

Plate-Based Assay Methods for the Assessment of Cellular Health

Andrew L. Niles, Senior Research Scientist

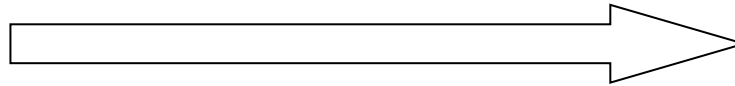


Biological Outcomes in Cell Culture

Treatment

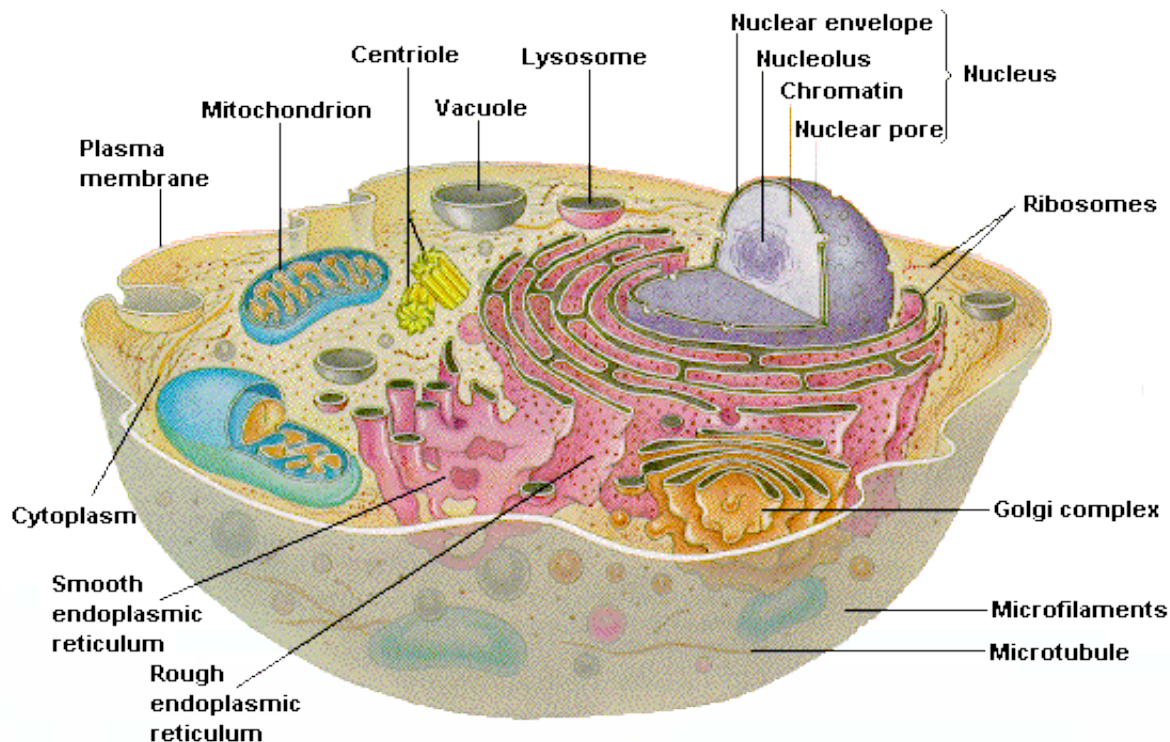
- Small molecule
- Bio-molecule
- Transgene
- Physical insult

Cause



Effect

- Normal Proliferation
- Enhanced Proliferation
- Cell-cycle Arrest
- Oxidative Stress
- 1° Necrosis
- Apoptosis
- 2nd Necrosis



Which response is it?

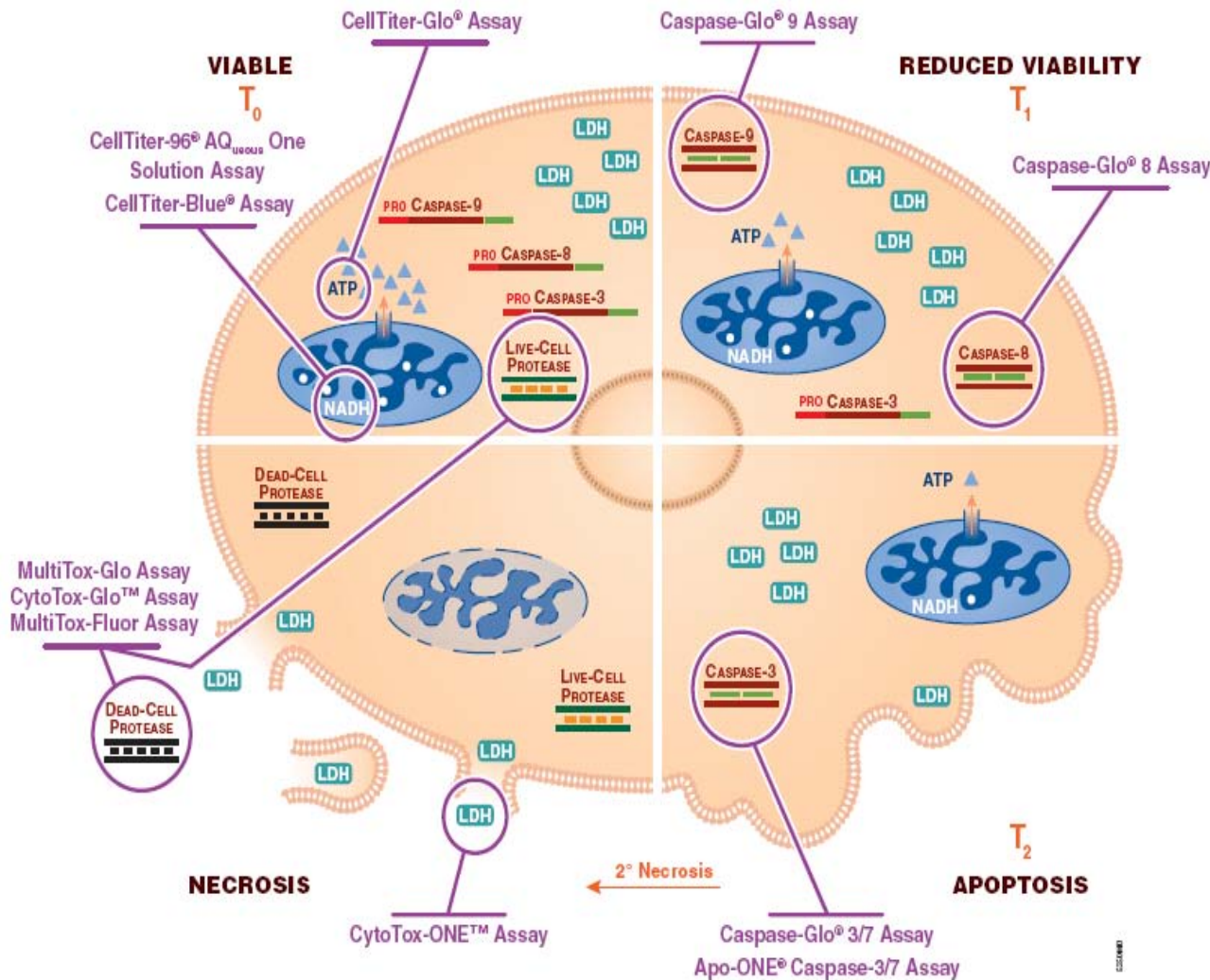


Rarely a simple answer. Often subject to experimental qualification.

The biological “**profile**” of any treatment is dependent upon:

1. Dosage
 - Addressed through serial dilution series
2. Exposure Time (cells with compound contact)
3. Mechanism of action of the test compound
4. Cell Type
 - specific target
 - off target

Biomarkers and the Cytotoxic Response



Biomarkers of cell health:

Can decrease due to cytotoxicity

Can increase due to cytotoxicity

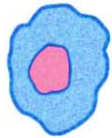
No single parameter assay can fully characterize cytotoxicity

Kinetics of Cell Death Affect Assay Results

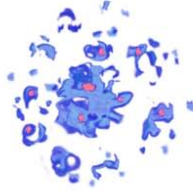
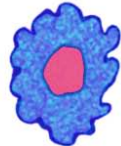


Time Zero 30 minutes to 4 hours 4 to 48 hours >48 hours

Apoptosis



Viable Cell



2° Necrosis

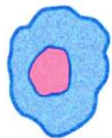
Enzyme Release	0	+	++++	+
Caspase	0	++	++++	0
ATP	++++	+++	+	0
Tetrazolium	++++	+++	+	0
Resazurin	++++	+++	+	0
Enzyme Retention	++++	+++	+	0

Choosing an appropriate biomarker and appropriate cell model is critical

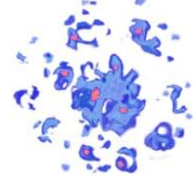
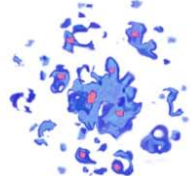
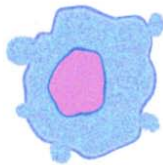
Primary Necrosis occurs quickly, apoptosis may take up to 48hrs to affect toxicity (secondary necrosis)

Cytotoxicity and caspase activities are transient but definitive

Necrosis



Viable Cell



Cell Debris

Enzyme Release	0	++++	++	0
Caspase	0	0	0	0
ATP	++++	0	0	0
Tetrazolium	++++	0	0	0
Resazurin	++++	0	0	0
Enzyme Retention	++++	0	0	0

Viability assays always report the number of viable cells remaining but offer little information about mechanism (anti-proliferative vs cytotoxic effects).

Tangible Examples of In Vitro Cytotoxicity



Typical Cell Health Work-Flow

Different mechanistic toxins

- serially diluted for potency calculations

+

Cell model(s)

- On target
- Off target

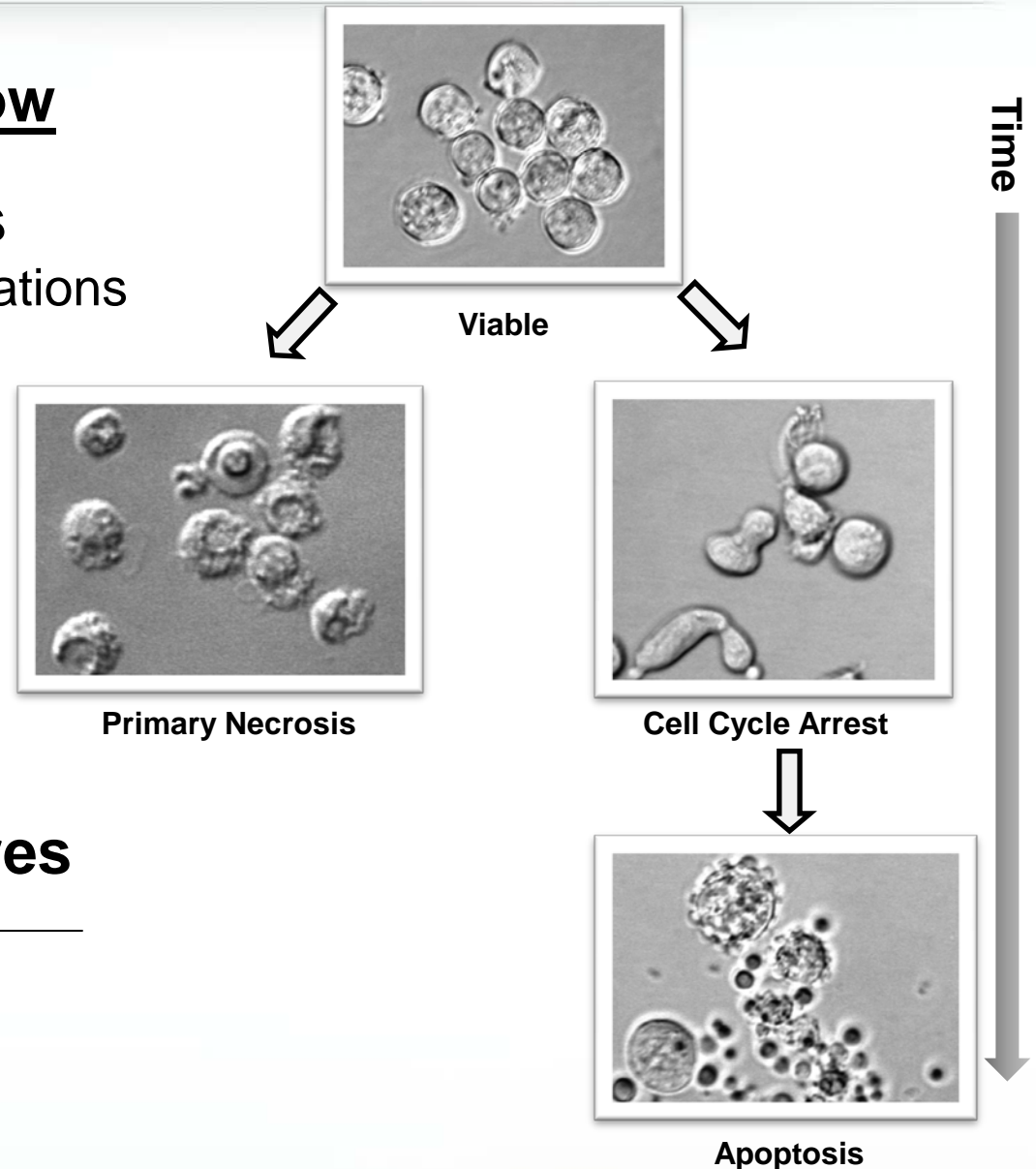
+

Time Course Exposures

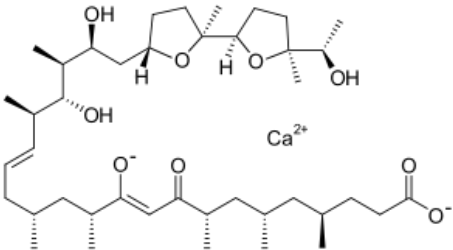
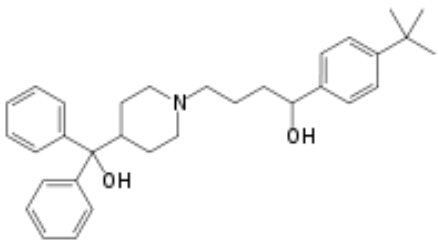
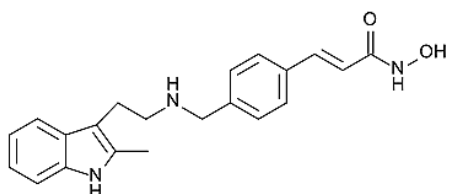
+

Different Cell Health Measures

= Data Interpretation(s)



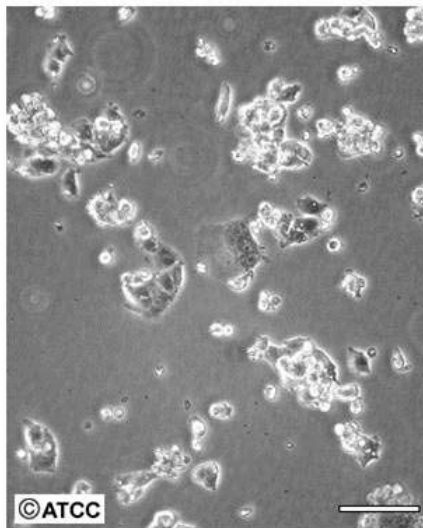
Three Model Toxins...

<u>Compound</u>	<u>Structure</u>	<u>Mode of Action</u>	<u>Cytotoxic Mechanism</u>
Ionomycin		Ionophore Ca ⁺⁺ flux	Primary necrosis
Terfenadine		Incompletely characterized pro-drug toxicity	Apoptosis (fast On-set)
Panobinostat		Histone deacetylase Inhibitor (HDACi)	Apoptosis (late On-set)

The Cell Model...

ATCC Number: **HB-8065**
Designation: **Hep G2**

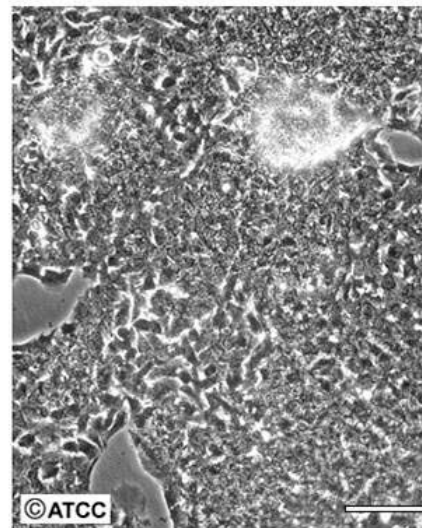
Untreated at “zero time”



Low Density

Scale Bar = 100µm

Untreated with
normal proliferation



High Density

Scale Bar = 100µm

HepG2 cells are a human hepatocellular carcinoma (epithelial morphology) commonly used as a model system for studies of liver metabolism and toxicity of xenobiotics.

The Experiment

Compounds diluted
in 10-fold dilutions of medium
and added in equal volumes
of sub-confluent HepG2 cells.
Vehicle control served as
“Untreated control”

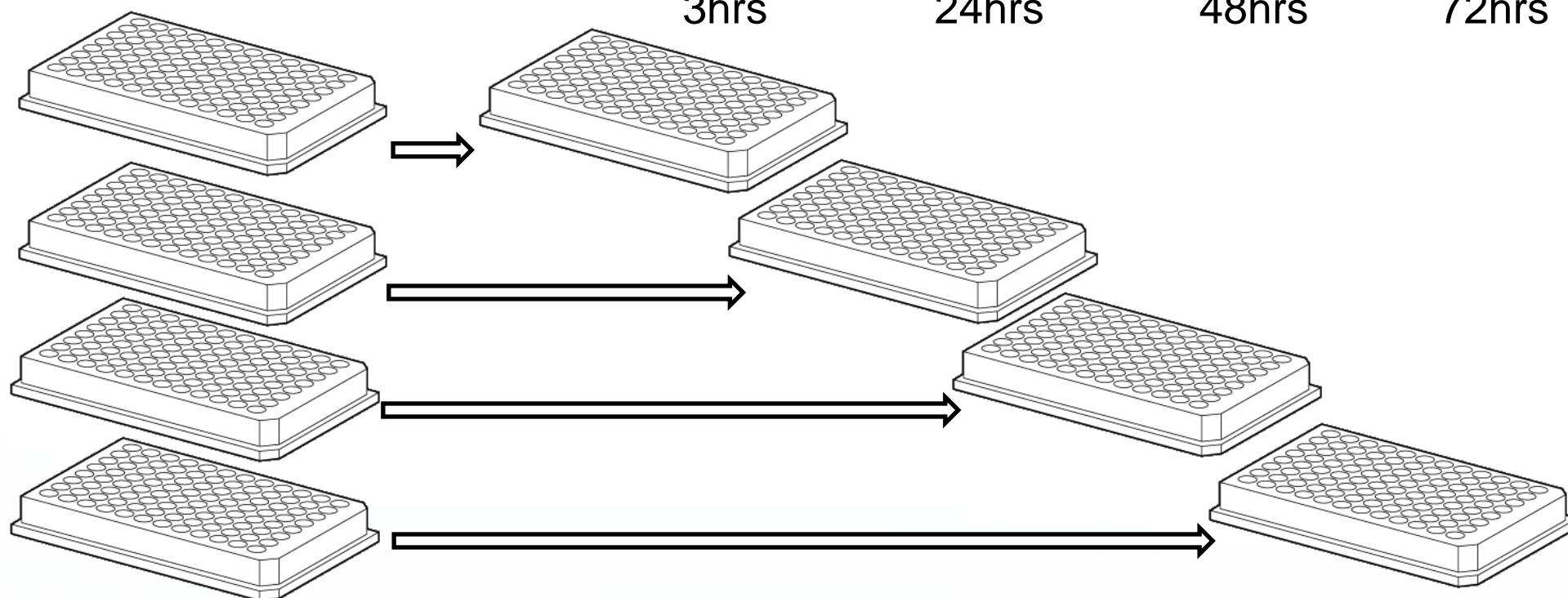
Hours of Compound Exposure with Cells

3hrs

24hrs

48hrs

72hrs



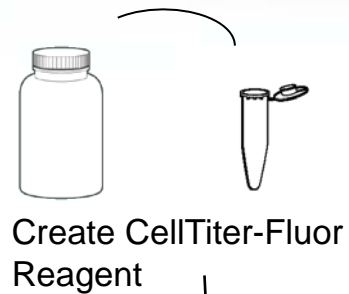
Three Model Cell Health Assays...



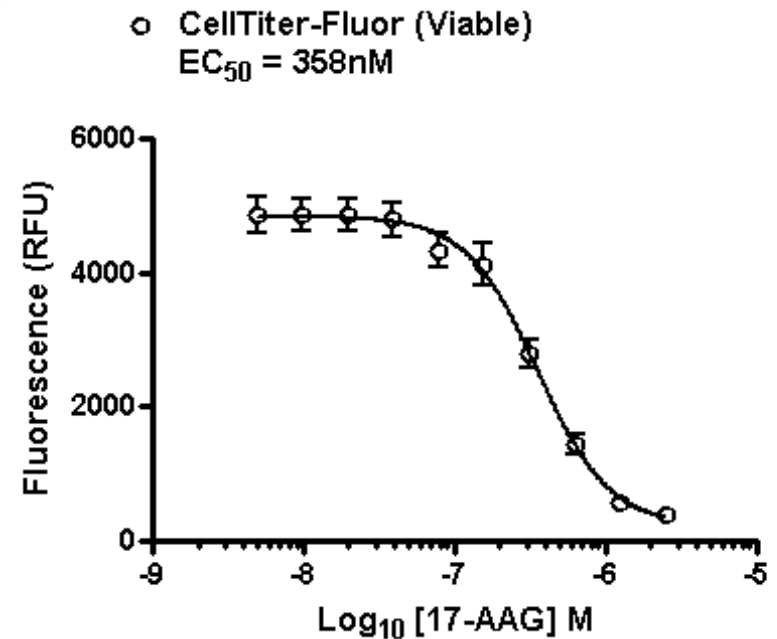
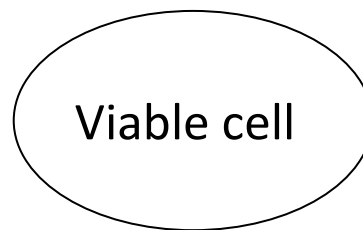
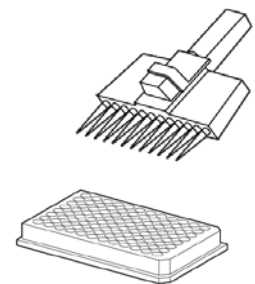
<u>Assay Name</u>	<u>Assay Type</u>	<u>Biomarker</u>	<u>Measurement</u>
CellTiter-Fluor™	Viability	Live Cell Protease	Fluorescence (AFC, 400 _{ex} /505 _{em})
CytoTox-Fluor™	Cytotoxicity	Dead Cell Protease	Fluorescence (R110, 485 _{ex} /530 _{em})
Caspase-Glo® 3/7	Cytotoxicity	Caspase Activity	Luminescence

Homogenous, “Add-Mix-Measure” Formats

CellTiter-Fluor™: A Protease Viability Assay

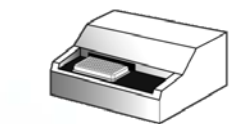


CellTiter-Fluor Reagent
(cell-permeable profluorogenic peptide substrate)



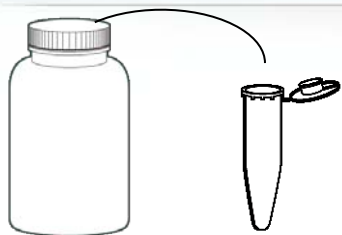
Advantages:

- Sensitive
- Scalable
- Compatible for multiplexing

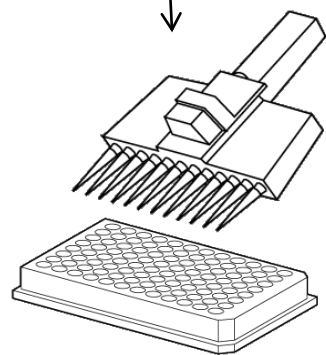


Constitutive
“live cell protease”
liberates AFC
fluorophore

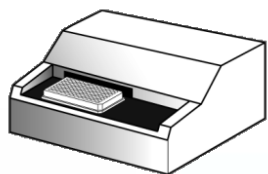
CytoTox-Fluor: Protease Cytotoxicity Assay



Create CytoTox-Fluor Reagent

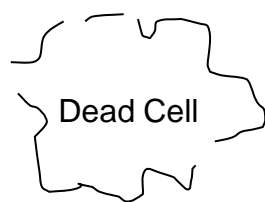


Incubate 30 min



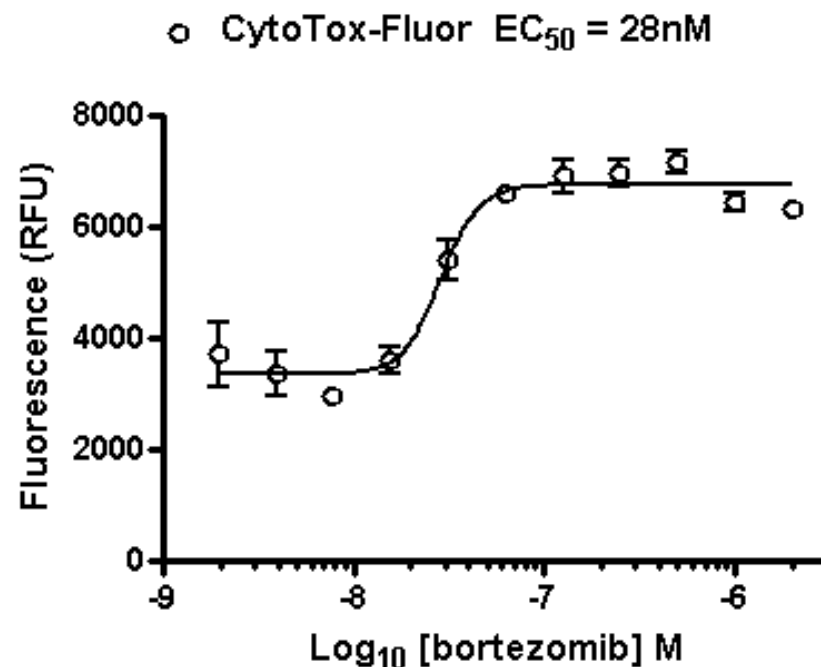
Fluorescence

CytoTox-Fluor Reagent
(non-cell permeable, profluorogenic Substrate)

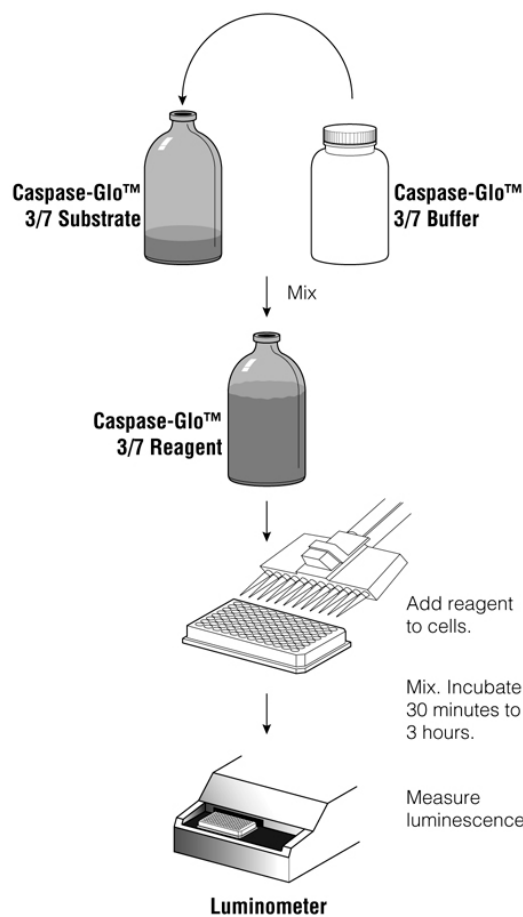


Leaked "Dead cell protease"

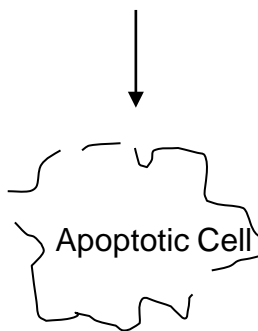
R110 is liberated
485ex/520em



Caspase Activity Assay

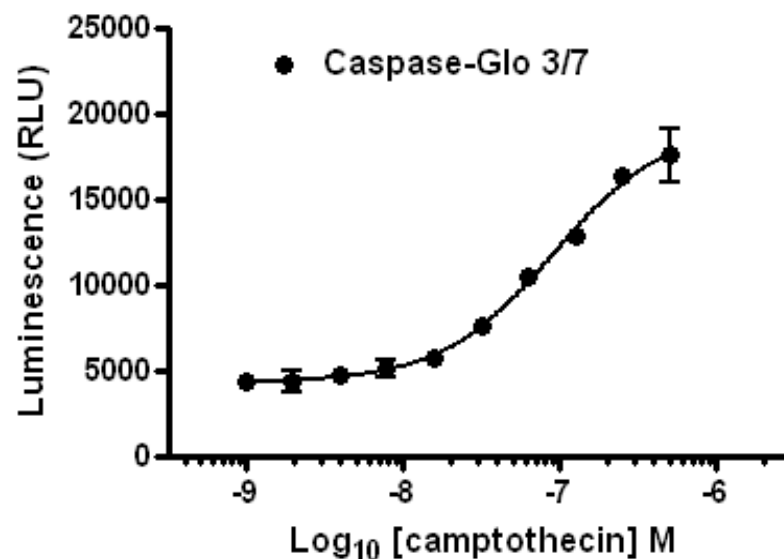


Caspase-Glo 3/7 Reagent
(Z-DEVD-luciferin + Ultra-Glo
+ ATP in lytic buffer)



Activated caspase
cleaves substrate

Aminoluciferin is liberated
and consumed by
luciferase reaction

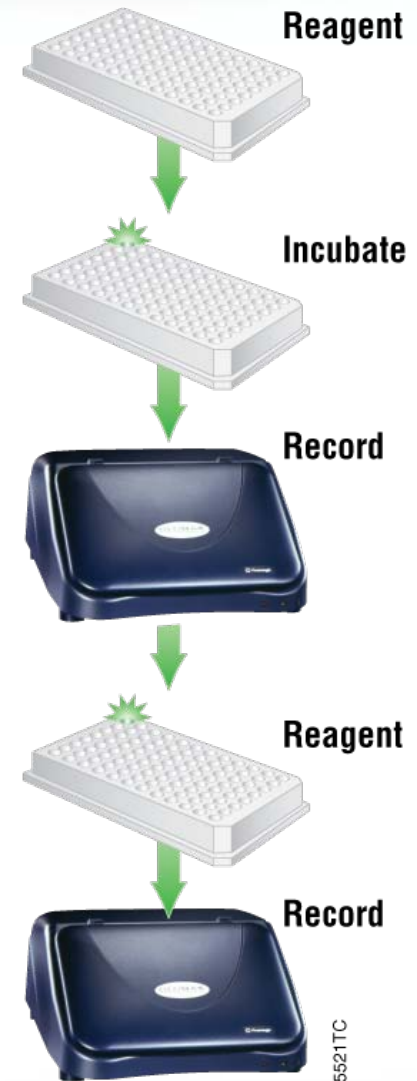
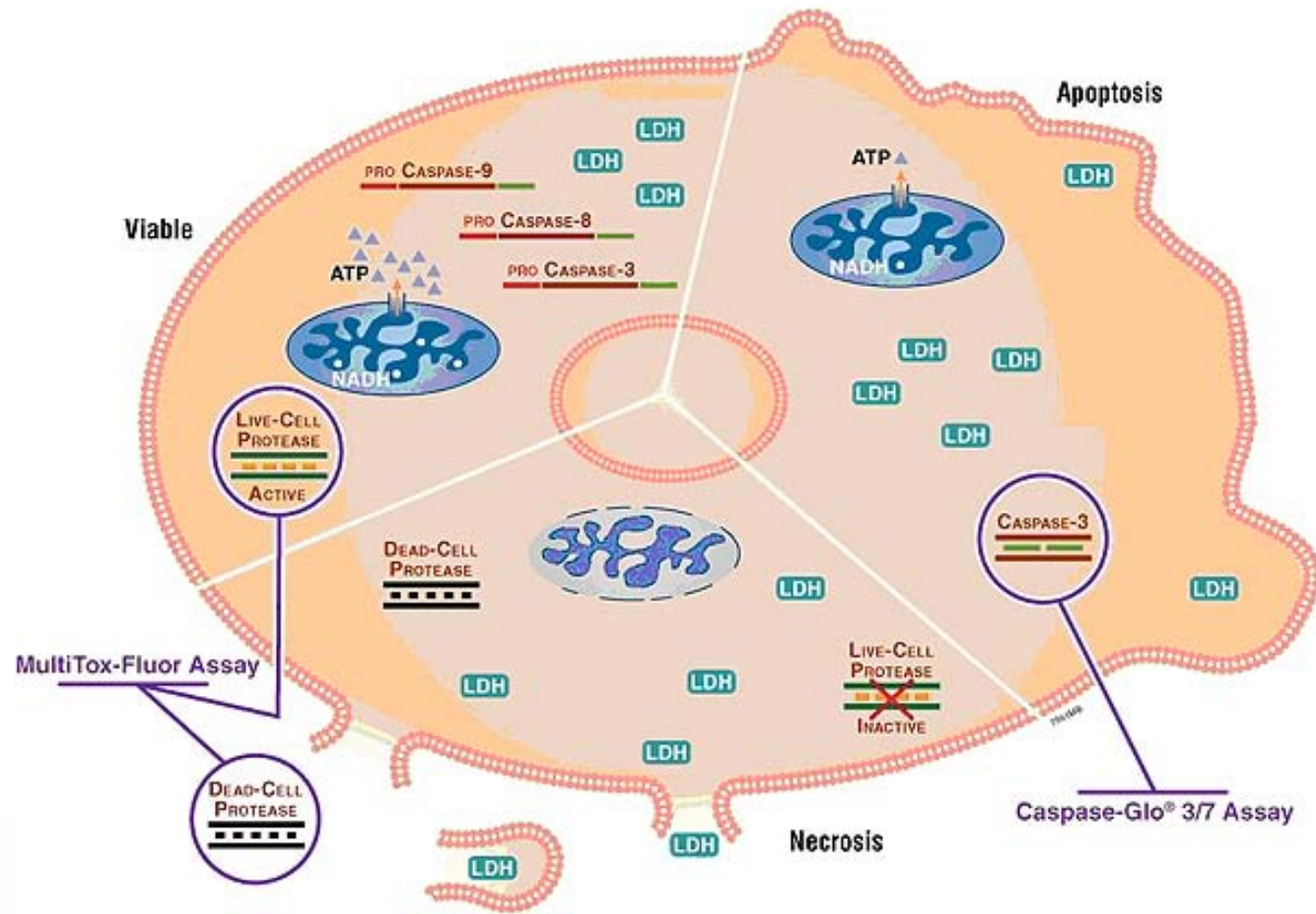


4055MA03_3A

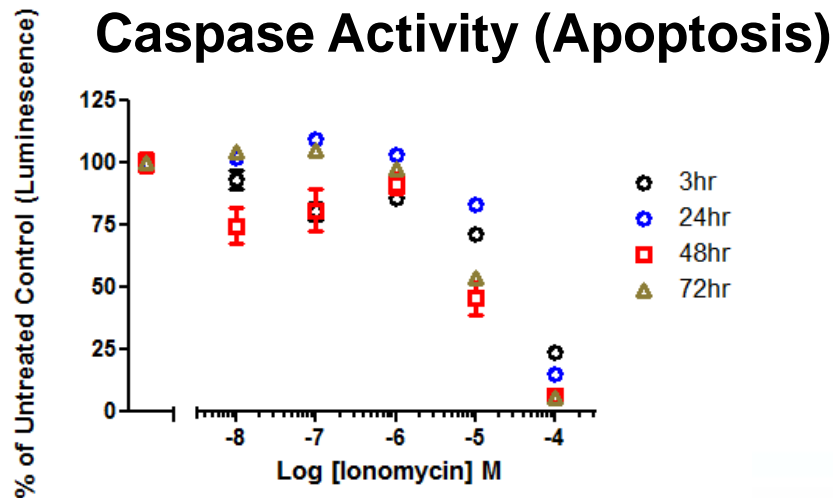
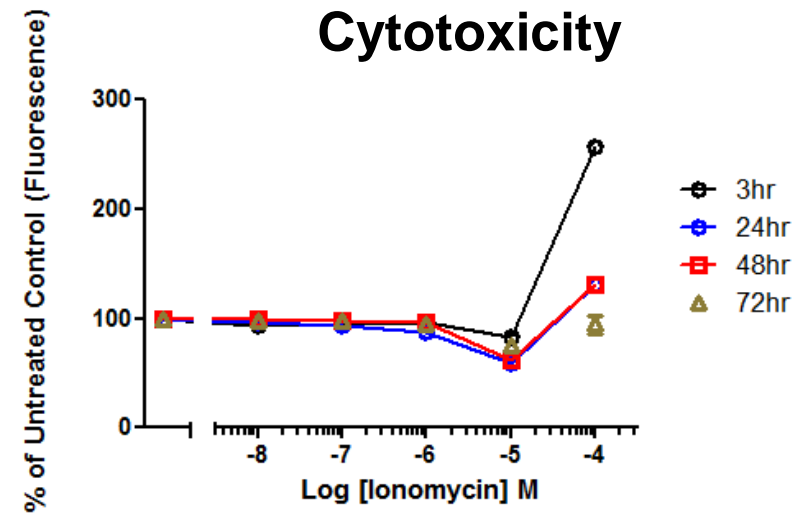
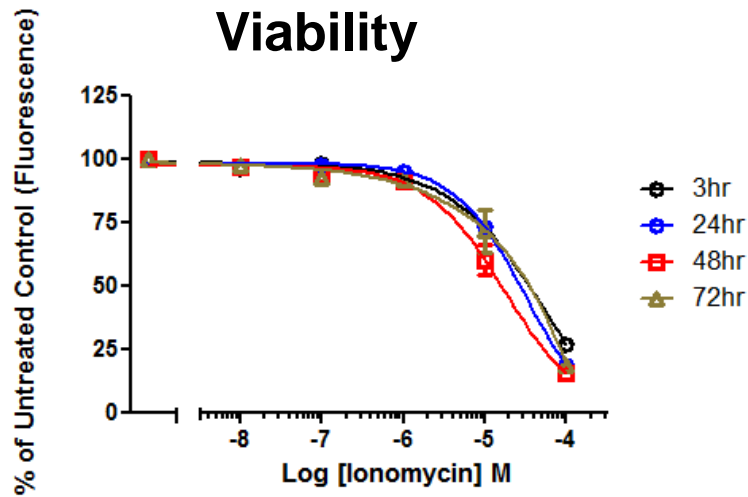
Three Assays Applied in Sequential, Same Well Multiplex = ApoTox-Glo™



Multi-parametric analysis facilitates dissection of the cytotoxic process



Ionomycin (Fast-acting, 1° Necrosis Inducer)

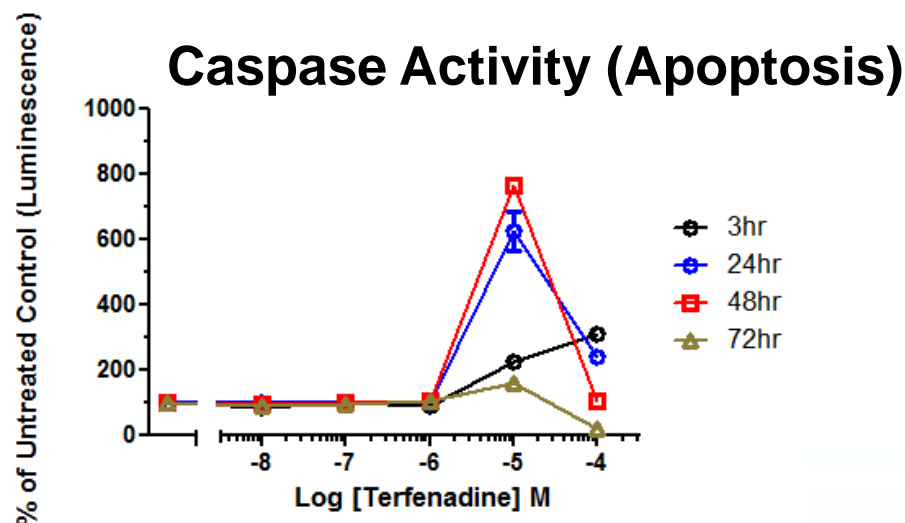
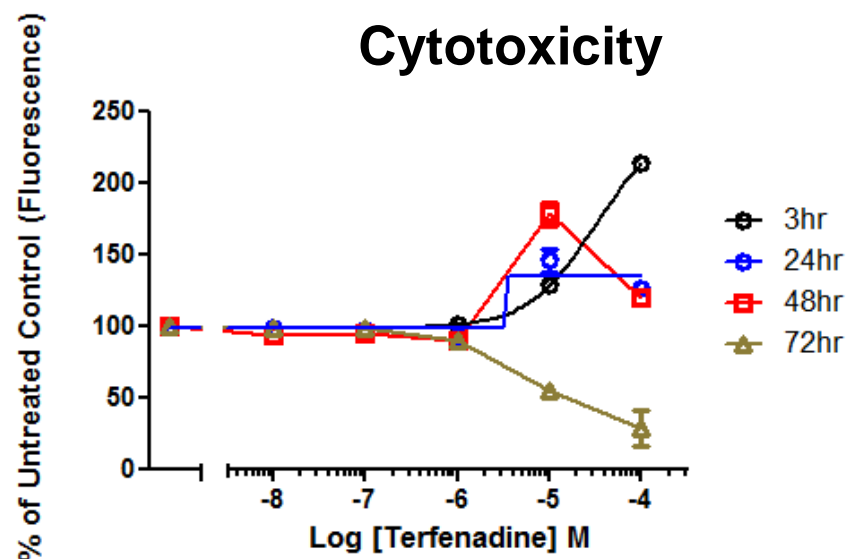
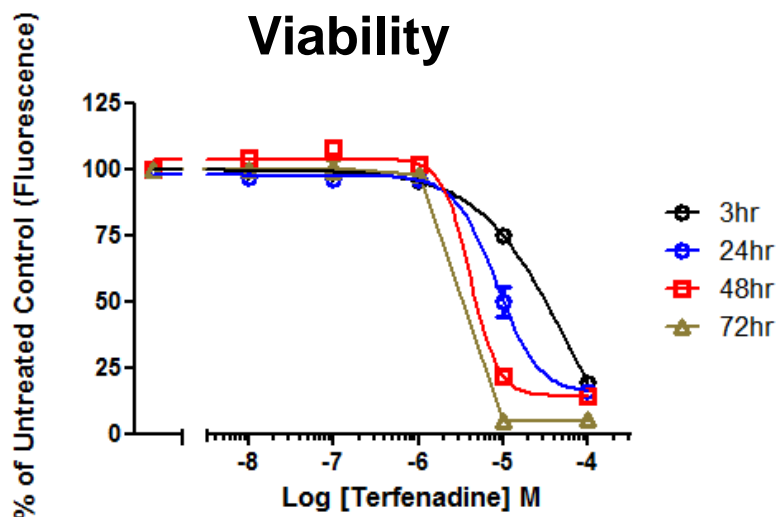


Viability assays always tell you the relative number of cells left after treatment.

Activity-based cytotoxicity markers are definitive for cell death, but subject to degradation as a function of time.

All cell populations have a basal caspase activity. In the absence of apoptosis, this activity declines with necrotic cell death.

Terfenadine (Fast-acting, Apoptosis Inducer)

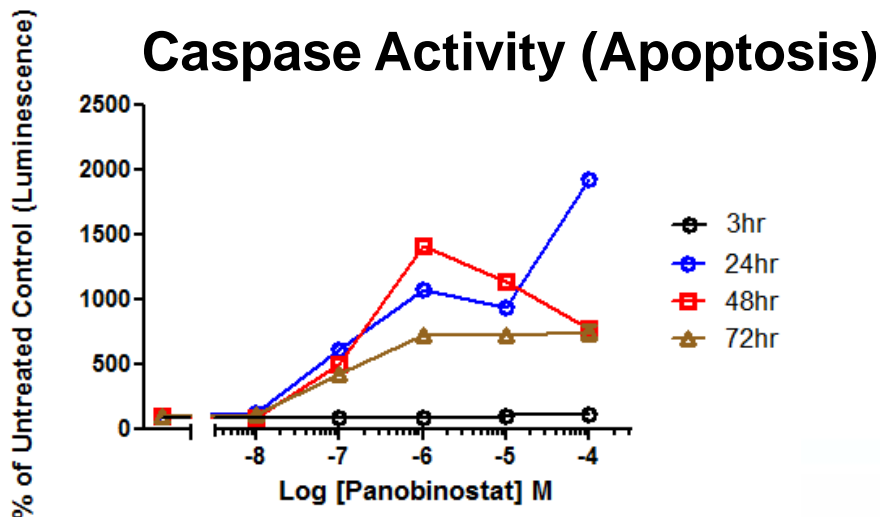
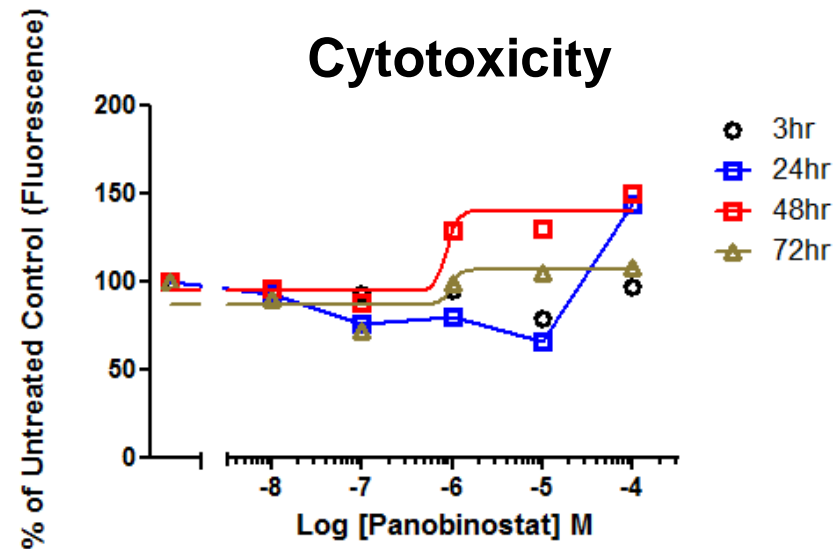
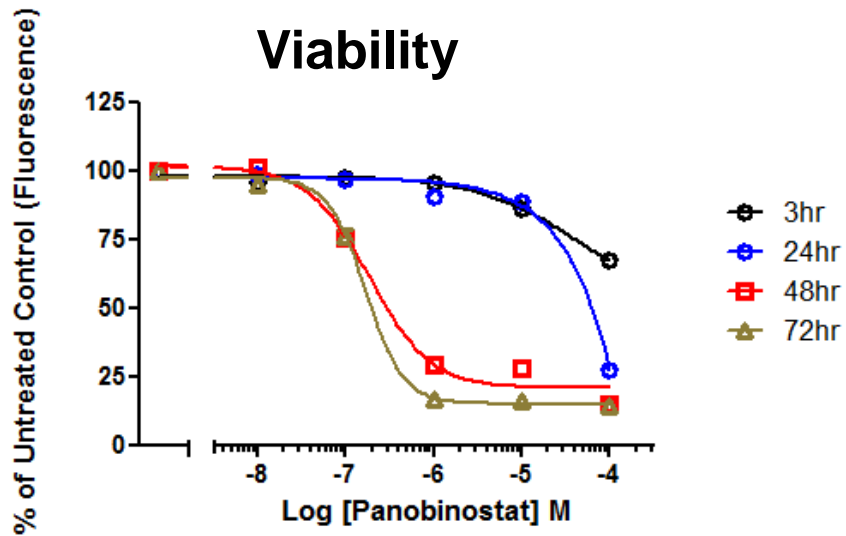


Viability assays always tell you the relative number of cells left after treatment.

Activity-based cytotoxicity markers are definitive for cell death, but subject to degradation as a function of time.

Caspase activity above basal levels is definitive for apoptosis. Caspase activity declines when cells reach secondary necrosis (natural enzyme degradation).

Panobinostat (Slow-acting, Apoptosis)

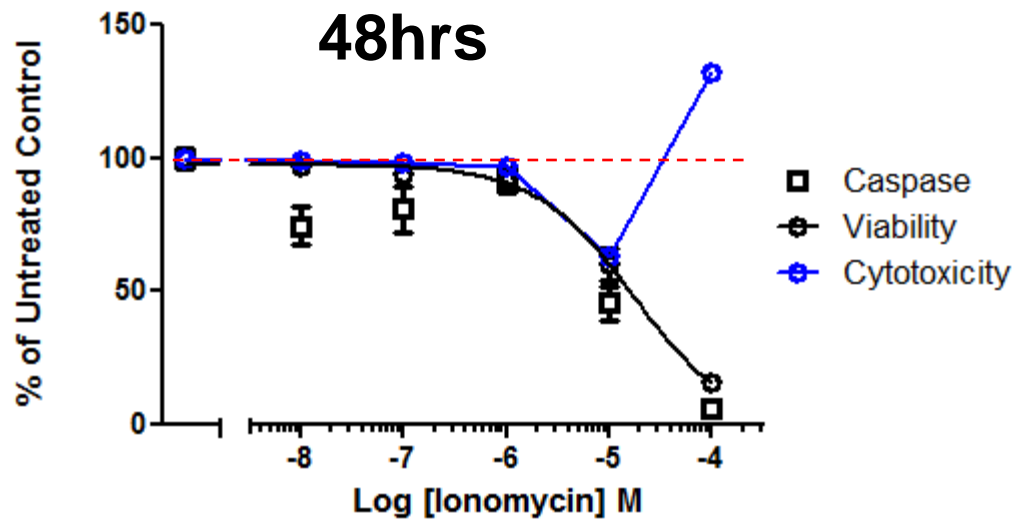


Viability assays always tell you the relative number of cells left after treatment.

Activity-based cytotoxicity markers are definitive for cell death, but subject to degradation as a function of time.

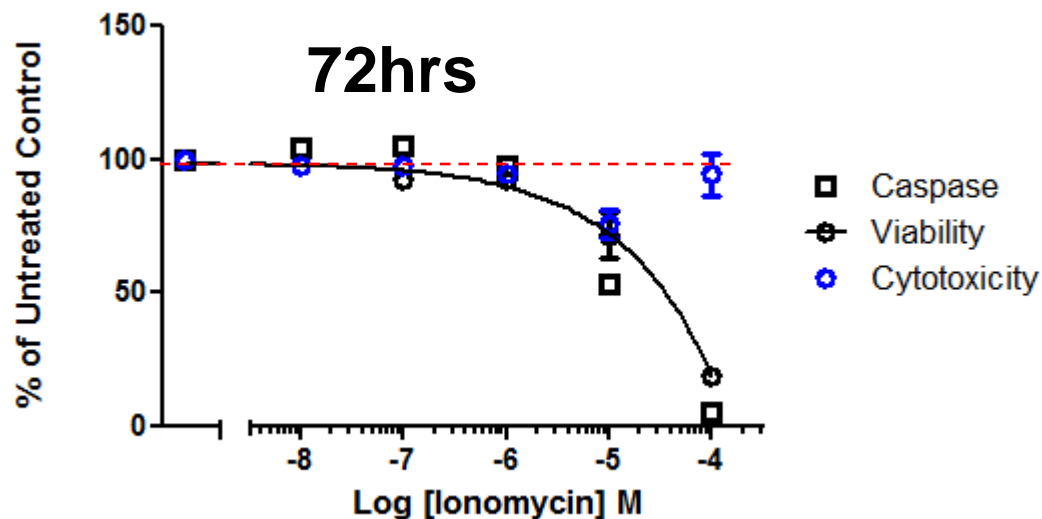
Caspase activity above basal levels is definitive for apoptosis. Caspase activity declines when cells reach secondary necrosis (natural enzyme degradation).

Cytotoxic Time Courses Preferred... Endpoints Can be Meaningful!



Caspase ↓
 Viability ↓
 Cytotoxicity ↑

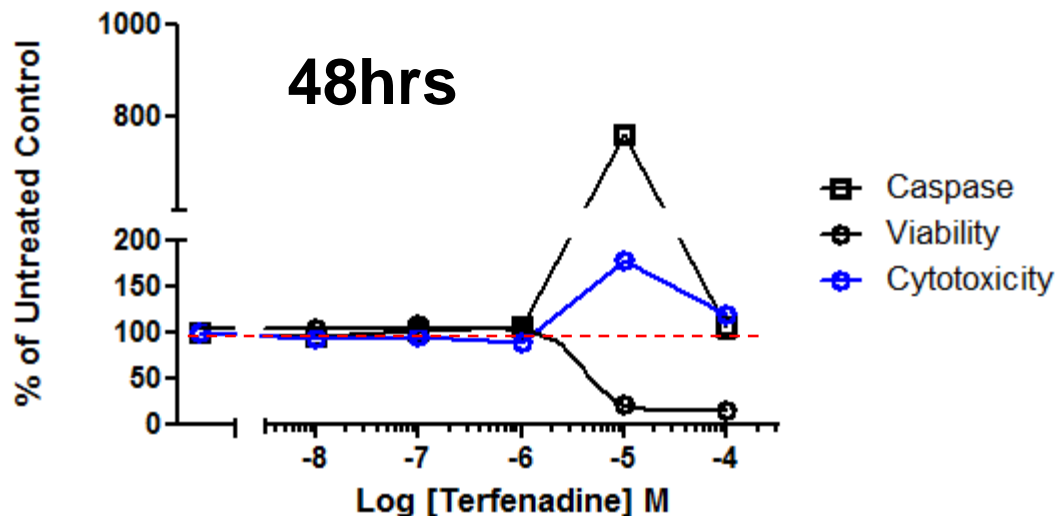
-Consistent with 1^o necrosis



Caspase ↓
 Viability ↓
 Cytotoxicity →

-Consistent with 1^o necrosis
 -Inconsistent with cell cycle arrest due to exposure period

Cytotoxic Time Courses Preferred... Endpoints Can be Meaningful!

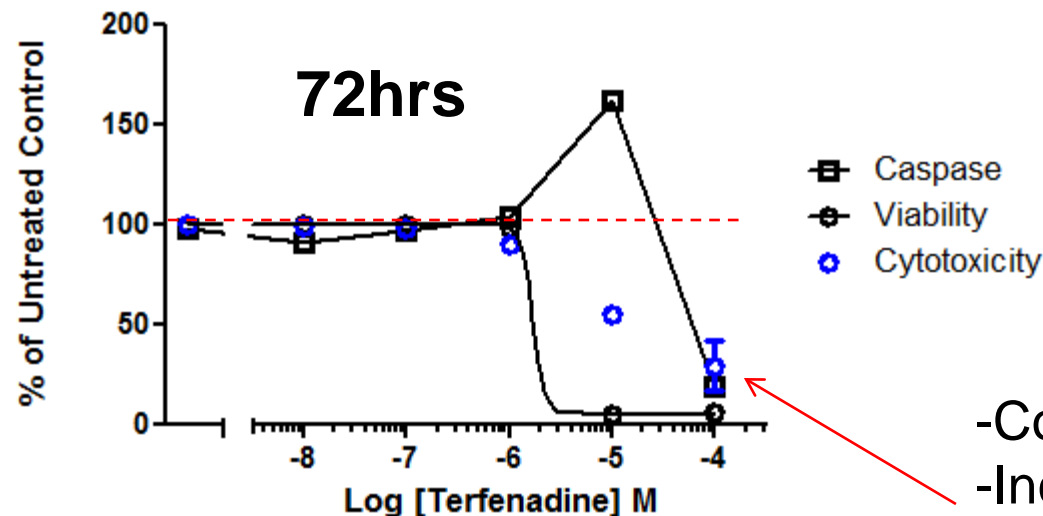


Caspase

Viability

Cytotoxicity

- Consistent with late apoptosis



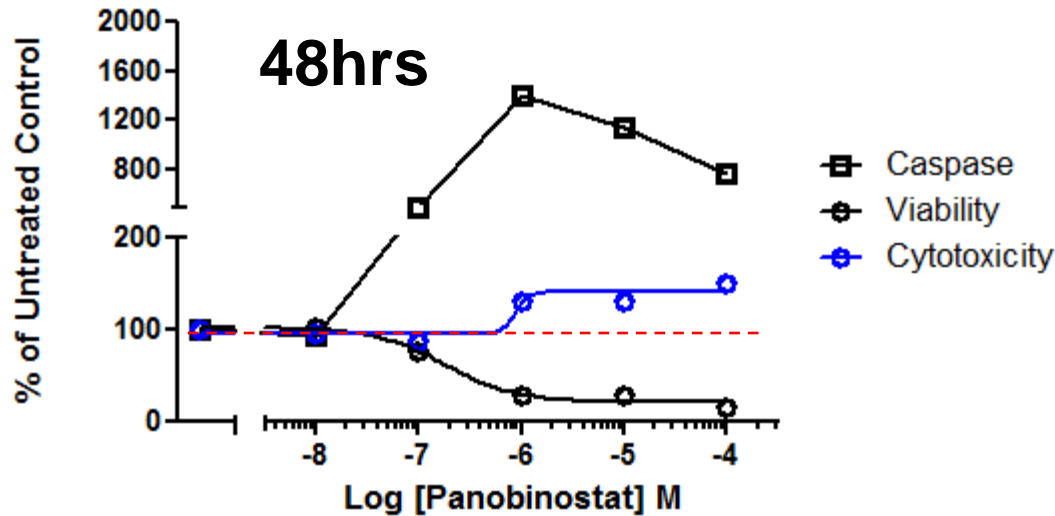
Caspase

Viability

Cytotoxicity

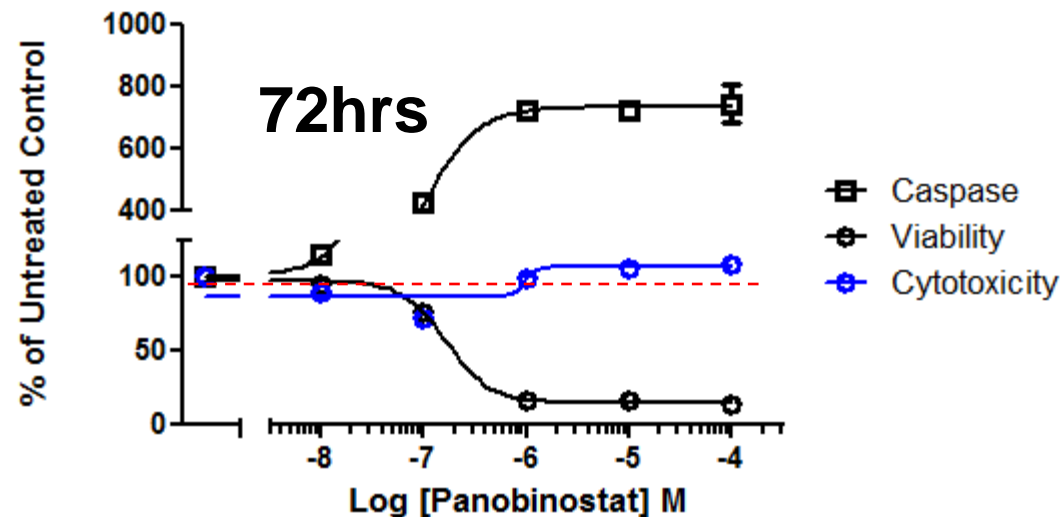
-Consistent with late(r) apoptosis
-Inconsistent with early cell cycle arrest (due to cytotoxicity and exposure)

Cytotoxic Time Courses Preferred... Endpoints Can be Meaningful!



Caspase ↑
 Viability ↓
 Cytotoxicity ↑

-Consistent with apoptosis



Caspase ↑
 Viability ↓
 Cytotoxicity ↑

-Consistent with apoptosis

ApoTox-Glo™ Multiplexed MOA Assay



Advantages

- Provides a detailed profile of the cytotoxic response
 - Primary Necrosis
 - Cell Cycle Arrest
 - Apoptosis
 - Secondary Necrosis
- Convenient, same-well multiplex
- Measures are inversely correlated, reducing optical interferences and providing “flagging” opportunities

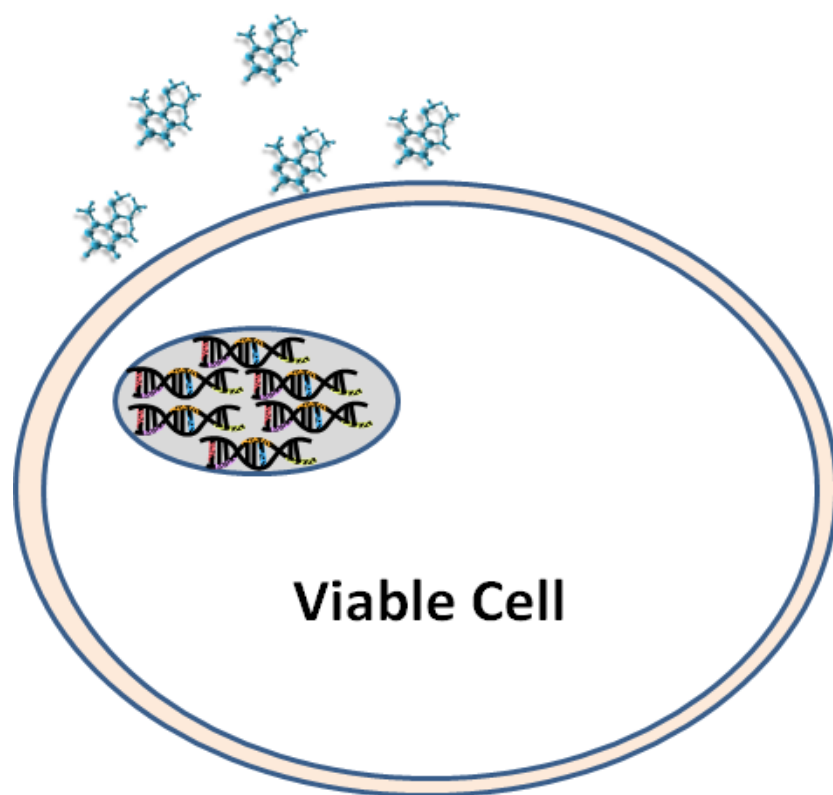
Disadvantages

- Best employed in time course experiments requiring multiple times and data points per test compound
- Can be applied at specified endpoints...but biomarker degradation may complicate data interpretation

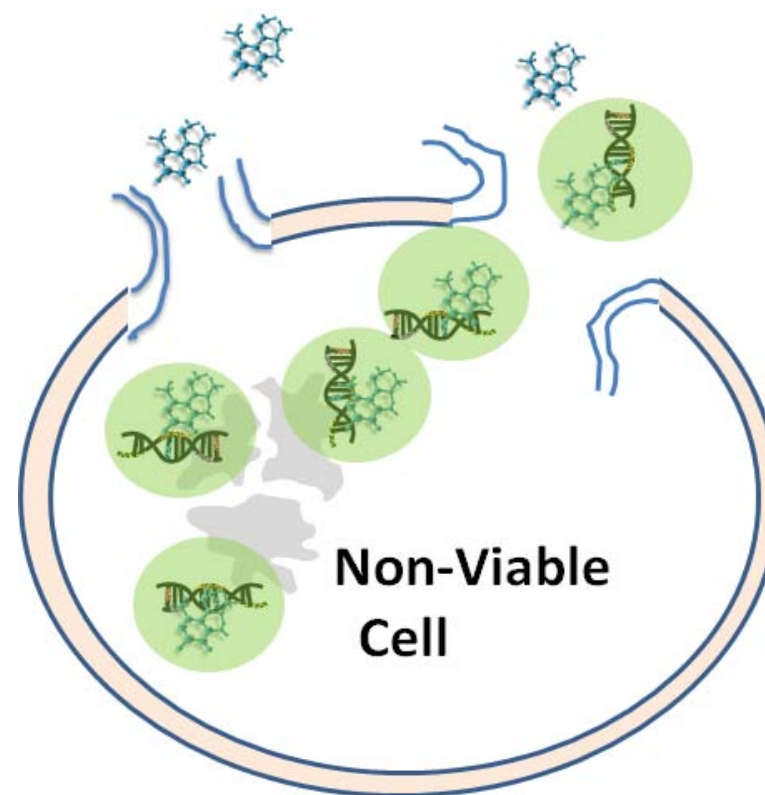
Other complimentary and orthogonal assay options for cytotoxicity determination?

Non-Activity Based Cytotoxicity Measures? Differential DNA Dye Excludability?

Membrane integrity “sensed” by environmental dye.



Excluded dye yields no increase in fluorescence.

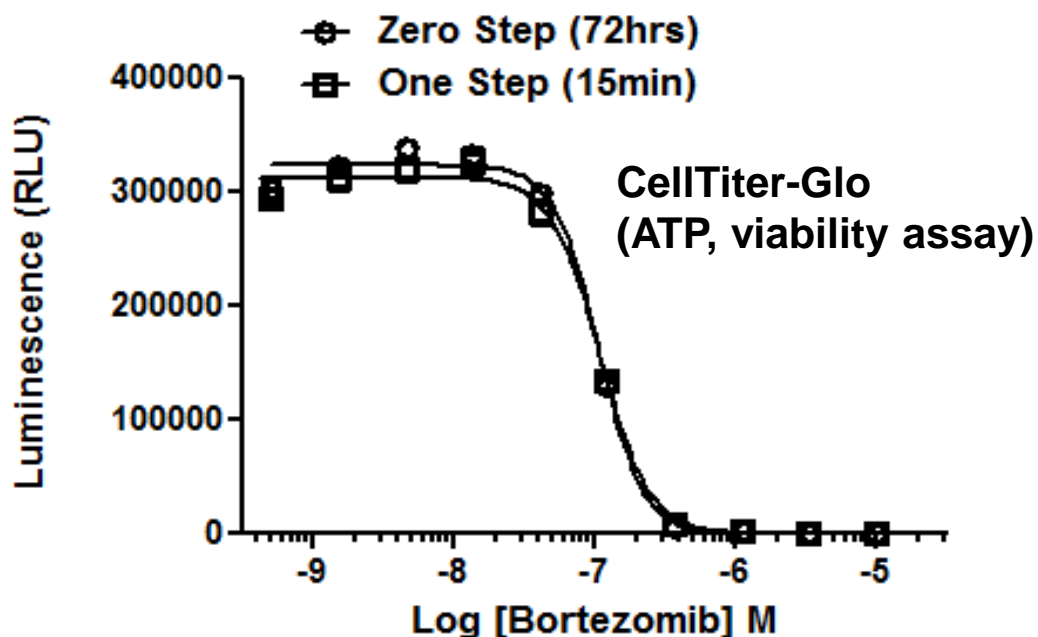


Non-excluded dye yields increase in fluorescence

Probe “Inertness” and “No Step* Format”



*No Step means adding the dye at cell dosing...
either in drug dilutions or with cells.



CellTox-Green™ does not alter viability or
impact dose-dependence of cytotoxic model
compounds in extended co-incubations.
HeLa at 72hrs shown.

Kinetic Assay Format

Add DNA probe at cell seeding



Add diluted test article



Measure fluorescence at 4hrs



Measure fluorescence at 24hrs



Measure fluorescence at 48hr

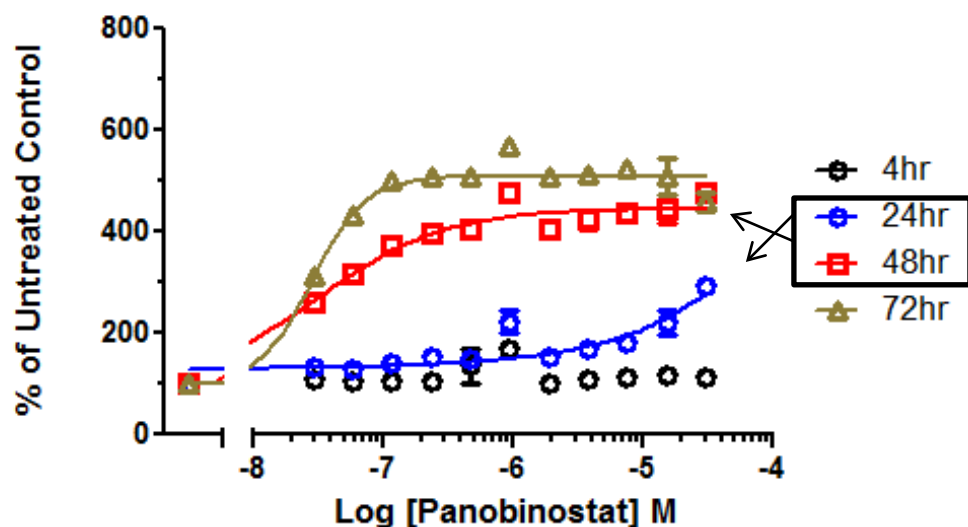
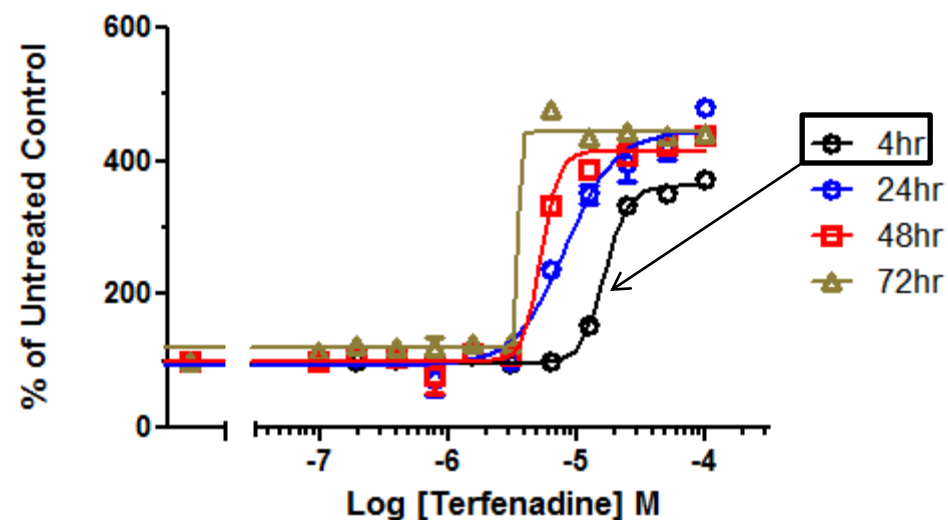
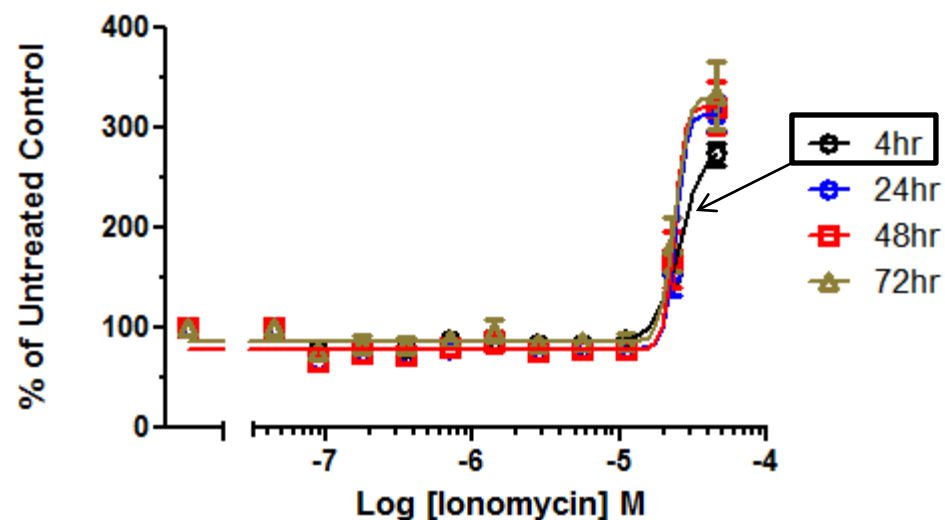


Measure fluorescence at 72hr



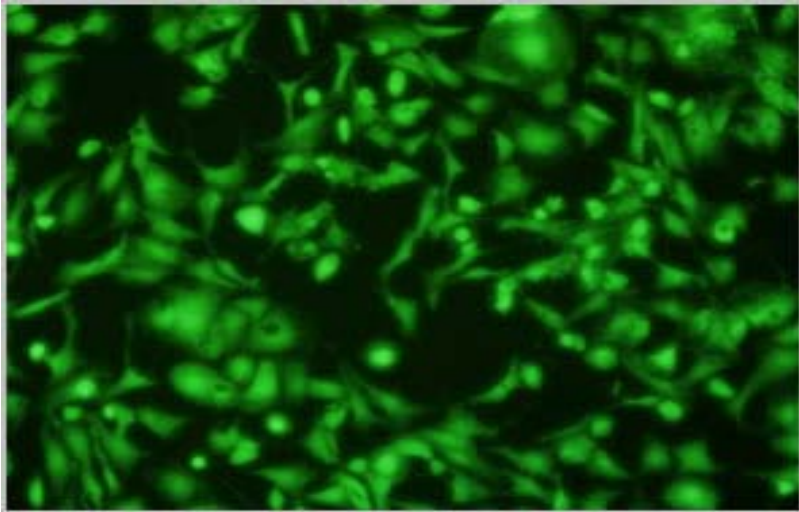
Optional sequential multiplex
with viability assay (CellTiter-Glo,
CellTiter-Blue or CellTiter-Fluor)

CellTox-Green™: A Kinetic Cytotox Assay



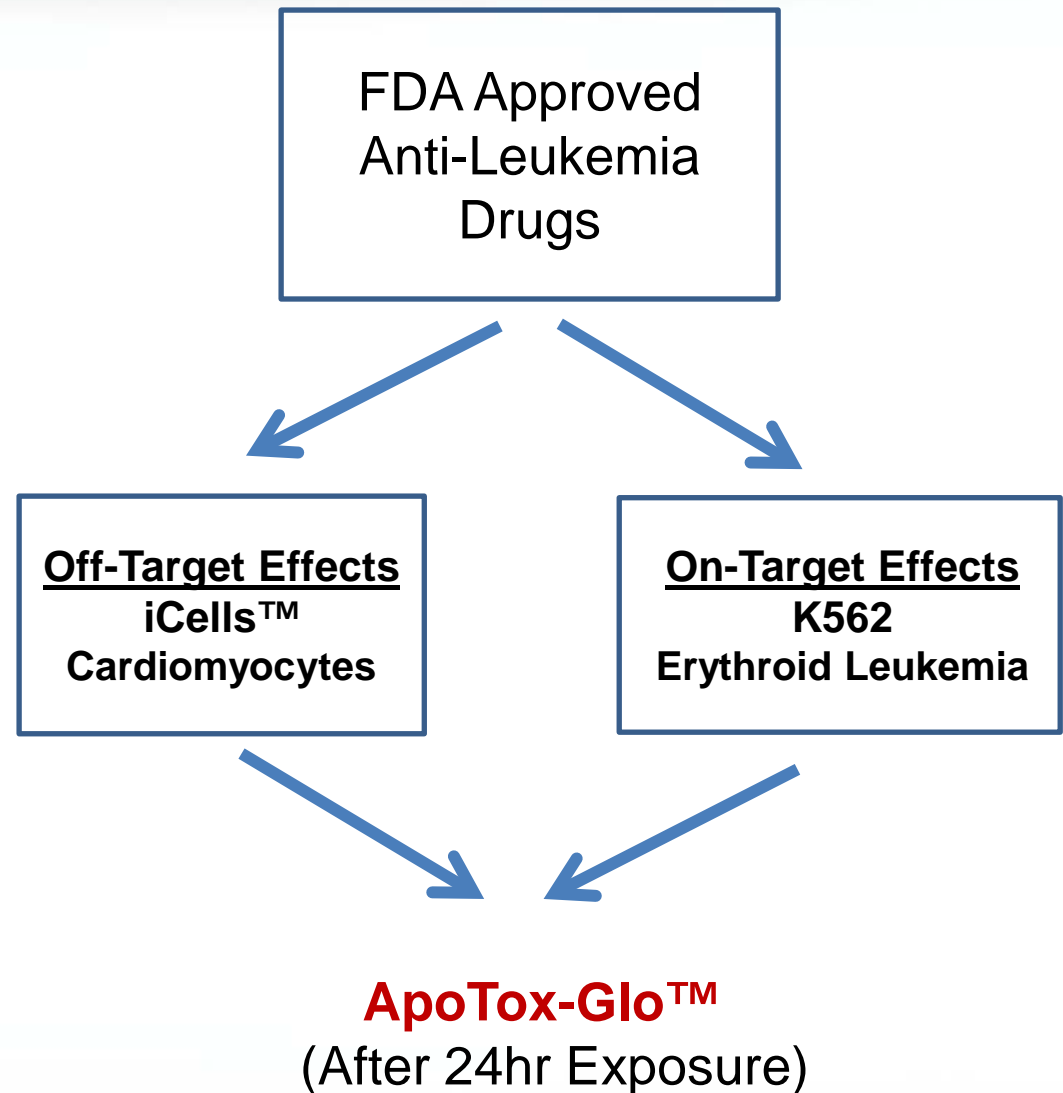
- Provides flexible, non-duplicative time course data for cytotoxicity
- Can be sequentially multiplexed at first emergence of cytotoxicity
- Data can be used to explore the complete cytotoxic response with a subsequent ApoTox-Glo assay

Potency and Safety Evaluation: Validation of ApoTox-Glo™ 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”.*

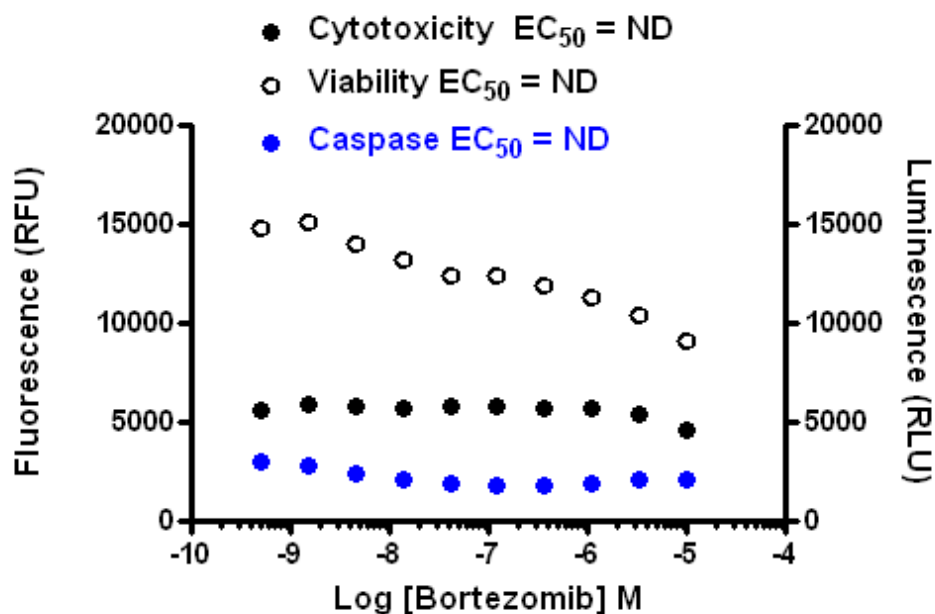
-Cellular Dynamics International



Proteasome Inhibitor (Velcade™)



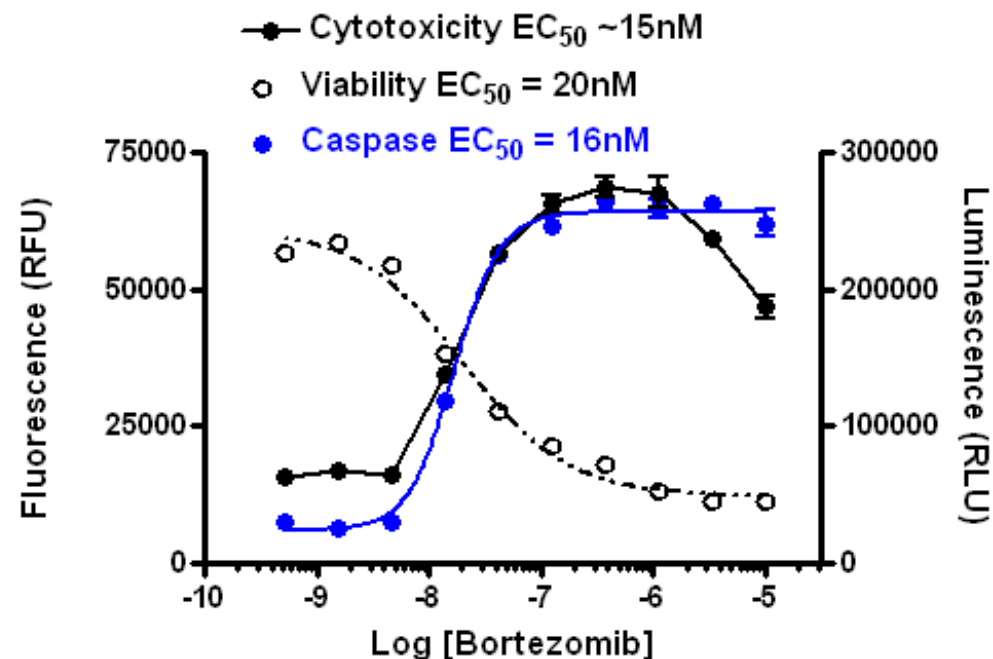
iCell™



**No apparent cytotoxicity*
or caspase activation**

*Bortezomib (and other proteasome inhibitors) are known to partially inhibit the viability assay protease biomarker at concentrations greater than 1µM

K562

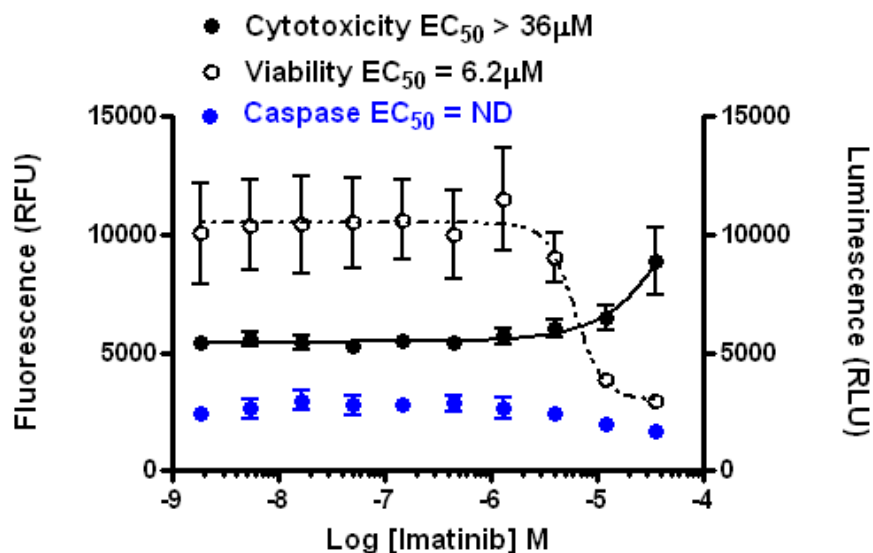


Cytotoxicity by apoptosis

Tyr Kinase Inhibitor (Gleevec™)

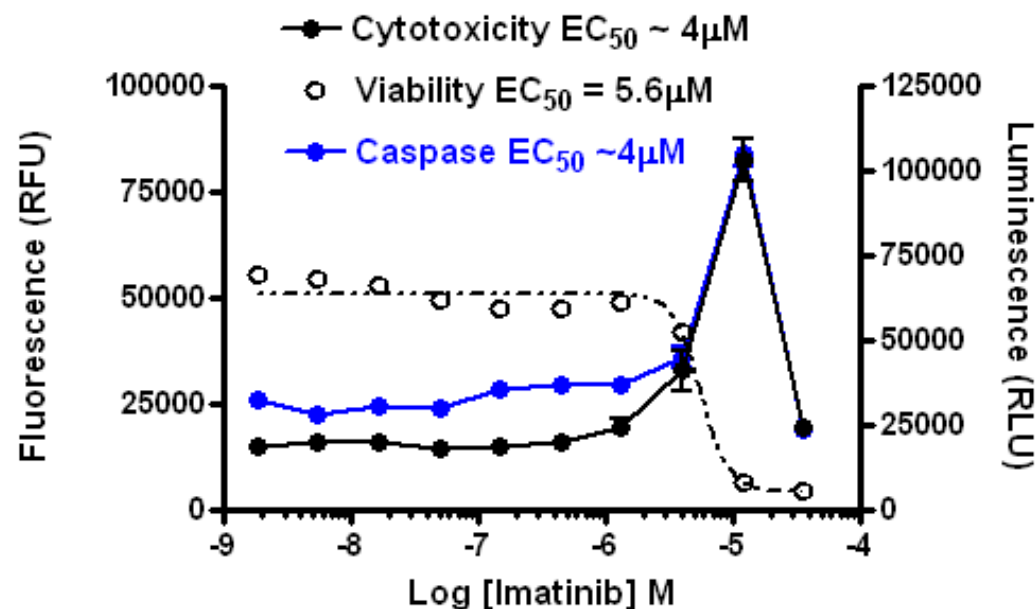


iCell™



**Caspase-independent
cytotoxicity at concentrations
greater than $1\mu\text{M}$**

K562



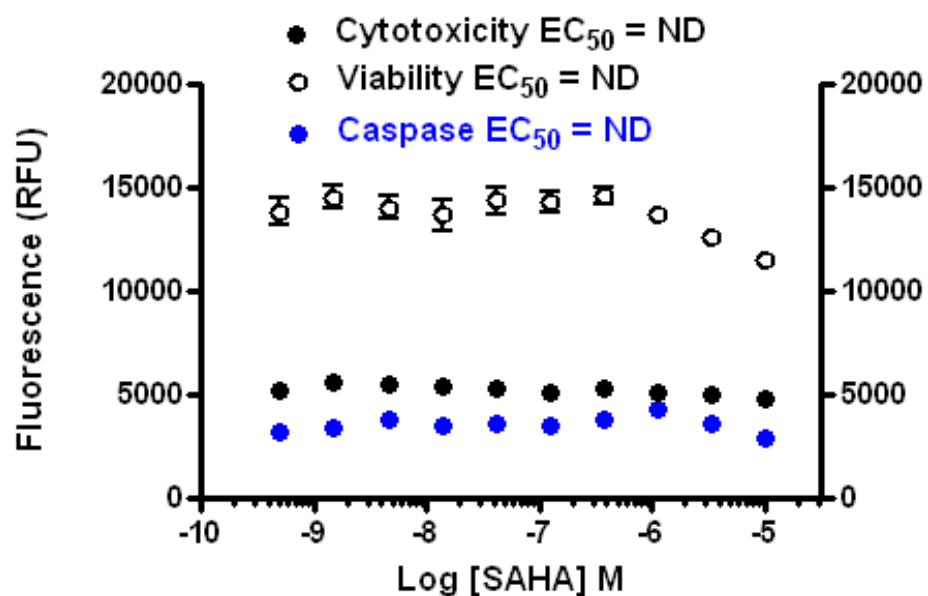
Cytotoxicity by apoptosis

Note: Diminution of caspase and cytotoxicity biomarker signals at highest doses of imatinib (with K562) are consistent with activation kinetics and time-dependent biomarker decay.

Histone Deacetylase Inhibitor (Vorinostat™)

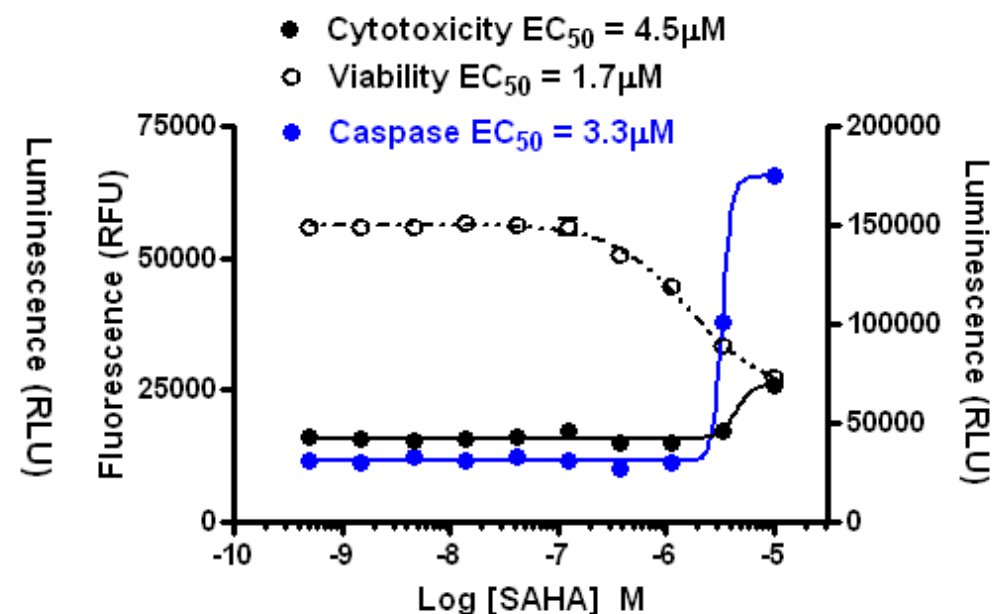


iCell™



No apparent cytotoxicity
or caspase activation.

K562

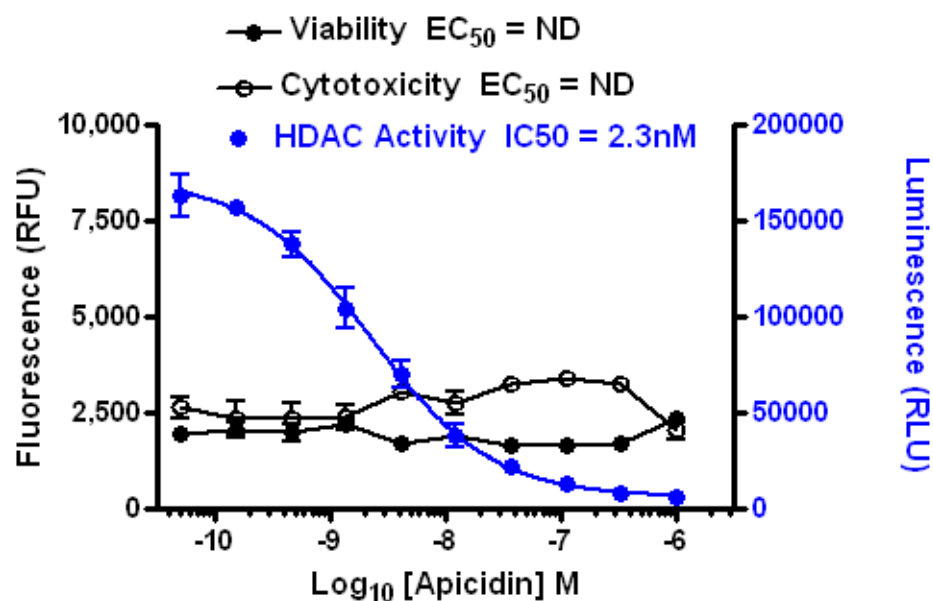


Cytotoxicity by apoptosis

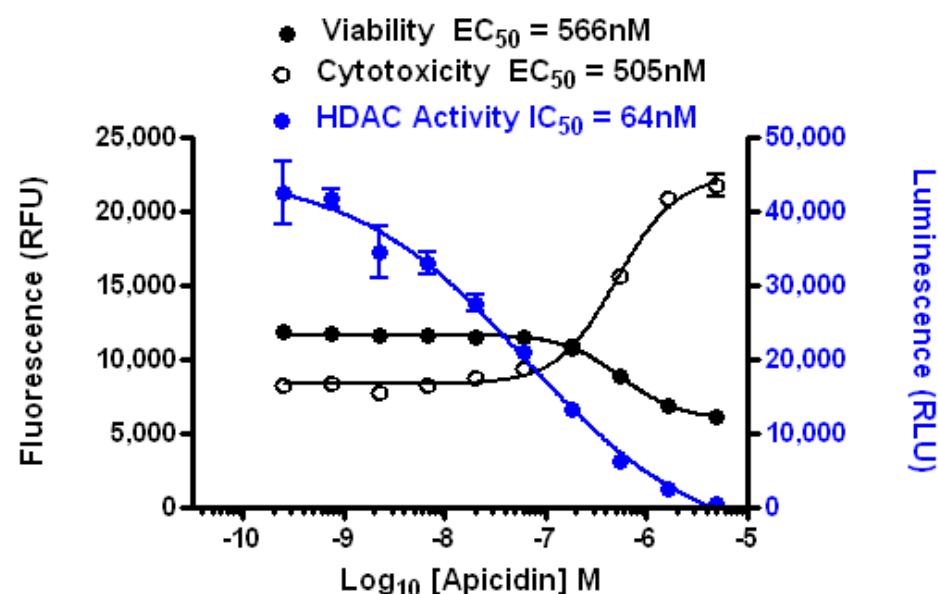
Targeted Activities...Differential Cytotoxicity



iCell™



U937 (Cancer Cells)



[24hr compound exposure]

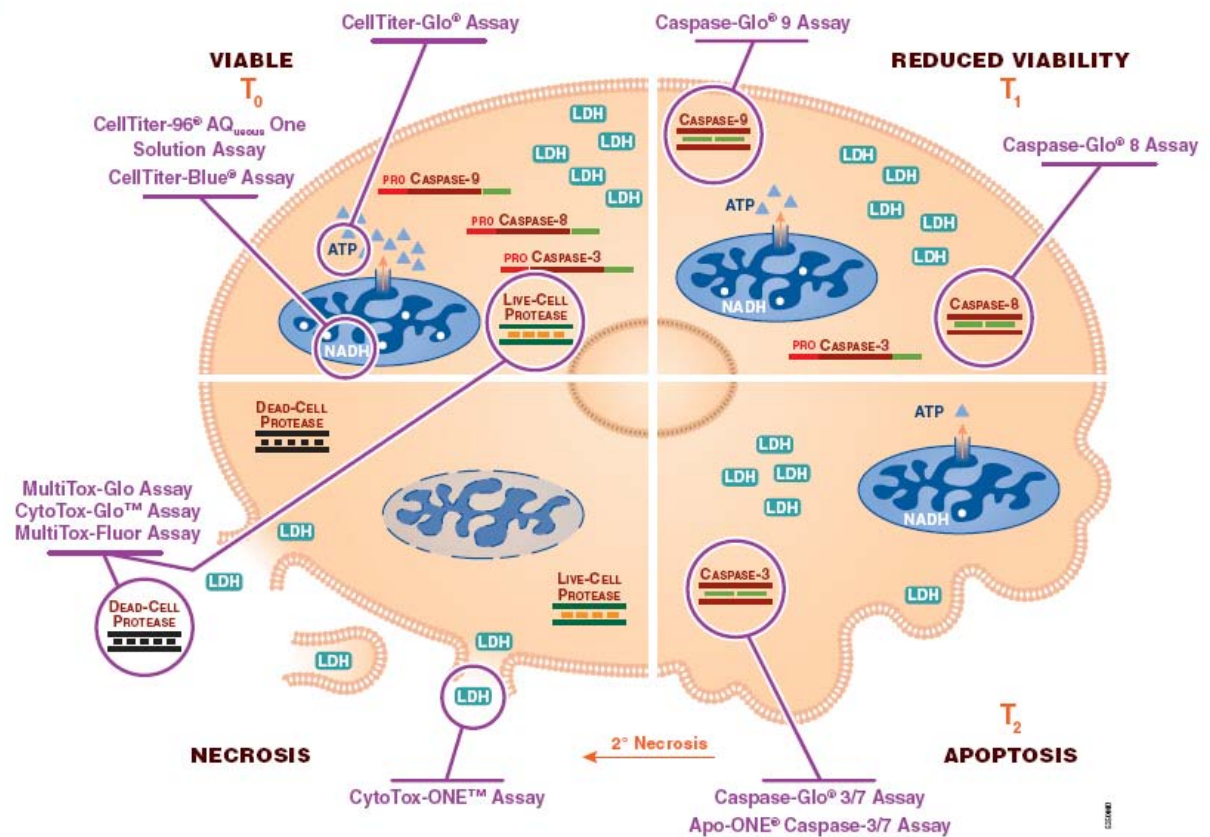
MultiTox-Fluor followed by HDAC-Glo™ I/II

Summary



In vitro cell health data is influenced by:

- Dosage
 - Addressed through serial dilution series
- Exposure Time (cells with compound contact)
- Mechanism of action of the test compound
- Cell Type
 - specific target
 - off target



...and ApoLive-Glo[™], ApoTox-Glo[™],
HDAC-Glo[™] I/II Assays!

For more information contact:
andrew.niles@promega.com

Acknowledgements



Promega Scientists and Support Staff Worldwide!

