



A Real-Time Cytotoxicity Assay That Delivers More Relevant Data

Andrew L. Niles

Senior Research Scientist

Biological Outcomes in Cell Culture

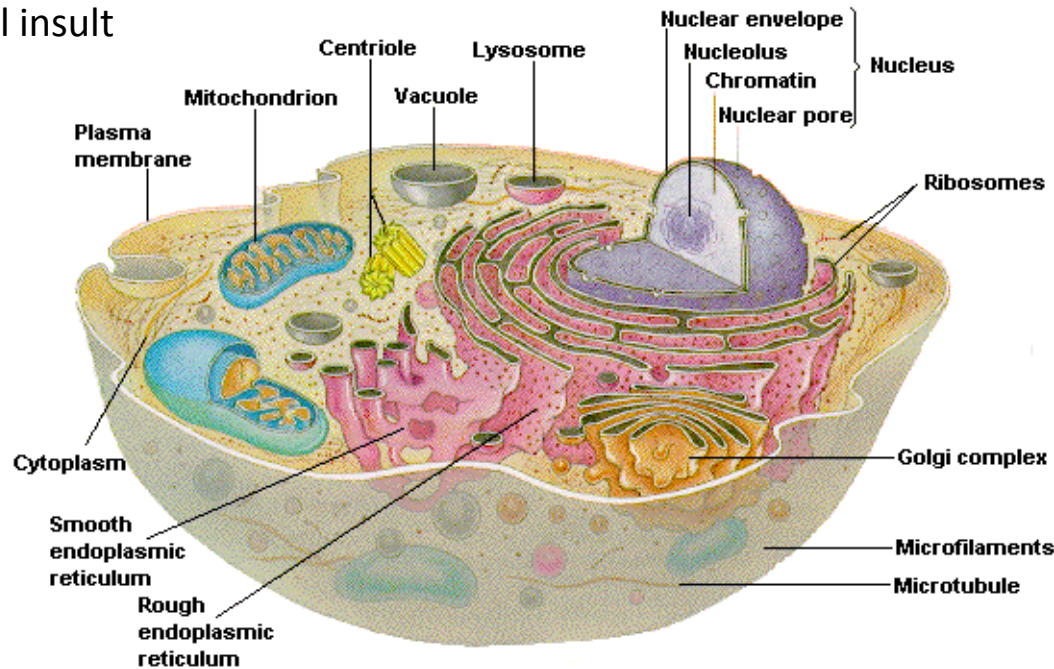
Treatment

- Small molecule
- Biological
- Industrial chemical
- Transgene
- Physical insult

Cause



Effect



Normal Proliferation

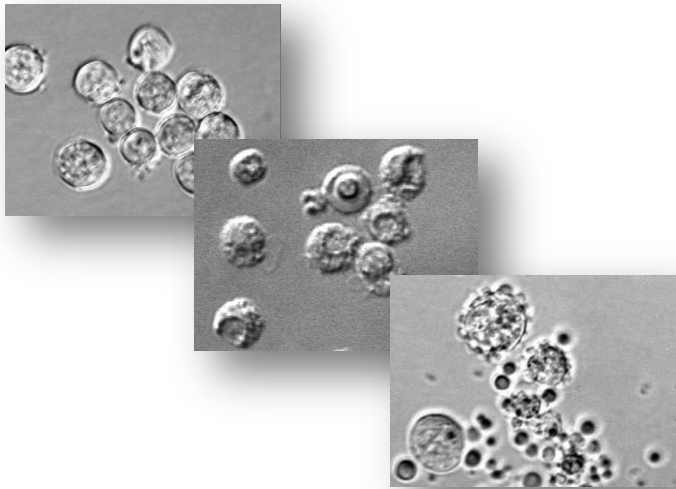
Enhanced Proliferation

- Cell-cycle Arrest
- Senescence/Quiescence
- Autophagy
- Primary Necrosis
- Apoptosis
- Necroptosis
- Secondary Necrosis

**“Cytotoxicity
Phenotype”**

Is My Compound or Treatment Toxic?

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

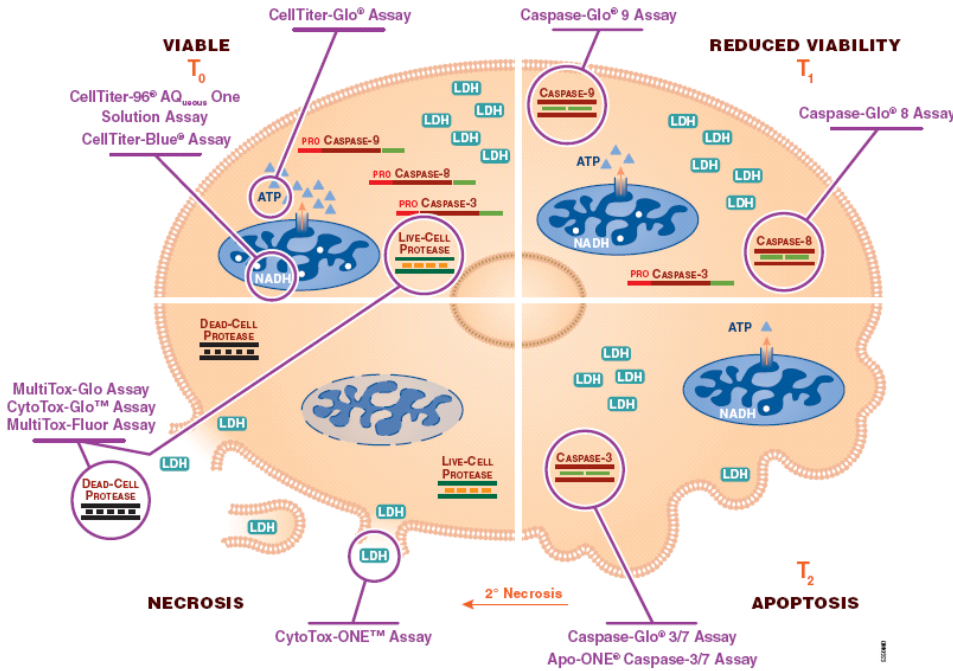
2. Exposure time (cells with compound contact)

3. Mechanism of action of the test compound

4. Cell type

- Specific Target
- Off Target

Biomarkers of a cytotoxic response



As cells lose vitality and health,

Viability biomarkers

- Metabolism biomarkers (ATP, reductase)
- Constitutive enzymatic biomarkers



Cytotoxicity biomarkers

- Constitutive enzymatic biomarkers
- Induceable biomarkers



Viable

“Stress” biomarkers

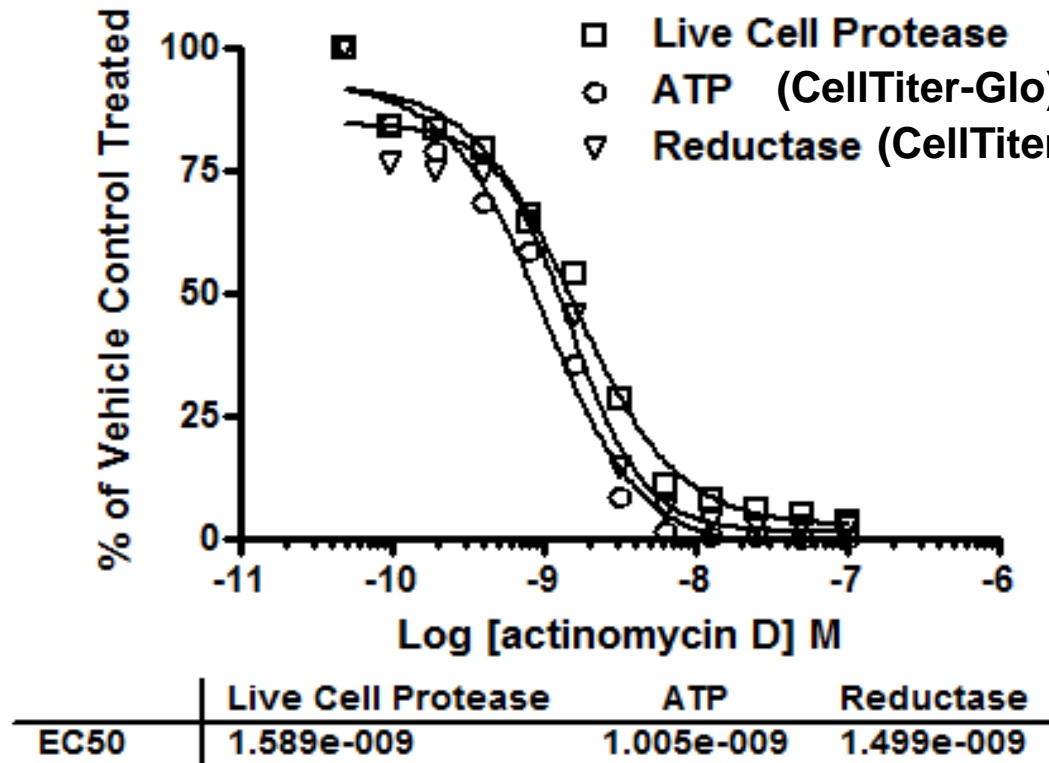
anti-proliferative effects

MOA death

Non-Viable

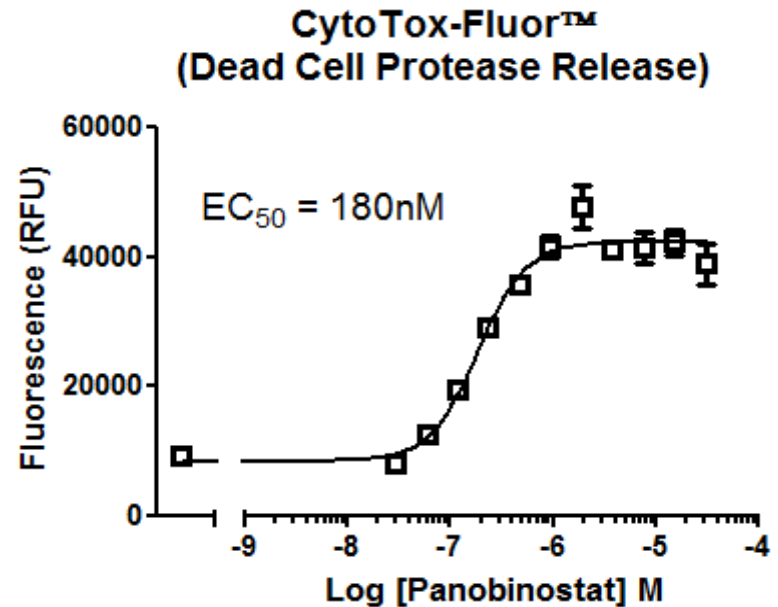
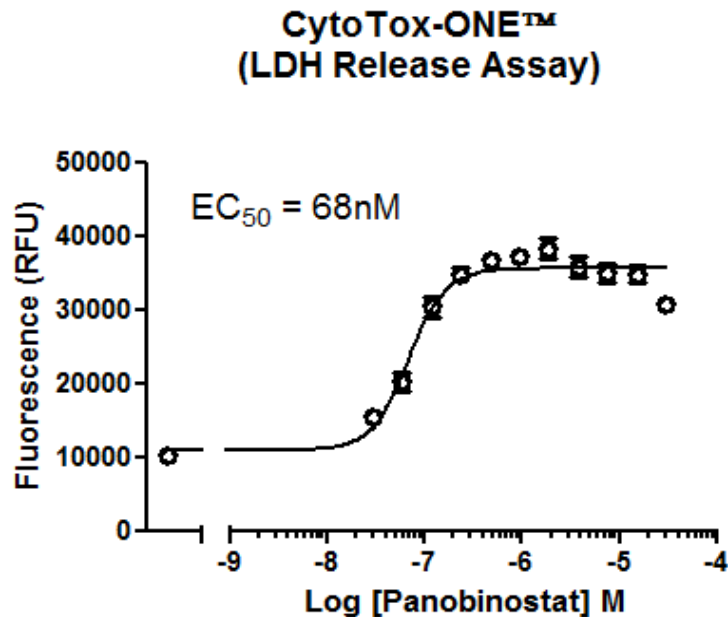
Elapsed Time

Viability Assays (in general) show good concordance with loss of membrane integrity



Reliably report the relative number of viable cells... but don't differentiate between cytotoxicity and cytostasis

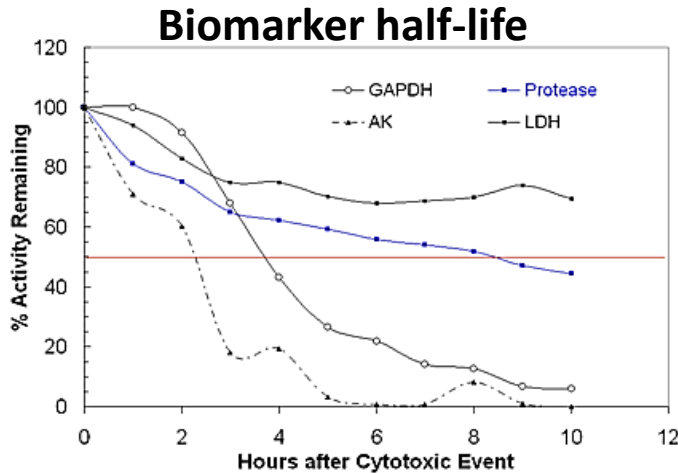
Conventional Cytotoxicity Assays



iCell™ Hepatocytes dosed for 72hr with panobinostat, endpoint chemistries added.

Definitive for cell death. Utility dependent upon kinetics of cytotoxic event...

Current Cytotoxicity Assay Limitations

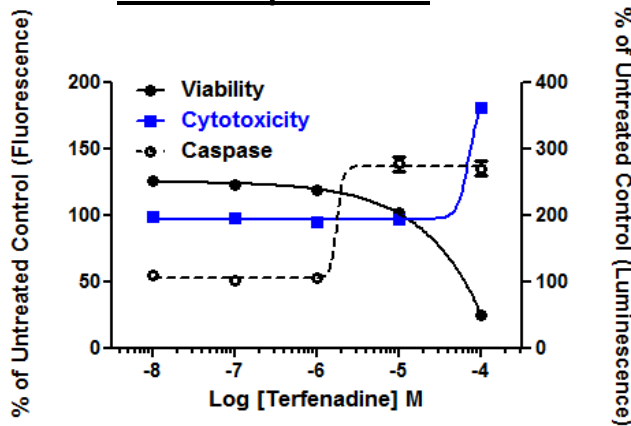


Activity-based cytotoxicity measures:

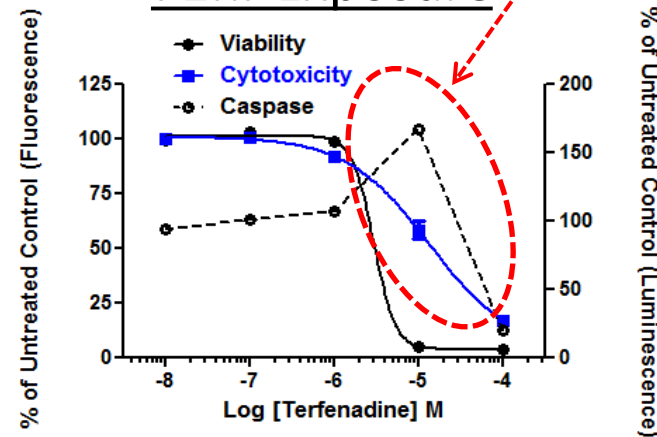
1. are susceptible to biomarker degradation
2. can underestimate cytotoxicity in long-term exposure models.

Caspases have a half-life, too.

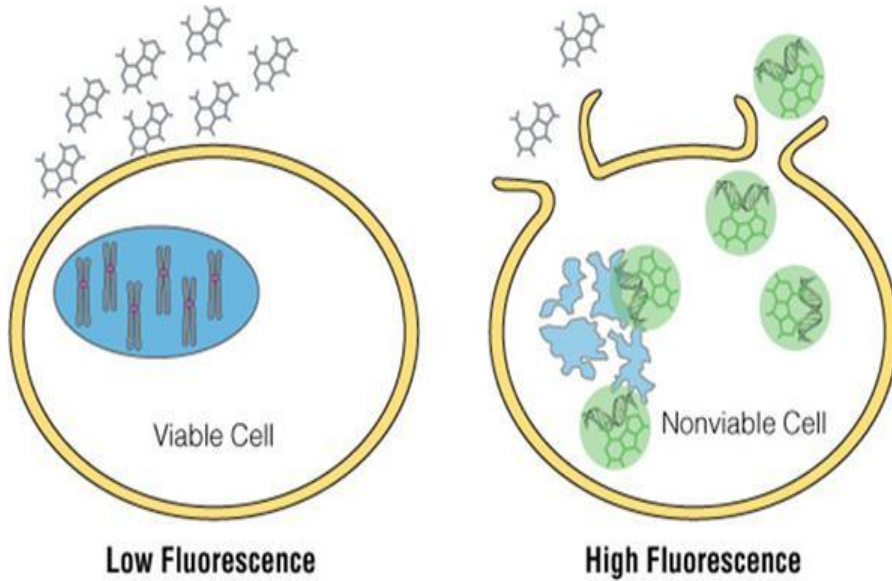
6hr Exposure



72hr Exposure



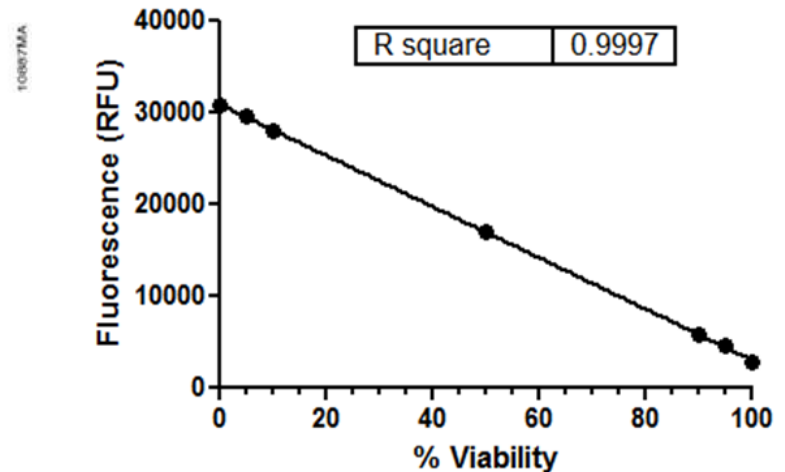
What About Dye Excludability?



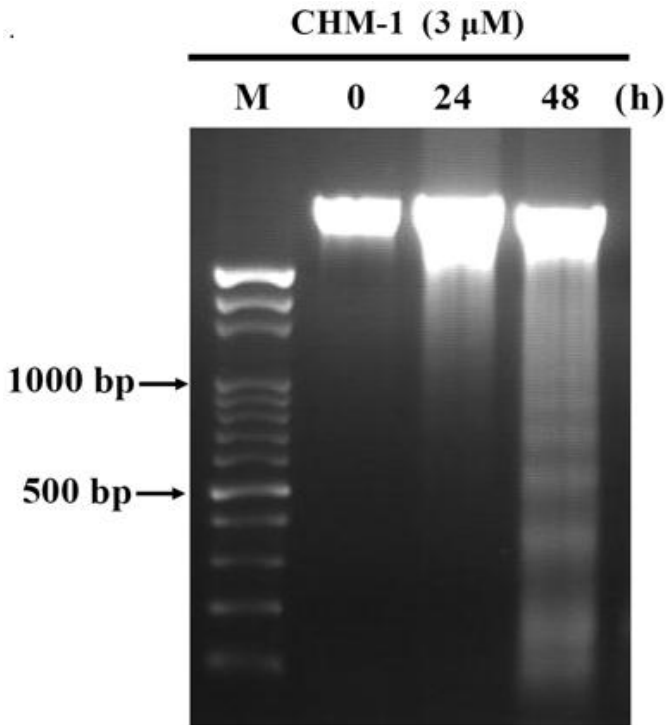
Signal intensity
is proportional to
cell death

CellTox™ Green Dye is a cell impermeant asymmetric cyanine dye with affinity for genomic DNA.

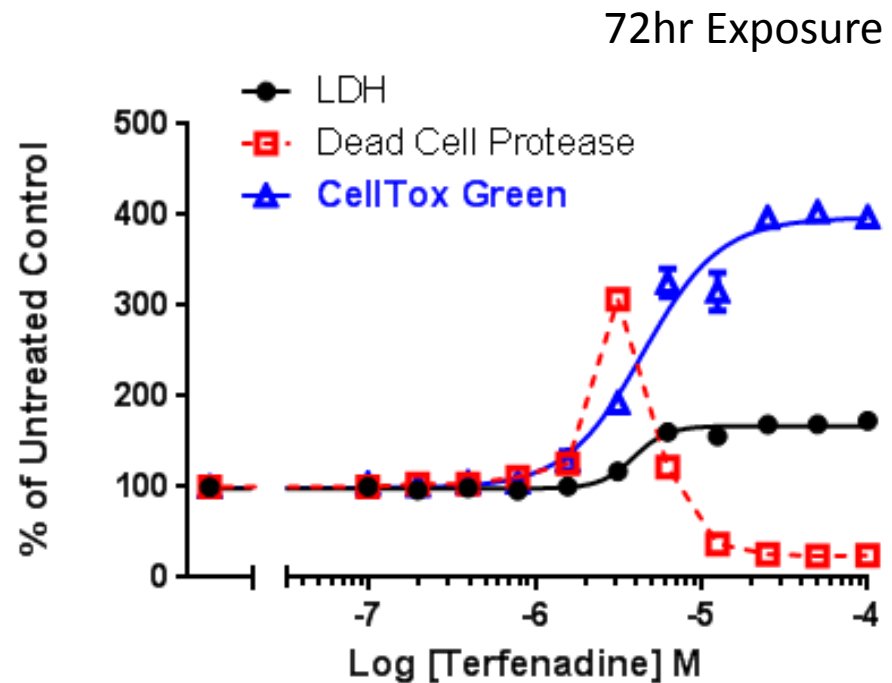
	4174	$\text{Ex} = 512 \text{ nm}$
		$\text{Em} = 532 \text{ nm}$
		$\epsilon = 88550$
		$\Phi_{\text{dsDNA}} = 0.28$
		$\Phi_{\text{oligo}} = 0.54$



DNA as a Biomarker of Cytotoxicity

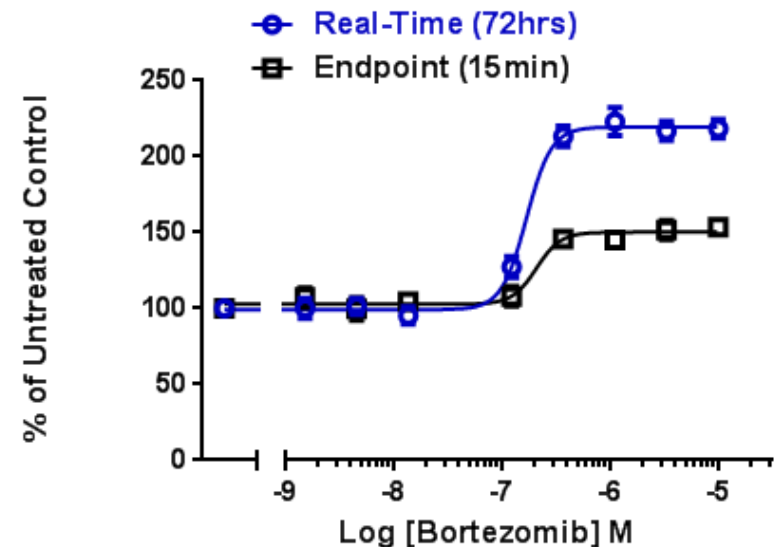
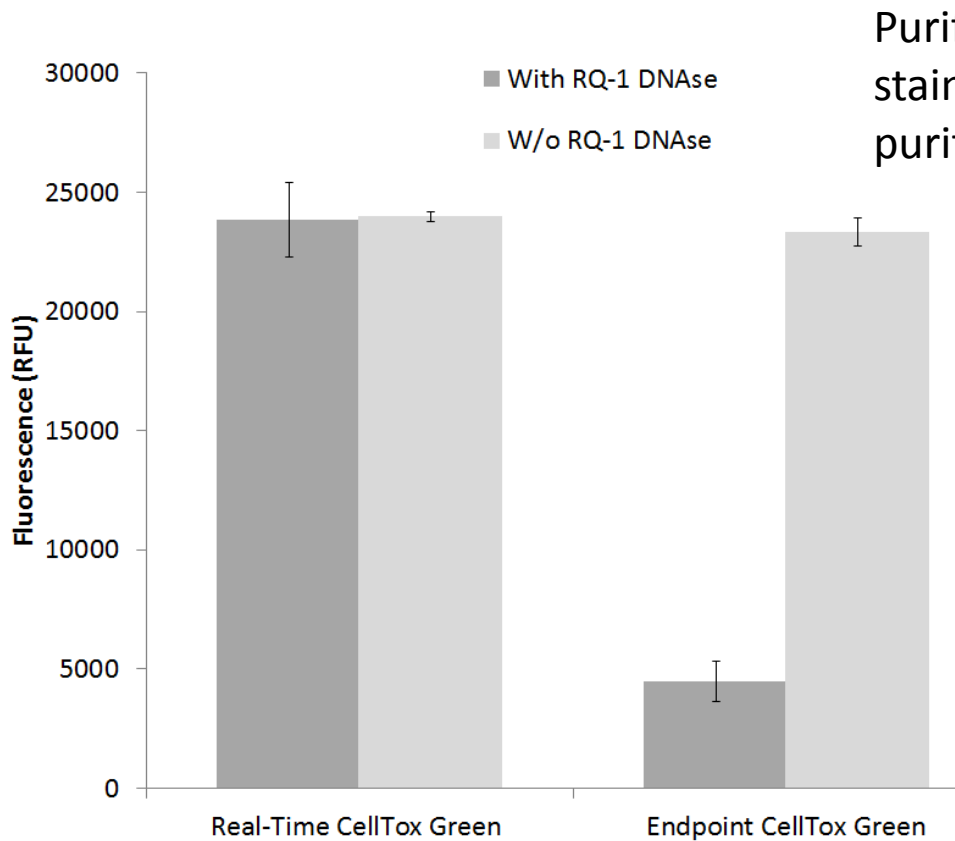


Chen et al (2010) Anti-Cancer Research
30:4187



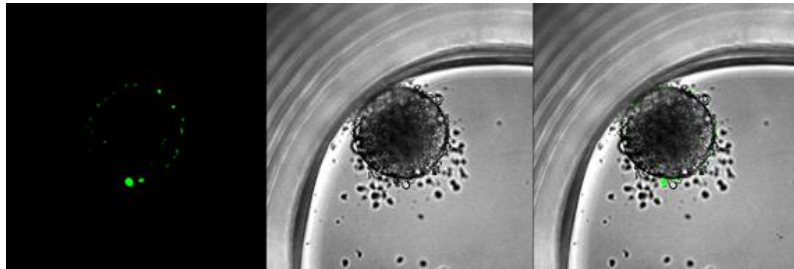
DNA degrades too, but retains sufficient base-pair pairing to facilitate staining

CellTox™ Green Staining Actually Protects DNA from Nuclease Activity



Non-Toxic, Allowing Real-Time Format

Untreated HCT-116 spheroid at 72hrs

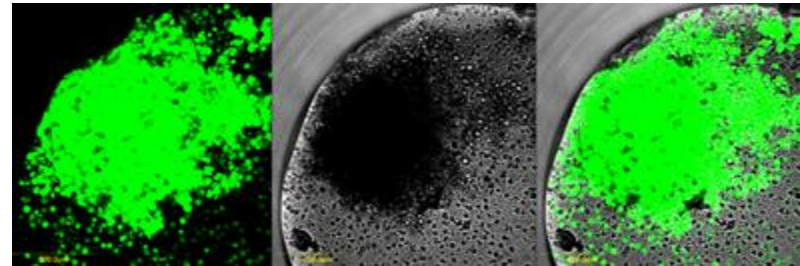


Fluorescence

Bright Field

Merge

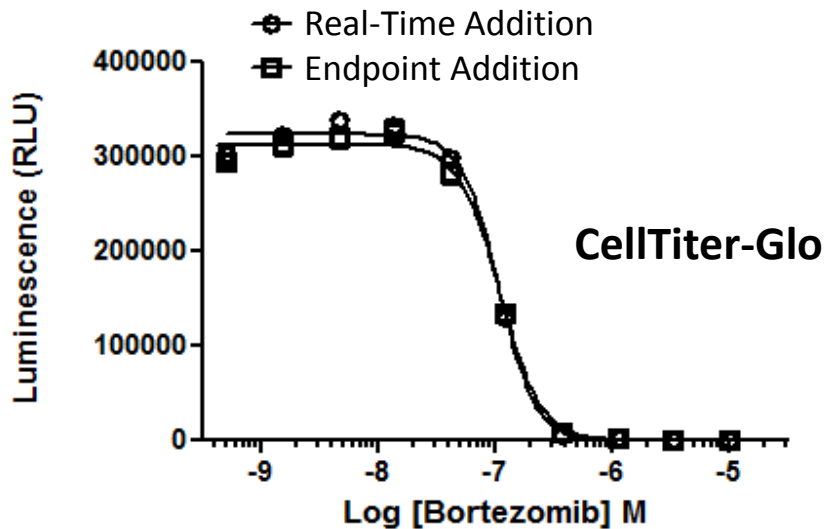
Panobinostat-treated HCT-116 at 72hr



Fluorescence

Bright Field

Merge



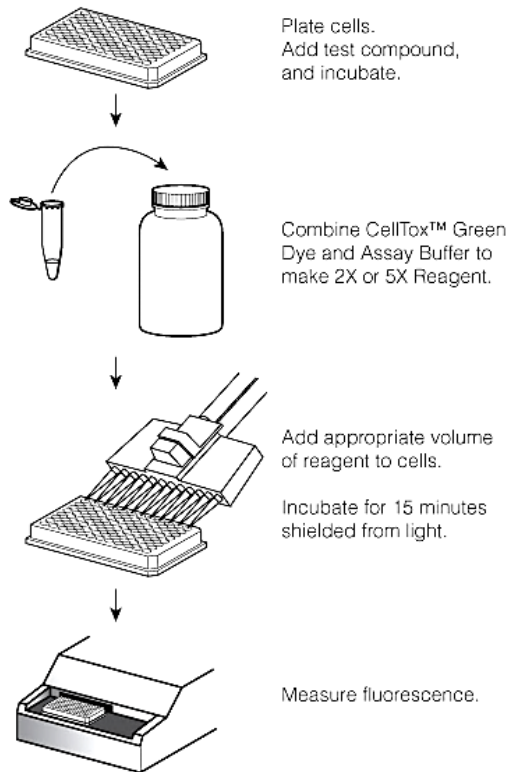
Spheroids were formed using “hanging drop method” courtesy of InSphero

CellTox™ Green is non-toxic and doesn't influence dose-dependent cytotoxicity measures

Three Reagent Delivery Formats

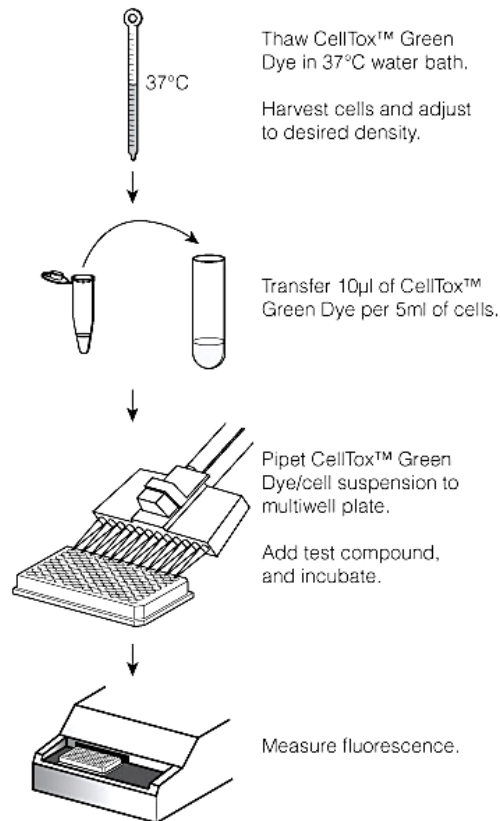
Endpoint

A. Endpoint Method, 2X or 5X Reagent Addition



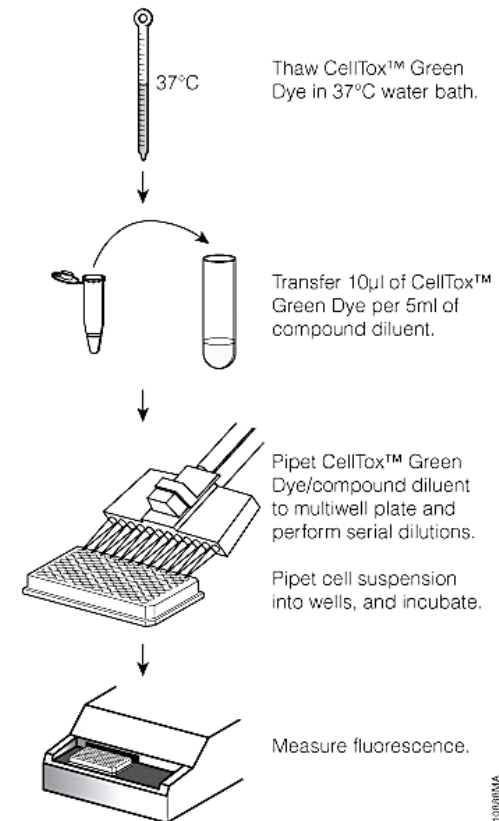
...at Seeding

B. Express, No-Step Addition at Seeding



...at Dosing

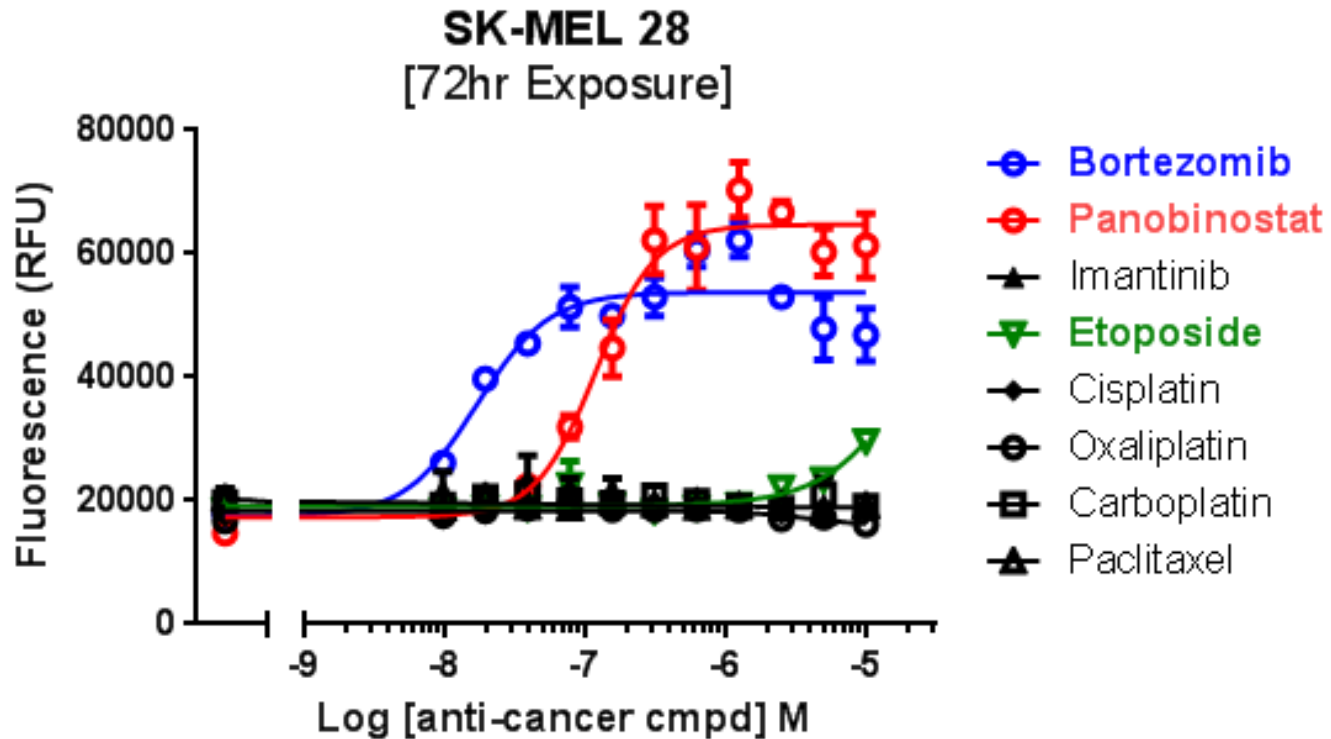
C. Express, No-Step Addition at Dosing



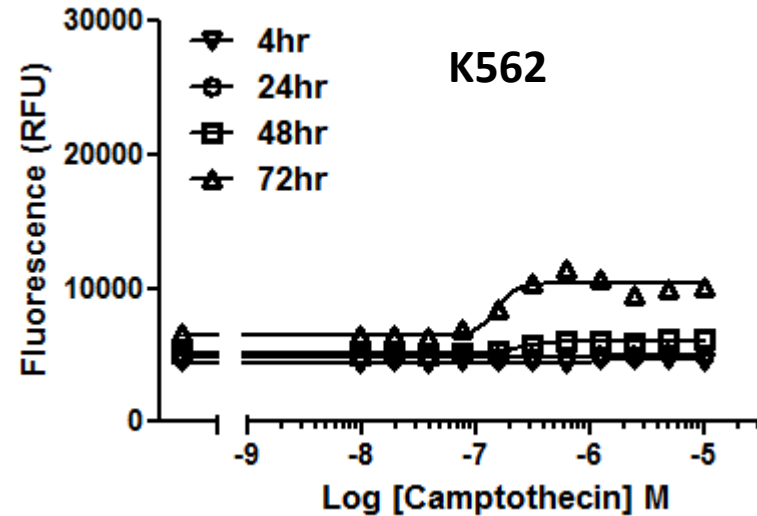
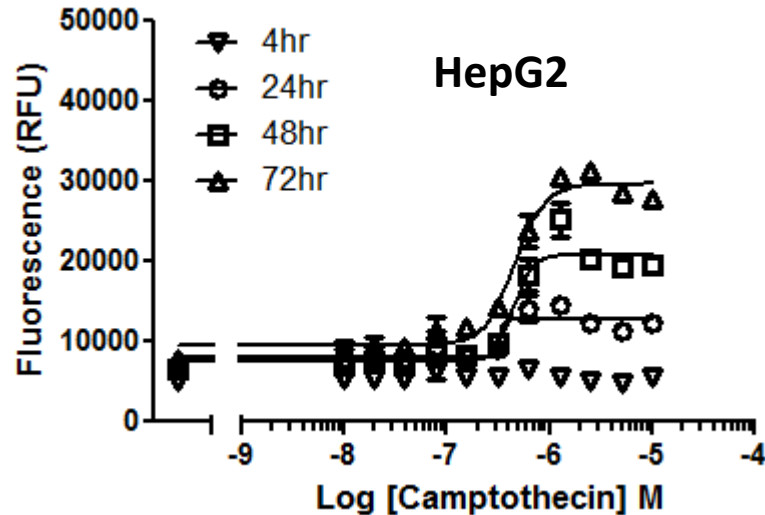
100881MA

Rank Order Cytotoxicity for Chemo-sensitivity Testing

No confounding results due to cell death detection chemistry used.



Real-Time Cytotoxicity Format Reduces Effort



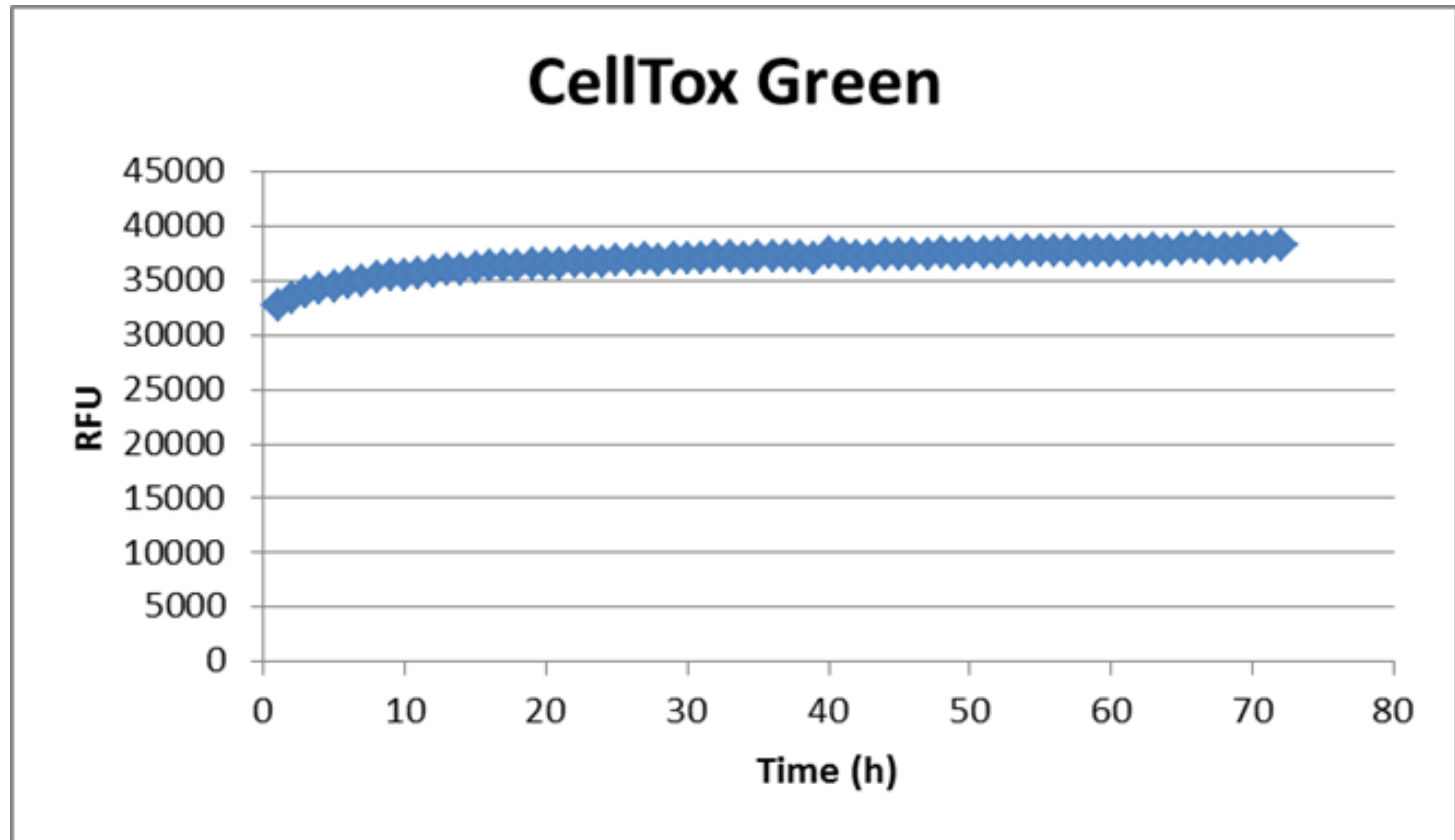
CellTox™ Green Dye added at plating or dosing:

1 plate measured/re-measured at 4-72 hours

Typical End-Point Assays:

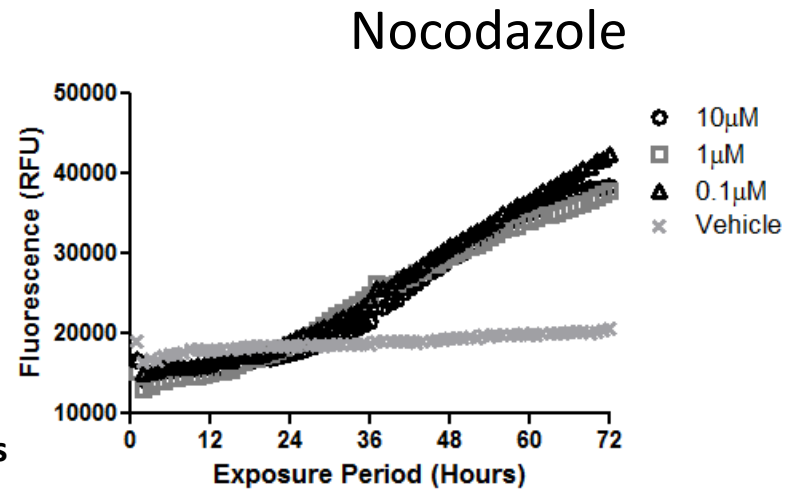
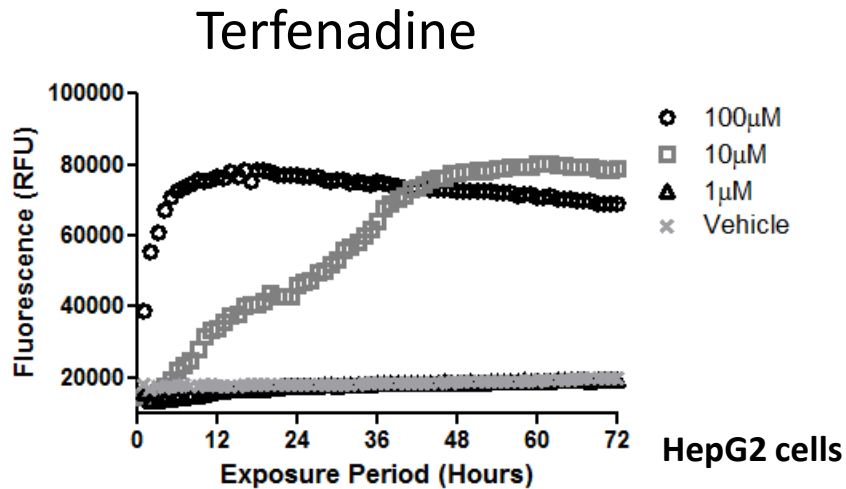
4 plates to measure at 4, 24, 48 and 72 hours

Photo-Stable Reagent Format



Fluorescence measured every hour for 72h

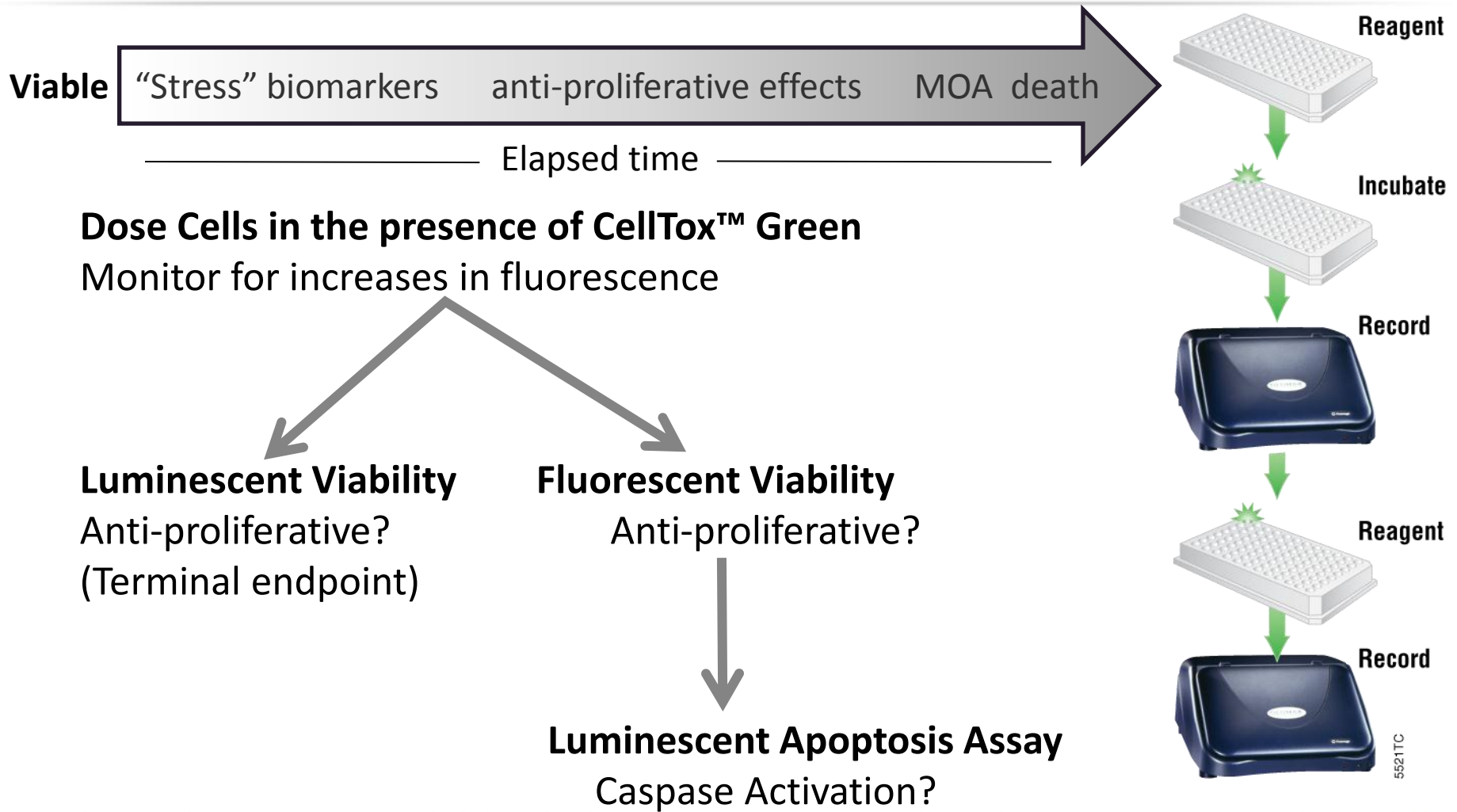
Real Time Kinetic Emergence



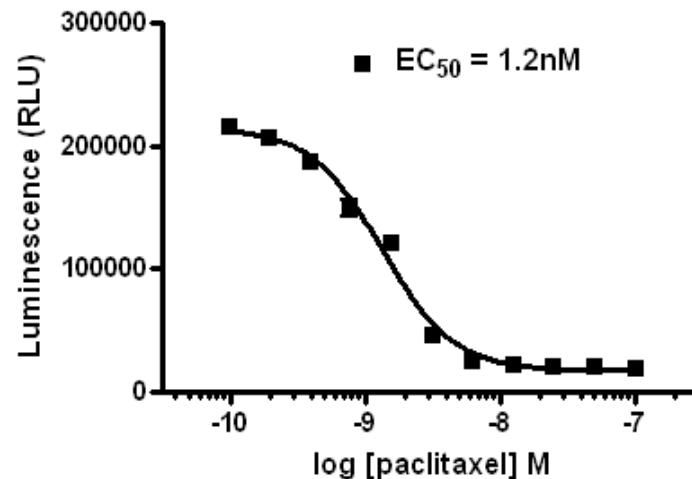
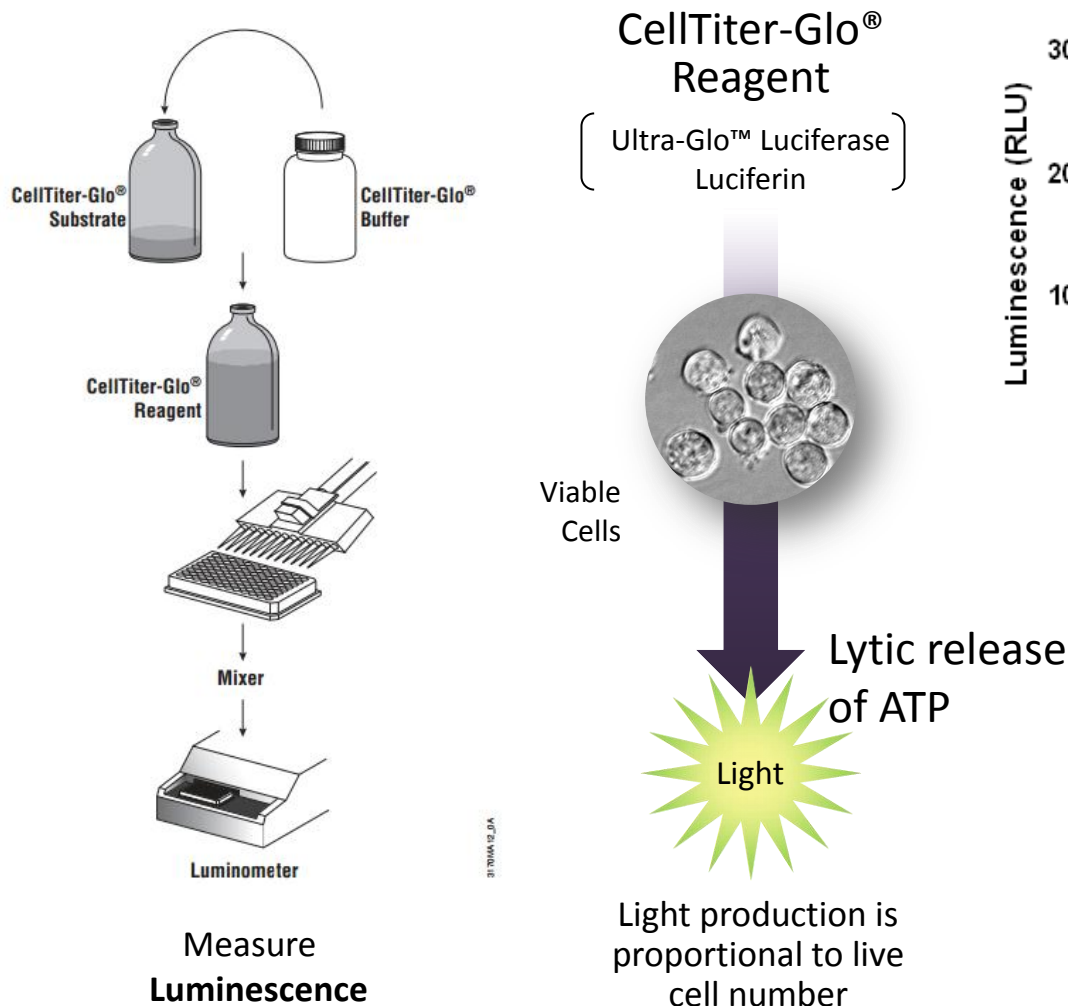
CellTox™ Green Dye added at dosing. Fluorescence measured every hour using Tecan Infinite™ 200 Pro with Gas Control Module™.



Examining Cytotoxic MOA by Multiplexing



CellTiter-Glo[®]: An ATP Viability Assay



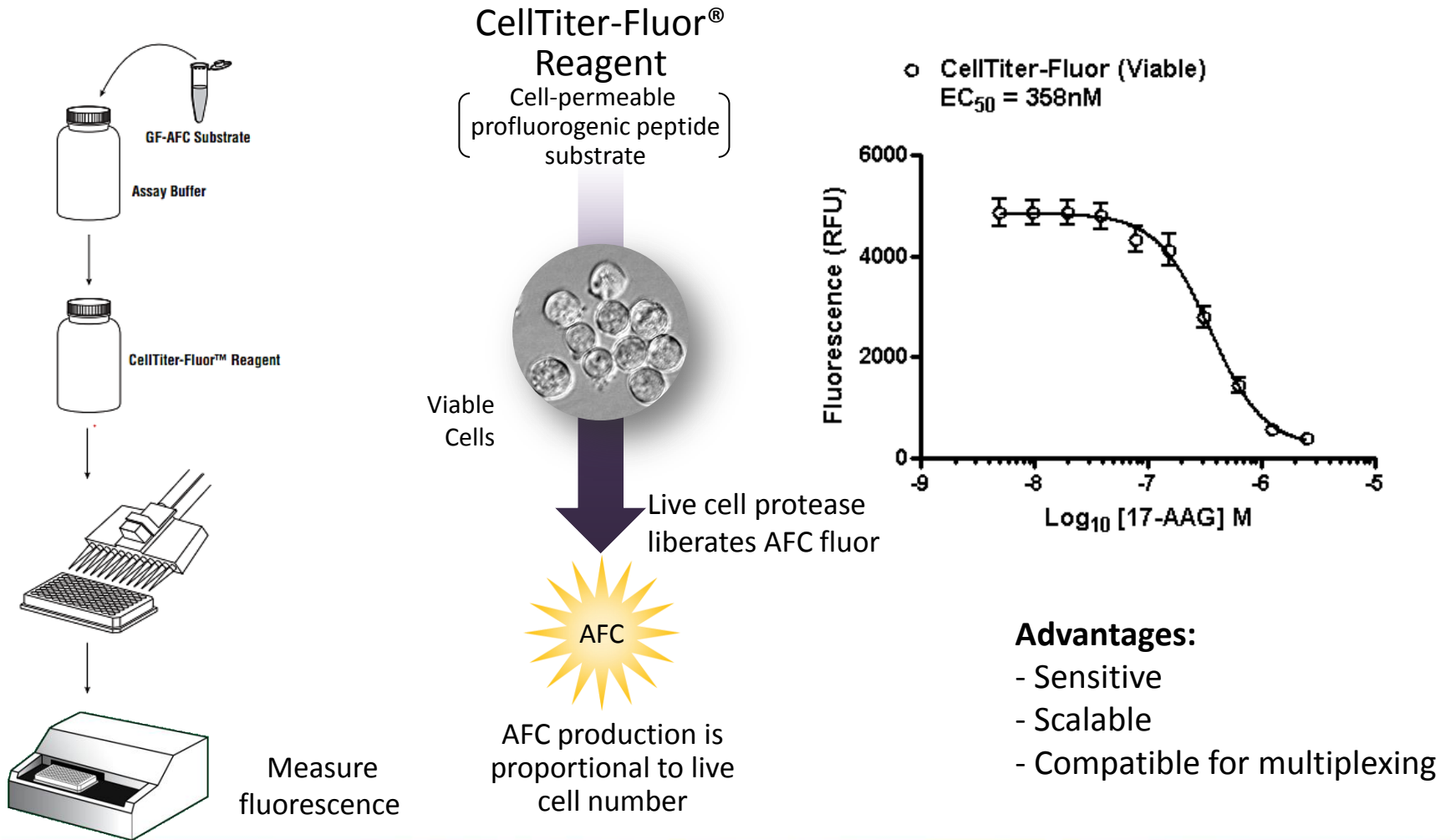
Advantages:

- Sensitivity
- Scalability
- Speed to first result

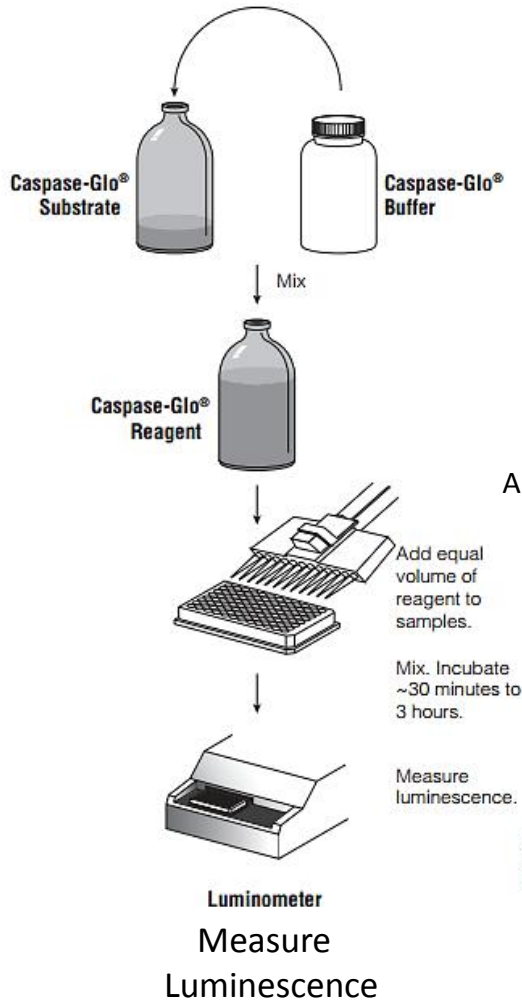
Disadvantages:

- Terminal assay
- Requires luminometer

CellTiter-Fluor™: A Protease Viability Assay

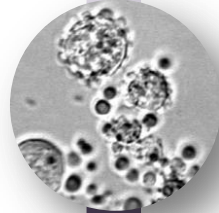


Caspase-Glo® 3/7: Caspase Activity Assay



Caspase-Glo® 3/7 Reagent

Z-DEVD-Luciferin
Ultra-Glo™ Luciferase
ATP



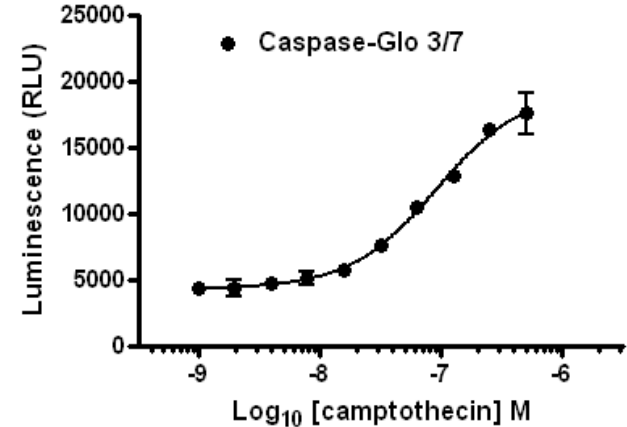
Apoptotic Cells

Activated caspase cleaves substrate liberating aminoluciferin



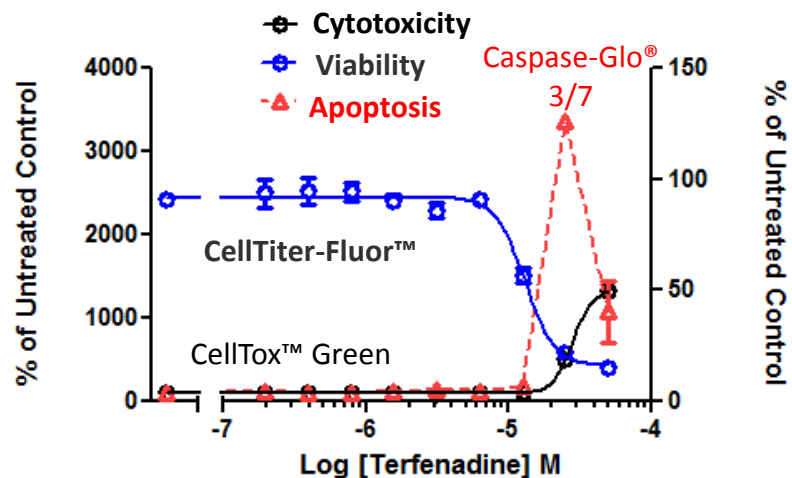
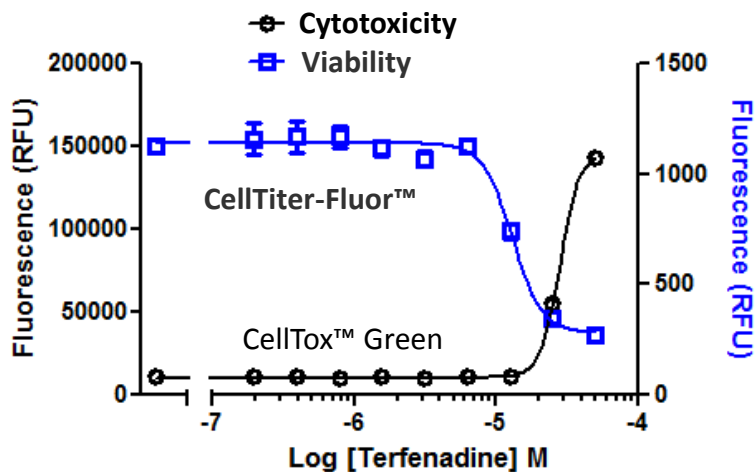
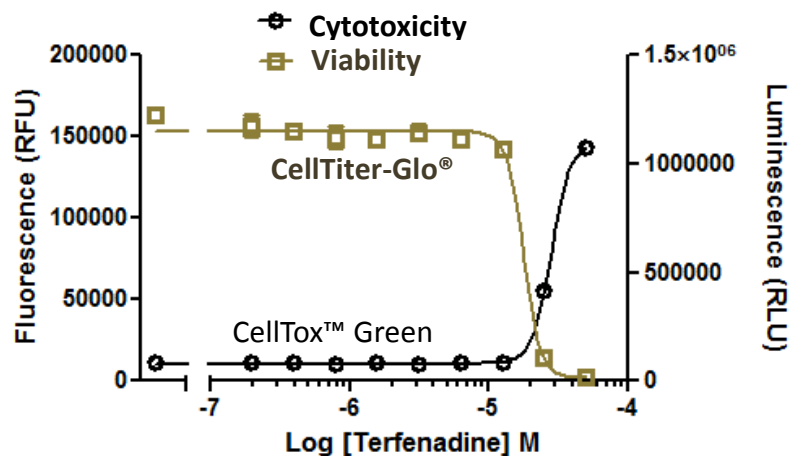
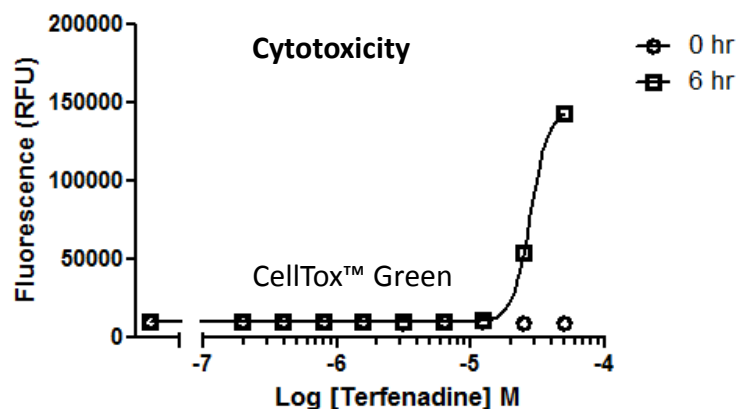
Light

Light production is proportional to caspase-3/7 activity

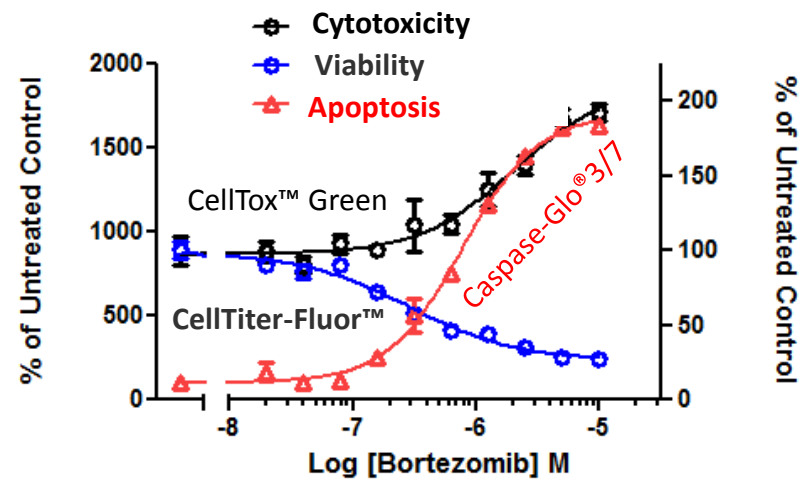
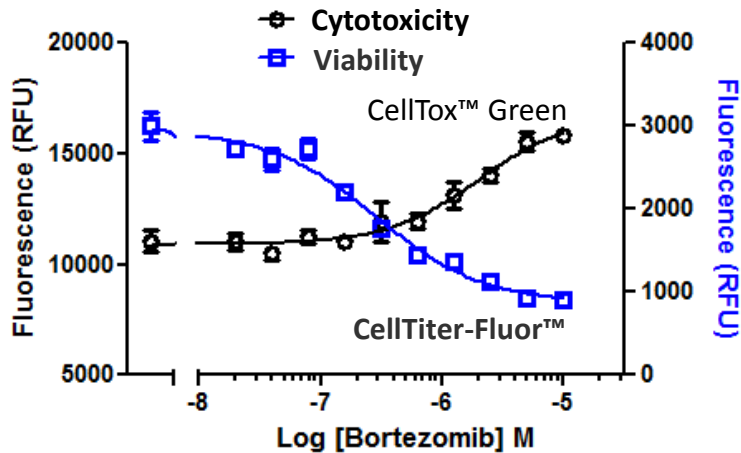
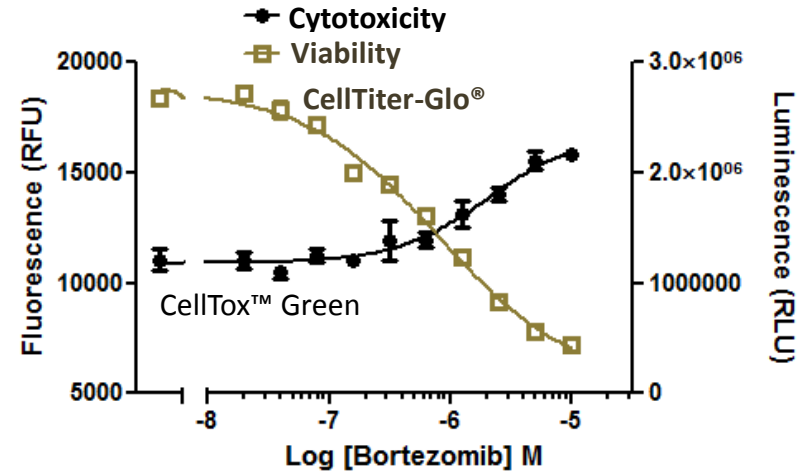
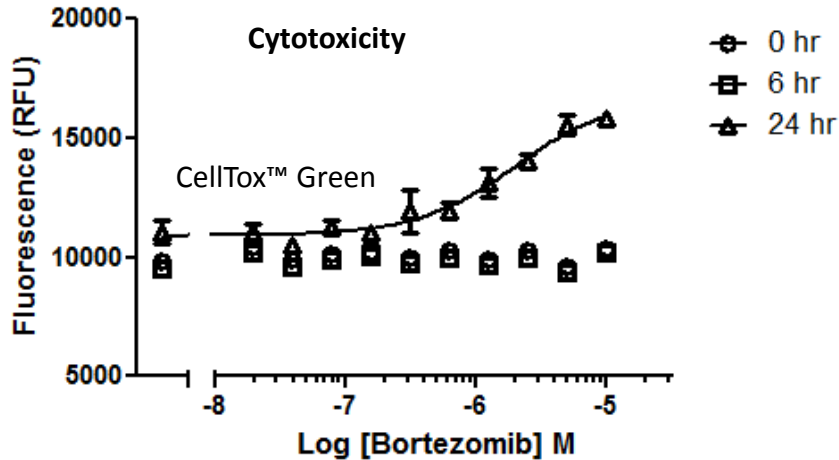


Caspase activity is definitive for apoptosis

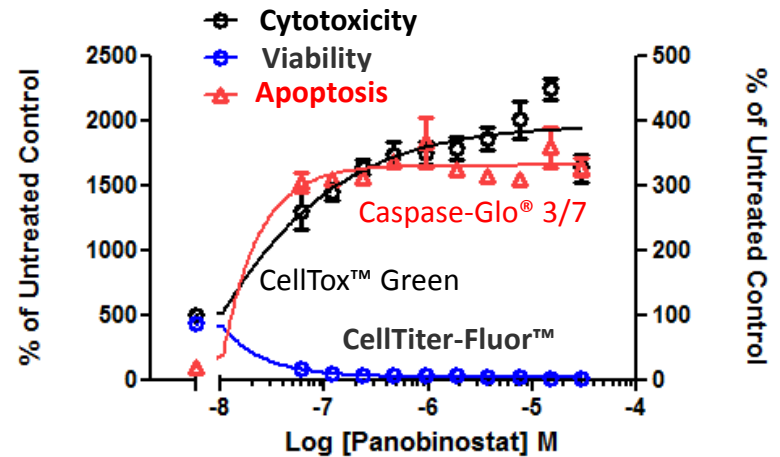
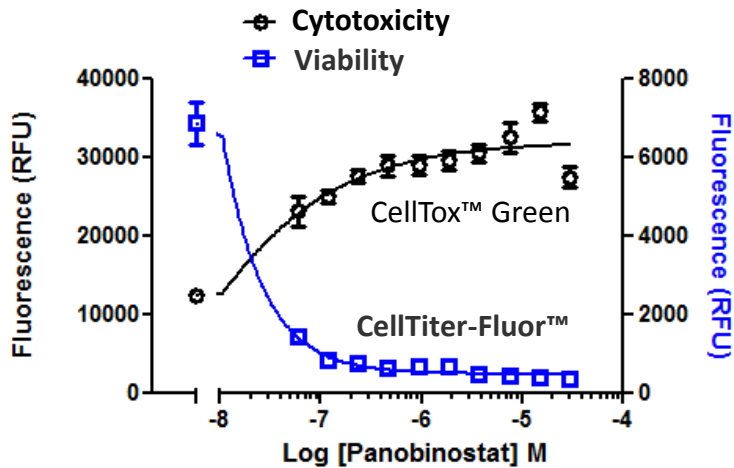
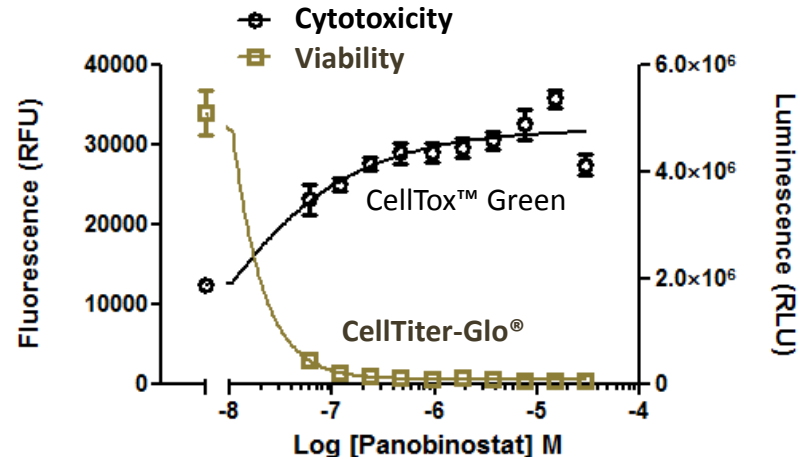
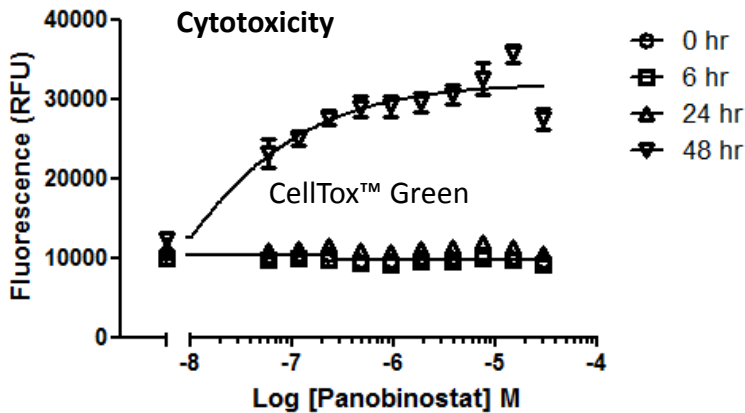
Terfenadine (Non-Sedating Antihistamine)



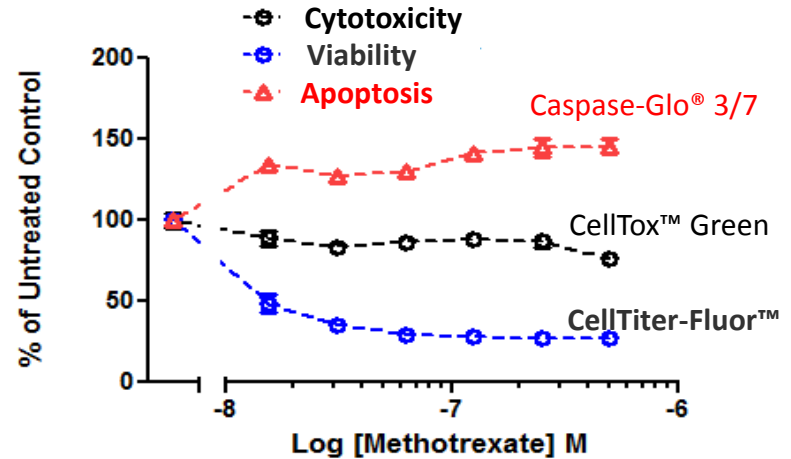
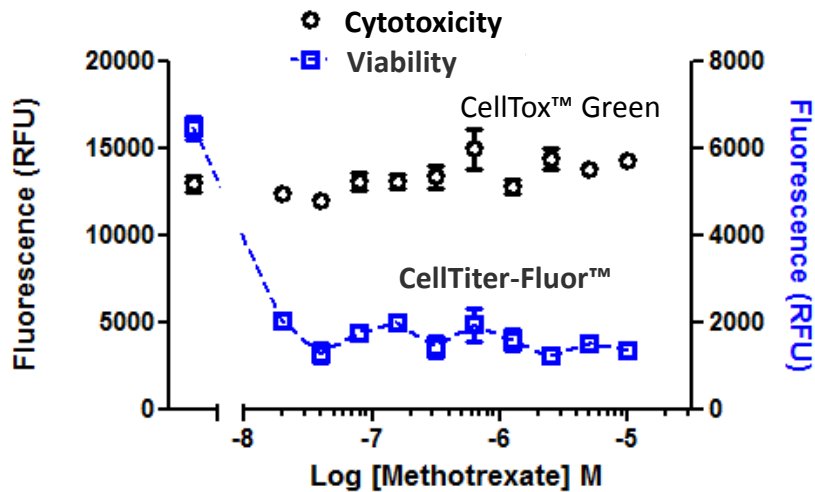
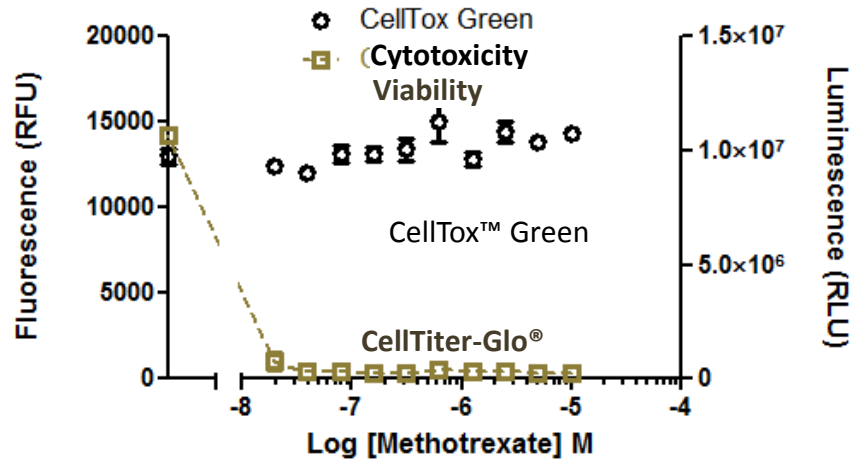
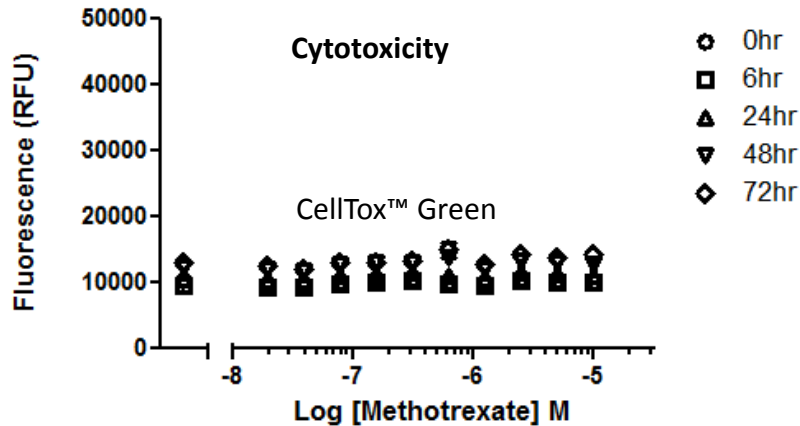
Bortezomib (Proteasome Inhibitor)



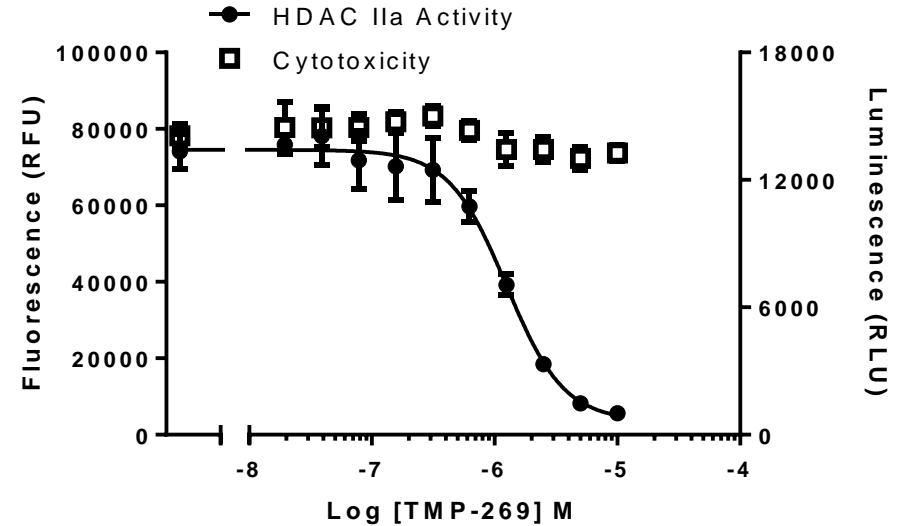
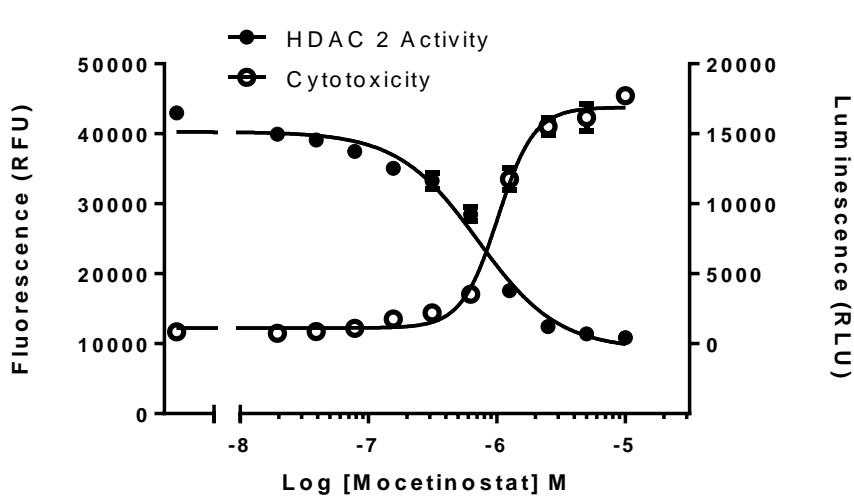
Panobinostat (HDAC Inhibitor)



Methotrexate (anti-cancer metabolite)

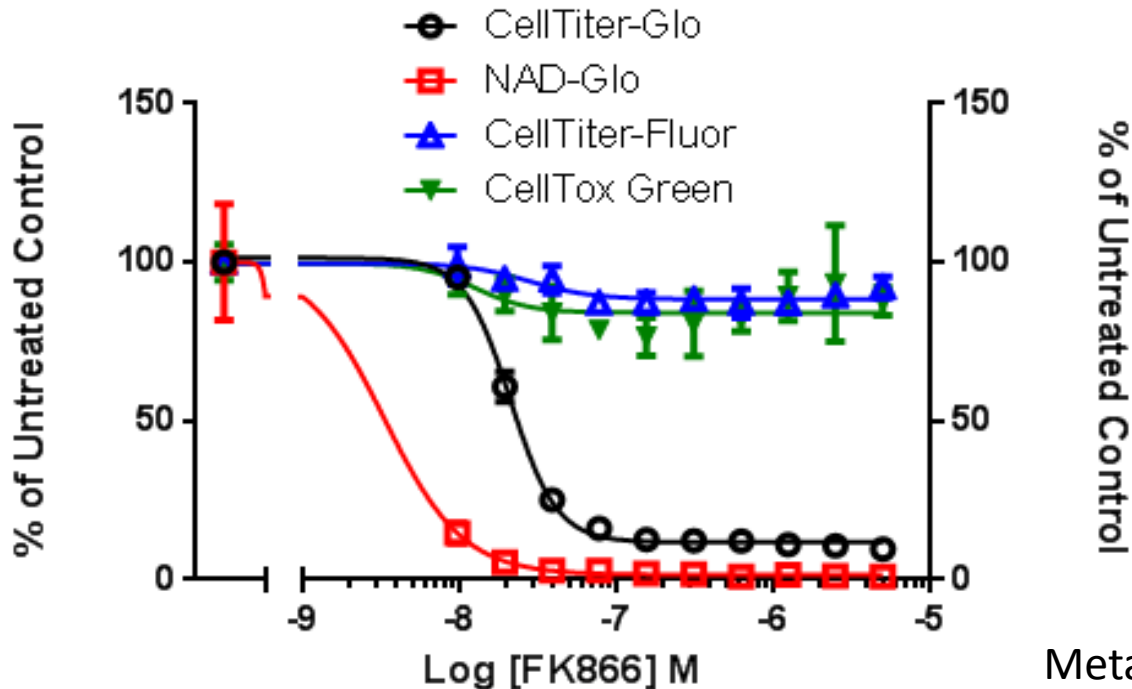


Consequences of Selective HDAC Class Inhibition



K562 were dosed with HDACi for 72hr with CellTox™ Green. Cytotoxicity fluorescence was gathered then HDAC 2 or IIa activity determined by HDAC-Glo 2 or IIa Assays in a multiplexed, same-well format.

General vs. Selective Pathway Effects



Metabolism biomarkers (NAD⁺ and ATP) can be reduced without loss of membrane integrity (CellTox Green) or reduction in non-metabolic biomarkers (CellTiter-Fluor)

Summary and Conclusions

Determination of cytotoxicity is greatly dependent upon:

- Dosage (“The dose makes the poison.”- Paracelsus in 1500’s)
- Exposure period with cells
- Mechanism of Action (primary vs. secondary necrosis, cytostasis)
- Intrinsic Cell Susceptibility (on- vs. off-target toxicity)

Not all cytotoxicity assays are created equal.

- Activity assays using endpoints are useful but imperfect
- CellTox™ Green allows for real-time measurement of cytotoxicity, and sequential multiplexing with viability and MOA assays for more complete characterization of the response

Acknowledgements

Promega Scientists and Support Staff Worldwide!



Thank You!