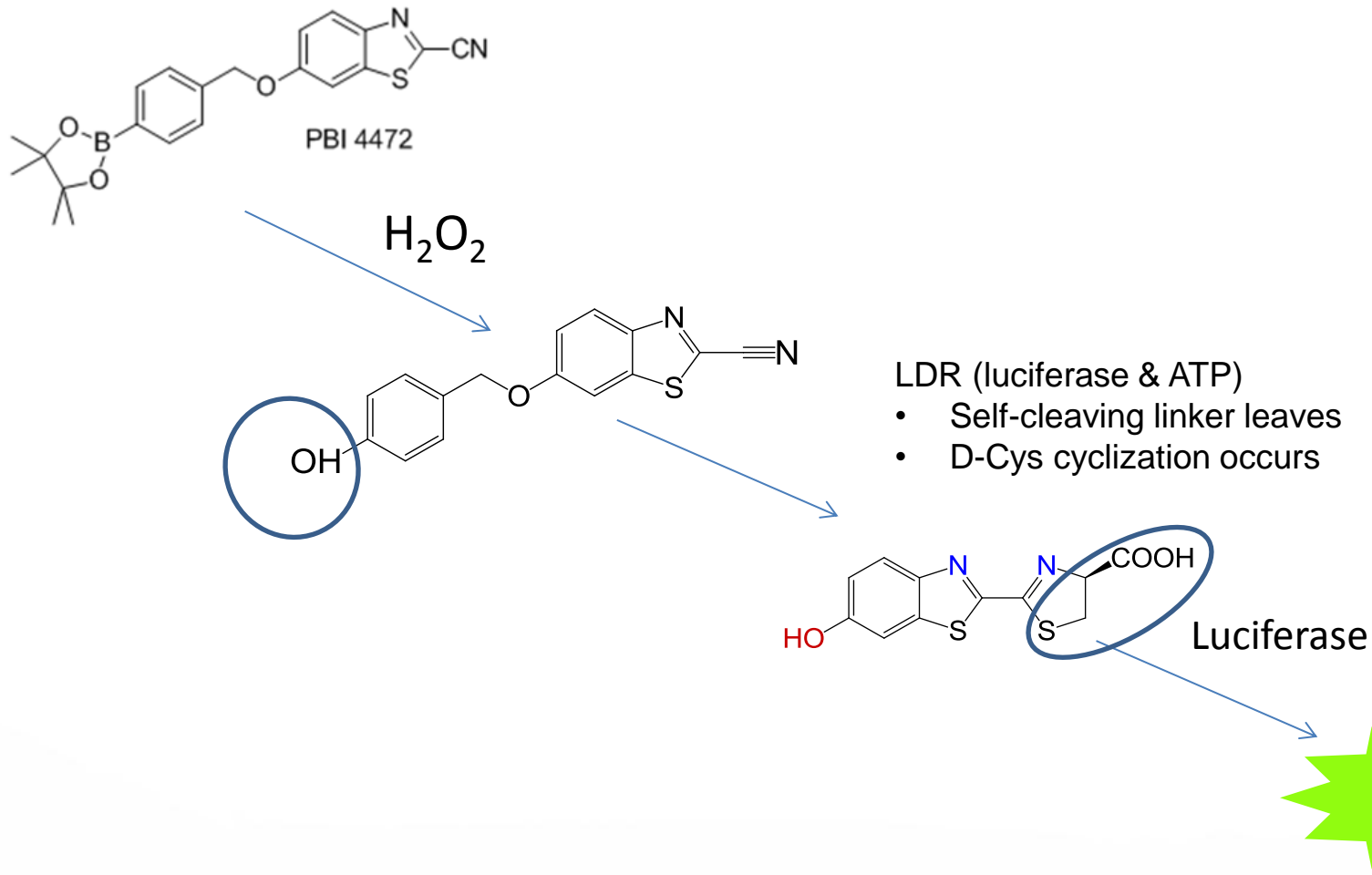
- Altered GSH:GSSG ratio (lowered GSH, increased GSSG)
- ROS (super oxide, hydroxyl radical, nitric oxide, hypochlorite convert to more stable  $\text{H}_2\text{O}_2$ )

## ***ROS-Glo $H_2O_2$ Assay***

- Direct  $H_2O_2$  detection without using Horseradish Peroxidase (HRP)
  - Mitigates HRP mediated false hits
- Homogeneous Bioluminescent Assay
  - Add-mix-read
  - No fluorescence interference
- Cell based assay
  - Detects  $H_2O_2$  content of culture wells

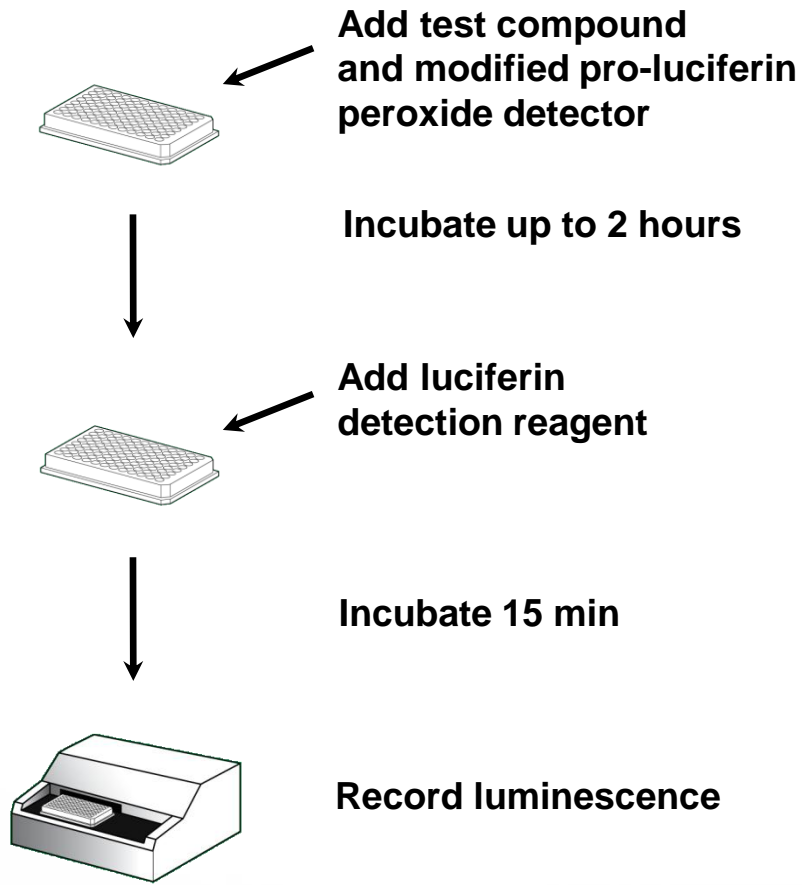
# ROS-Glo Assay Chemistry Based on Pro-Luciferin

## Modified Pro-luciferin (Peroxide Sensor)

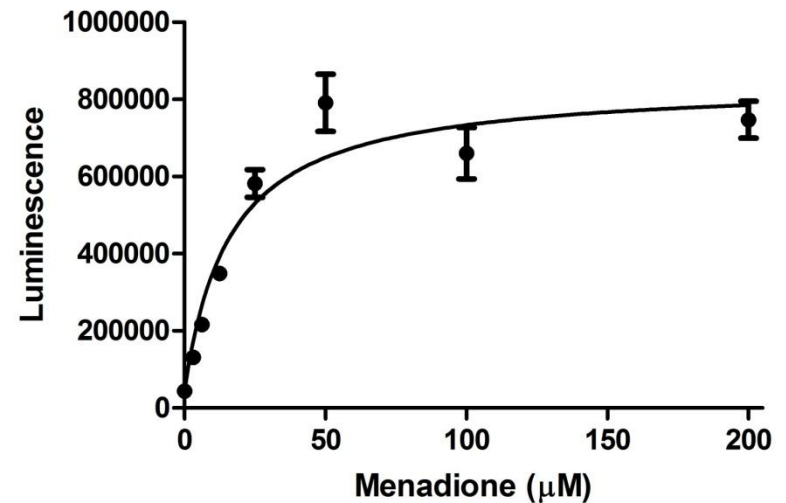




# ROS-Glo™ $H_2O_2$ Assay Protocol



## ROS-Glo™ $H_2O_2$ Assay of Hep G2 Cells Treated with Menadione



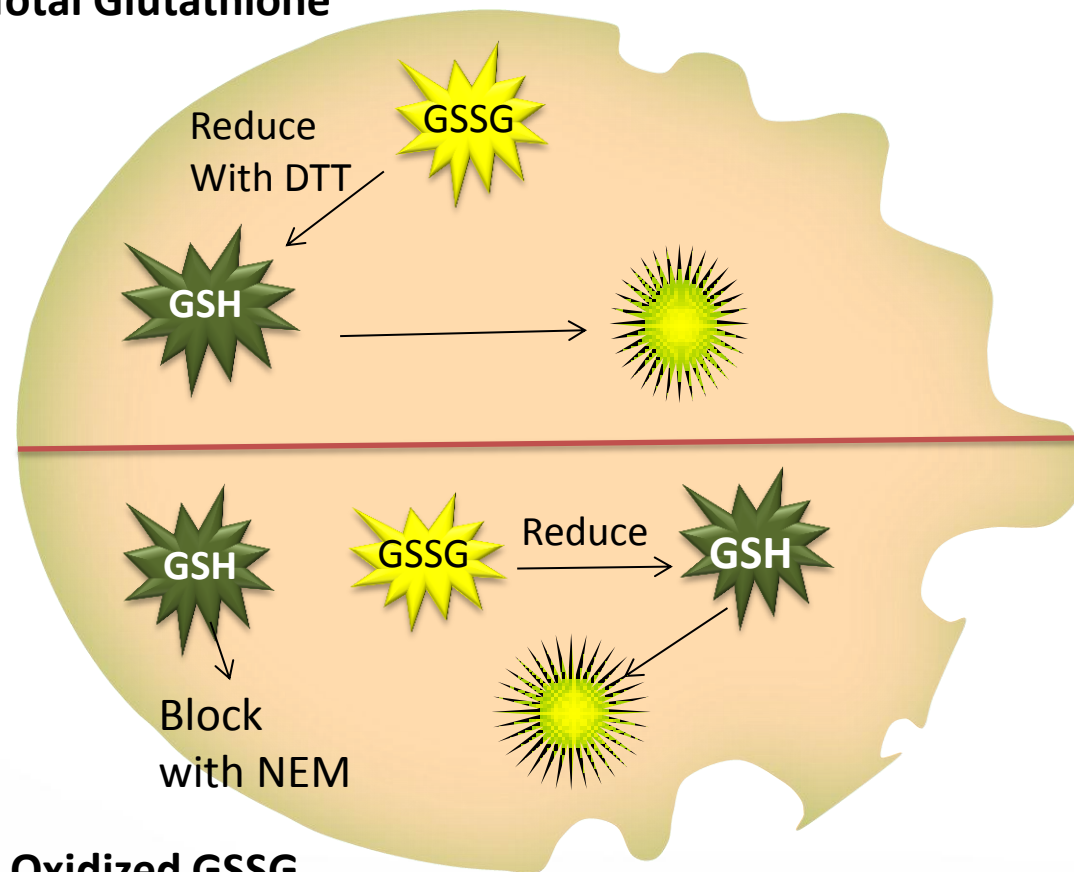
## ***GSH Assay as Marker for Oxidative Stress***

- Reduced form of glutathione (GSH) serves as an antioxidant in cells
- Decreased levels of GSH are associated with oxidative stress
- GSH and GSSG can be measured separately with a luminescent assay using Glutathione S Transferase (GST) and luciferase
- A fluorescent cell viability assay can be sequentially multiplexed with the luminescent GSH assay

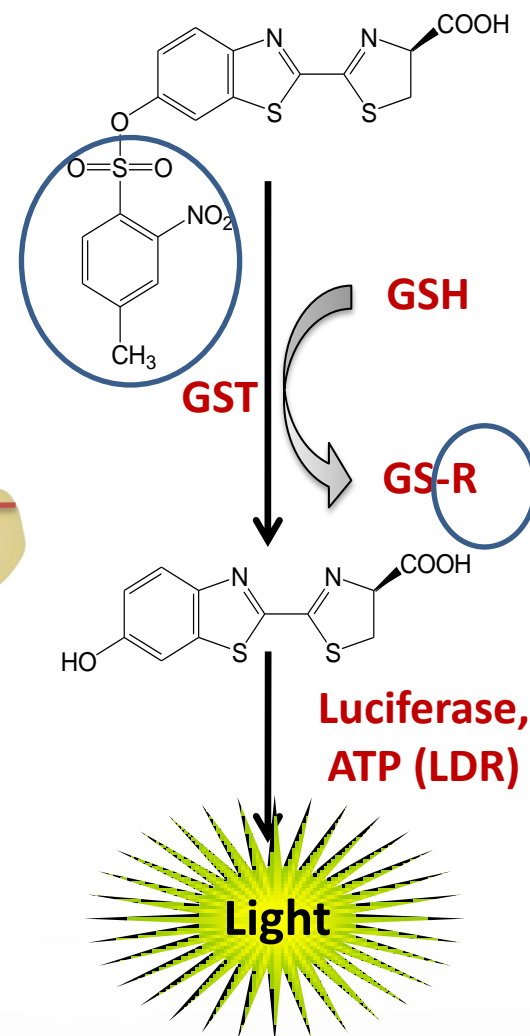
# Principal of GSH:GSSG Ratio Assay

*(Assays must be run in parallel in separate wells.)*

## Total Glutathione



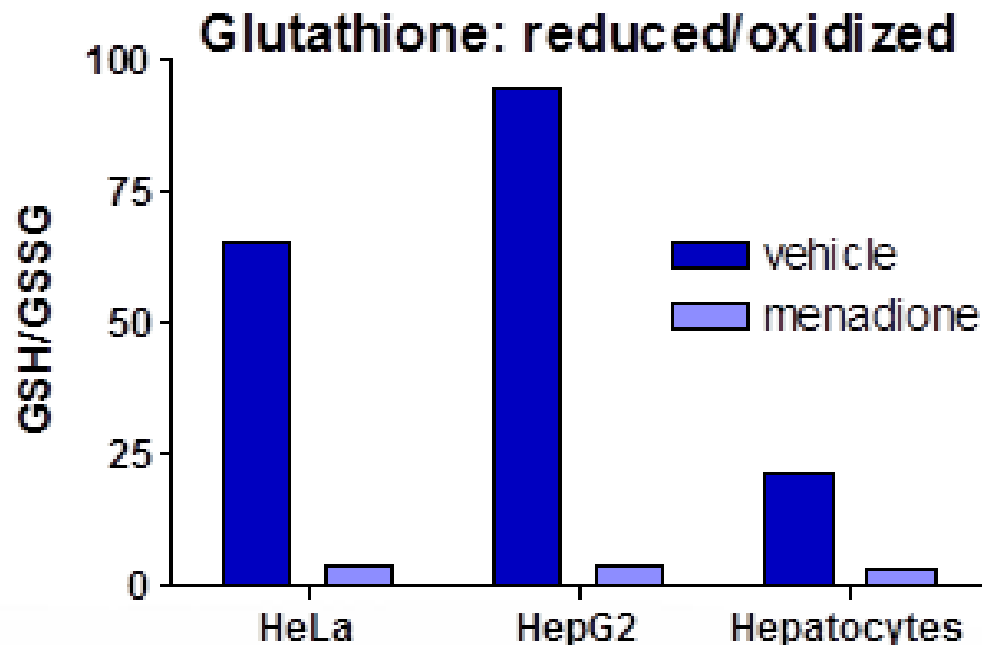
## Oxidized GSSG



## ***Menadione Treatment Drops GSH:GSSG Ratio***

GSH:GSSG changes indicate:

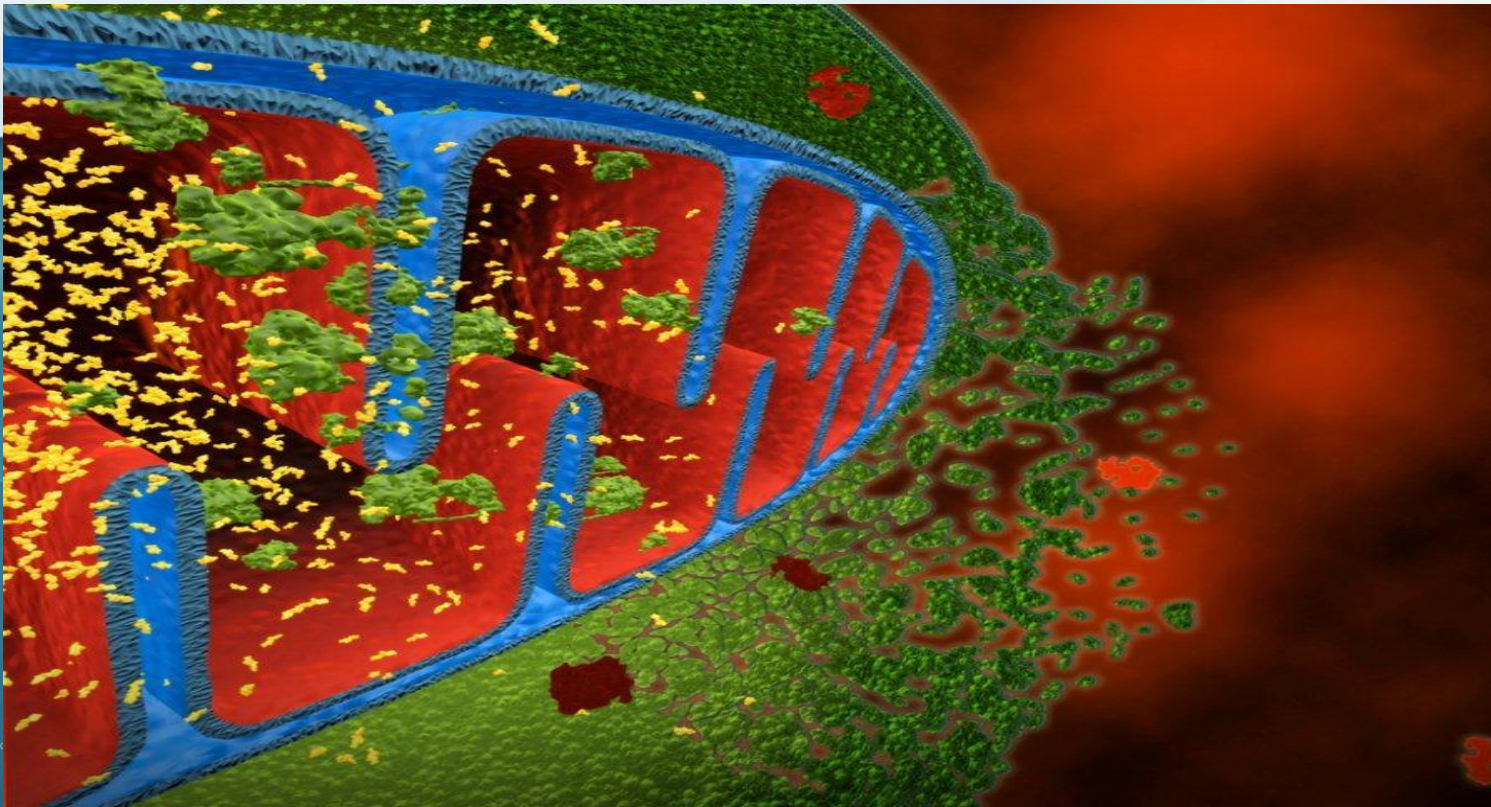
- Oxidative Stress
- Compound toxicity
- Reactive metabolite formation





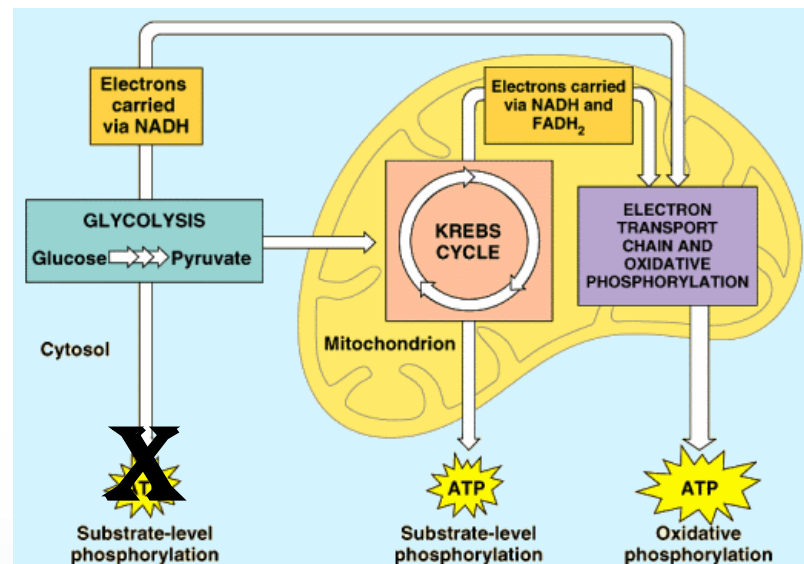
# ***Detecting Mitochondrial Toxicity***

***Rapid multiplexed cell-based assays using plate reader***



## Detecting Mitochondrial Toxicity

- ATP can be used as a marker of functioning mitochondria
- Net ATP production from glycolysis can be blocked by using glucose-free medium\* (without serum reduces background)
- Decrease in ATP marker (without general necrosis) during 1-4hr incubation suggests mitochondrial toxicity
- ATP and membrane integrity assays can be multiplexed



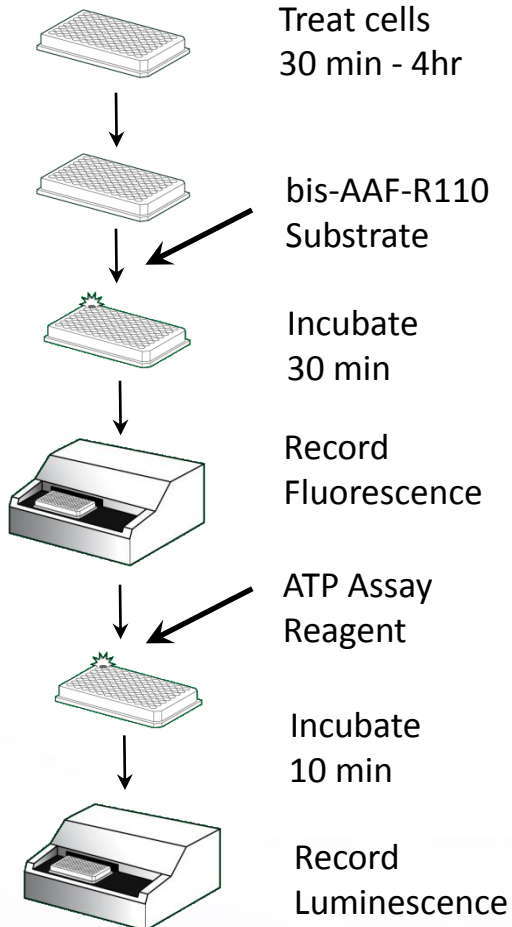
\* Lisa D. Marroquin, James Hynes, James A. Dykens, Joseph D. Jamieson, and Yvonne Will. Circumventing the crabtree effect: Replacing media glucose with galactose increases susceptibility of HepG2 cells to mitochondrial toxicants. *Toxicol. Sci.* 97:539 – 547, 2007.



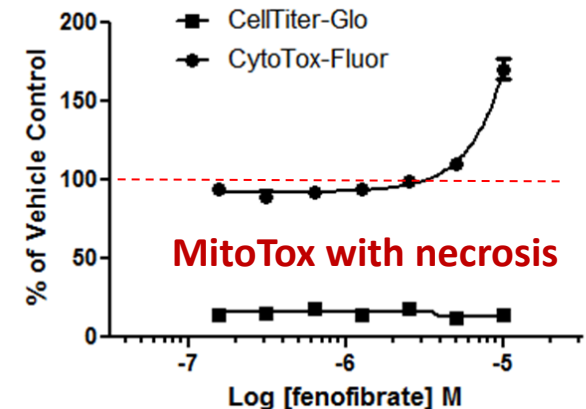
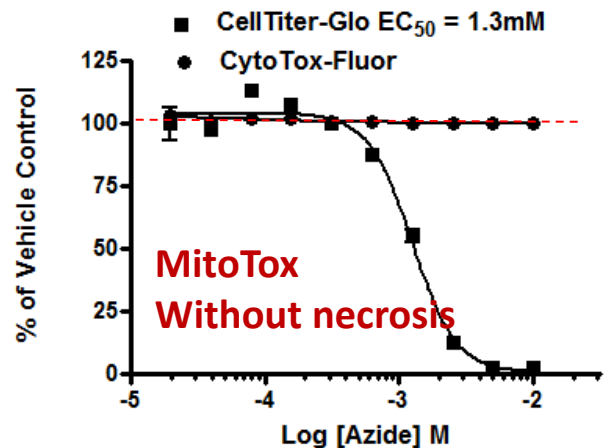
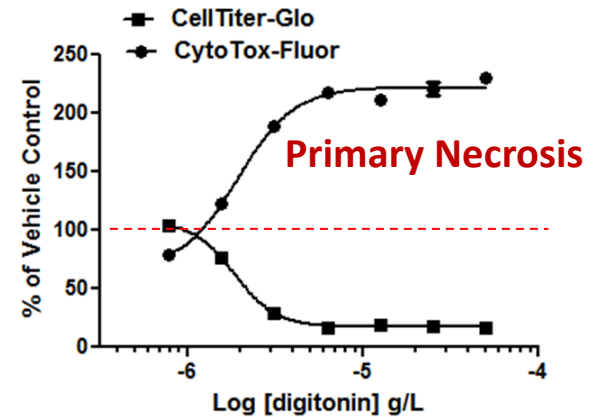
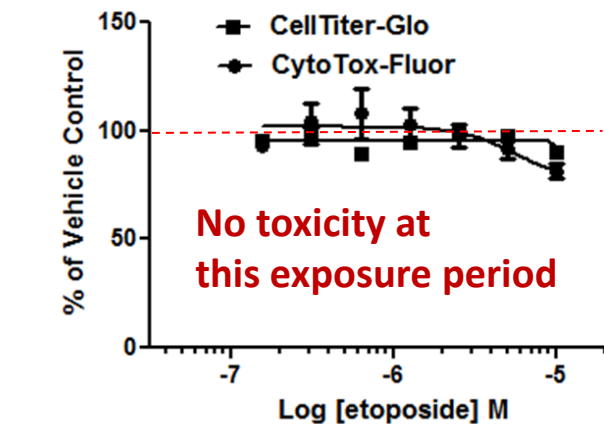
# Mitochondrial ToxGlo™ Assay

## Multiplex membrane integrity and ATP content

Change to glucose-free serum-free medium + galactose



### Expected Assay Profiles



# ***Cell Stress Response Pathway Reporters***

## ***Stress Response Pathways Leading to Cytotoxicity***

- Stress response pathways are toxin activated signal transduction events that modulate transcription factors to trigger expression of cytoprotective genes to enable the cell to attempt to restore homeostasis.\*
- Triggering cell response pathways occurs at lower toxin doses or exposure times than what is needed to trigger necrosis or apoptosis.
- If stress cannot be overcome to re-establish homeostasis, the result is induction of apoptosis and removal of the cell.

\*Simmons, S.O. et al., Cellular stress response pathway system as a sentinel ensemble in toxicological screening. Tox. Sci. 111(2): 202-225, 2009.

# Stress and Toxicity Pathway Vectors

<http://www.promega.com/resources/articles/pubhub/novel-pgl4-reporter-vector-panel-for-profiling-cellular-stress-and-chemical-toxicity/>

MAPK/JNK	Activation Protein 1 (AP1)	AP1 Response Element (AP1 RE)	pGL4.44[ <i>luc2P</i> /AP1 RE/Hygro]	E4111
Inflammation	Nuclear Factor $\kappa$ B (NF- $\kappa$ B)	NF- $\kappa$ B Response Element	pGL4.32[ <i>luc2P</i> /NF- $\kappa$ B-RE/Hygro]	E8491
Oxidative Stress	NF-E2-related factor 2 (Nrf2)	Antioxidant Response Element (ARE)	pGL4.37[ <i>luc2P</i> /ARE/Hygro]	E3641
DNA Damage	p53	p53 Response Element (p53 RE)	pGL4.38[ <i>luc2P</i> /p53 RE/Hygro]	E3651
Endoplasmic Reticulum Stress	Activating Transcription Factor 6 (ATF6)	ATF6 Response Element (ATF6 ERSE)	pGL4.39[ <i>luc2P</i> /ATF6 RE/Hygro]	E3661
Heavy Metal Stress	Metal-Responsive Transcription Factor-1 (MTF1)	Metal Regulatory Element (MRE)	pGL4.40[ <i>luc2P</i> /MRE/Hygro]	E4131
Heat Shock	Heat Shock Factor 1 (HSF1)	Heat Shock Element (HSE)	pGL4.41[ <i>luc2P</i> /HSE/Hygro]	E3751
Hypoxia	Hypoxia-Inducible Factor 1 $\alpha$ (Hif1 $\alpha$ )	Hypoxia Response Element (HRE)	pGL4.42[ <i>luc2P</i> /HRE/Hygro]	E4001
Xenobiotic Stress	Aryl hydrocarbon receptor (AhR)	Xenobiotic Responsive Element (XRE)	pGL4.43[ <i>luc2P</i> /XRE/Hygro]	E4121

# Cell lines available as custom products

<http://www.promega.com/a/forms/custom-assays/custom-assay-services.html>

## Cell Lines for Toxicity Pathway Analysis

		Pathway	Response Element	Cell line background
<input type="checkbox"/>	1	Antioxidant	ARE	HEK293
<input type="checkbox"/>	2	Hypoxia	HIF1	HEK293
<input type="checkbox"/>	3	Hypoxia	HIF1	HEPG2
<input type="checkbox"/>	4	MAPK	AP1	HEK293
<input type="checkbox"/>	5	Ras/MEK-1	SRE	HEK293
<input type="checkbox"/>	6	RhoA (Gα12/13)	SRF	HEK293
<input type="checkbox"/>	I am interested in a different toxicity pathway cell line			



# ***Validating Performance of Cytotoxicity Assays Applied to 3D Culture Models***



# One of many Examples to Determine Predictive Value of 3D Culture Models

## Activity of Anticancer Agents in a Three-Dimensional Cell Culture Model

Victor Sanjit Nirmalanandhan,<sup>1,2</sup> Alicia Duren,<sup>1</sup> Peter Hendricks,<sup>1</sup> George Vielhauer,<sup>1,3</sup> and Gurusingham Sitta Sittampalam<sup>1,2</sup>

<sup>1</sup>The Institute for Advancing Medical Innovation, <sup>2</sup>Department of Pharmacology, Toxicology, and Therapeutics, and <sup>3</sup>Department of Urology, The University of Kansas Medical Center, Kansas City, Kansas.

### ABSTRACT

Cell-monolayer-based assays for chemotherapeutic drug discovery have proven to be highly artificial compared with physiological systems. The objective of this study was to culture cancer cells in a simple 3-dimensional (3D) collagen gel model to study the antiproliferative activity of known lung cancer drugs. The validity of our 3D model was tested by measuring the activity of 10 lung cancer drugs (Paclitaxel, Alimta, Zactima, Doxorubicin, Vinorelbine, Gemcitabine, 17AAG, Cisplatin, and 2 experimental drugs from the University of Kansas [KU174 and KU363]) in 2 lung cancer cell lines (A549 and H358) and comparing the activity in a traditional 2-dimensional (2D) *in vitro* cellular assay. Both potency and efficacy of these drugs were calculated to evaluate the activity of the drugs. Our results demonstrate that the activity of these drugs showed significant differences when tested in 3D cultures, which varied with individual drugs and the cell line used for testing. For example, the cytotoxicity of Paclitaxel, KU174, Alimta, Zactima, Doxorubicin, Vinorelbine, KU363, and 17AAG was significantly changed when tested in the 3D model, whereas the potency of Cisplatin and Gemcitabine in H358 cell line remained unaffected. A similar pattern, with some differences, was observed in A549 cells and is discussed in detail in this article. The observed differences in potency and efficacy of the cancer drugs in 3D models suggest that the biological implications

of screening configurations should be taken into account to select superior cancer drug candidates in preclinical studies.

### INTRODUCTION

Cancer is the second leading cause of death in the United States, with lung cancer having the highest mortality rate.<sup>1</sup> With this high rate of death, it is becoming more imperative to find new molecular entities (NME) that can be rapidly advanced to the clinic by using improved high-throughput screening (HTS) and lead optimization strategies. Currently, NME for biological and pharmacological activities are screened by growing cells as a 2-dimensional (2D) monolayer culture. Cell monolayers are physiologically different compared with 3-dimensional (3D) human tissues or tumors.<sup>2,3</sup> Monolayer cell cultures are not a physiologically relevant environment for NME screening, and therefore may lead to the low predictability of these 2D models in preclinical trials.<sup>4</sup> Although the conventional focus of chemotherapeutic drug development has been on tumor cells, recent studies show that interactions with the stroma (Fig. 1) might influence drug sensitivity and development of drug resistance.<sup>5</sup> The extracellular matrix (ECM), which is considered part of the stroma and has been shown to activate signaling pathways, might not have been present when the drugs were tested in 2D cell-based assays, rendering this approach less representative of a human tissue or tumor<sup>6</sup> and therefore could contribute to less success in animal studies and clinical trials.

To closely mimic the *in vivo* behavior of normal or cancerous human tissue, a 3D model needs to be used.<sup>4,6–10</sup> Although 3D cultures of lung cancer cells have been used to study various aspects of tumorigenesis and tumor progression,<sup>6,11</sup> to the best of our knowledge, these models have yet to be used to screen new lung cancer drugs. Using a 3D tumor model as a strategy to test NMEs is more physiologically relevant than the traditional 2D cell culture

*“The observed differences in potency and efficacy of the cancer drugs in 3D models suggest that the biological implications of screening configurations should be taken into account to select superior cancer drug candidates in preclinical studies.”*

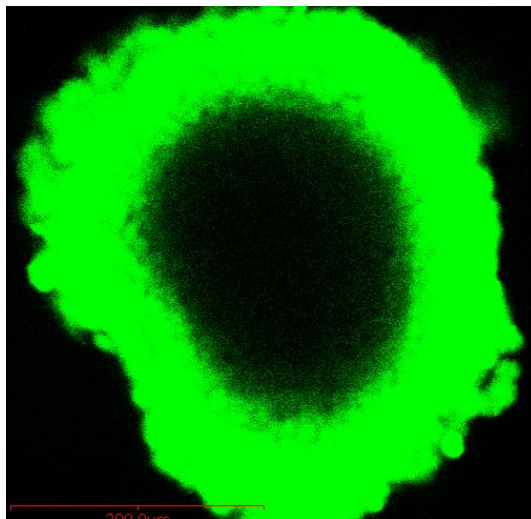
**ABBREVIATIONS:** ANOVA, analysis of variance; AUC, area under the curve; DMSO, dimethyl sulfoxide; ECM, extracellular matrix; EGFR, epidermal growth factor receptor; FBS, fetal bovine serum; HTS, high-throughput screening; NME, new molecular entities; PBS, phosphate-buffered saline; 3D, 3-dimensional; 2D, 2-dimensional; VEGFR2, vascular endothelial growth factor receptor 2.

# ***Critical Experiment that Prompted Much of our Effort to Begin Validating Promega Assays***

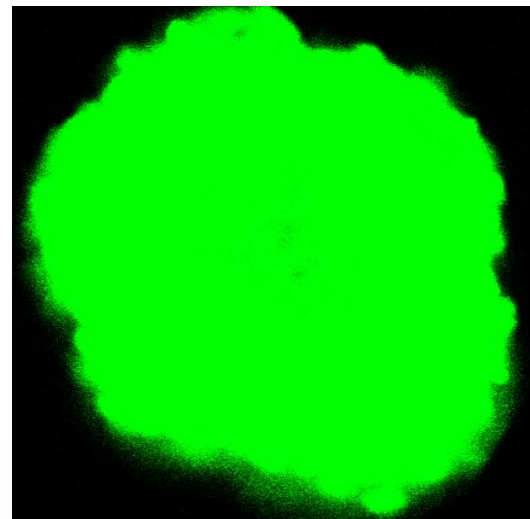
## ***Observation of Lytic Efficiency of ATP Assay Reagents***

- Microspheres grown to  $\sim 350\mu\text{m}$  using hanging drop method
- Add ATP assay reagents + DNA dye to indicate lytic effectiveness
- Photograph using laser confocal microscopy

ATPLite 1-Step Reagent



CellTiter-Glo<sup>®</sup> Reagent



$\sim 350\mu\text{m}$   
spheroids

# *Improved ATP Assay Reagent Protocol*

## **CellTiter-Glo® Assay**

- Equilibrate plate to ambient temperature
- Add equal volume of reagent
- Wait 10 min for signal to equilibrate
- Record luminescence

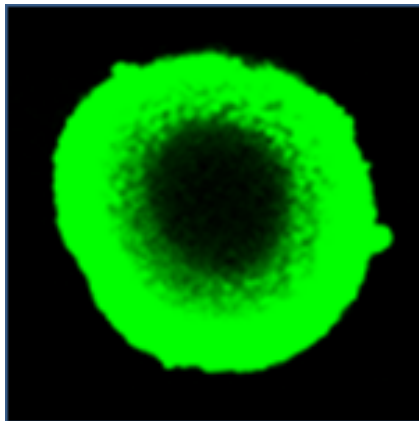
## **CellTiter-Glo® 3D Assay**

- Equilibrate plate to ambient temperature
- Add equal volume of reagent
- **Shake 5 min 450 rpm**
- Wait **25 min** for signal to equilibrate
- Record luminescence

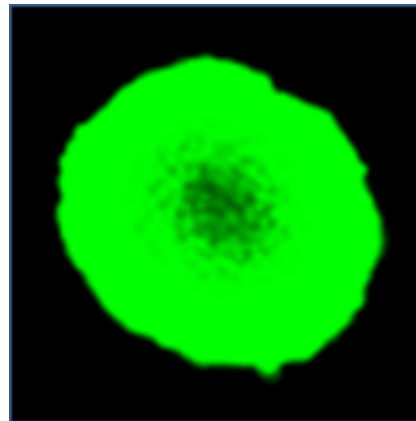
# ***Caspase-Glo® 3/7 Assay: Protocol Optimization to Facilitate Enhanced Microtissue Cell Lysis***

- HCT116 cell spheroids grown to ~330µm using hanging drop method
- Add Caspase-Glo® 3/7 assay reagent + DNA dye to indicate lytic effectiveness
- Shake with assay reagent for 5 or 30 min
- Image with confocal after a total of 30 min incubation with reagent

5 min shake



30 min shake



- Increased shake time results in near complete spheroid cell lysis
- ***Note: Human liver microtissues undergo complete lysis and disruption using these conditions (data not shown)***

## *Take Home Message*

- Most cell-based assays were designed for 2D culture models
- Assay effectiveness depends on size of microtissue
- Validation should be done for each assay and 3D culture model
- Promega is in the process of validating many of our cell health and genetic reporter
- Collaborations to validate assays on model systems are welcome

# *Summary*

- Bioluminescent assays
- Live and Dead Cell Assays
- “Real Time” Assays
- Apoptosis
- Stress Events Leading to Cytotoxicity
- Validation of assays for 3D culture models



תודה  
 Dankie Gracias  
 Спасибо شكراً  
 Merci Takk  
 Köszönjük Terima kasih  
 Grazie Dziękujemy Děkojame  
 Ďakujeme Vielen Dank Paldies  
 Kiitos Tänname teid 谢谢  
**Thank You** Tak  
 感謝您 Obrigado Teşekkür Ederiz  
 Σας ευχαριστούμε 감사합니다  
 ขอบคุณ  
 Bedankt Děkuje vám  
 ありがとうございます  
 Tack