

Tips for Multiplexing Cell-Based Assays: Plan for success

Fall 2010



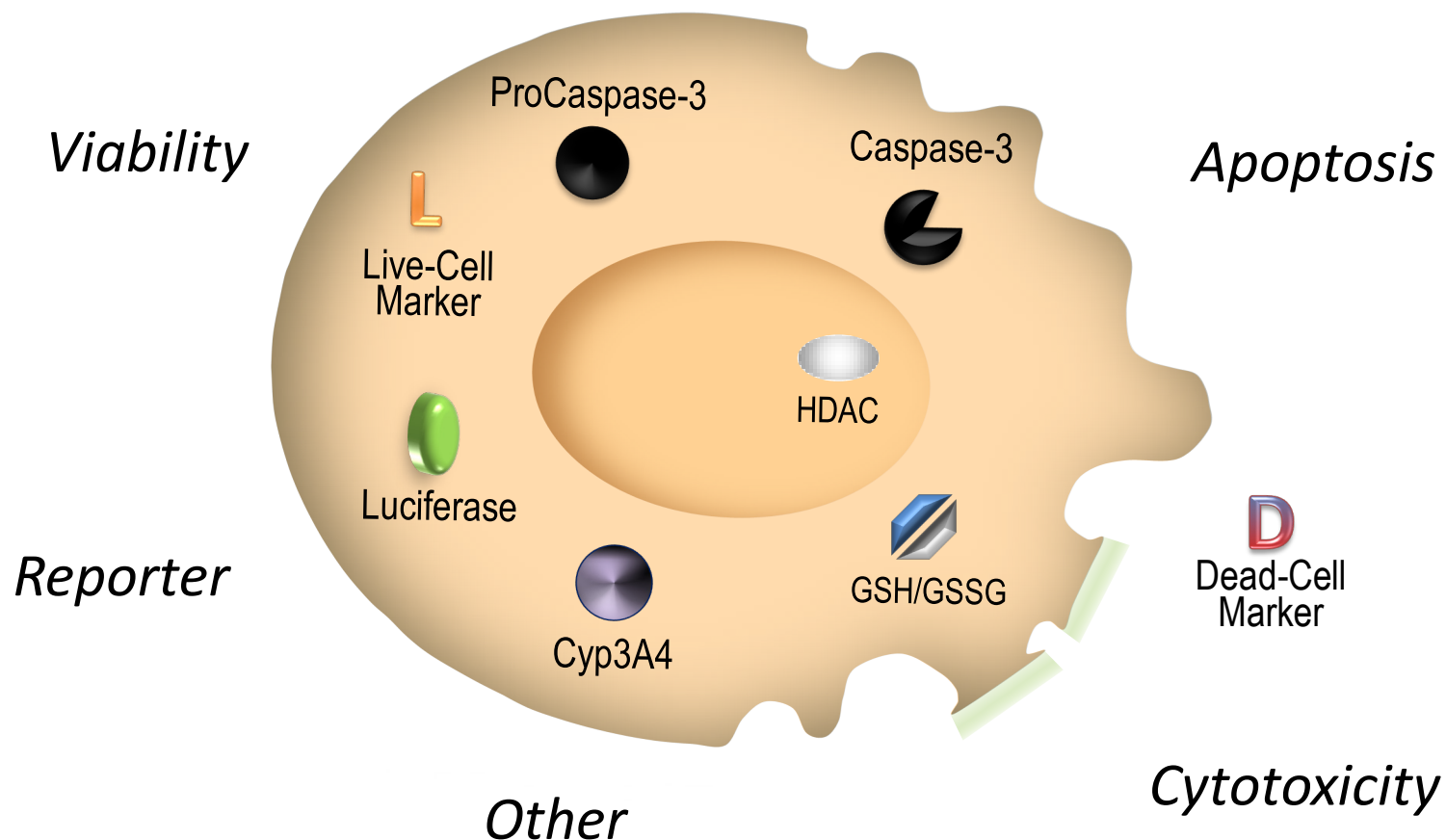
Click this icon to view
speakers notes for
each slide.



Considerations for Successful Cell-Based Assays



Assay Choices

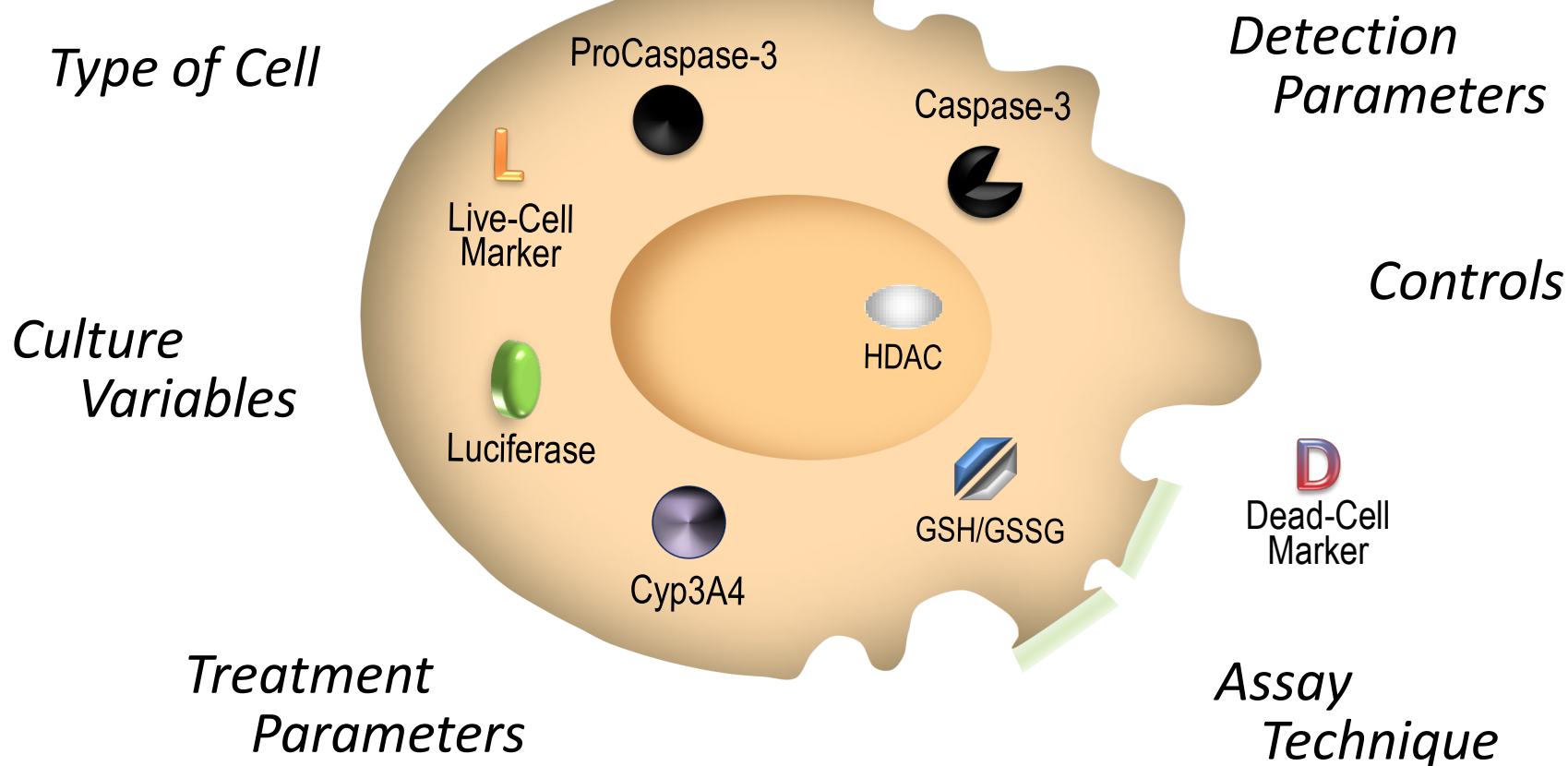




Considerations for Successful Cell-Based Assays



Assay Design/Process

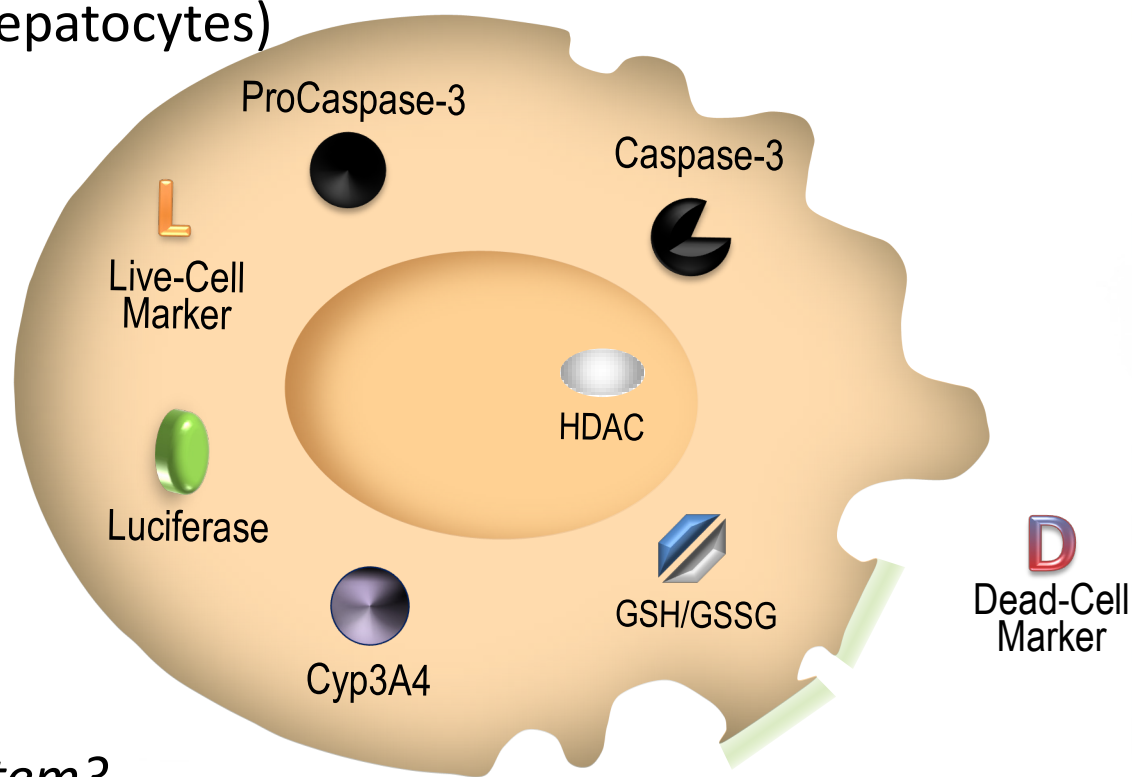




Considerations for Successful Cell-Based Assays

- Fibroblasts (HEK293, Cos)
- Cancer cell lines (HepG2, PC-3)
- Primary cells (HUVECs, hepatocytes)

Type of Cell



- *Amenable to assay?*
- *Faithfully represent system?*
- *Express factors, signaling intermediates?*



Choice of cell type influences assay design

- **Viability** - metabolism higher in cell lines, lower in primary cells
- **Cytotoxicity** – background higher in primary cells
- **Apoptosis** - differences in susceptibility to inducers

Assay parameters that may need to be altered:

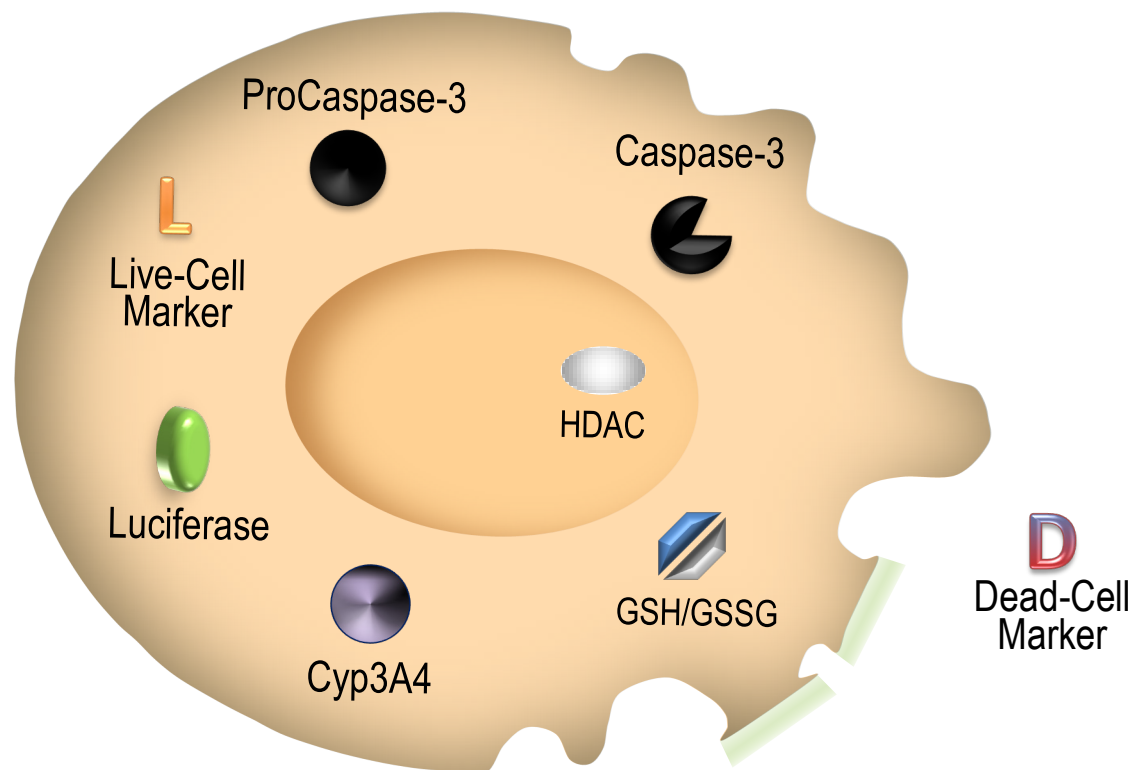
- Seeding density
- Treatment choice, concentration, duration
- Assay incubation time



Considerations for Successful Cell-Based Assays



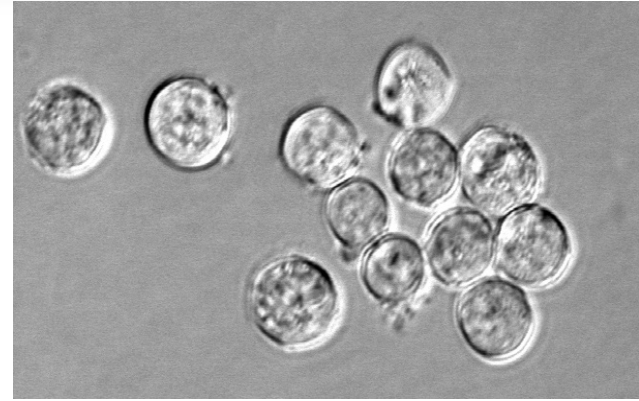
Culture Variables





Basic Cell Culture Considerations

- Cell health
 - Medium
 - *High quality, fresh*
 - *Consistent source*
 - Handling
 - *Trypsinization*
 - *Pipeting*
 - *Counting (viable cells)*
 - Contamination
- Cell line identity
 - *Verification required by NIH, many journals*



*Consistent, careful handling
of the cells is crucial
for reliable, repeatable results*



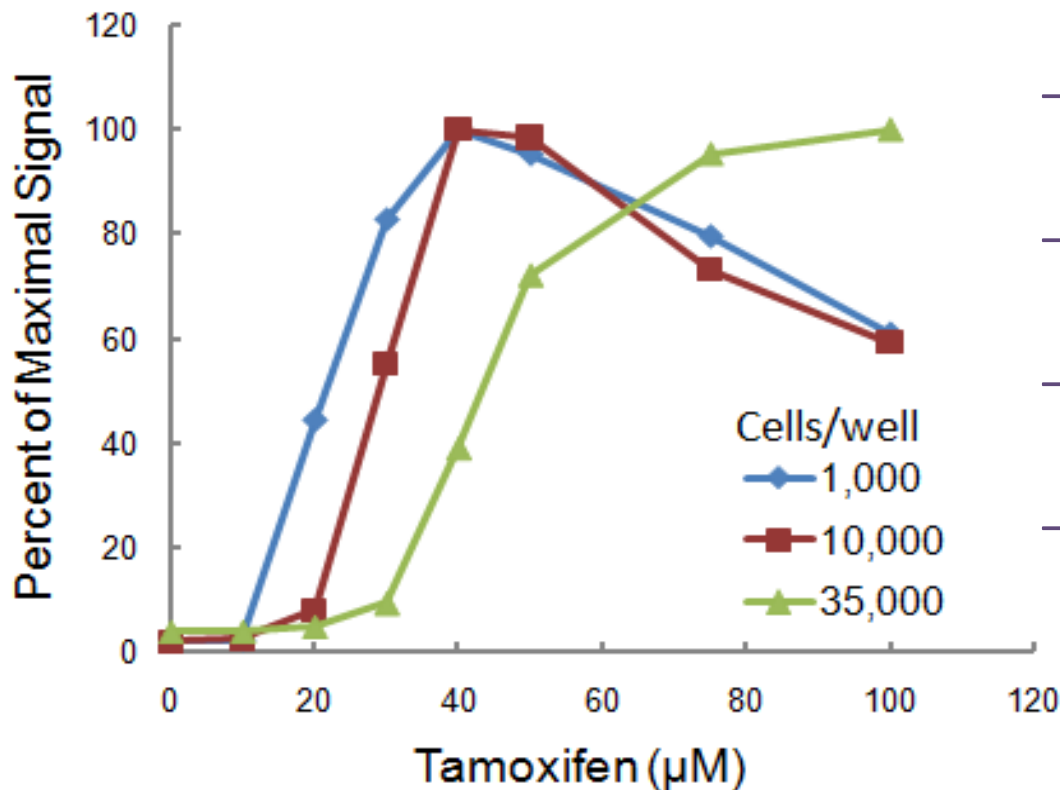
More Cell Culture Considerations

- Cell line passage
 - As passage number increases, cells may change character:
 - *Change in response to treatments?*
 - *Line may change character if allowed to reach confluence*
 - *Use low passage number – keep passage number consistent*
 - *Keep track of cultures when bulking up & during experiment*
- Cell density
 - Seeding density & cell density at time of treatment
 - Will the cells become confluent by treatment?
 - *Will cell density influence response to treatment?*
 - *Will the culture be transfected? (% confluence important parameter)*
 - *Do cells need to be equilibrated, or assay begun immediately?*



Cell number/density may alter dose response

- Induction of apoptosis as measured by Caspase-Glo® 3/7 Assay
- Replicate Dose Response Curves (DRC) from different seeding densities



- *Sparser cultures show increased sensitivity*
- *Higher densities may provide protective effect...*
- *Higher density cultures may need longer exposure*
- *Overgrown cultures will have a higher background*



Cell number optimization for viability and/or cytotoxicity assay



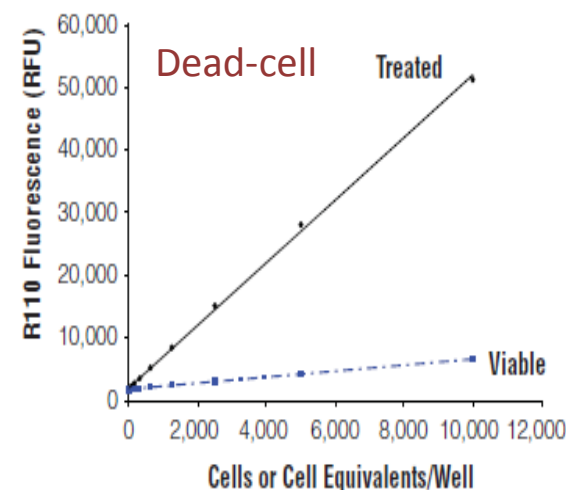
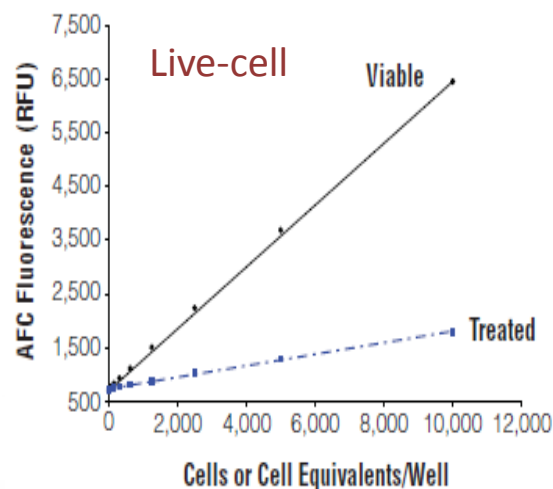
Control assay with serial dilution of cells

- *Determines sensitivity & range of assay*
- *Guides decision on seeding density, timing between plating, treatment, assay*

For live/dead multiplex, duplicate sets, one treated & one untreated

Table 1. Schematic of 96-well plate layout.

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|-----------|-------------------|---|---|---|---|---|---------|---|---|----|----|----|
| A | 10,000 Cells/Well | | | | | | | | | | | |
| B | 5,000 Cells/Well | | | | | | | | | | | |
| C | 2,500 Cells/Well | | | | | | | | | | | |
| D | 1,250 Cells/Well | | | | | | | | | | | |
| E | 625 Cells/Well | | | | | | | | | | | |
| F | 313 Cells/Well | | | | | | | | | | | |
| G | 156 Cells/Well | | | | | | | | | | | |
| H | 0 Cells/Well | | | | | | | | | | | |
| untreated | | | | | | | treated | | | | | |

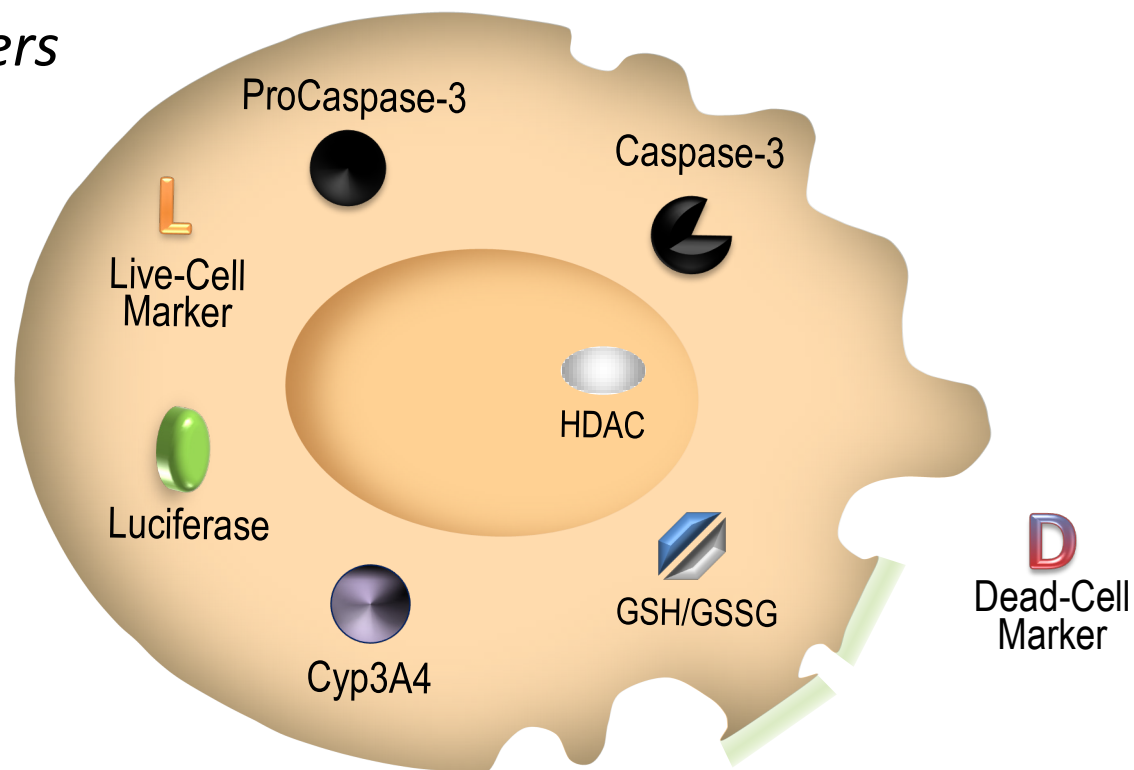




Considerations for Successful Cell-Based Assays

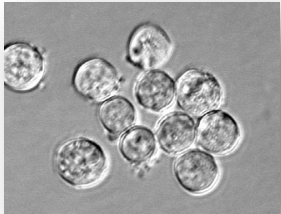


Treatment Parameters



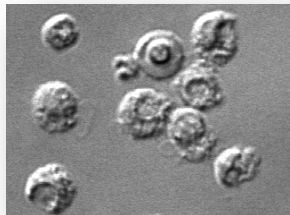


Treatment Parameters



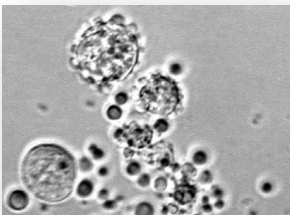
Dose

- Affects the synchrony & timing of response
- May even determine *type* of response
 - *Some drugs which induce apoptosis may be acutely cytotoxic at higher doses*



Timing

- Treatment → Assay
 - *Optimal timing affected by type of cell, compound, dose*

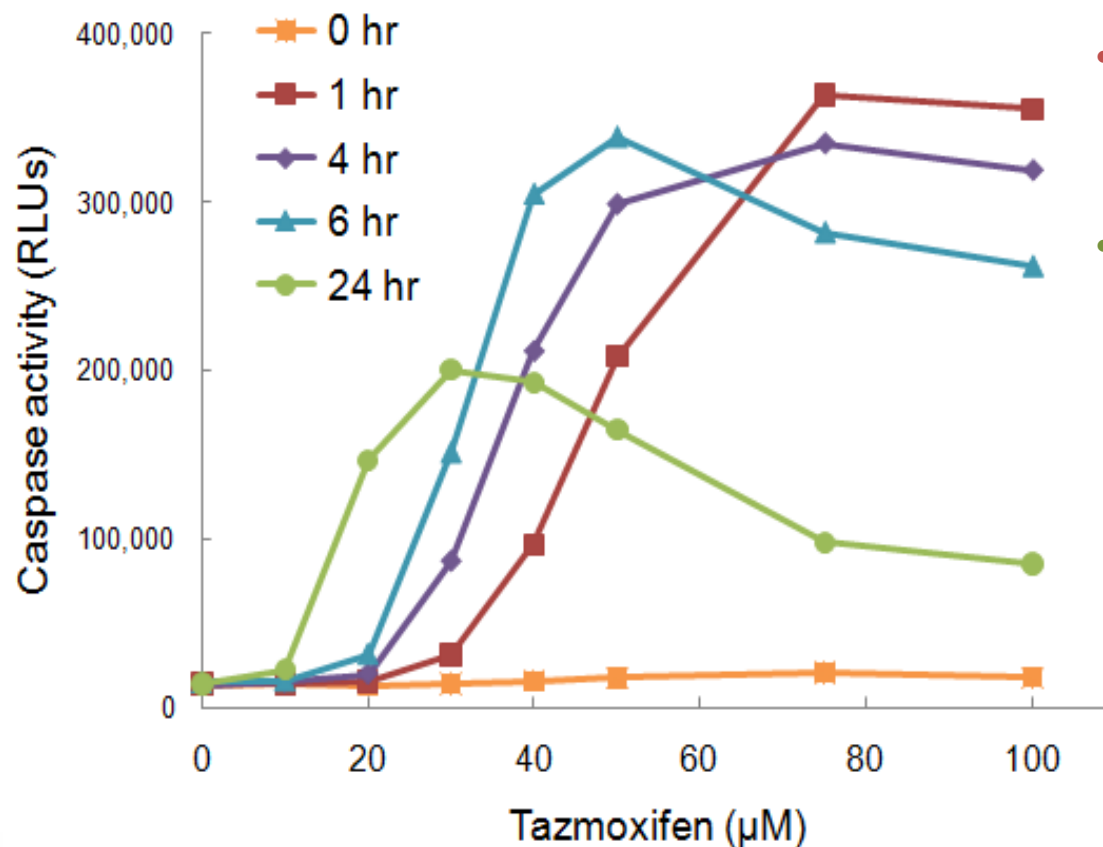


All assay parameters may interact, but dose timing & cell type are the most important & intertwined



Timing and dose interact to shape response

- Induction of apoptosis as measured by Caspase-Glo® 3/7 Assay
- Replicate Dose Response Curves (DRC) measured at different times

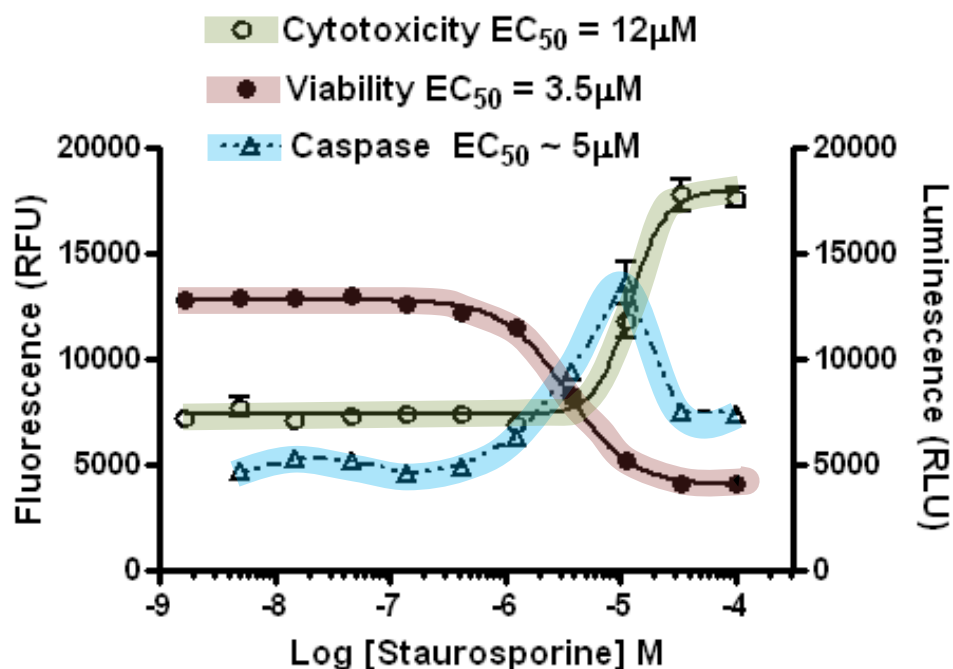


- *At 1 hour max response at higher doses; no response seen at lower doses*
- *At 24hr max response at lower doses; at higher doses response missed, analyte has greatly diminished*

DRC & timecourse control assays are often worth the investment

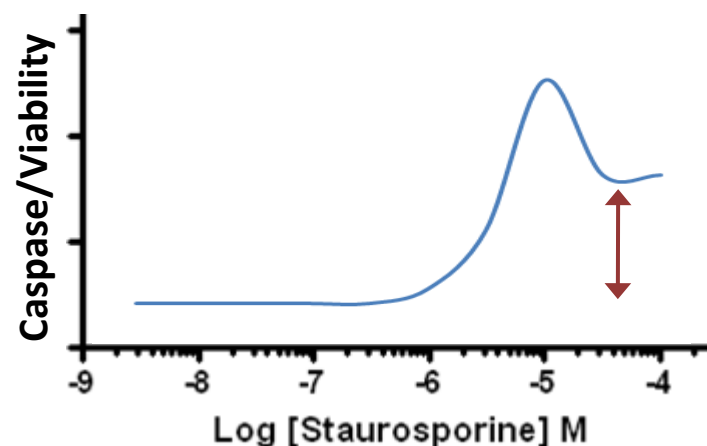


Multiplexing makes more robust assays: Caspase data compared to live/dead data



- Induction of apoptosis is asynchronous in a population of cells
- Cells undergoing apoptosis are still "alive"

- At highest doses caspase signal is low ...
- But it is because most cells have already completed 2° necrosis...
- Normalizing caspase to viability score shows there is still a strong induction at those doses!

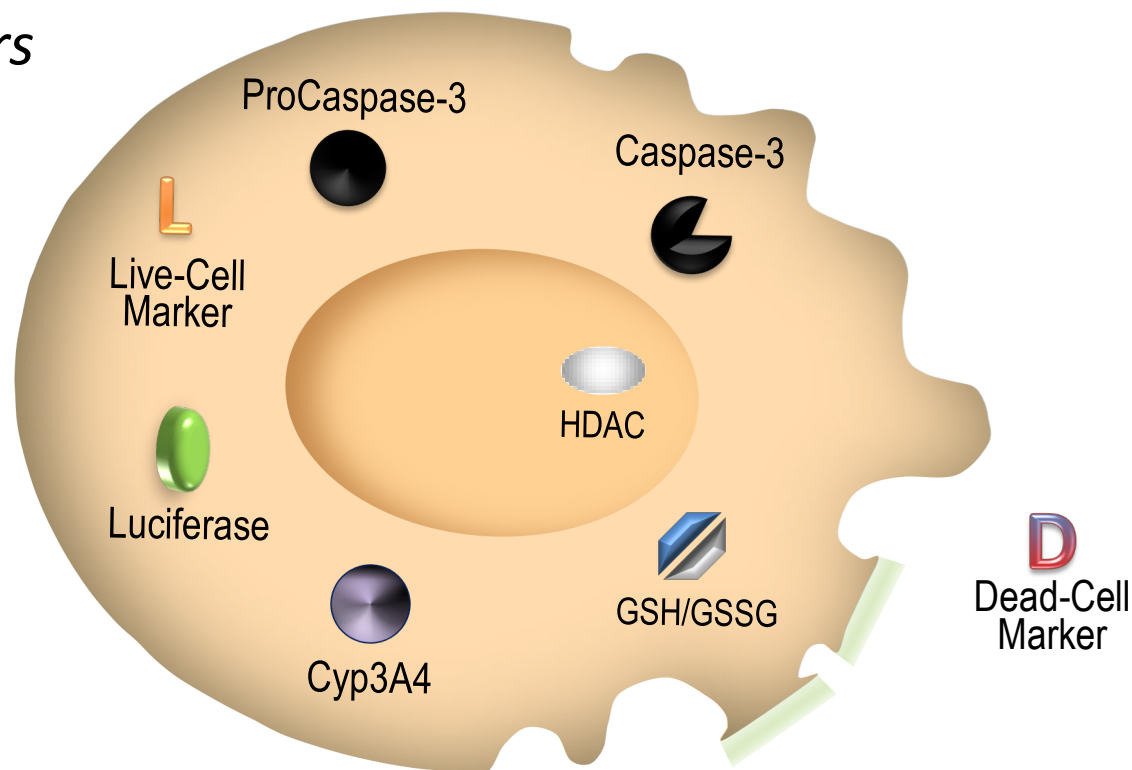




Considerations for Successful Cell-Based Assays



Detection Parameters

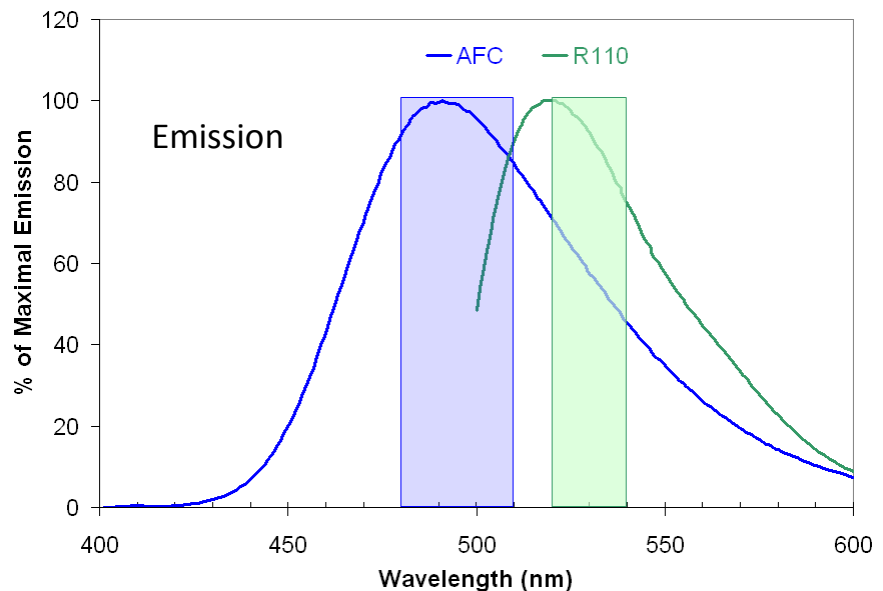
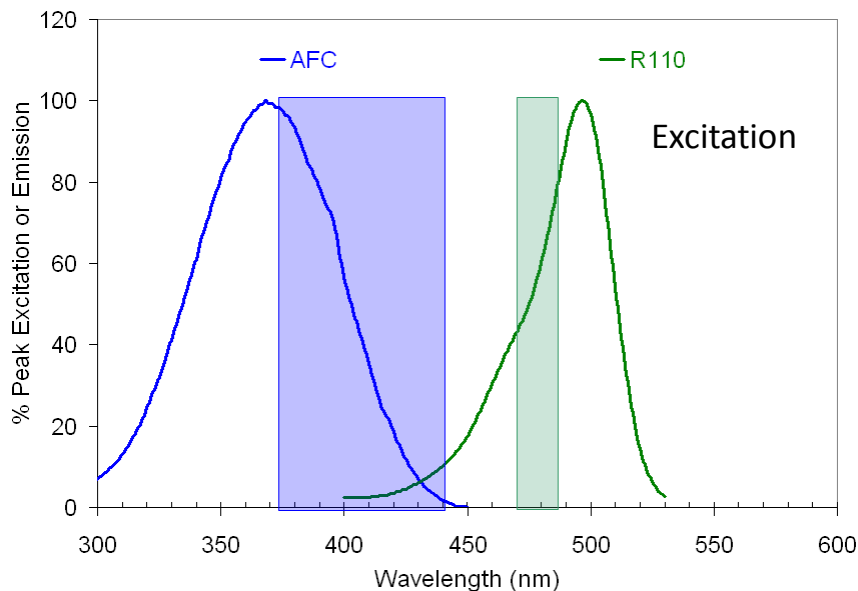




Detection Parameters – Filters & Fluorophores

The best filter settings for your multiplex may not equal the peak ex/em of the fluorophores...

- Consider the spectral overlap of the multiplexed fluorophores
- Consider the capabilities of your detection platform

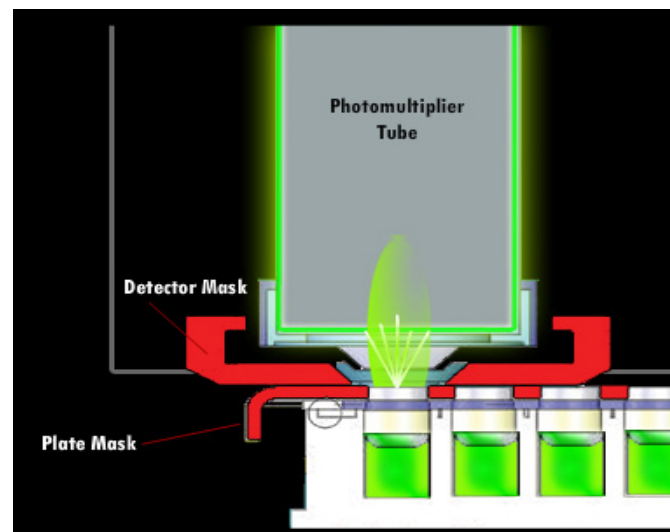
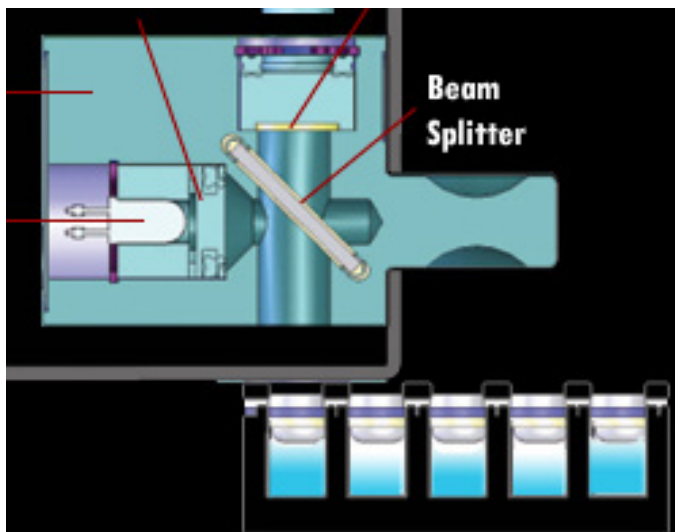


MutliTox-Fluor AFC & R110 dye spectra and example of compatible filter sets

Detection Parameters – Gain & Integration

- *Gain* – sensitivity setting
- *Integration* – time of signal collection

- *User adjustable on many instruments*
- *Settings too high or too low can cause problems*





Detection parameters affect assay performance

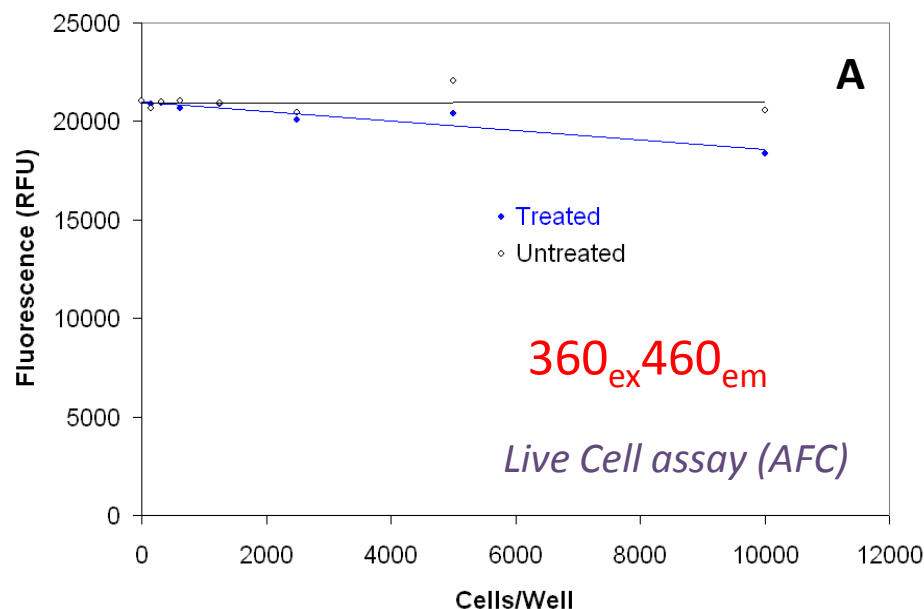


Symptoms:

*No signal \uparrow with cell # \uparrow ;
no Δ with treatment*

Cause:

Incorrect filters used!

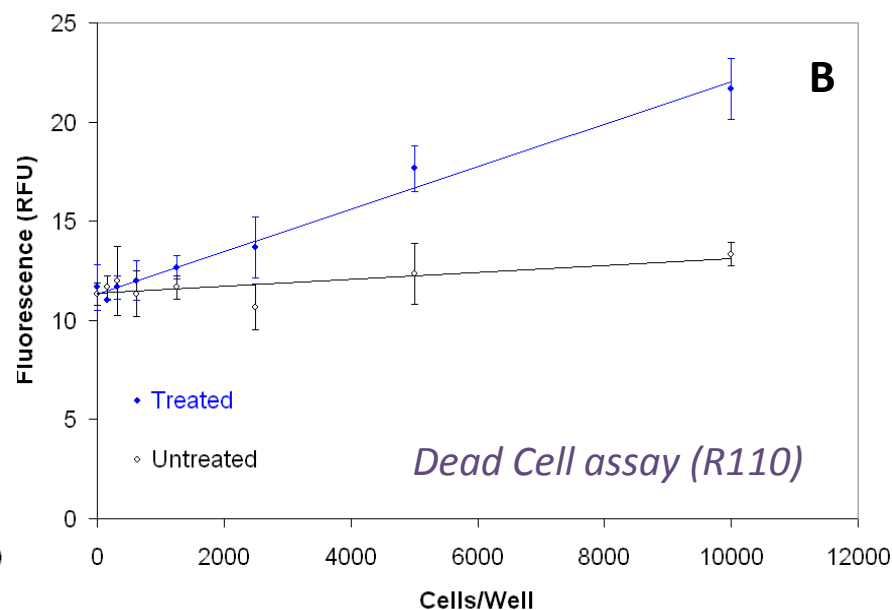


Symptoms:

*Low signal, high variation;
little change with treatment*

Cause:

Gain is too low!

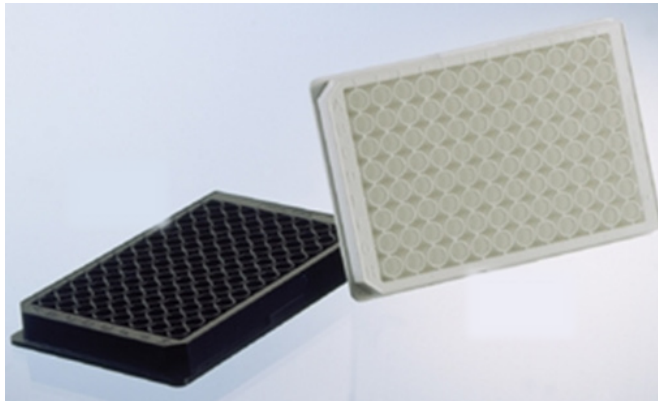


MultiTox-Fluor assay on serial dilutions of cells, +/- toxic treatment



Detection Parameters – plate choice

- *White or Black?*

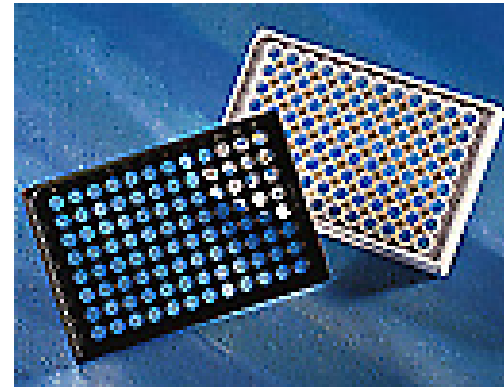


General recommendation:

- Black for fluorescence
- White for luminescence

- In reality, either plate can be used for both assays
- For multiplex we tend to use white

- *Solid or Clear Bottom?*



- Clear bottom plates can be used for viewing cultures prior to assay
- Opaque bottom gives better signal, less cross-talk

- Generally small effect on signal and cross-talk
- Opaque tape can be applied



Detection Parameters – Software/analysis issues

- The instrument software may be automatically...
 - ...*subtracting background*
 - ...*assigning & using a control sample for normalization*

If so, are the correct wells being used?

What automatic functions are being applied?

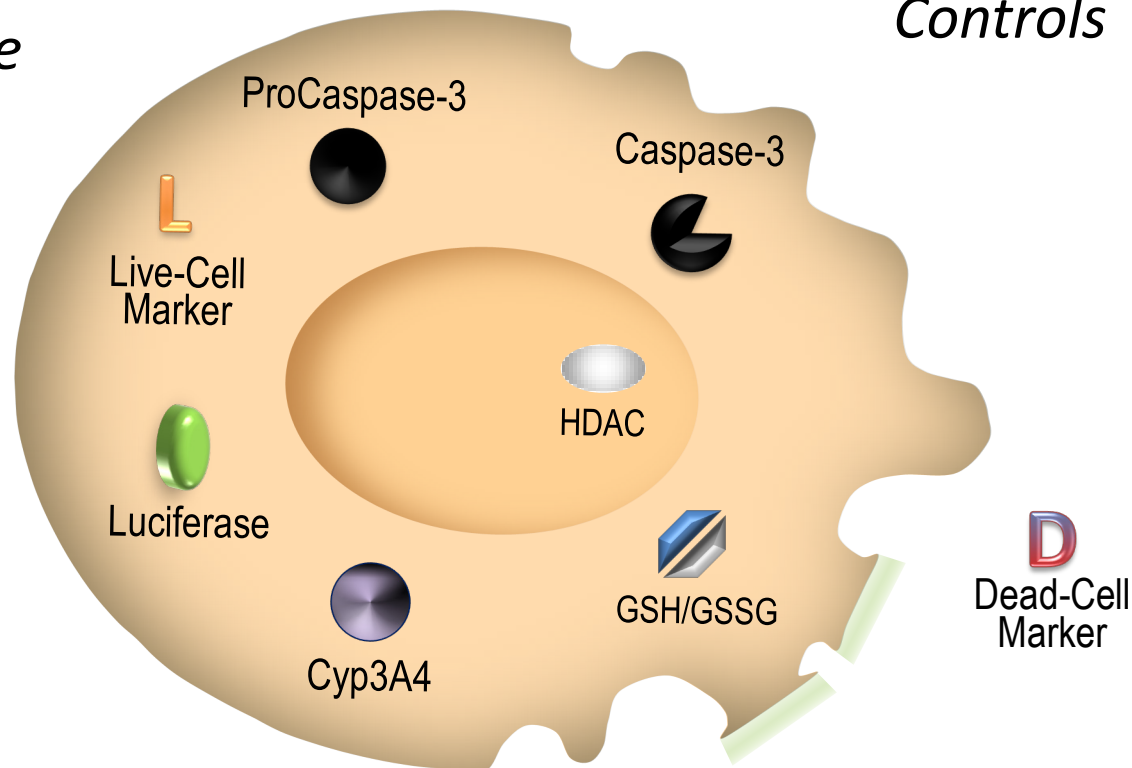
| Read 1 | 1 | 2 | 3 | 4 | 5 |
|--------|---------|---------|-----------|----------|---------|
| A | 1022620 | 976023 | | | |
| B | 745541 | 804604 | 12391.6 | | |
| C | 1011300 | 1440090 | 578110 | 11656.7 | 81.0004 |
| D | 86737.9 | 101260 | 580505 | 505791 | 94.0005 |
| E | 89070 | 101812 | 549744 | 495521 | 119.001 |
| F | 83276.4 | 92886 | 703031 | 602741 | 181.002 |
| G | 12396.1 | 17421.6 | 940443 | 680742 X | |
| H | 14852.7 | 15333.5 | 481679 | 699395 X | |
| | | 2418.31 | 471340 X | | |
| | | | 2346.29 X | | |

Considerations for Successful Cell-Based Assays



*Assay
Technique*

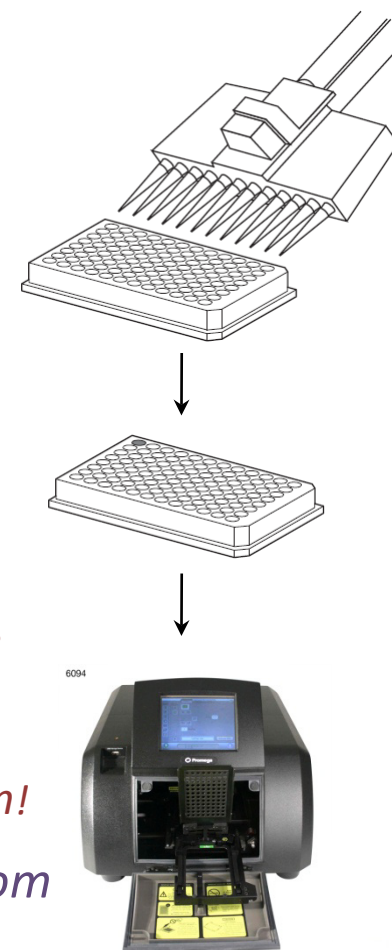
Controls



Common technical mistakes in cell-based assays



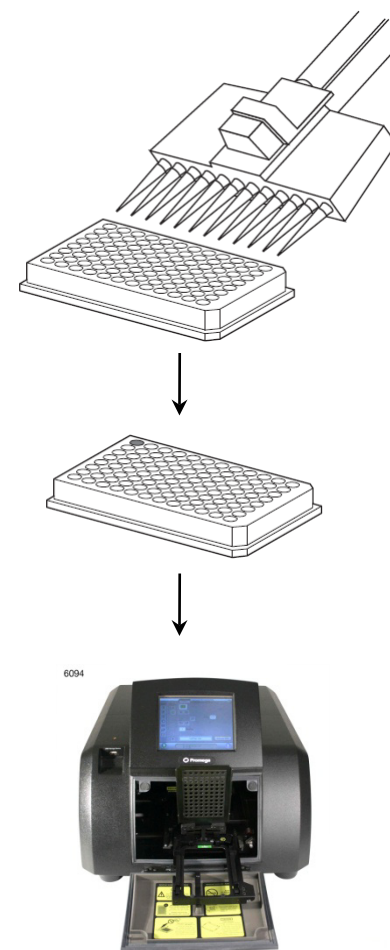
- Problems with reagent addition
 - *Introducing bubbles*
 - *Droplets of reagent on upper surfaces well*
 - *Insufficient mixing or inconsistent mixing*
- Poor temperature control during reagent incubation
 - *Enzyme assays are temperature dependent!*
 - Fluorogenic assays may be incubated at room temp or 37°C...
 - *either way, temperature should be consistent across plate*
 - *use of water bath is ideal*
 - *Beware cell culture incubators for 37°C reagent incubation!*
 - *Luminescence assays should always be incubated at room temperature (equilibrate reagent before addition)*



Common technical mistakes in cell-based assays



- Improper timing of reagent incubation
 - *Reading too soon may*
 - \downarrow *signal:noise*
 - \downarrow *sensitivity*
 - \downarrow *range*
 - \downarrow *discrimination*
 - *Incubating too long may*
 - \downarrow *range* – *higher levels of analyte may saturate assay or detector*
 - *Optimal incubation period will vary with temperature*
 - *e.g. MultiTox-Fluor assays incubated at 37°C require ~30 minutes; incubation at room temperature may require up to 3 hours*



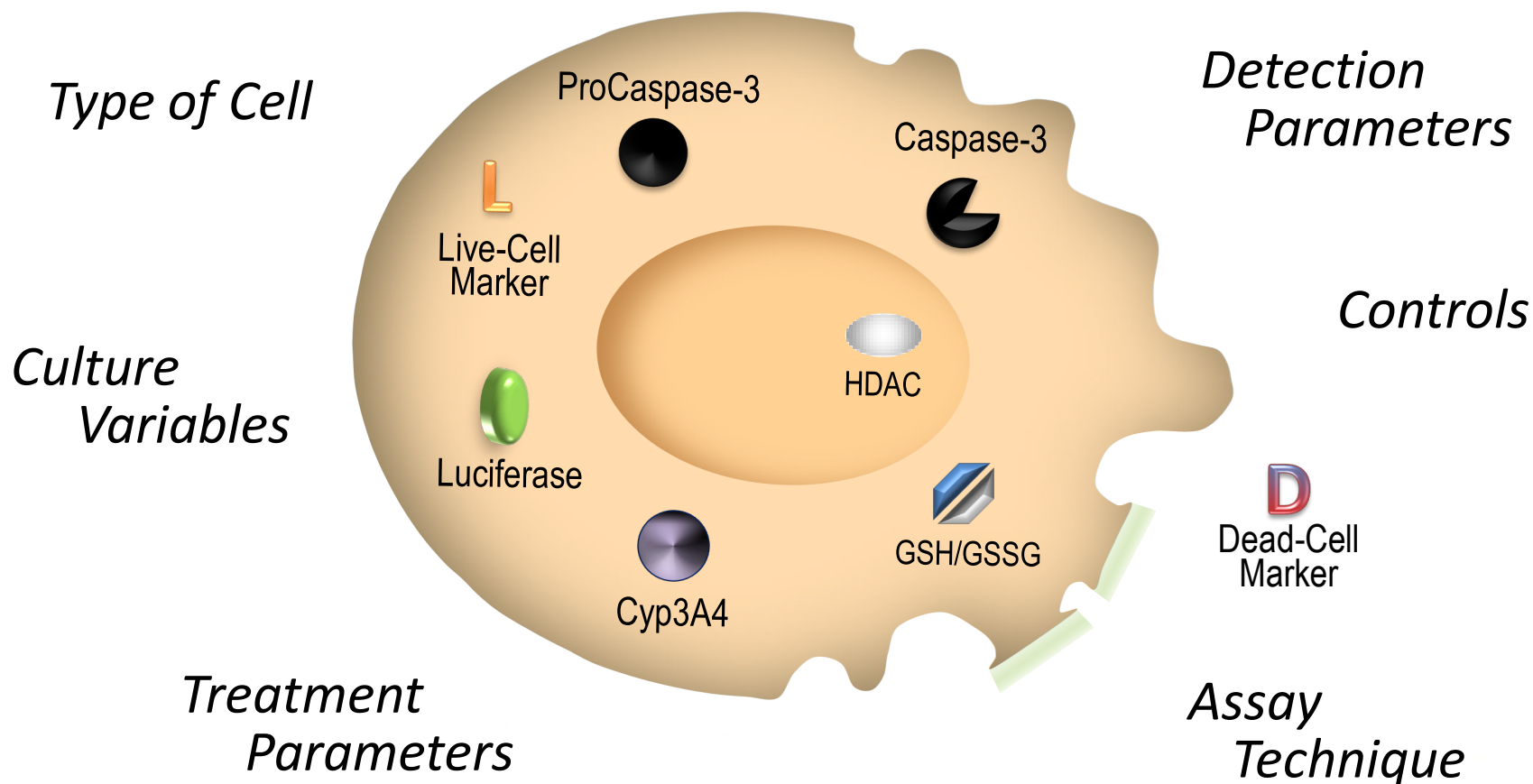
Assay controls

- No-Cell control
 - *Determines background of the assay (media/plate/reagent/detector)*
 - *Used in data processing, background subtraction*
- Untreated Cells Control
 - *Vehicle only*
 - *Also used in analysis - determining fold response, 100% viability &/or 0% cytotoxicity for relative response ratio*
- Positive Controls
 - *For cytotoxicity/apoptosis/pathway effect*
 - *Also used in analysis - determining fold response 0% viability and/or 100% cytotoxicity for relative response ratio*
- Multiplate assays
 - *Replicates of untreated control and a positive control*
 - *Controls for possible variation in processing, temp, timing, etc.*

Considerations for Successful Cell-Based Assays



Assay Design/Process



Questions?



Rely on Promega Technical Services

- Experienced & highly trained scientists
*>150 years cumulative bench experience,
>10 yrs average*
- Varied technical expertise
reporters, cell culture, HTS, etc.
- Varied scientific expertise
model systems, genetics, development, etc.
- ***Easy!*** – phone, chat, e-mail

