



# GloMax<sup>®</sup> Discover

An instrument for your assays

**Dr. Harald Winteler**  
Product Manager, Promega AG, Switzerland

Promega Corporation

©2013 Promega Corporation.

# Getting the Most From Your Plate-Based Assays

---

- Scientific discoveries involve not only finding the right assay, but also the right instrument.
- *Assay sensitivity, dynamic range, sample-to-sample cross-talk, and ease-of-use* are critical considerations when choosing an assay and instrument.
- Promega develops and QC tests many of its luminescence assays using GloMax<sup>®</sup> instruments, because of their:
  - Superior luminescence performance
  - World-wide acceptance and publications in peer-reviewed research journals

# Introducing ...

- Multi-mode detection (Luminescence, Fluorescence, UV-Vis Absorbance, BRET, FRET, filtered luminescence)
- Easy to use: Plug-n-play setup
- Tablet PC touchscreen control and wireless connectivity to network and Promega.com
- 6 to 384-well plate formats
- Heating & shaking
- Optional: Dual injector
- Automation-friendly
- IQ / OQ Service

**GloMax**  
DISCOVER



➤ Integrated with Promega assays

# Integrated with Promega Assays



## The Perfect Partner for Promega Assays

Preloaded Promega protocols  
or customize your own

### Cell Signaling & Metabolism Assays:

Including:

- ADP-Glo™
- Kinase-Glo®
- P450-Glo™
- cAMP-Glo™

### Cell Health Assays:

Including:

- CellTiter-Glo®
- CellTox™ Green
- Caspase-Glo®
- BacTiter-Glo®

### Luciferase Reporter Assays: Including:

- Nano-Glo®
- ONE-Glo™
- Dual-Glo® & DLR
- Bright-Glo™

### BRET and FRET Assays: Including:

- NanoBRET™
- Renilla/YFP
- Commercial and Homebrew assays

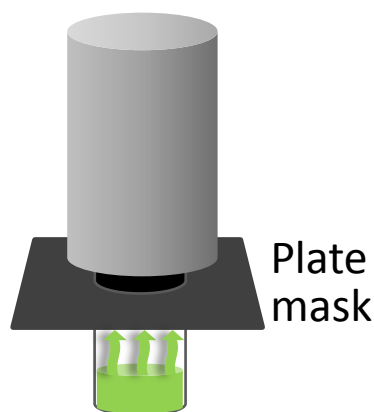
*...plus many, many more*



# Built-In Detectors for each Module

## -behaves like three separate instruments

Head-on PMT for  
photon counting

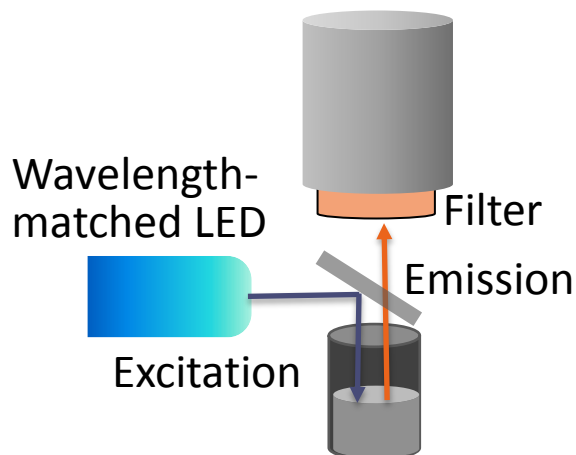


### Luminometer

Plate mask switching for 96/384-well measurements

- $3 \times 10^{-21}$  moles Luciferase sensitivity
- 9 logs dynamic range

PiN-Photodiode

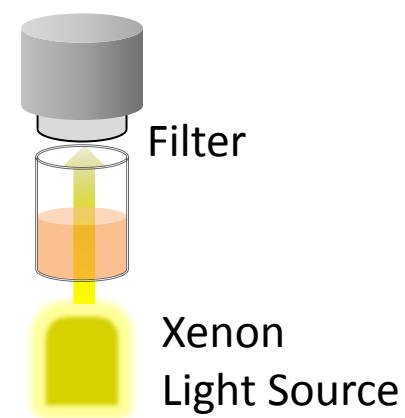


### Fluorometer

Filter slides for automatic filter switching

- 2fmol/200ul fluorescein sensitivity
- 6 logs dynamic range

PMT



### Photometer

Filter Wheel with UV-Vis Abs filters

- 0.01 OD sensitivity
- 0-4 OD range

# Luminescence Filters

## Luminescence Filters (standard)



Position	Filter wavelength	Assay
5	Empty (user configurable)	
4	530nm LP	Click Beetle Luciferase; BRET1
3	540nm SP	Firefly Luciferase; ChromaGlo
2	600nm LP	HaloTag
1	495nm SP	NanoLuc (NanoBRET <sub>618</sub> ), Renilla Luciferase

- All of the filters needed for NanoBRET™<sub>618</sub> are included
- Custom Luminescence Emission filters are available

# Fluorescence Filters

5 Standard Excitation filters and  
2 custom positions.

Empty Emission filter paddles for  
6 customized positions are  
available as accessory

	Excitation Peak Wave length	Emission wave length	Assays
UV	365	415-445	Hoechst dye, 4-MU
Blue	475	500-550	EGFP, or hMGFP, DNA, RNA or protein quantitation dyes, QuantiFluor™, Fluorescein, Rhodamine-110
Green	525	580-640	Rhodamine, Cy®3, resorufin
Red	605	660-720	Cy®5, RNA quantitation dyes
AFC	405	495-505	Aminofluorocoumarin

Custom Excitation and Emission filters  
are available

# Absorbance Filters

9 UV-Vis Absorbance filters are provided with the instrument

The wavelength just needs to be close, not exact. A 450nm filter is actually 445-455nm

**i.e.** Pierce 660 Reagent calls for a 660nm filter, but 600nm works great, that's 60nm off-peak

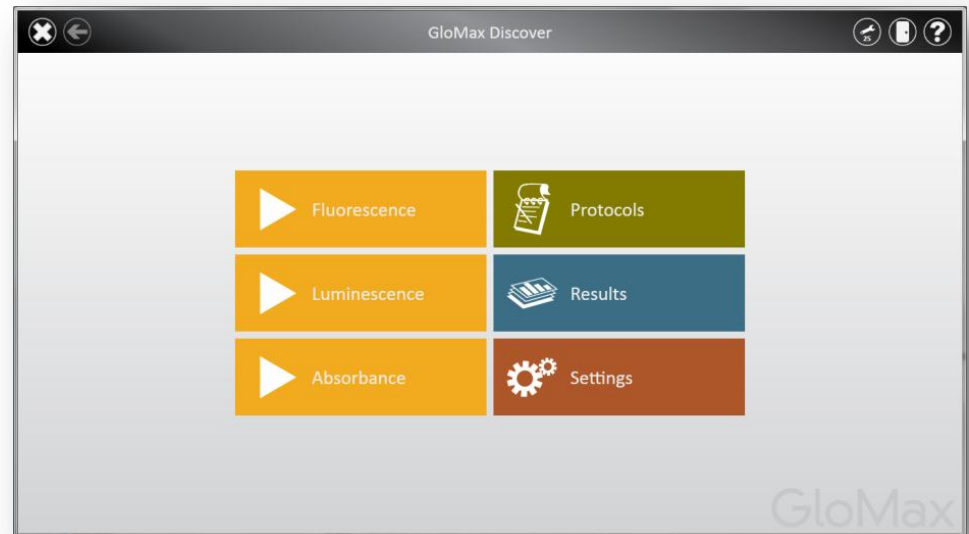
Filters (10nm bandpass)	Assays
230nm	Contaminants (Guanidine, Phenol, carbohydrates)
260nm	Nucleic Acid Quantitation
280nm	Protein Quantitation (Nucleic Acid Purity)
320nm	Background subtraction for Nucleic Acid/Protein Quantitation
405nm	Colorimetric Promega Assays (CaspACE)
450nm	ELISA Assays
490nm	Colorimetric Promega Assays (CellTiter 96/Aqueous/Aqueous ONE / CellTiter Blue)
560nm	BCA Protein Assays
600nm	Bradford Protein Assays Coomassie® Blue Protein Assays



# Intuitive Software Makes it Easy

## Simple GUI Interface

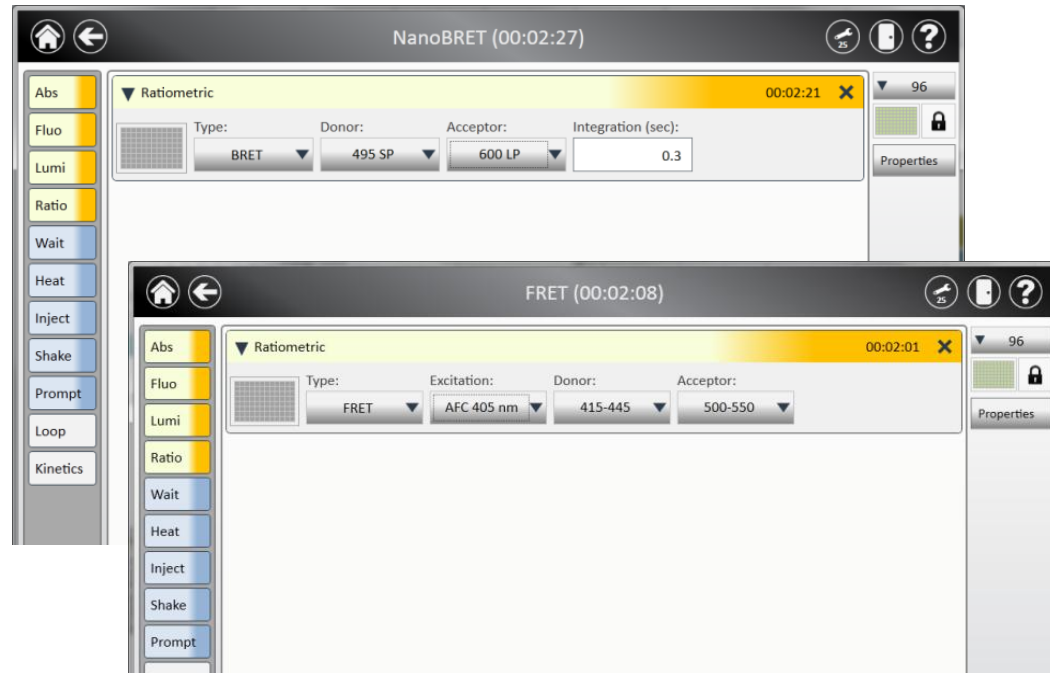
- Quick Start reads
- Select / Create Protocols
- View / Export Results
- User Settings



# Intuitive Software Makes it Easy

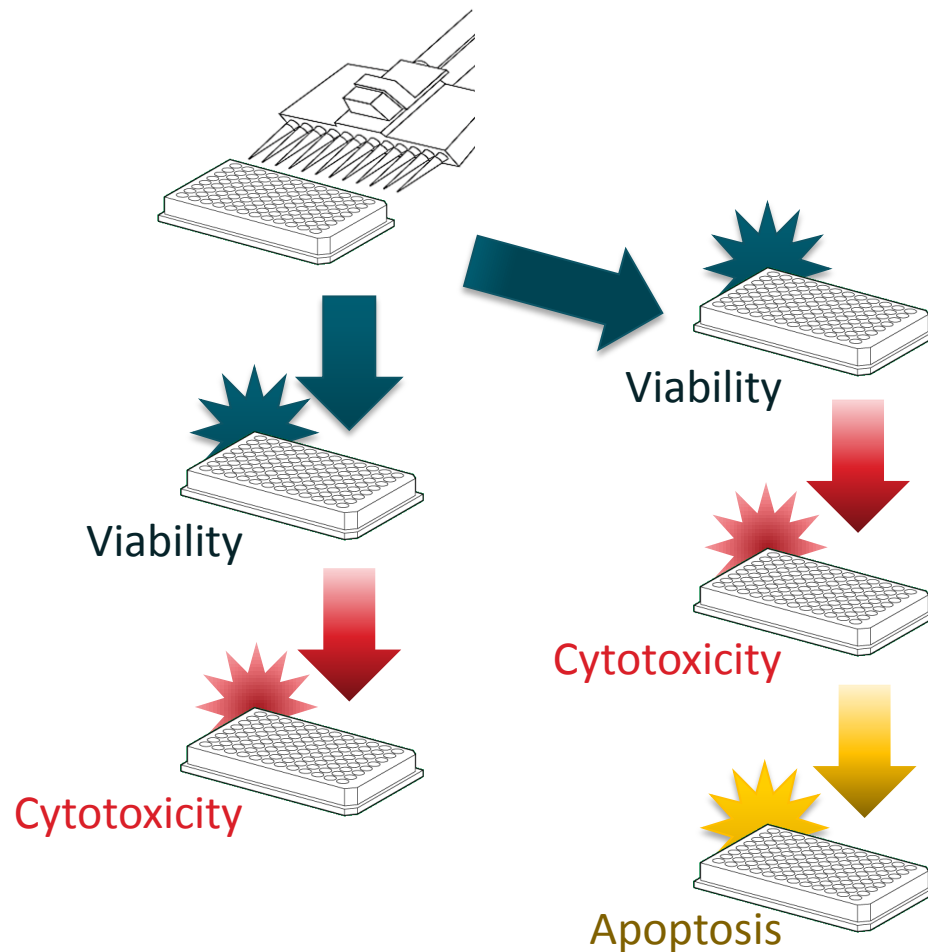
## Flexible protocol-builder

- User customizable
- State-of-the-art touchscreen response
- Drag and drop navigation



# Multiplex Assays for a More Complete Picture

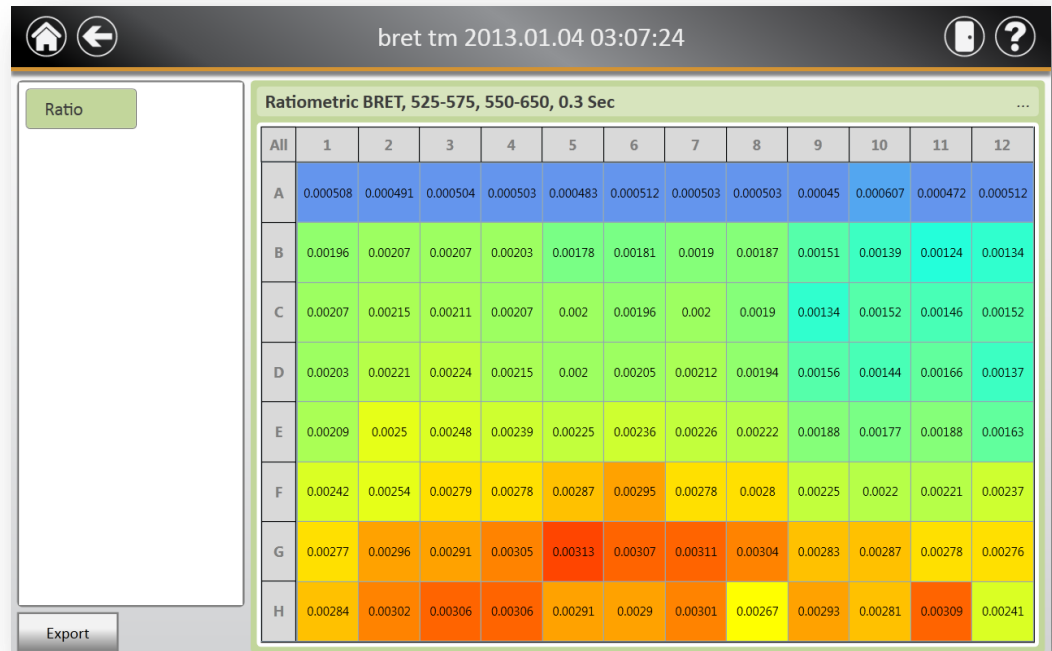
- Automatic filter slides enable multiple reads without user intervention
- Intuitive software provides flexible and easy protocol design
- Easily perform BRET and FRET studies
- Customize filters for your needs



# Intuitive Software Makes it Easy

## Data Portability

- Network-ready
- Heat-map display
- Multi-touch pinch/zoom gestures
- Export to USB, microSD, network, LIMS, Cloud

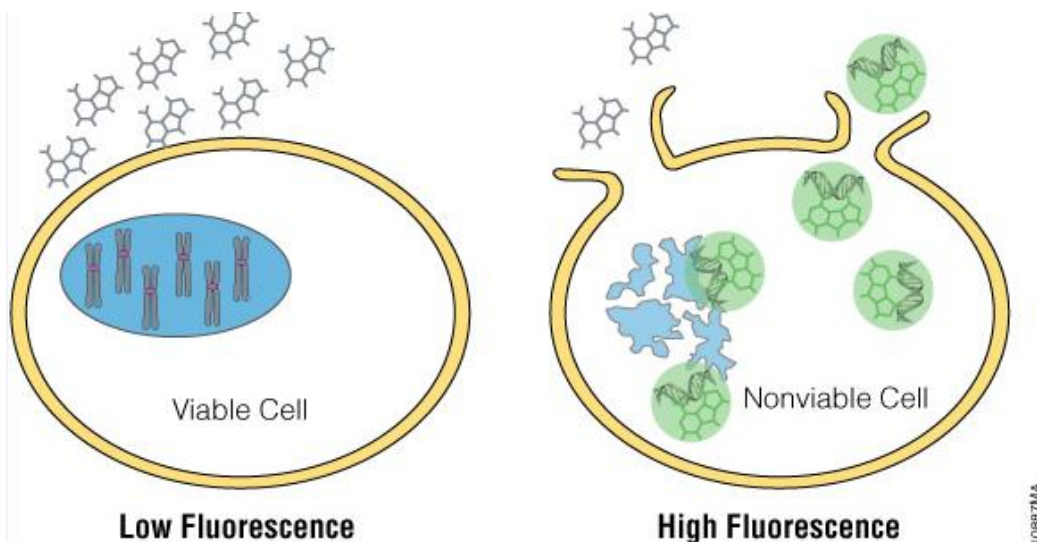




# Performance

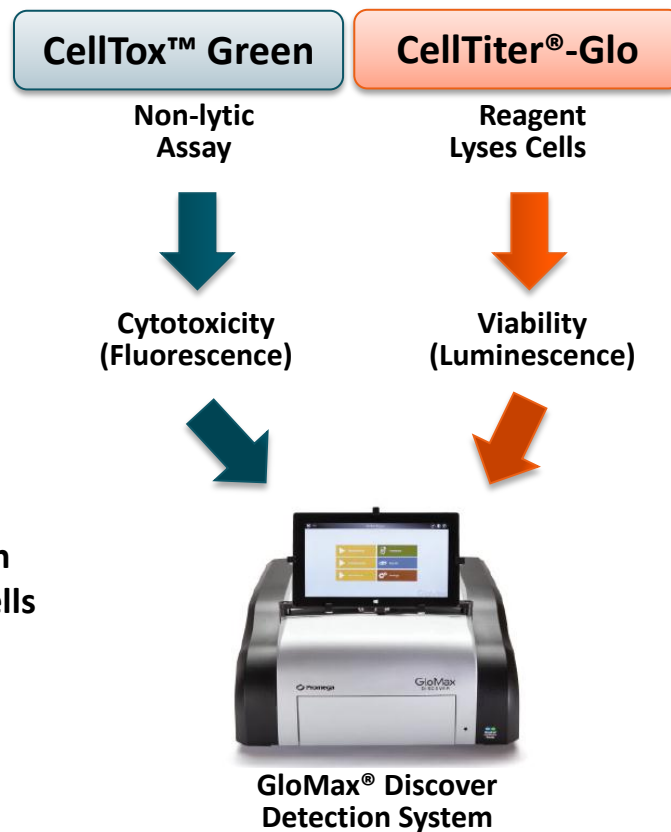
Multiplexing: CellTox™ Green and CellTiter®-Glo  
NanoBRET™ Technology for Protein Interactions

# Multiplexing: CellTox™ Green and CellTiter® Glo



Excluded dye yields **no increase** in fluorescence with viable cells

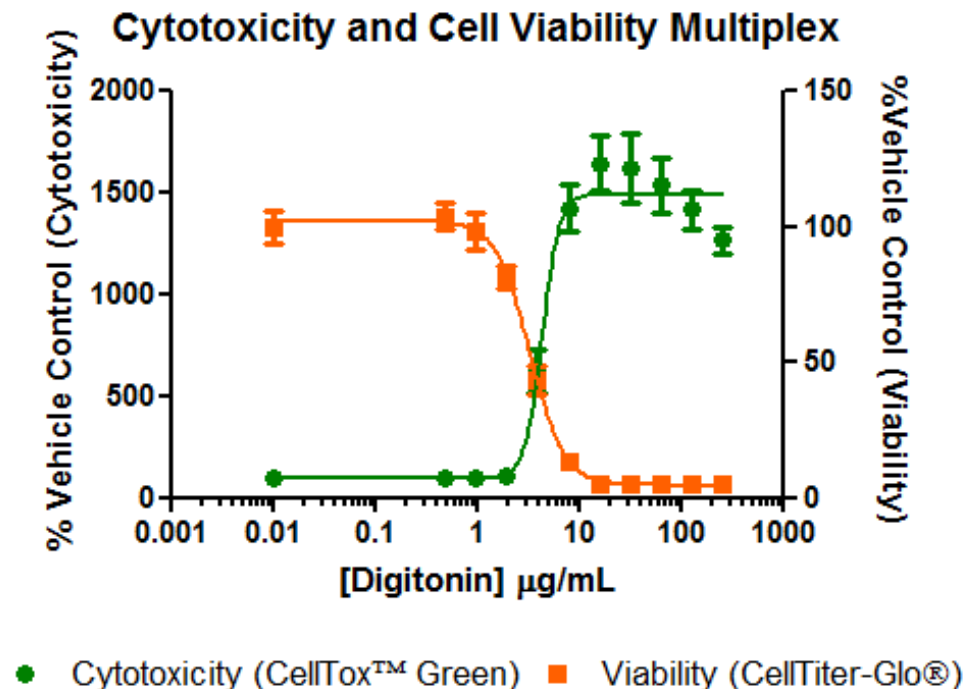
Non-excluded dye yields **increase** in Fluorescence with compromised cells



# Multiplexing: CellTox™ Green and CellTiter® Glo

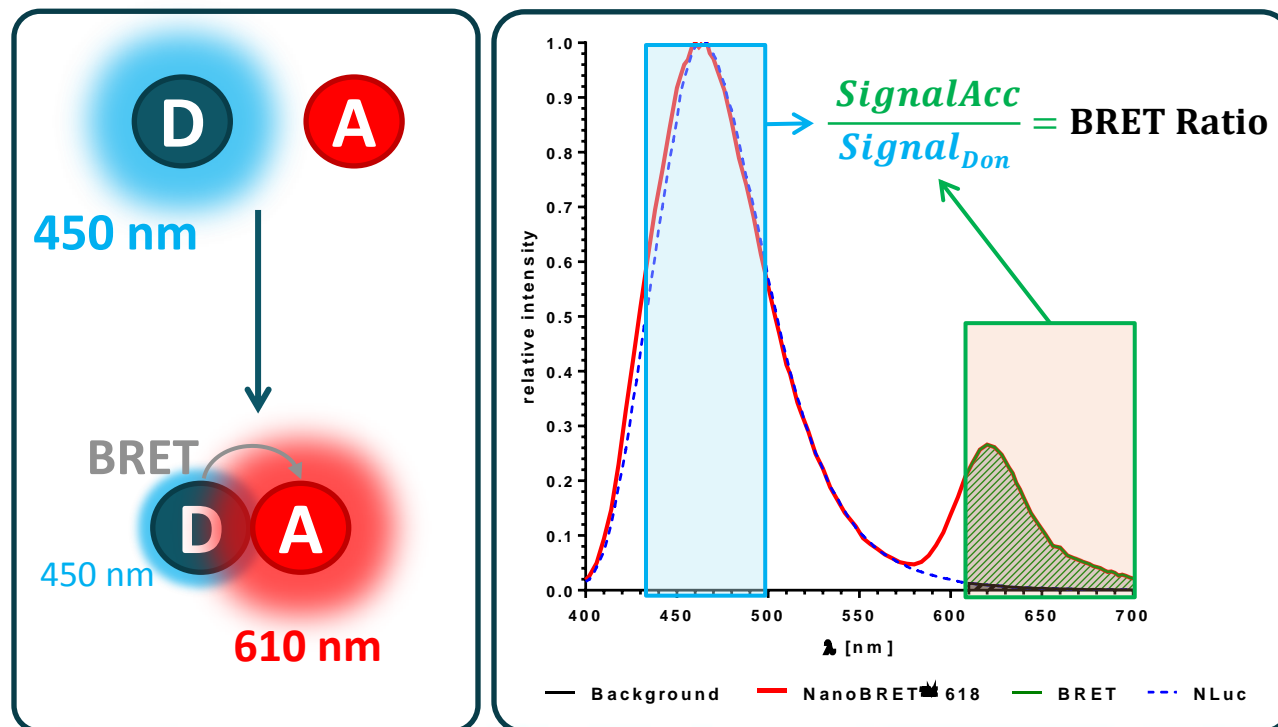
## Expected Results:

- A dose-dependent effect on K562 cell viability. As digitonin concentration increase, so does cytotoxicity
- Luminescent ATP detection decreases due to decreased cell viability
- GloMax® Discover provides easy setup to multiplex Promega assays
- GloMax® Discover results provide expected biology



# NanoBRET™ Technology for Protein Interactions

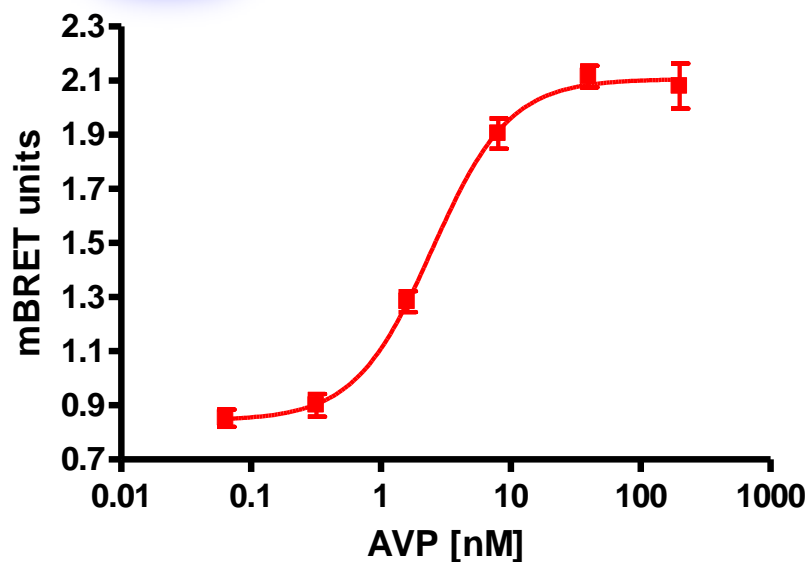
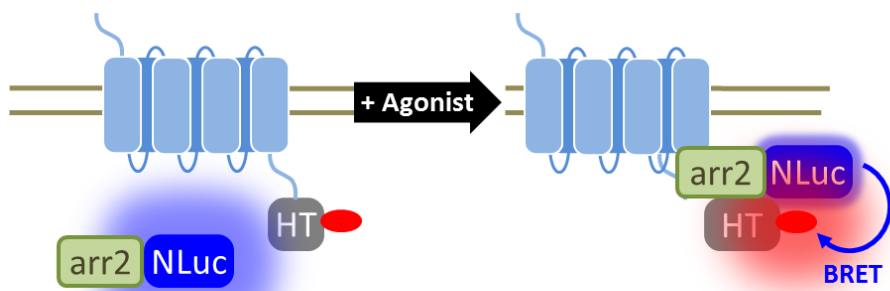
- BRET provides real time measurement *in living cells*
- Superior luminescent Donor signal from NanoLuc™
- Flexible choice of Donor/Acceptor Separation
- Low Donor/Acceptor ratios provides best dynamic range





# NanoBRET™ Application

## AVPR2 / $\beta$ -arr2 Model



**Donor:**

b-arrestin-NanoLuc

**Acceptor:**

Arginin-Vasopressin  
receptor-HaloTag fusion

**AVP** = Arginin Vasopressin  
(Agonist)

In HeLa cells

**NanoBRET™**

Express Donor and  
Acceptor protein fusions

Label Cells with HaloTag  
(Acceptor fusion)

Induce interaction



**GloMax® Discover**  
Detection System

# Automation Integration

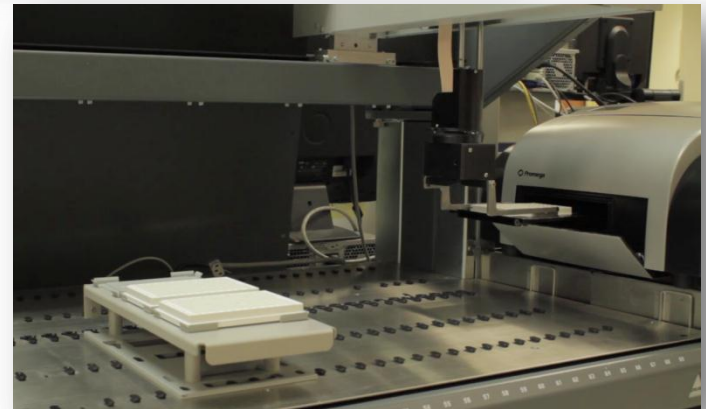
## Demonstrated Integration with:

- **Tecan Freedom EVO®** liquid handler
- **Hudson Robotics Solo™** liquid handler and PlateCraneEX™ robotic arm
- Additional platforms planned



**3<sup>rd</sup> party software control** of Discover

**Integrator's Kit** (PDF command-line instructions)



# GloMax<sup>®</sup> Discover

<b>Integrated</b>	Seamless workflow with Promega Cell and Reporter assays.
<b>Performance</b>	Broader dynamic range, better sensitivity, and lower well-to-well cross-talk for more usable data from your experiment.
<b>Easy-to-use</b>	Simple Tablet PC touchscreen navigation with full PC capabilities and a state-of-the-art Graphical User Interface makes the workflow simple, smooth, yet flexibly.
<b>Connected to your Workflow</b>	<p>Stand-alone instrument or integrate with automation.</p> <p>Export data to your laboratory network, LIMS, or Cloud.</p> <p>Enable 21CFR Part 11 electronic signature compliance.</p>

# Promega Supports Your Entire Workflow for Reporter and Cell-Based Assays

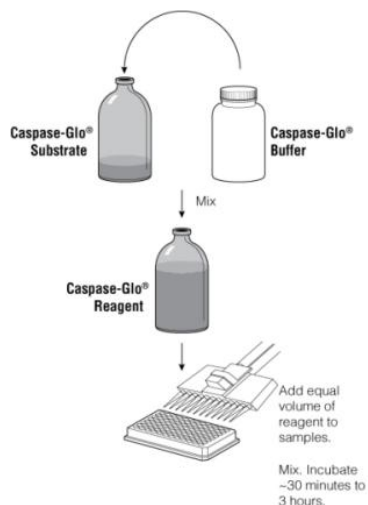
Investigate

Quantitate

Analyze

## Add – Mix – Read :

Cell Viability, Apoptosis,  
Gene Reporter assays,  
Drug Discovery assays



**GloMax<sup>®</sup> Discover**

**Request a free demo at  
[www.promega.com/discover](http://www.promega.com/discover)**

## One Call Supports it All



Scientists to help you:

- Interpret results
- Troubleshoot issues
- Design experiments

# Hands-On: Find the dirtiest place

## Promega's BacTiter-Glo™ Microbial Cell Viability Assay (reagent) and GloMax® Discover (Instrument)

- 1) Swab with a cotton swab a surface that you think is dirty
- 2) Add 400ul water (sterile) in Eppendorf Tube (2ml)
- 3) Put cotton swab into the Ependorftube that contains 400ul water
- 4) Remove the cotton swab
- 5) Take 100ul of the water and put into a well of a white 96-well plate
- 6) Add 100ul BacTiter-Glo™
- 7) Incubate 5 mins
- 8) Measure with Glomax® Discover

**Request a free demo at  
[www.promega.com/discover](http://www.promega.com/discover)**



# Using GloMax® Discover

1) Open door



2) Add plate  
(note: A1)



3) Select  
protocol



4) Select wells



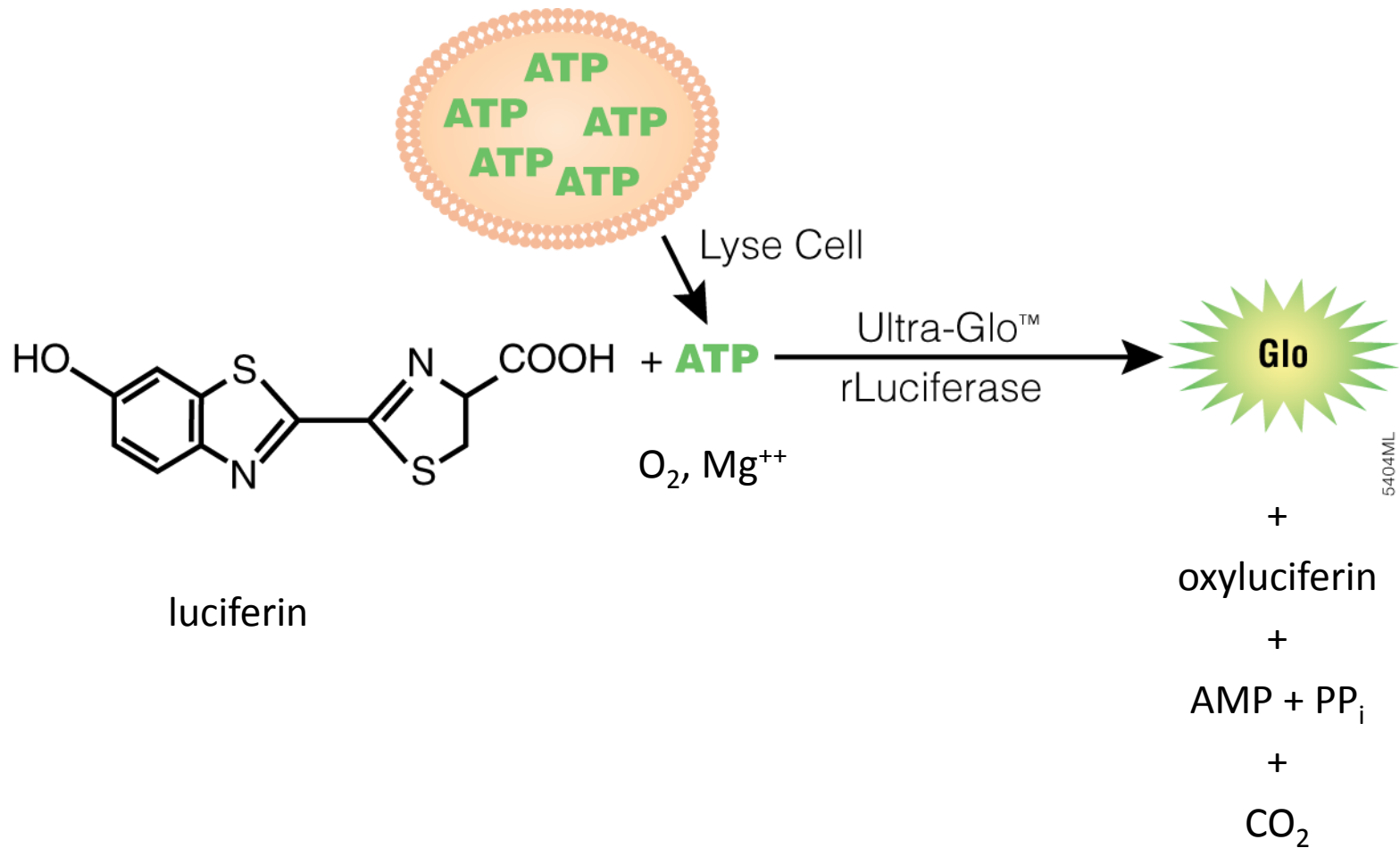
5) Start read



6) Export  
results



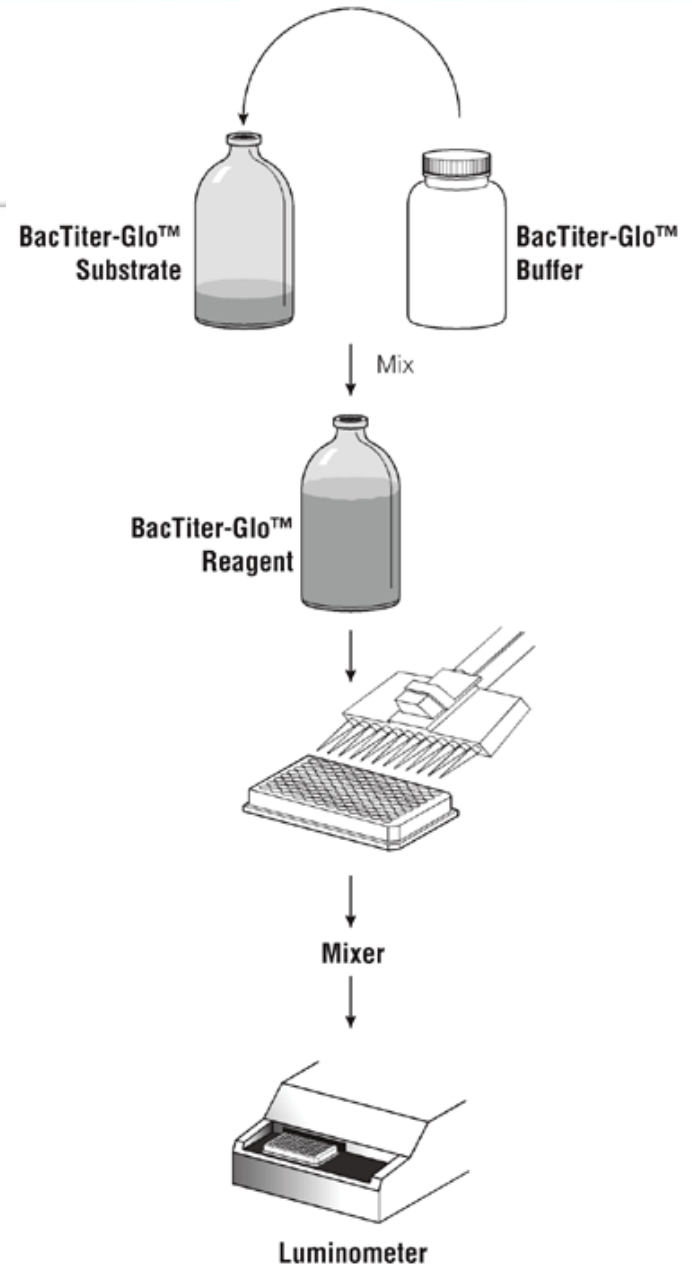
# Bactiter-Glo™ Experiment





# Find the dirtiest spot.....

- Use a cotton swab to sample  
Hands, pen, toilet, money, cell phone
- Return the cotton swab with the form
- Specify your sample on the form
- Incubate the cotton swab in water
- Transfer sup to white 96-wells plate
- Add 1:1 Bactiter-Glo™ reagent
- Incubate 5 min
- Read luminescence in the Discover
- Win an orange labcoat





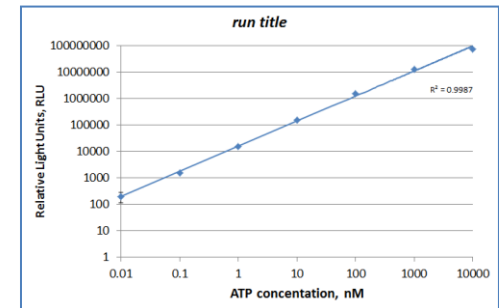
# Hands-On: Smart Protocols

1. Select Smart Protocol from menu



2. Discover loads the Kit protocol & guides FAS / customer through assay setup

3. Discover reads the plate & automatically analyzes data



- 1) GloMax<sup>®</sup> Luminescence Functional Test Kit: 7 point serial dilution of ATP and Promega's CellTiter<sup>®</sup>-Glo assay.
- 2) GloMax<sup>®</sup> Fluorescence Functional Test Kit: 7 point serial dilution of DNA and the QuantiFluor<sup>™</sup> dsDNA System.
- 3) GloMax<sup>®</sup> Absorbance Functional Test Kit: 7 point BSA protein dilution and Pierce 660nm Protein assay