

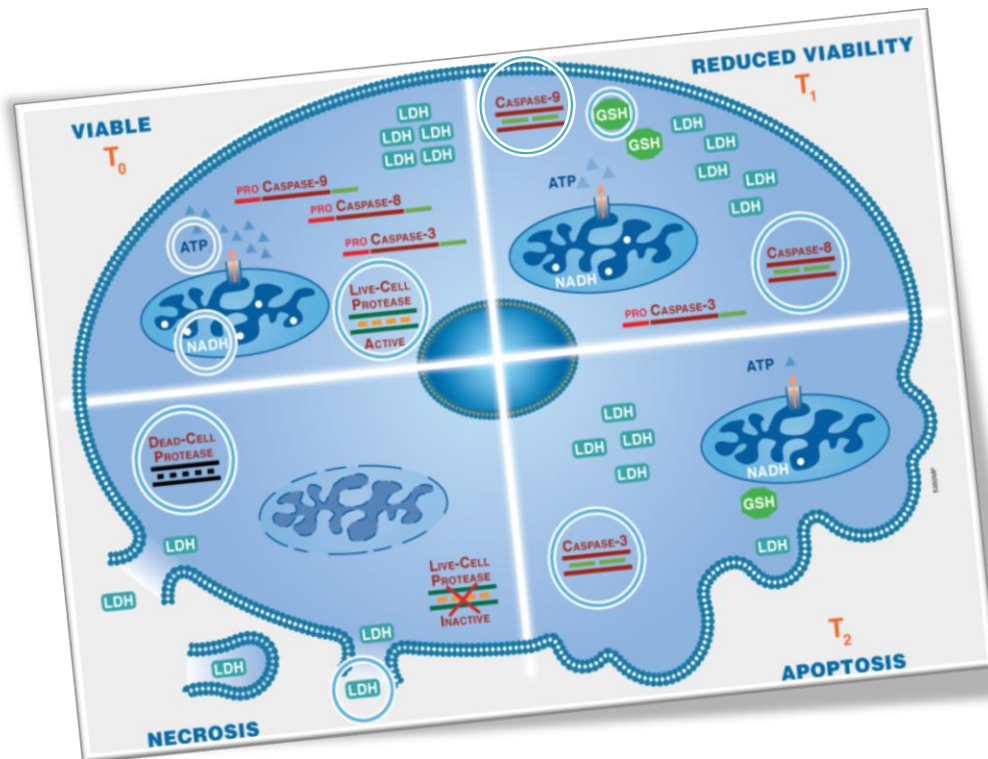
Multiplexing Cell-Based Assays: Get More Biologically Relevant Data

Fall 2010



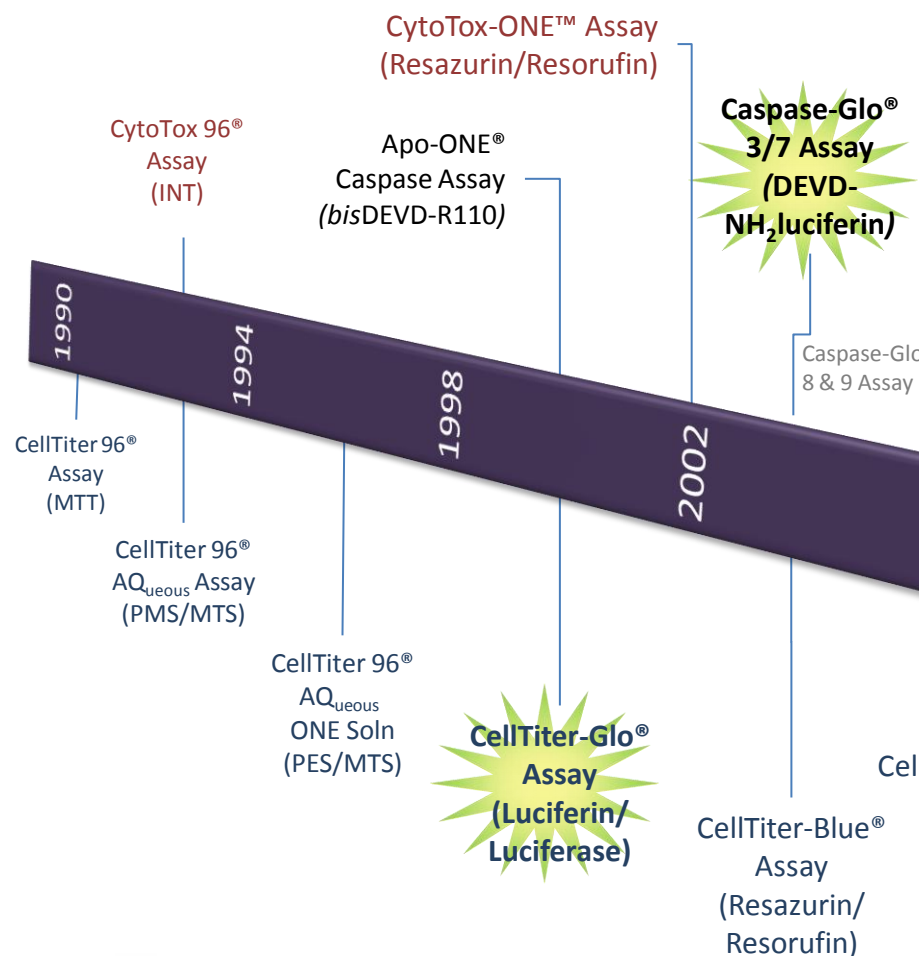
Click this icon to view
speakers notes for
each slide.

Multiplexing assays for more informative data



- Plate-based assays for viability, cytotoxicity and apoptosis measurement
- Using multiplex assays to understand cell death mechanism
- Monitoring cell response in multiple applications

Development timeline for cell-based viability, cytotoxicity and apoptosis assays

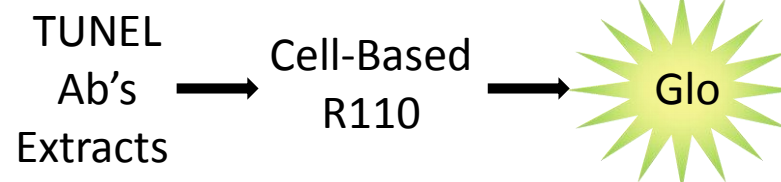


Ease-of-Use & Sensitivity

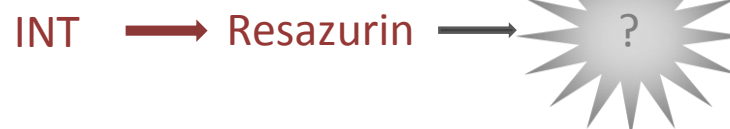
Viability Assays



Apoptosis Assays

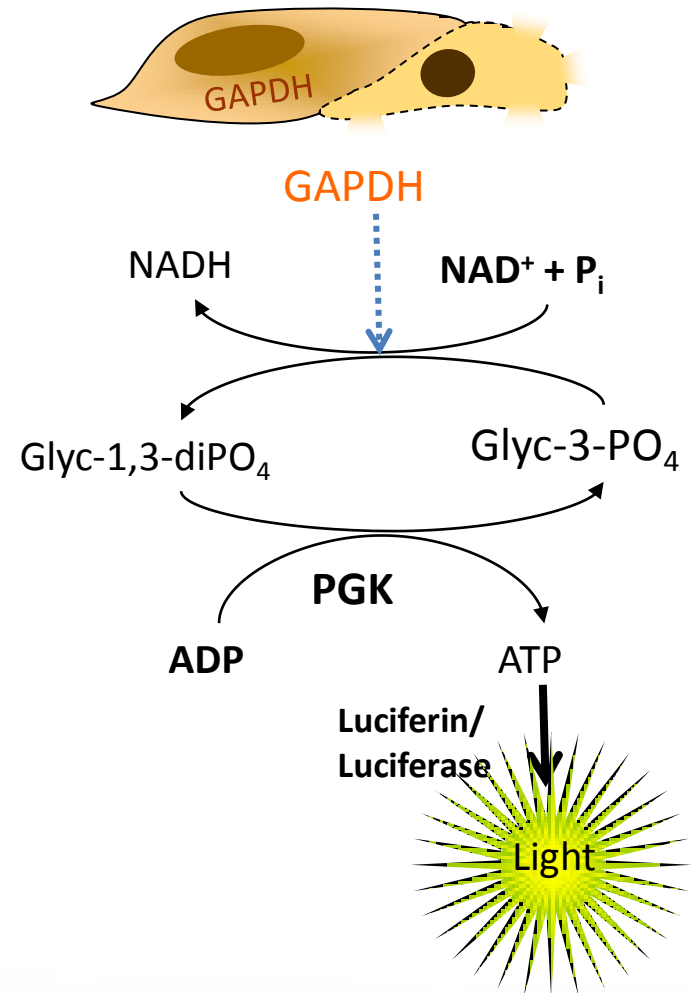
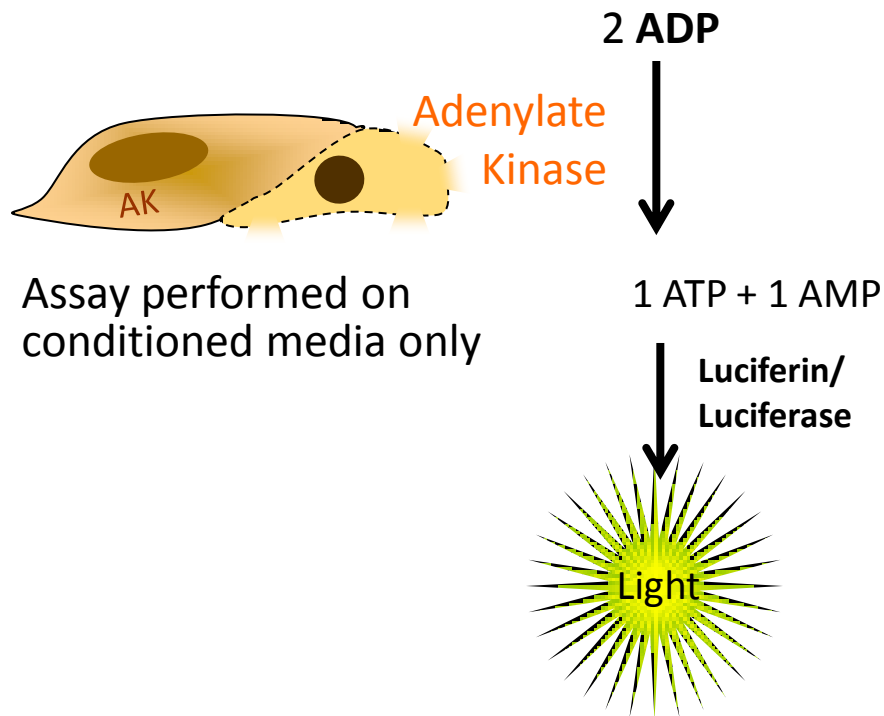


Cytotoxicity Assays



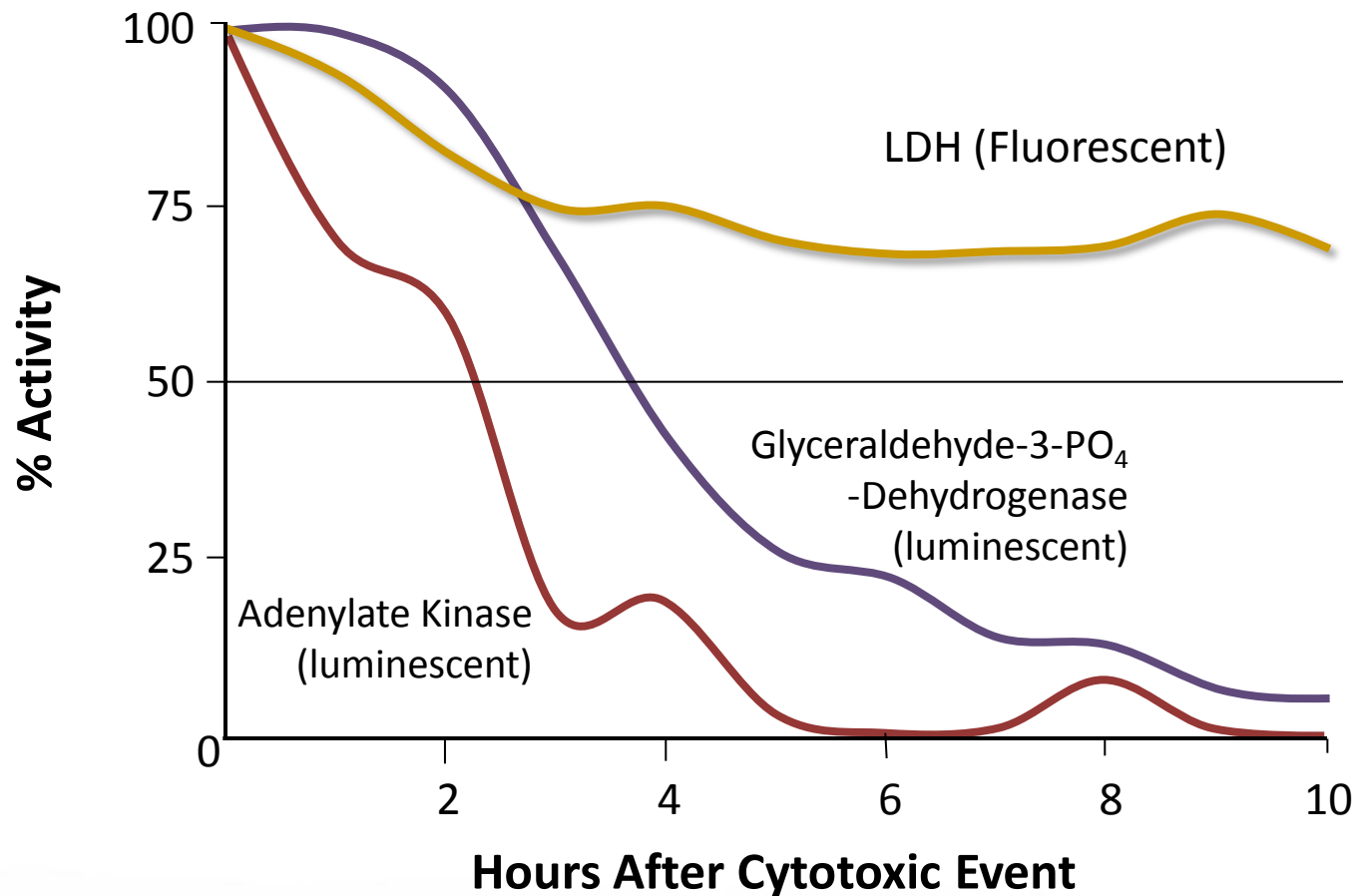


Other enzymatic markers of cytotoxicity





Luminescent methods where not ideal...



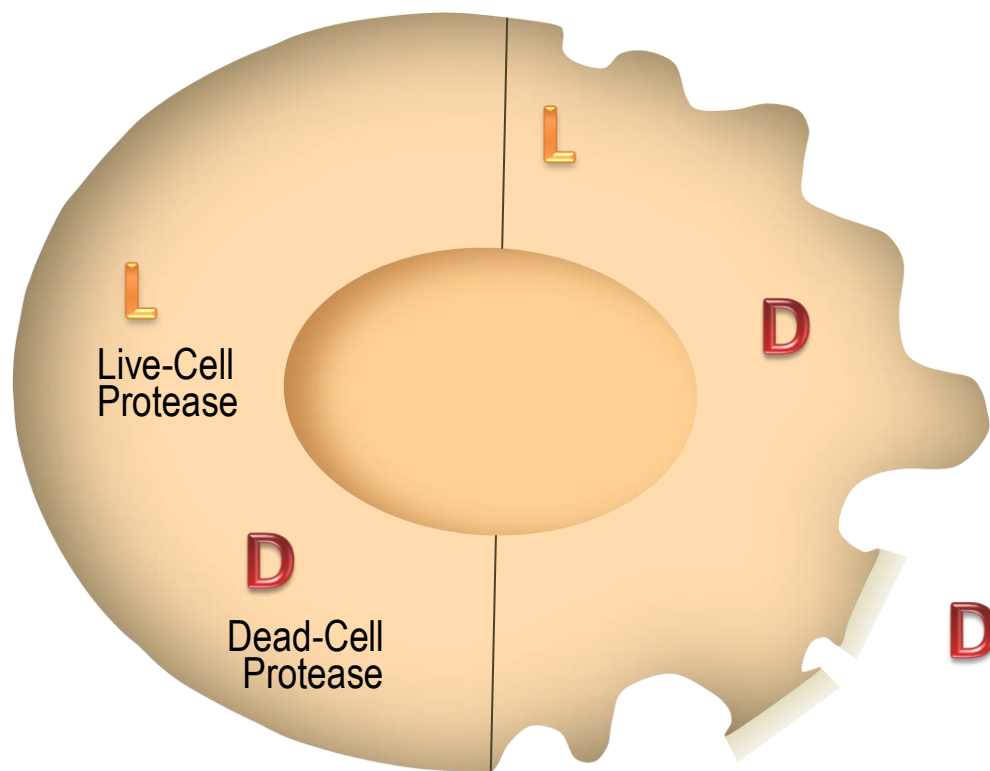


Two Protease activities = live/dead cell assay



Are my cells living?

Are my cells dying?



Live-Cell Protease quickly inactivated outside the cell.



Available online at www.sciencedirect.com

ScienceDirect

Analytical Biochemistry 366 (2007) 197–206

ANALYTICAL
BIOCHEMISTRY

www.elsevier.com/locate/yabio

A homogeneous assay to measure live and dead cells in the same sample by detecting different protease markers

Andrew L. Niles ^{a,*}, Richard A. Moravec ^a, P. Eric Hesselberth ^b,
Michael A. Scurrey ^b, William J. Daily ^b, Terry L. Riss ^a

^a Promega Corp., Madison, WI 53711, USA

^b Promega Bioscience, San Luis Obispo, CA 93401, USA

Received 15 January 2007

Available online 12 April 2007

Abstract

A method to simultaneously determine the relative numbers of live and dead cells in culture by introducing a combination of two fluorogenic substrates or a fluorogenic and a luminescent protease substrate into the sample is described. The method is based on detection of differential ubiquitous proteolytic activities associated with intact viable cells and cells that have lost membrane integrity. A cell-permeable peptide amine-fluorogenic substrate detects protease activity restricted to intact viable cells. Upon cell death, the viable cell protease marker becomes inactive. An impermeable peptide (chondroitinase 110 (or aminoluciferin) conjugated substrate detects protease activity from nonviable cells that have lost membrane integrity. The multiplex assay can detect 200 dead cells in a population of 10,000 viable cells. The protease substrate reagents do not damage viable cells over the course of the assay; thus the method can be multiplexed further with other assays in a homogeneous format. Ratios of measurement of viable and dead cells in the same sample provides an internal control that can be used to normalize data from other cell-based assays.

© 2007 Elsevier Inc. All rights reserved.

Keywords: Cell-based; Cytotoxicity; Viability; Multiplex; Fluorescence; Luminescence; Protease; Apoptosis; Homogeneous; High throughput

It is an important and necessary experimental practice to determine the viability of cells in culture after chemical, biological, or physical treatment and manipulation. Maintenance of membrane integrity is a common criterion for cell viability. Measurable changes in membrane permeability include trypan blue exclusion, nucleic acid staining, and ³Cr or lactate dehydrogenase release [1,2]. Conversely, measures of viability by metabolic capacity include tritiated thymidine incorporation, ATP content, tetrazolium dye reduction, and fluorescein diacetate labeling [3].

These existing techniques have a number of technical or practical drawbacks which limit their utility in multiplexed or high-throughput formats. For instance, cellular ³Cr release assays require significant prelabeling preparation, and all assays utilizing radiological tracers or mutagens/

teratogenic dyes impose significant exposure, handling, and disposal issues. In addition, tetrazolium or resazurin chemistries can significantly complicate additional downstream applications by color quenching of fluorescence or luminescence.

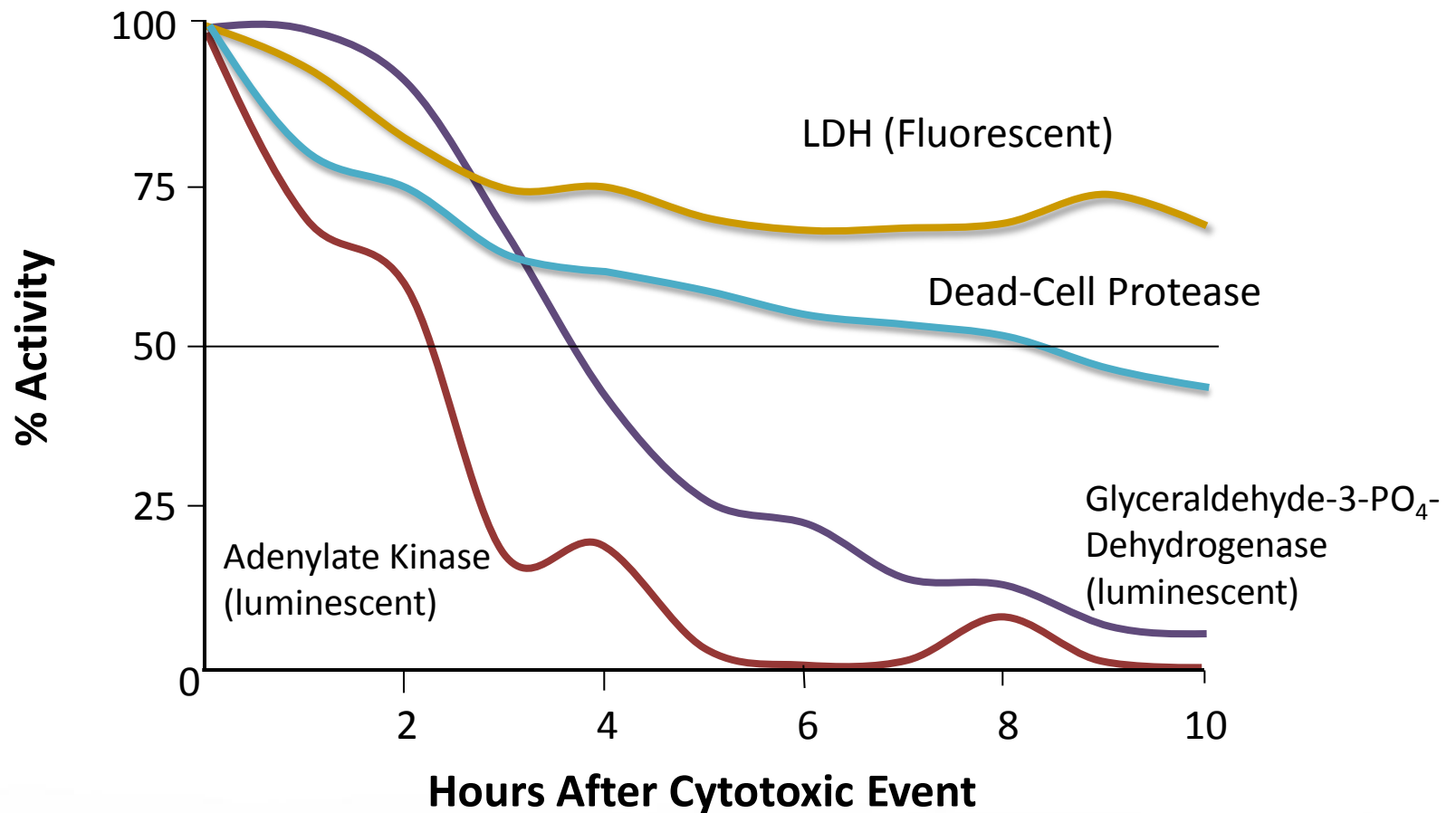
Cultured mammalian cells contain a rich milieu of protease, esterase, lipase, and nuclease activities which contribute to homeostatic maintenance. In particular, cytosolic, lysosomal, and transmembrane-bound proteases are involved in intracellular protein degradation, generation of immunogenic peptides, posttranslational modification, and cell division [4–6]. The activity of these enzymes is regulated by various mechanisms including specialized compartmentalization [7]. In response to extreme stress, environmental adversity, or committed progression of the apoptotic program, a loss of subcellular structure or membrane integrity is observed [8,9]. Therefore, we hypothesized that the release of stable protease activity into the

* Corresponding author. Fax: +1 608 286 4818.
E-mail address: andrew.niles@promega.com (A.L. Niles).

Niles, A.L., et al. (2007)
Analytical Biochemistry
366, 197–206.

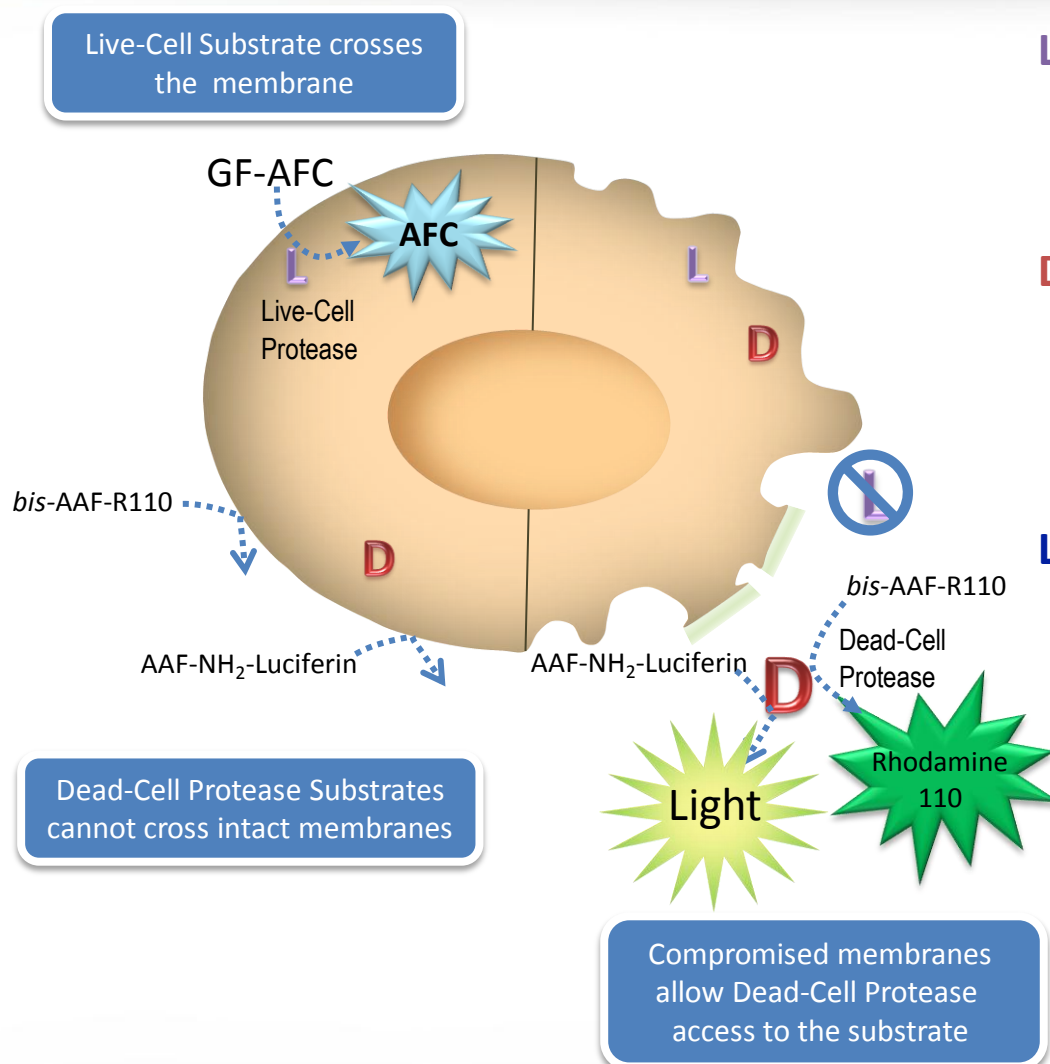


Dead-Cell Protease more like LDH





Measure Live, Dead or Both



Live-Cell Protease Assay

CellTiter-Fluor™ Cell Viability Assay

Dead-Cell Protease Assay

CytoTox-Fluor™ Cytotoxicity Assay

CytoTox-Glo™ Cytotoxicity Assay

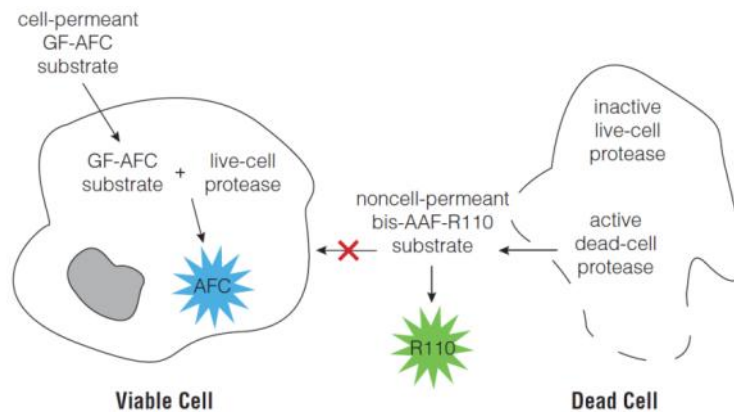
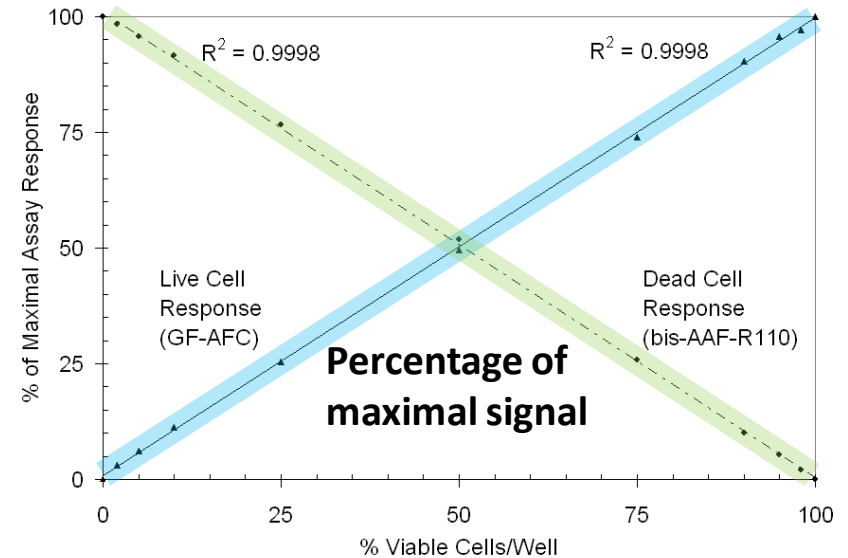
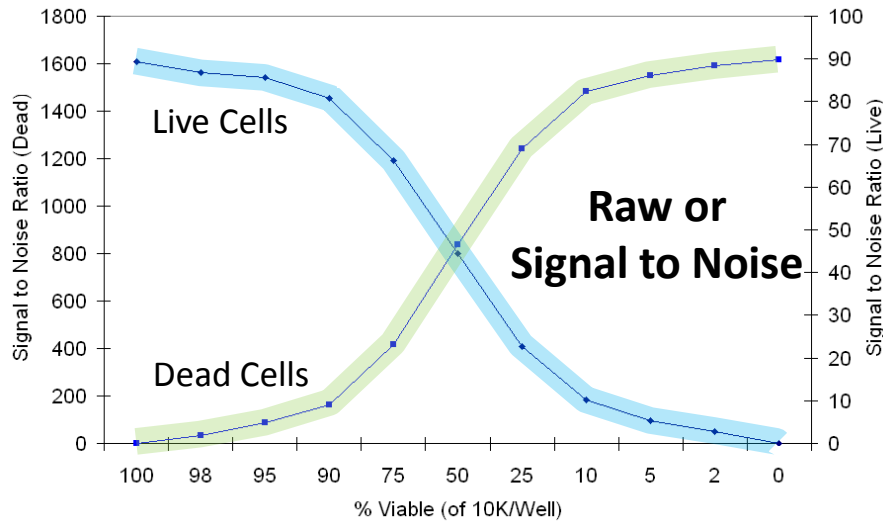
Live- & Dead-Cell Assay

MultiTox-Fluor Multiplex Cytotoxicity Assay

MultiTox-Glo Multiplex Cytotoxicity Assay



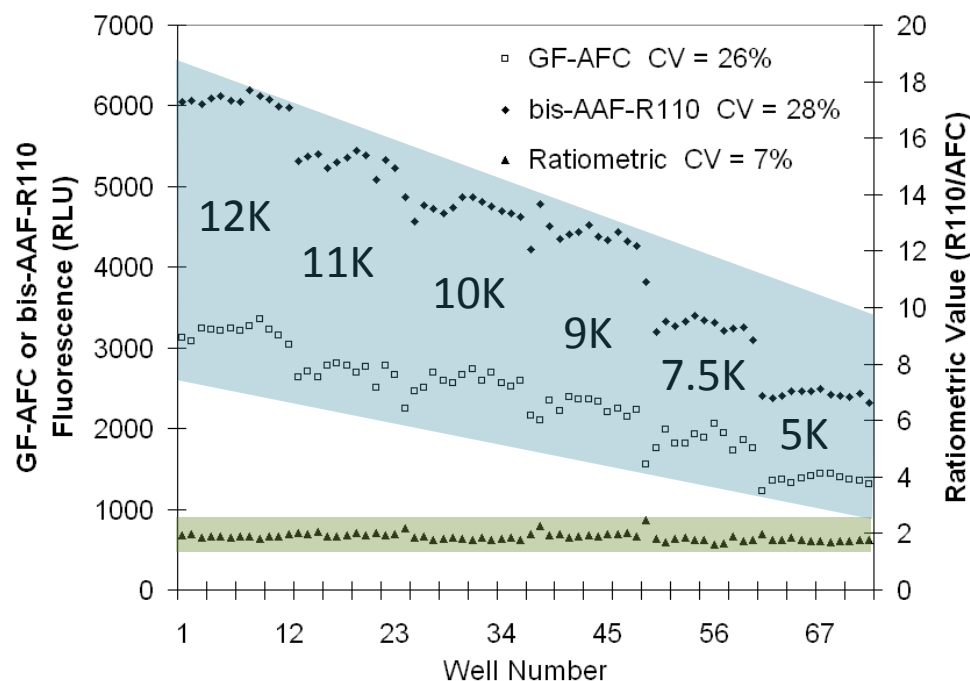
Inverse relationship between live & dead cell signals





Ratiometric measures address variability

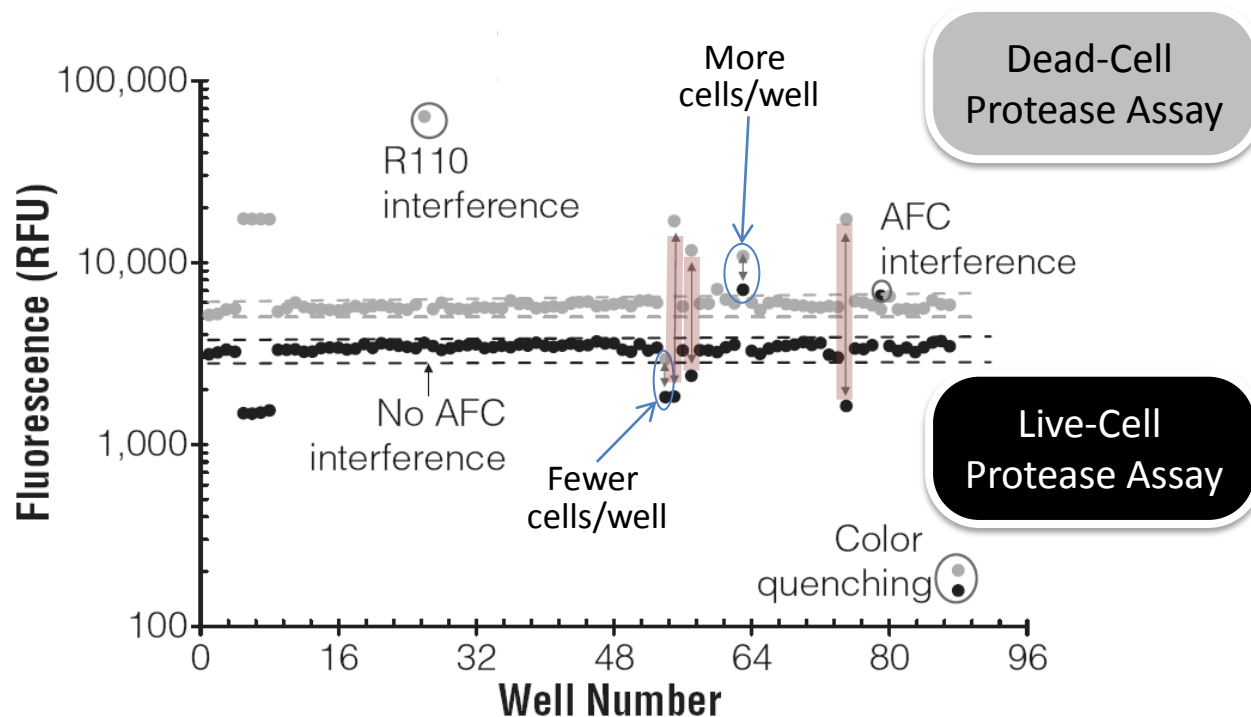
MultiTox-Fluor Assay Data



Variable number of 50% viability cells
plated per well.

- Single parameter responses are partially dependent on cell number
- Subtle clumping or pipetting error can make screens difficult to interpret
- Ratiometric measures decrease variation by normalizing the data

MultiTox-Fluor assay improves data confidence for cytotoxicity screens



A cytotoxic event must yield an **increase** in dead-cell protease activity and a **decrease** in live-cell protease activity



MultiTox-Fluor can be the perfect multiplexing partner



Assays must be chemically & biologically compatible

- Signals must be spectrally distinct (Fluorescence or Luminescence)
- Assay chemistries must be compatible
- The assays must fit in the available volume of the well or be separable.

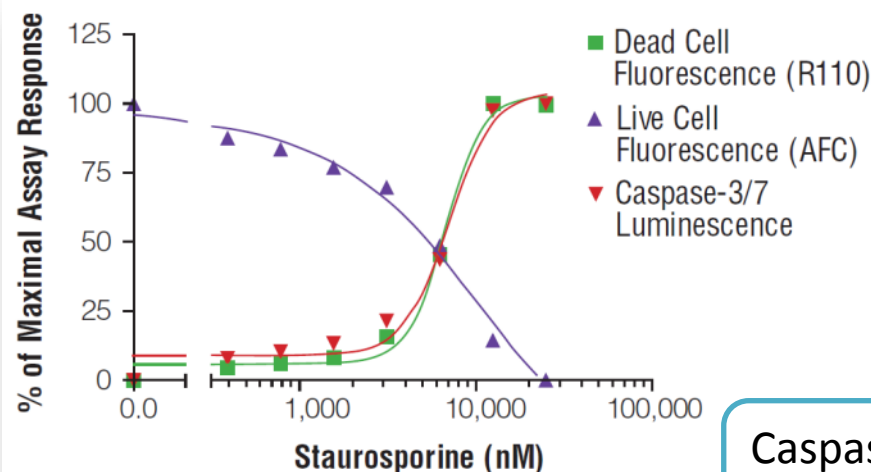
Multi-Tox Fluor
Non-Lytic
2 data points

Lytic Luminescent Assay





Multiplexing with Caspase-Glo[®] 3/7 Assay



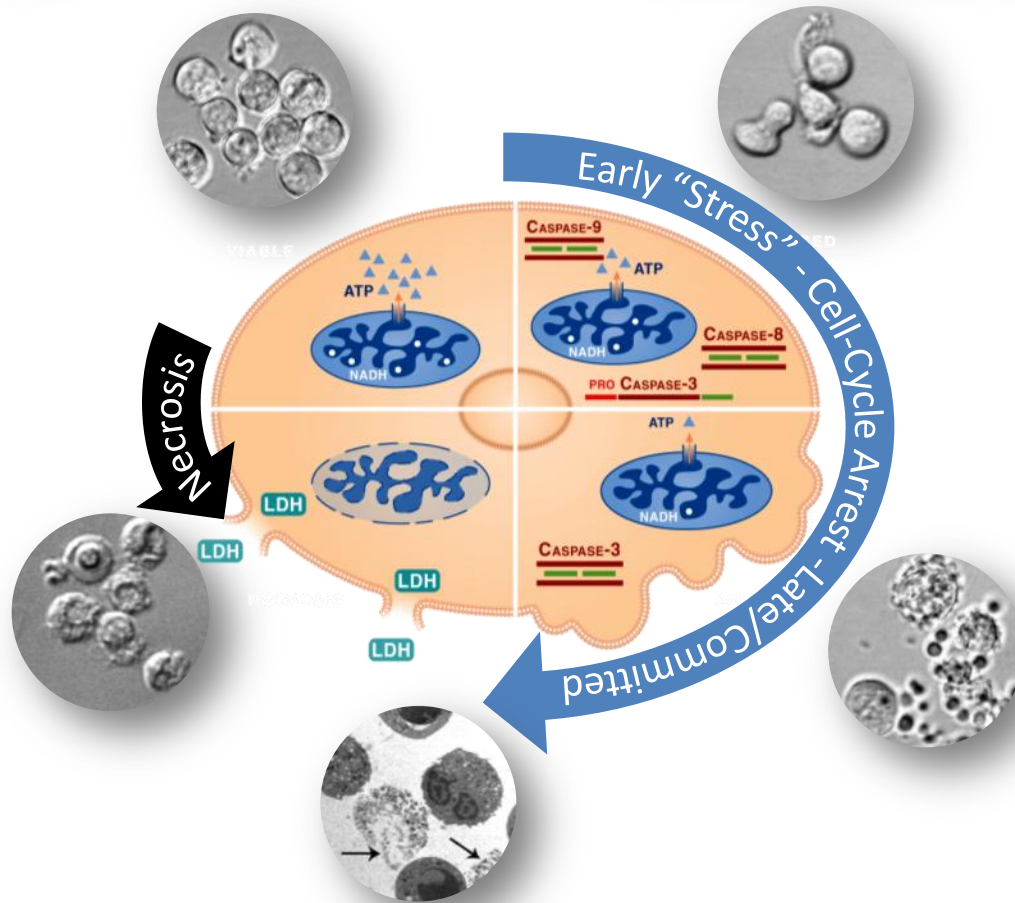
MultiTox-Fluor Multiplex Cytotoxicity Assay matches well with the Caspase-Glo[®] 3/7 Assay

- Used as the multiplexing example in the MultiTox-Fluor manual

Caspase-dependent apoptotic cell death

This combination can do so much more...

The Cytotoxicity Paradox: A Simple Concept with Inherent Biological Complexity



Did the treatment affect cell viability
-Yes/No?
-How?
-When?

How potent was the treatment?

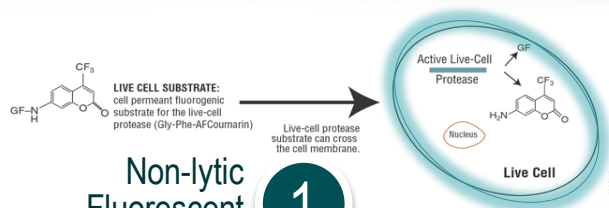
Is the treatment selective?

The cytotoxic phenotype is shaped by multiple factors:

1. Dosage
2. Exposure Time
3. Cellular susceptibility

No single parameter assay can fully characterize cytotoxicity

Deciphering a complicated process: Multiplexed, cytotoxicity signatures



Non-lytic
Fluorescent
Live Cell Assay

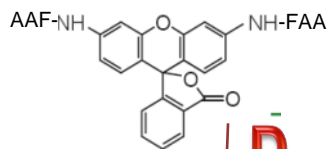
1

VIALE

REDUCED
VIABILITY

NECROSIS

APOPTOSIS



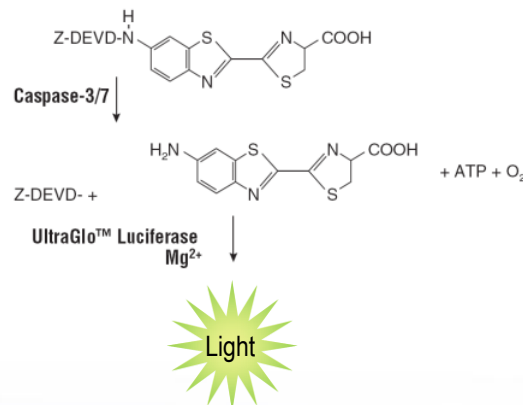
2

Non-Lytic
Fluorescent
Dead Cell Assay

Rhodamine
110

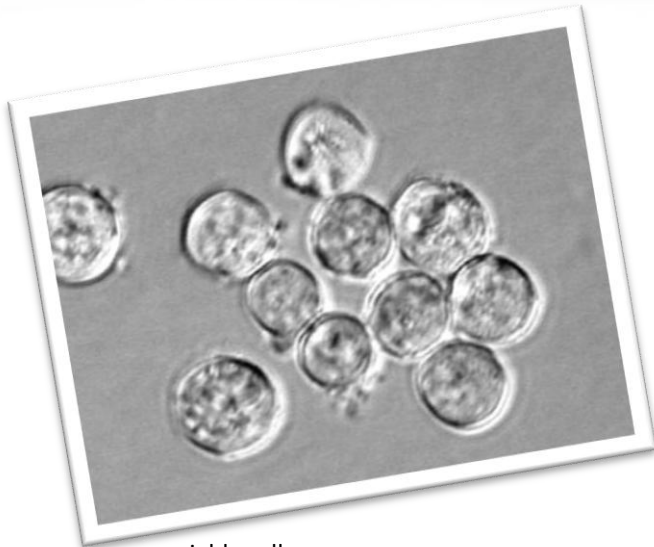
3

Lytic
Bioluminescent
Caspase-3/7
Assay

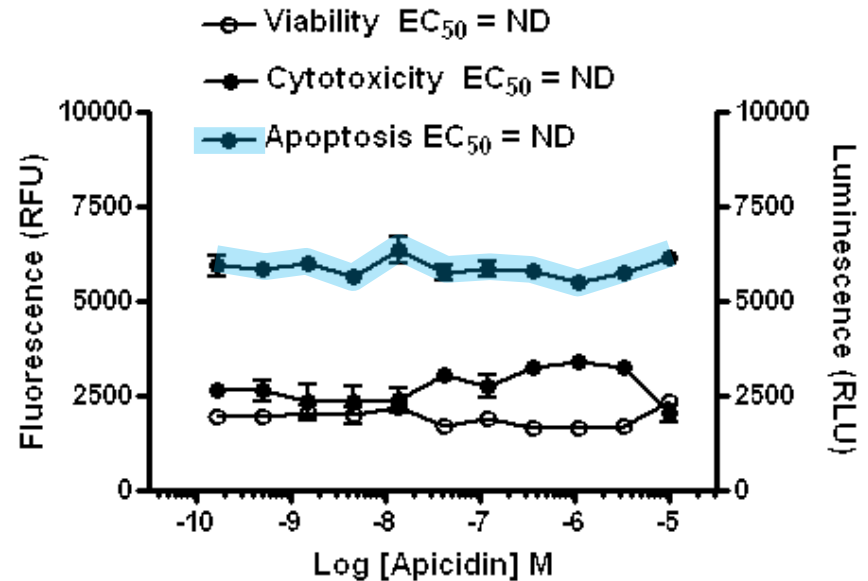




Signature #1. No Cytotoxic Effect



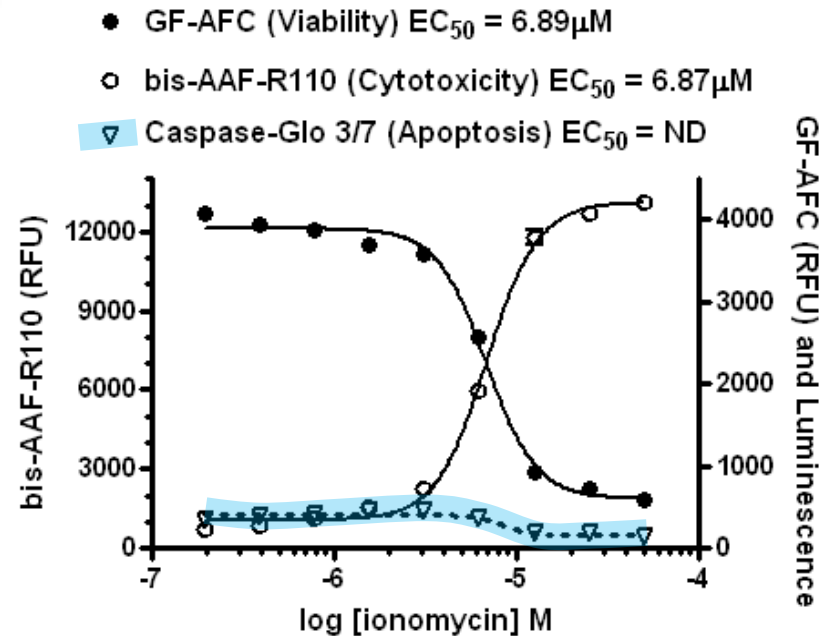
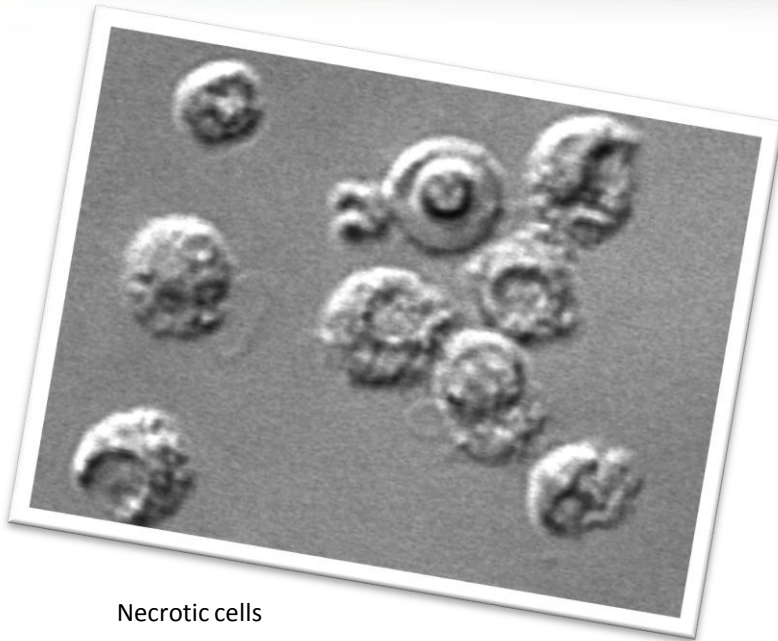
viable cells



Compound **exposure period** and **cell type** are critical parameter for establishing cellular inertness.

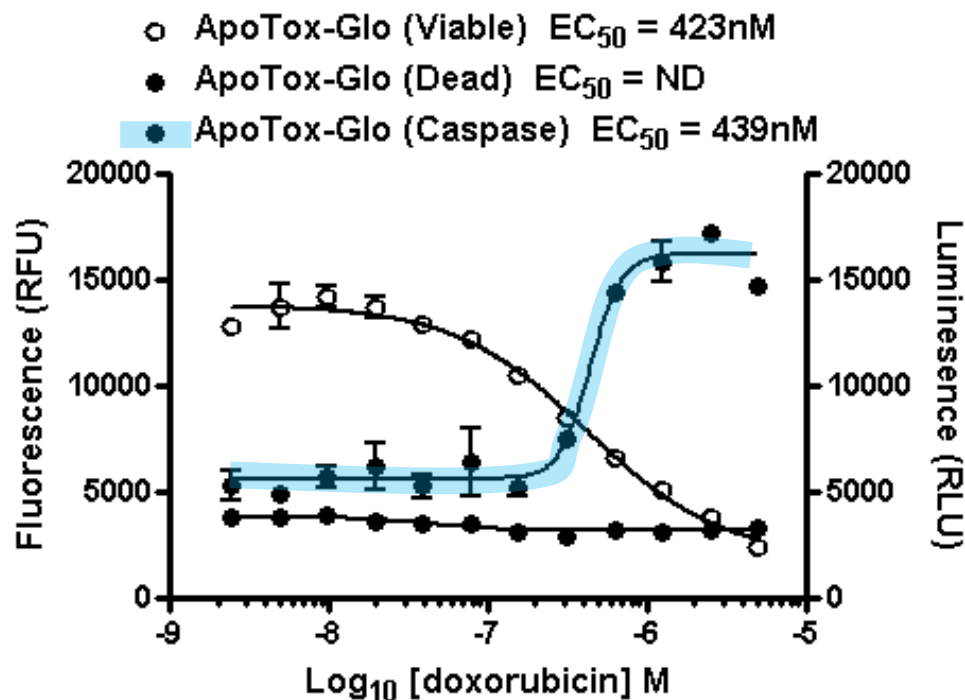
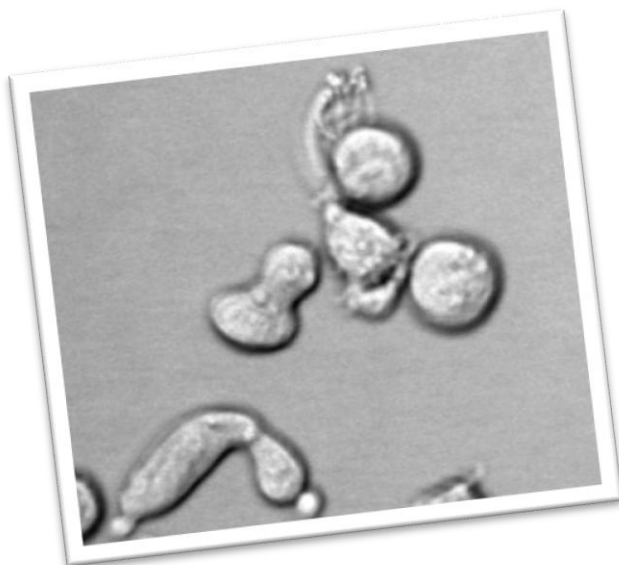


Signature #2. Primary Necrosis



Rapid loss of membrane integrity (<4hrs) **without caspase activation** is strongly indicative of primary necrosis.

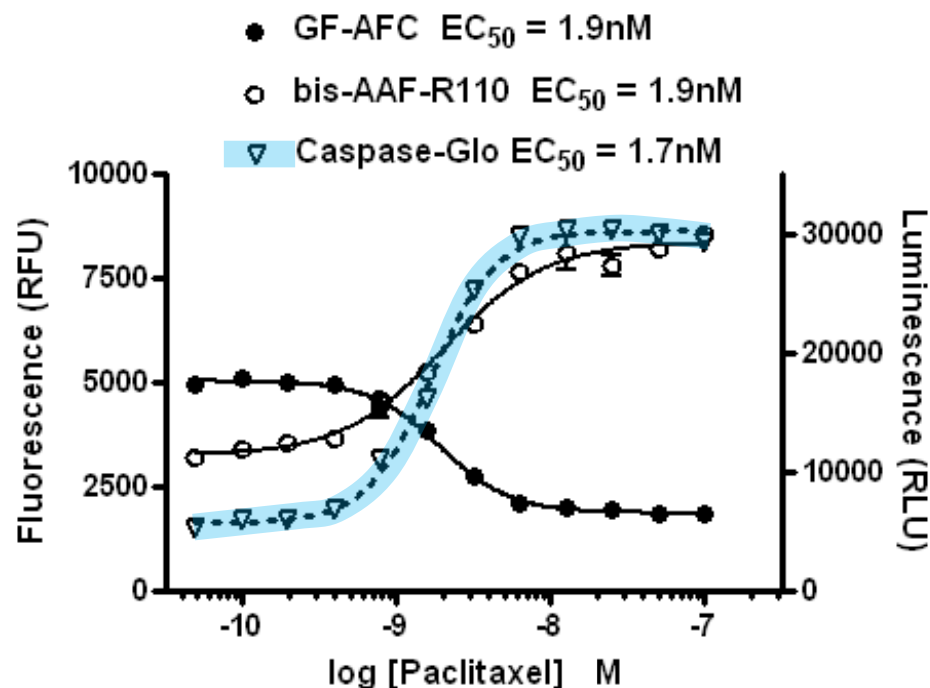
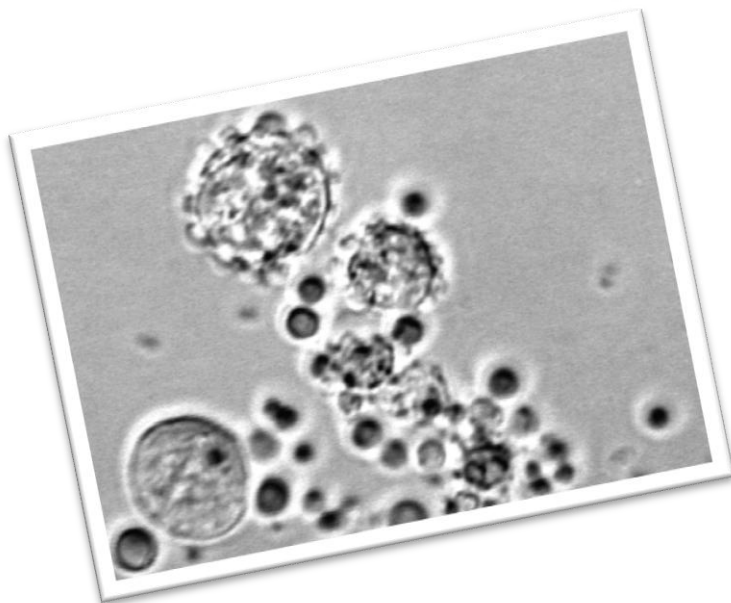
Signature #3 Cell Cycle Arrest...and early apoptosis



Decreases in apparent viability (viable cell number) with **increases** in caspase activation are consistent with cell-cycle arrest.



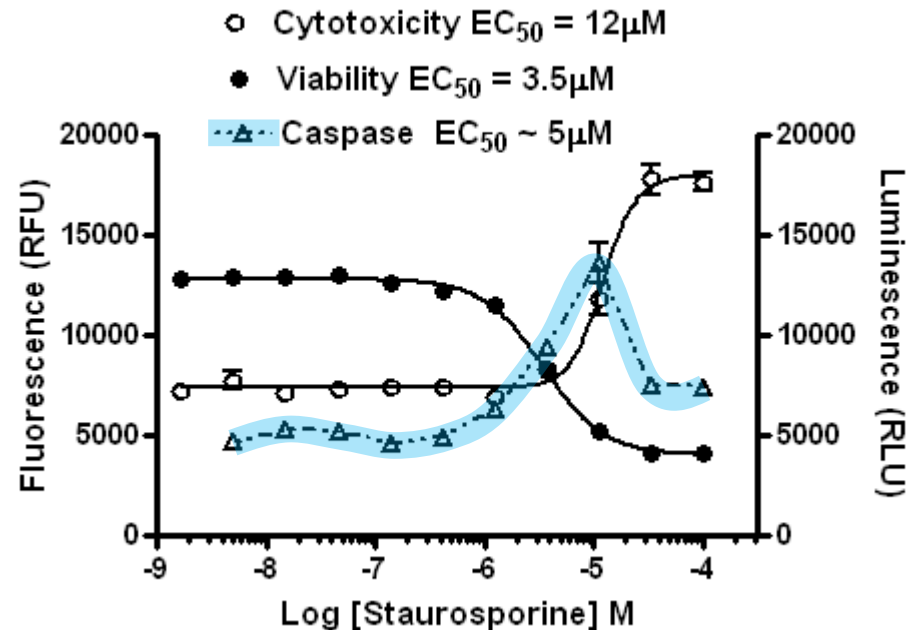
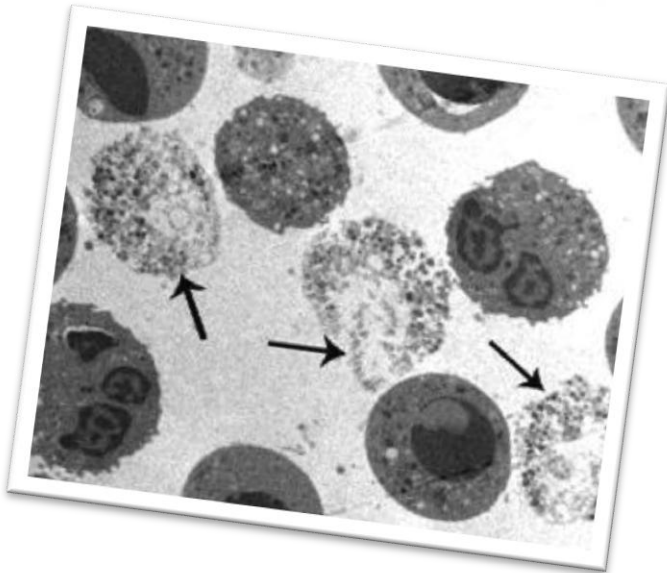
Signature #4: Apoptosis



Decreases in viability with a commensurate **increase** in cytotoxicity with caspase activation are consistent with **apoptosis and secondary necrosis**



Signature #5: Late State Apoptosis



Dose-dependent decrease in viability, increase in cytotoxicity with **caspase biomarker degradation** at highest concentrations is consistent with late stage apoptosis.



Does “Biological Relevance” Equate into Translational Relevance?



Translational Problem:

“Patients with [various cancers] experience poor outcomes, especially in metastasized disease, and treatment of all stages is associated with **strong side effects** [off-target] resulting in **impaired quality** of life. **Specific therapies** for such high-risk patients are therefore urgently needed to resolve this unsatisfactory situation.”

– Milde, T., et al. (2010) *Clinical Cancer Research* **16**, 3240-52.

Imaging, Diagnosis, Prognosis

**Clinical
Cancer
Research**

HDAC5 and HDAC9 in Medulloblastoma: Novel Markers for Risk Stratification and Role in Tumor Cell Growth

Till Milde¹, Ina Oehme¹, Andrey Korshunov^{2,5}, Annette Kopp-Schneider³, Marc Remke^{4,6}, Paul Northcott⁷, Hedwig E. Deubzer^{1,6}, Marco Lodrini^{1,6}, Michael D. Taylor⁷, Andreas von Deimling^{2,5}, Stefan Pfister^{4,6}, and Olaf Witt^{1,6}

Abstract

Purpose: Medulloblastomas are the most common malignant brain tumors in childhood. Survivors suffer from high morbidity because of therapy-related side effects. Thus, therapies targeting tumors in a specific manner with small molecules such as histone deacetylase (HDAC) inhibitors are urgently warranted. This study investigated the expression levels of individual human HDAC family members in primary medulloblastoma samples, their potential as risk stratification markers, and their roles in tumor cell growth.

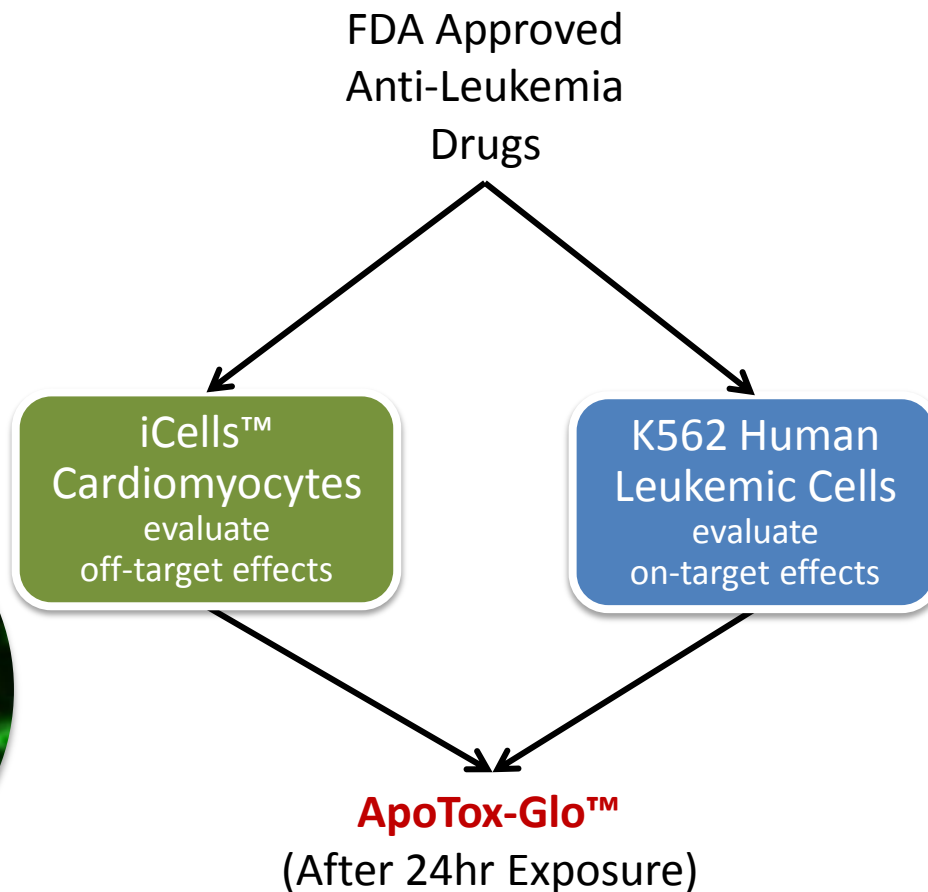
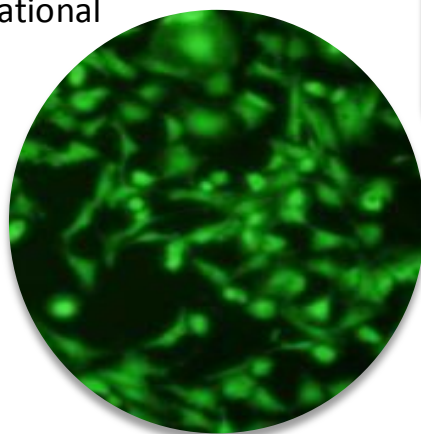


Potency & Safety Evaluation: Validation of ApoTox-Glo™ Assay with Clinical Cancer Therapeutics



"iCells™ are specifically designed to aid drug discovery and **improve the predictability of drug efficacy and toxicity** screens, weeding out ineffective and potentially toxic compounds early in the pharmaceutical pipeline process before significant time and resources have been invested."

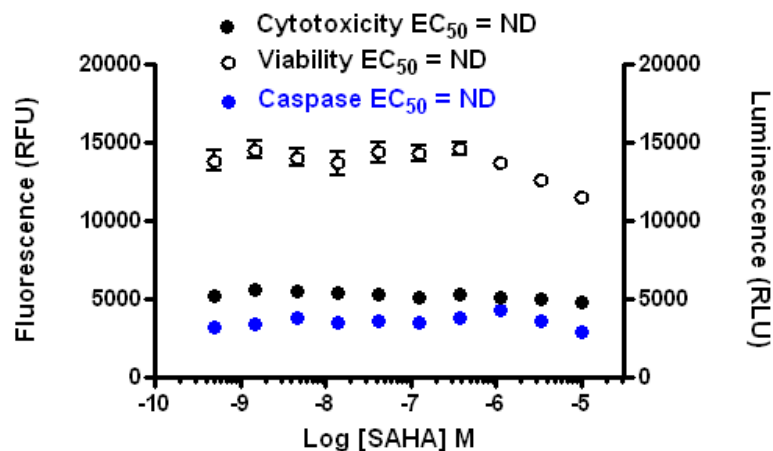
-Cellular Dynamics International





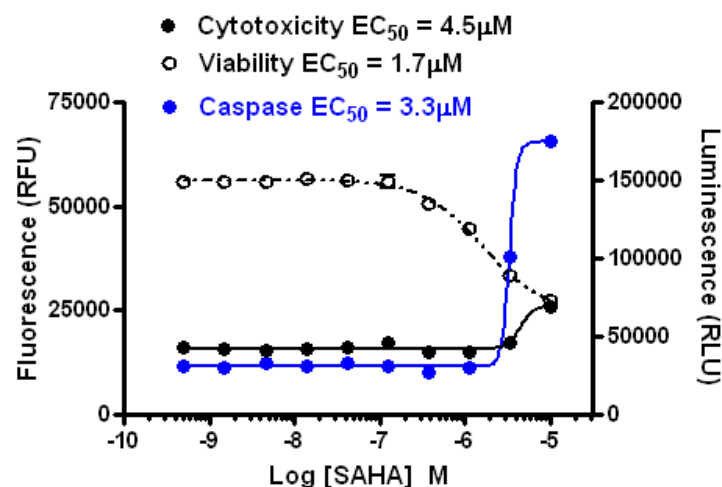
HDAC inhibition shows target specificity

iCell™



No apparent cytotoxicity or caspase activation.

K562



Cytotoxicity by apoptosis

Histone Deacetylase Inhibitor
SuberoylAnilide Hydroxamic Acid
(Vorinostat™)

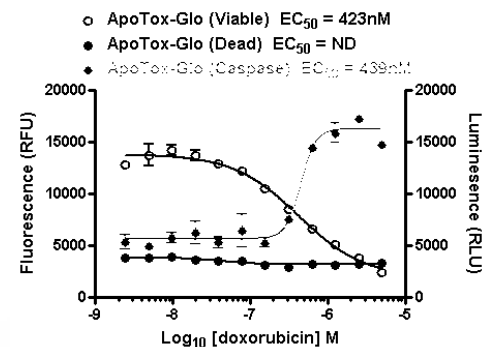
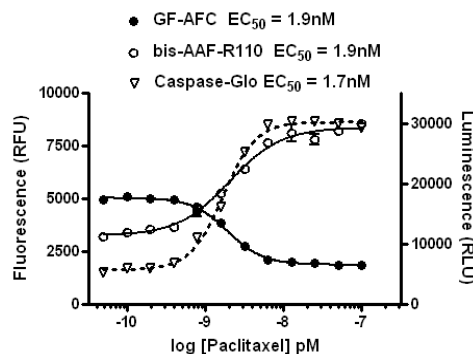
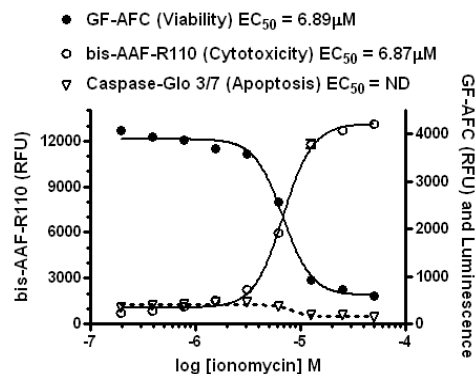
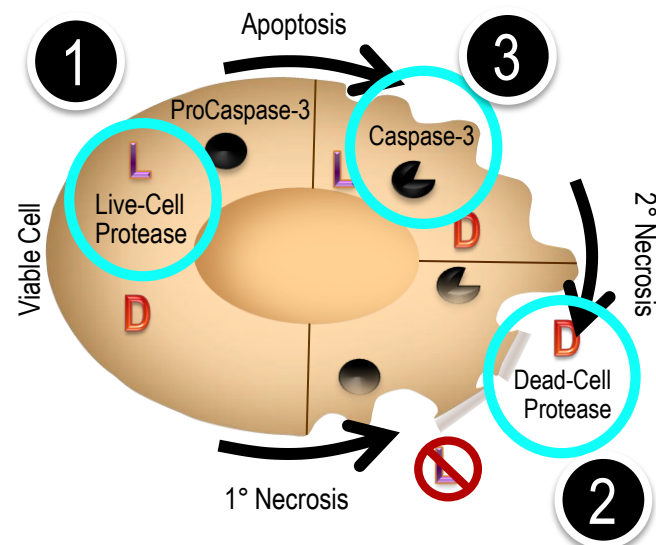
Determine Death Mechanism with ApoTox-Glo™ Triplex Assay



ApoTox-Glo Triplex Assay measures:

- Live Cells
- Dead Cells
- Apoptotic Cells

...and gives profile signatures



Need help?



Promega Technical Services



>150 years of combined lab experience

- Experienced & highly trained scientists
- Varied technical expertise
- Varied scientific expertise



Chat



eMail

techserv@promega.com



Phone

Promega Corporation - USA

Tel: 608-274-4330

Toll-Free: 800-356-9526

7:00am – 6:00pm CST

