

GloMax® Discover – Ideal Partnership of Assays and Reagents

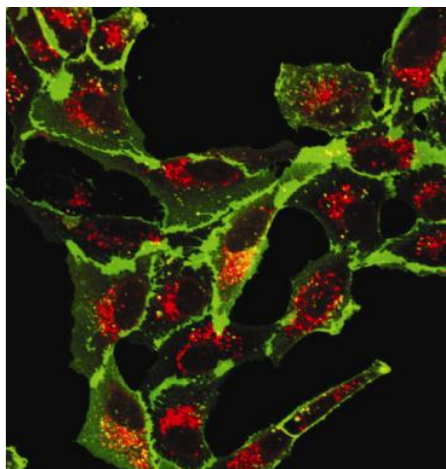
Craig Malcolm: Product Manager Cell Analysis & Proteomics

Cell-Based Assay Tour - March 2014

Outline

- Overview of GloMax[®] Discover
- Assay Performance
- Automation Integration

Promega capabilities



Cellular & Biochemical Technologies

- Assay Design
- Integrated Cellular Biology
- Macromolecular Design
- Protein Analysis
- Organic Chemistry



Nucleic Acid Technologies

- Purification
- Amplification
- Detection



Instrument & Reagent Technologies

- Instrumentation
- Reagents
- Software
- Services

Introducing ...

- Integrated with Promega assays
- Multi-mode detection (Luminescence, Fluorescence, UV-Vis Absorbance, BRET, FRET, filtered luminescence)
- Easy to use: Plug-n-play setup
- Superior sensitivity, dynamic range, and cross-talk performance
- Integrated filter paddles for assay multiplexing
- 6 to 384-well plate formats
- Automation-friendly
- Tablet PC touchscreen control and wireless connectivity to network and Promega.com
- Electronic signature control for 21CFR Part 11
- IQ / OQ Service



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™

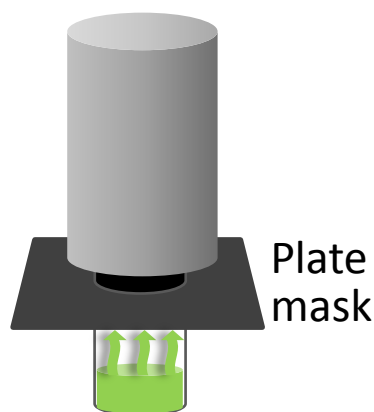
BRET and FRET Assays: Including:

- 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

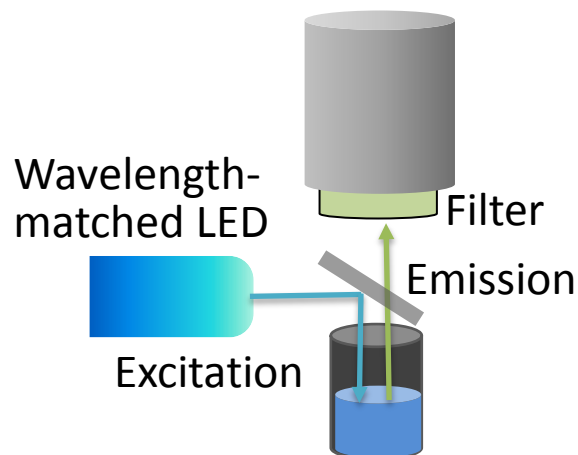


Luminometer

Plate mask switching for 96/384-well measurements

- 3×10^{-21} moles Luciferase sensitivity
- 9 logs dynamic range

PiN-Photodiode

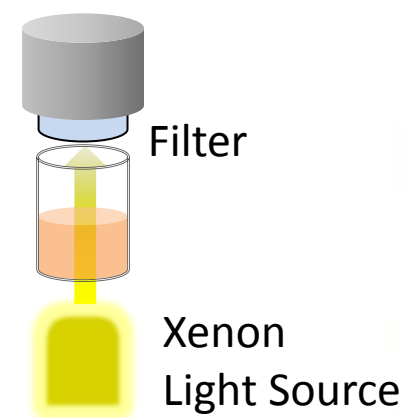


Fluorometer

Filter slides for automatic filter switching

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

PMT



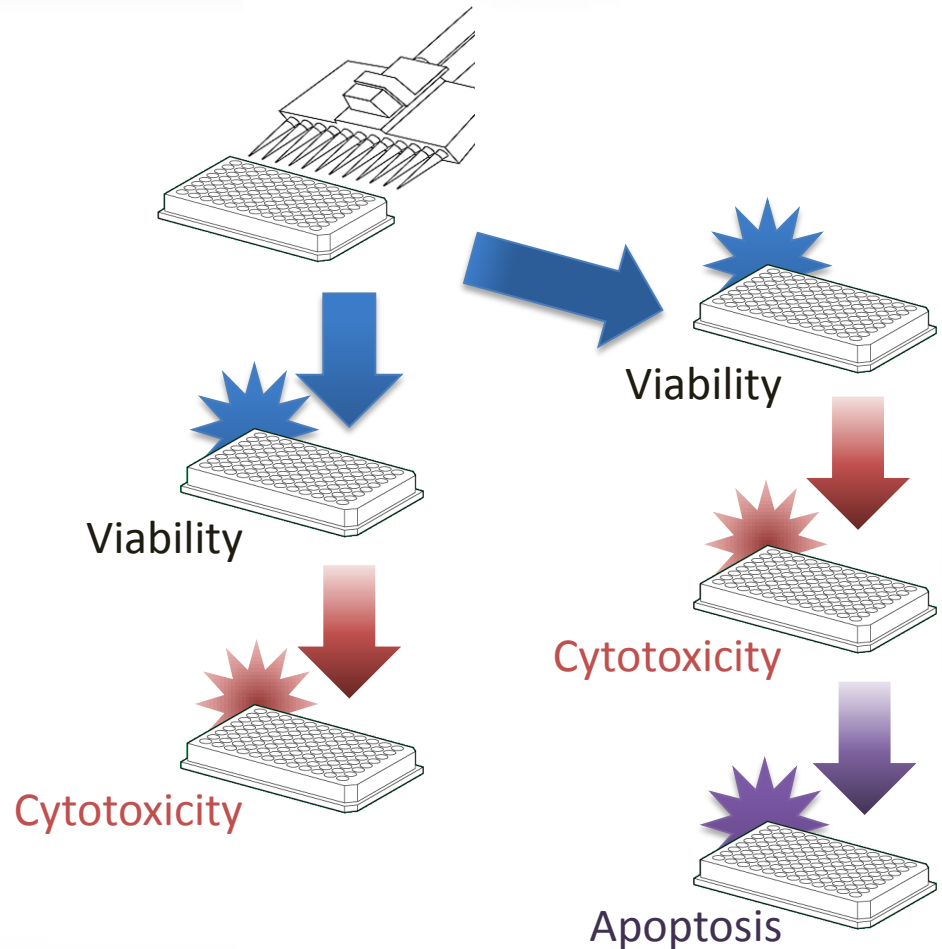
Photometer

Filter Wheel with UV-Vis Abs filters

- 0.01 OD sensitivity
- 0-4 OD range

Multiplex Assays for a More Complete Picture

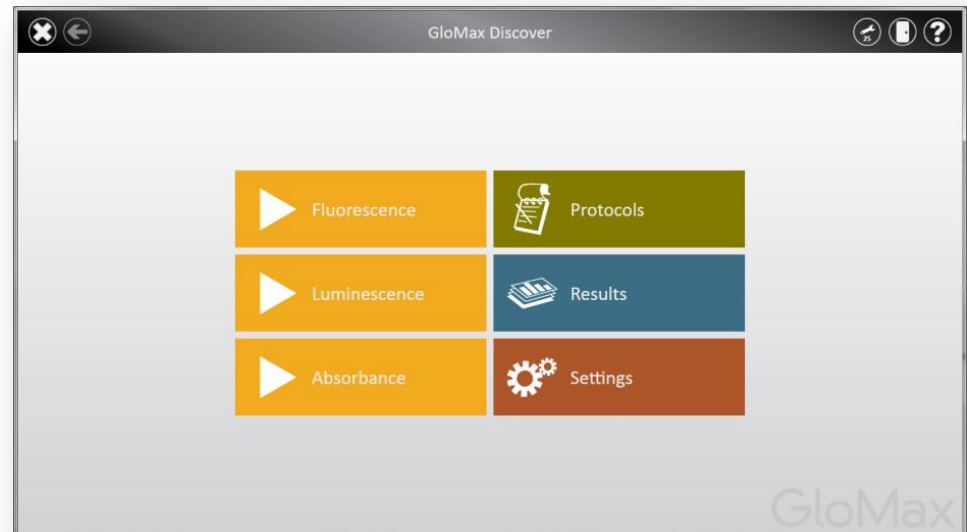
- Automatic filter slides enable multiple reads without user intervention
- Intuitive software provides flexible and easy protocol design
- Easily perform BRET and FRET studies
- Customize filters for your needs



Intuitive Software Makes it Easy

Simple GUI Interface

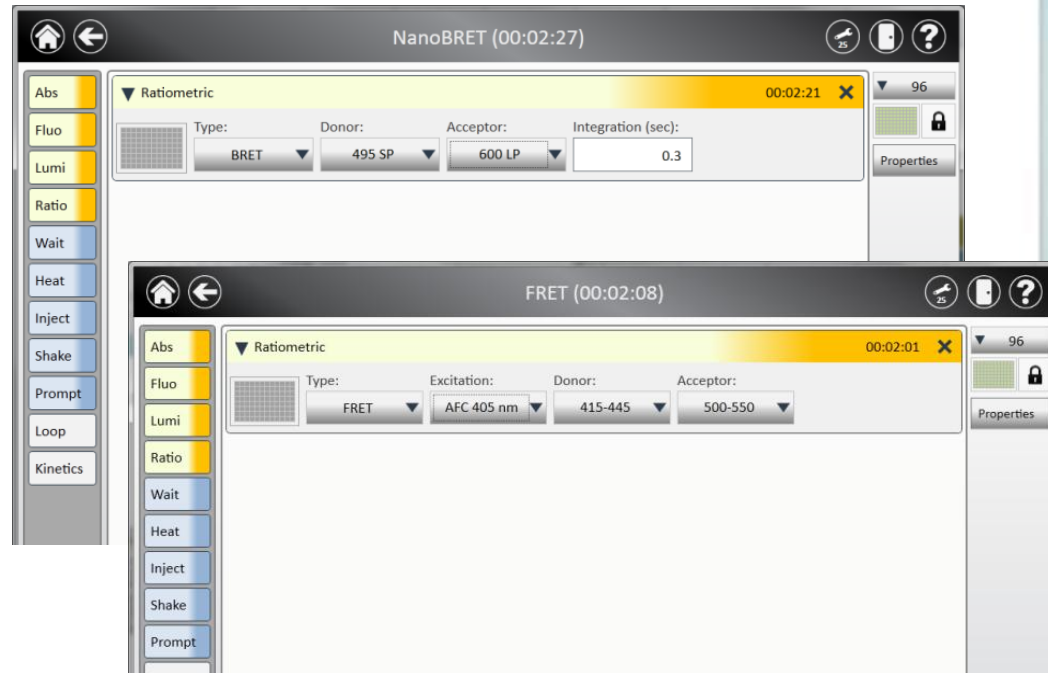
- Quick Start reads
- Select / Create Protocols
- View / Export Results
- User Settings



Intuitive Software Makes it Easy

Flexible protocol-builder

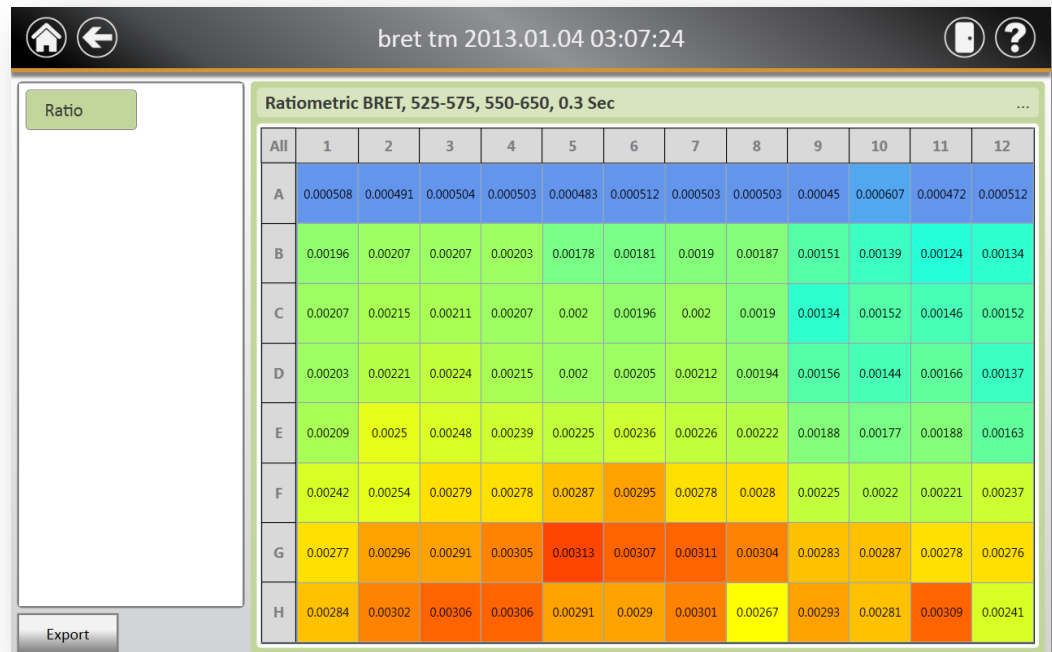
- User customizable
- State-of-the-art touchscreen response
- Drag and drop navigation



Intuitive Software Makes it Easy

Data Portability

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



An Extensive List of Applications on Promega.com

Application Notes: GloMax® Discover System



GloMax® Discover System

Bioassays

Measuring the ADCC Reporter Bioassay Complete Kit (WIL2-S) Signal on the GloMax® Discover System

Cell Health and Metabolism

Measuring the Output of the CytoTox-Fluor™ Cytotoxicity Assay on the GloMax® Discover System

Measuring the ONE-Glo + Tox Luciferase Reporter and Cell Viability Assay on the GloMax® Discover System

Measuring Cell Viability Using the CellTiter-Glo® Cell Viability Assay and GloMax® Discover System

Measuring P450-Glo™ Assays on the GloMax® Discover System

Measuring Bacterial Cell Viability Using the BacTiter-Glo™ Assay and GloMax® Discover System

Measuring Fluorescence Using the Apo-ONE® Homogeneous Caspase-3/7 Assay with the GloMax® Discover System

Measuring Fluorescence Using the CellTiter-Blue® Cell Viability Assay with the GloMax® Discover System

Measuring Fluorescence Using the ApoTox-Glo™ Toxin Assay With the GloMax® Discover System

www.promega.com/discover

Using GloMax[®] Discover is Simple

1) Open door



2) Add plate
(note: A1)



3) Select
protocol



4) Select wells



5) Start read



6) Export
results



What's Inside

To access the interior of the instrument:

- 1) Hold the door open
- 2) Use the Phillips screwdriver to unscrew the 2 screws (one at the left side, and one at the right side).

Note: Only a $\frac{1}{4}$ turn is necessary

- 3) Pull the front access panel off

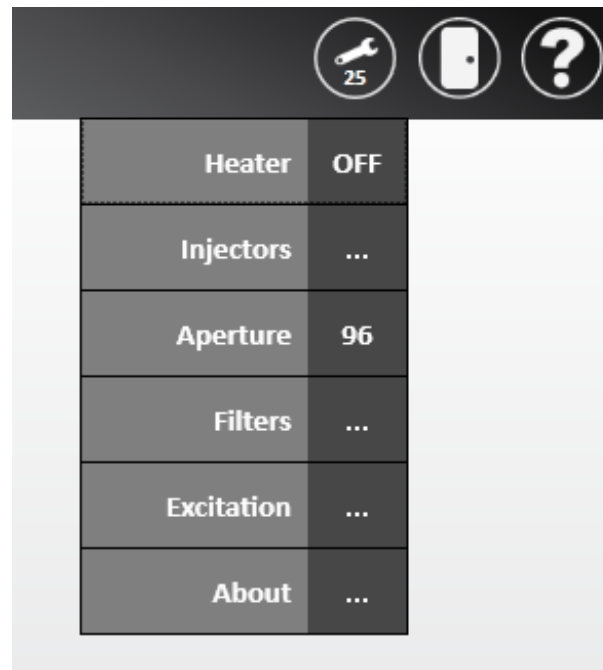


What's Inside

To change excitation filters, emission filters, or apertures:

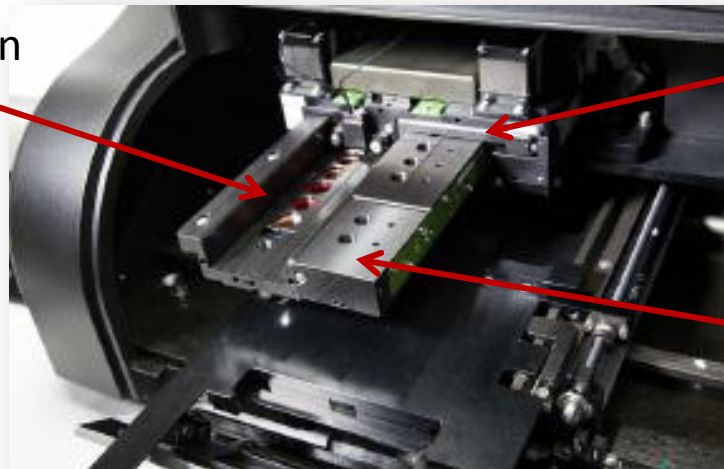
Touch the Tools icon
at the top of the
window and then
the desired function

Follow the on-screen
instructions



What's Inside

Luminescence Emission
filters



Fluorescence
Emission filters

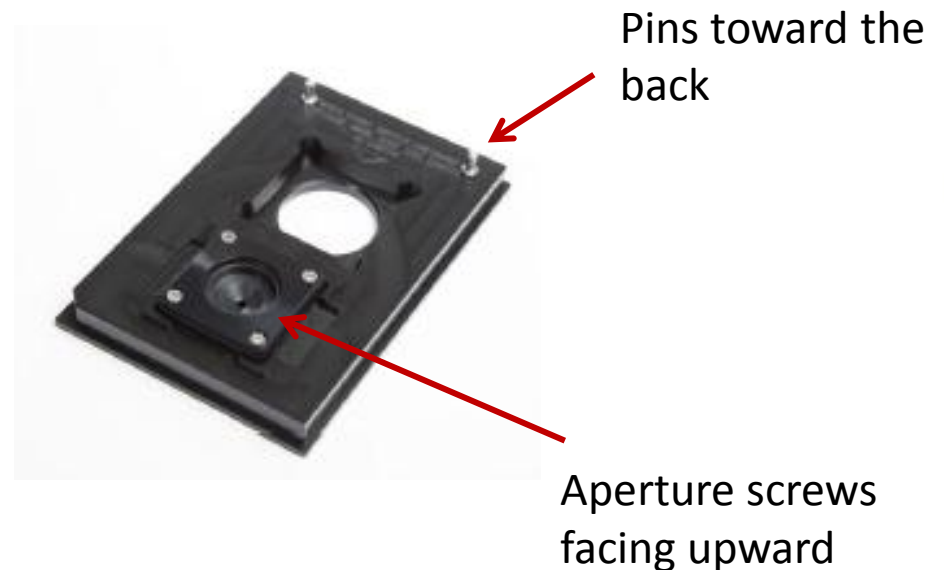
Fluorescence Excitation
modules



What's Inside

Semi-automated method for changing the internal aperture.

1. Follow the on-screen guide to change the aperture
2. Make sure to load the Aperture Plate correctly



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

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 30nm 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

Luminescence Filters

Luminescence Filters (standard)



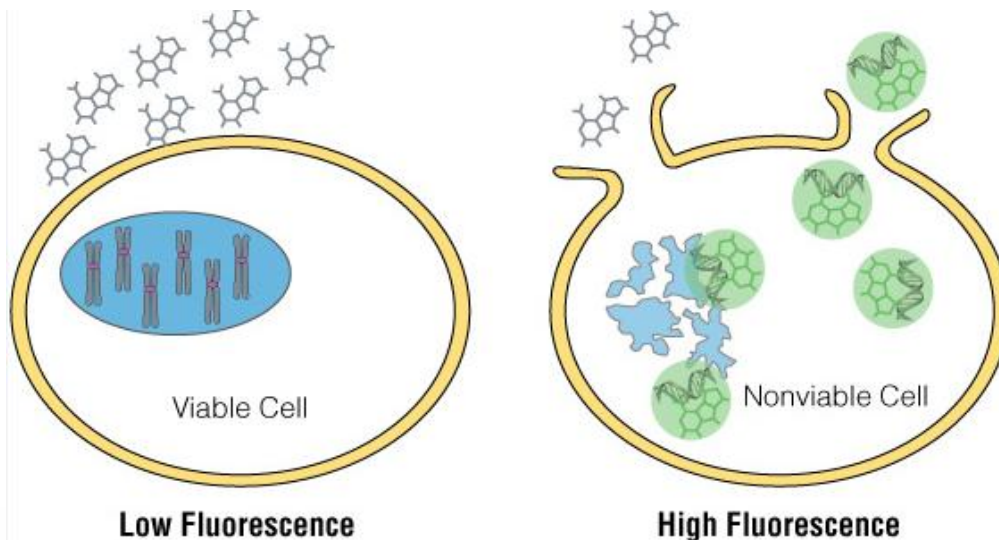
Position	Filter wavelength	Assay
5	Empty (user configurable)	
4	530nm LP	Click Beetle Luciferase; BRET1
3	540nm SP	Firefly Luciferase; ChromaGlo
2	600nm LP	HaloTag
1	495nm SP	NanoLuc (NanoBRET), Renilla Luciferase

- All of the filters needed for NanoBRET™ are included
- No additional custom filters are needed for NanoBRET™

Assay performance

CellTox™ Green for Real-Time Cytotoxicity

- Measures changes in membrane integrity as a result of cell death
- Dye preferentially stains dead cell DNA
- Seamless multiplex with viability assay to provide mechanistic information related to cytotoxicity



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

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

CellTox™ Green

Non-lytic Assay

Cytotoxicity (Fluorescence)

CellTiter®-Glo

Reagent Lyses Cells

Viability (Luminescence)

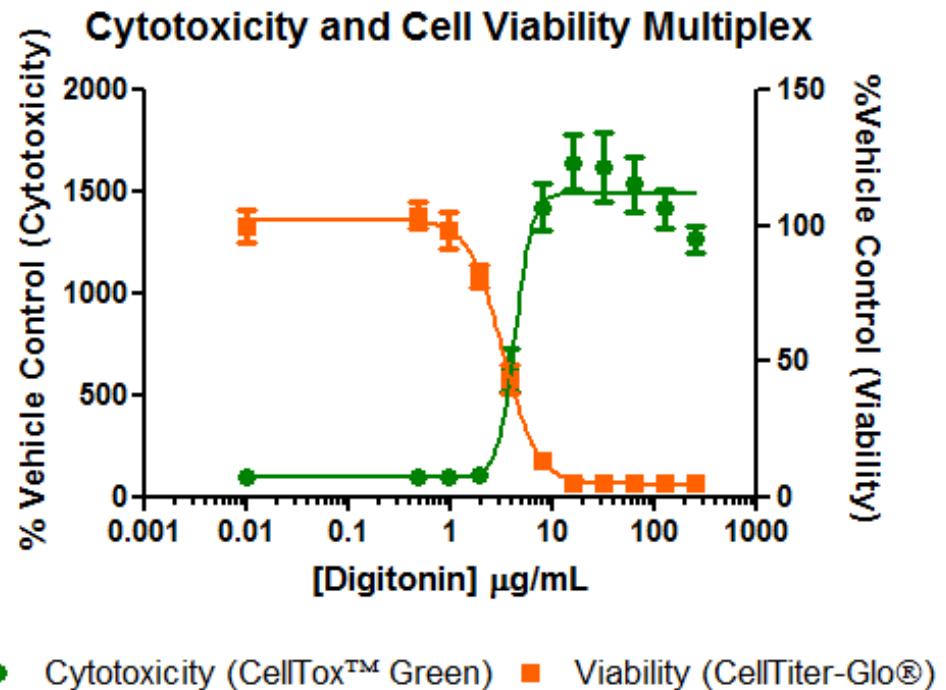


GloMax® Discover Detection System

CellTox™ Green for Real-Time Cytotoxicity

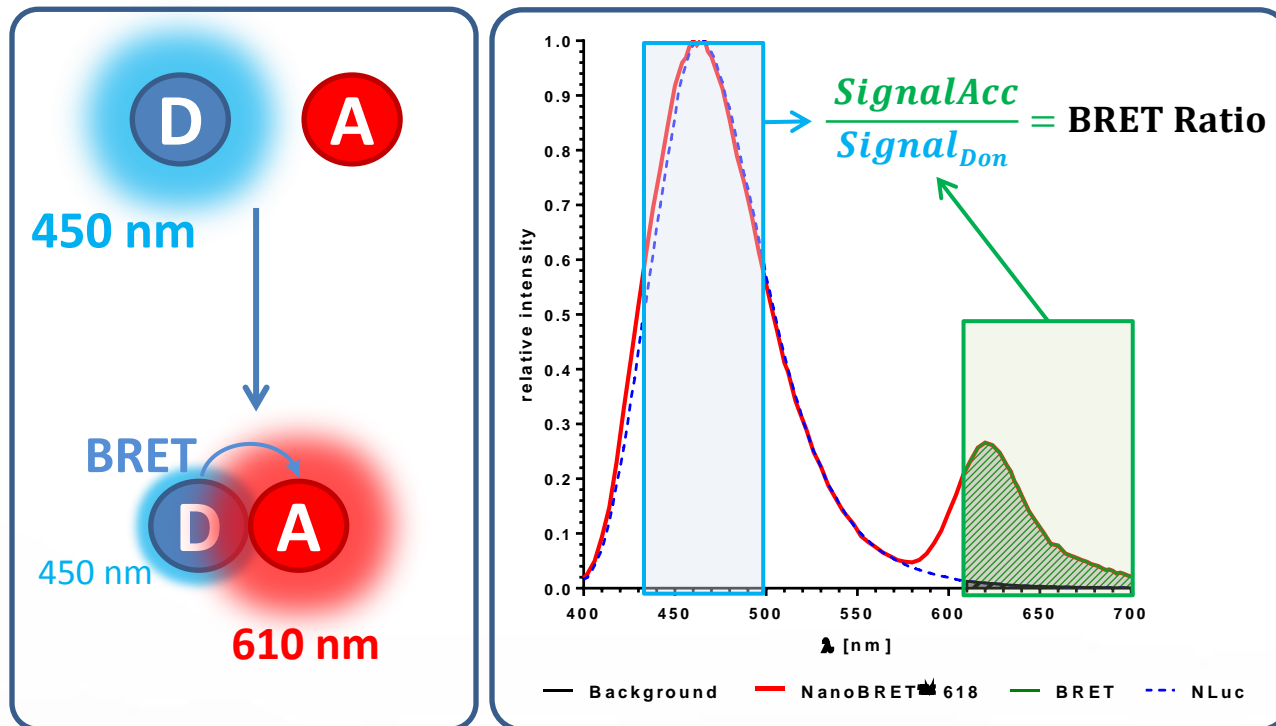
Expected Results:

- 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™

Express Donor and
Acceptor protein fusions

Label Cells with HaloTag
(Acceptor fusion)

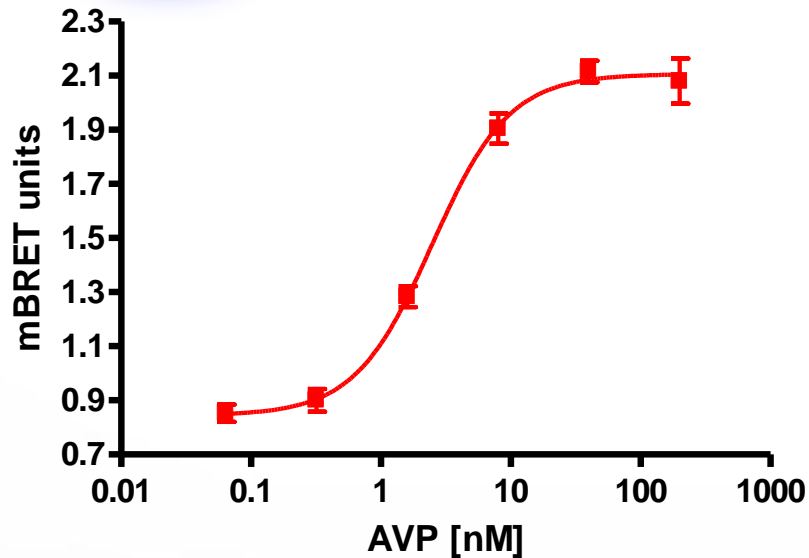
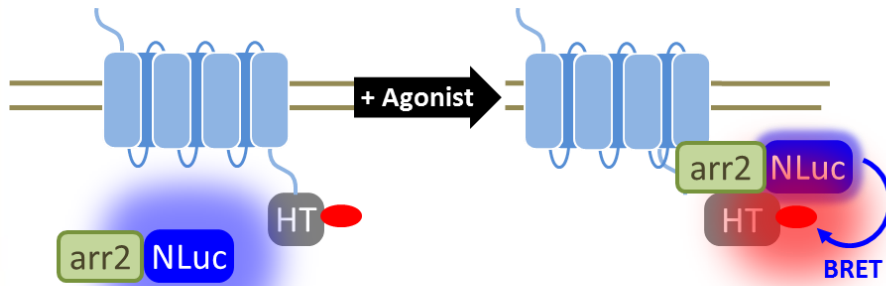
Induce interaction



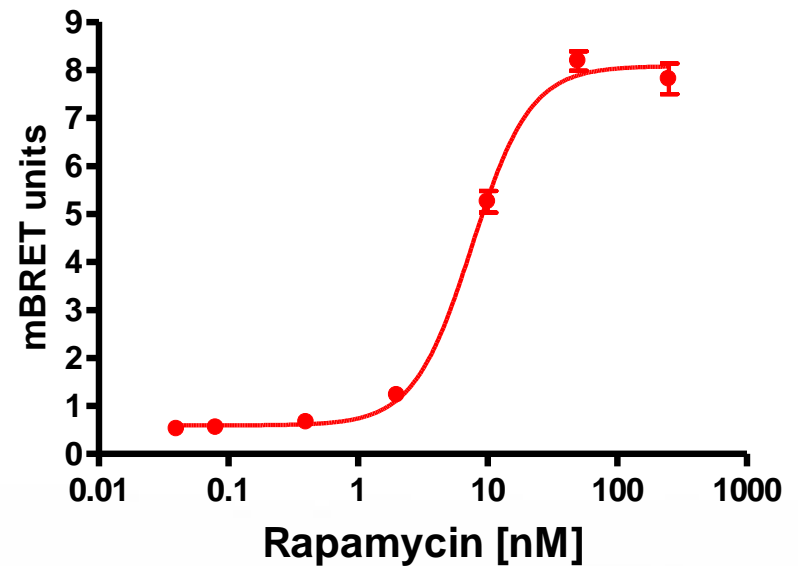
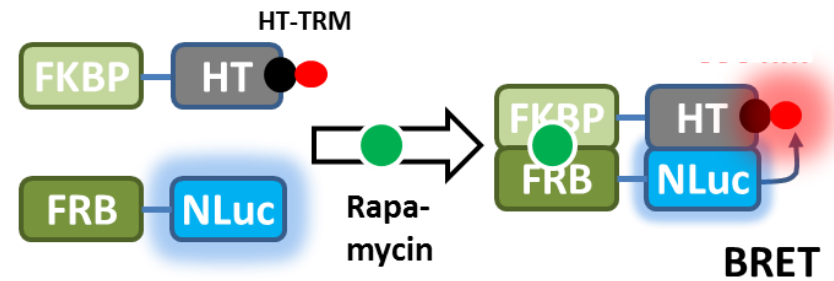
GloMax® Discover
Detection System

NanoBRET™ Application Examples

AVPR2 / β -arr2 Model



FKBP/Frb/Rapamycin Model



Other BRET and Filtered Luminescence Assays

Renilla / YFP

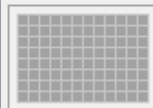
Donor Filter

495nm SP

**Acceptor
Filter**

530nm LP

▼ Ratiometric



Type:

BRET ▼

Donor:

495 SP ▼

Acceptor:

530 LP ▼

Integration (sec):

0.5

Chroma-Glo

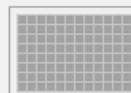
1st reading

600nm LP

2nd reading

540nm SP

▼ Luminescence



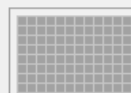
Filter:

600 LP ▼

Integration (sec):

0.3

▼ Luminescence



Filter:

540 SP ▼

Integration (sec):

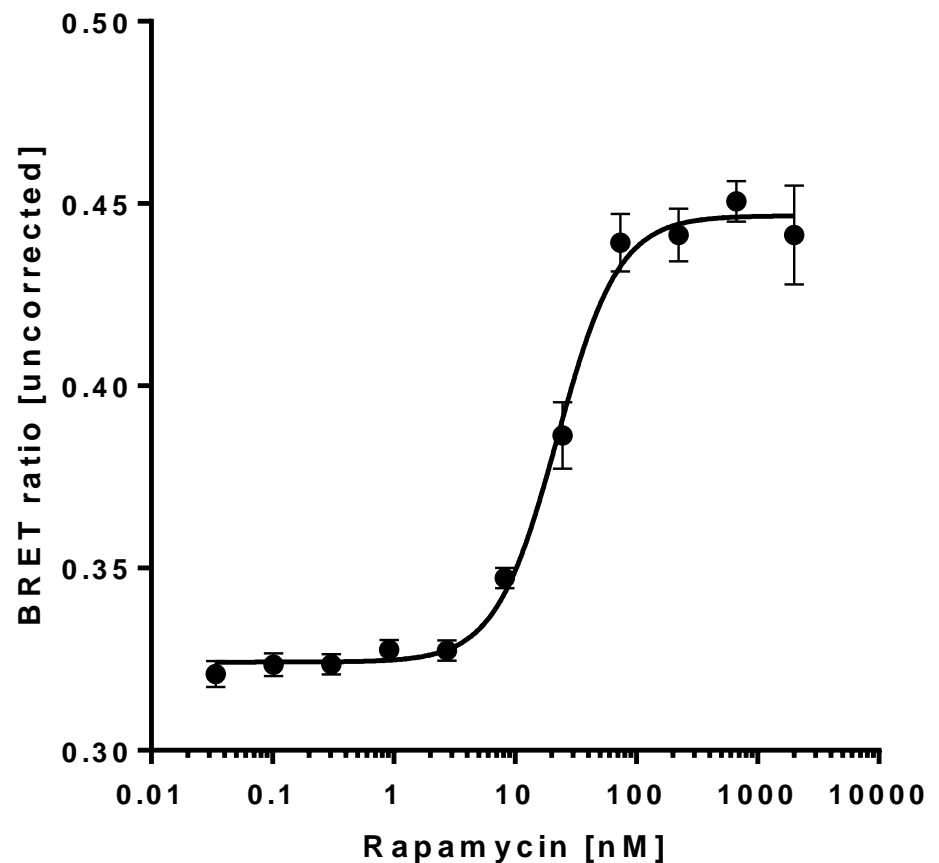
0.3

BRET1: Frb-RLuc8 / FKBP-YFP

Experiment

- HEK293 cells reverse transfected with expression constructs for Frb-RLuc8 and FKBP-YFP (ratio 1:4) using Fugene HD
- Cells were plated in white 96-well plates in DMEM.
- After 24 h medium was replaced with Optimem.
- Cells were treated with a serial dilution of Rapamycin for 15 minutes at 37C before Coelenterazine h was added to a final concentration of 20 μ M.
- Plate was read in GloMax Discover using the following settings
 - Donor filter 495 shortpass filter
 - Acceptor filter 530 nm longpass filter
 - 0.5 seconds integration time

HEK293 Frb-RLuc8 / FKBP-YFP
GloMax Discover

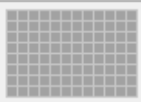


FRET Assays

GFP / RFP FRET

Fluorescein / Rhodamine FRET

▼ Ratiometric



Type:

Excitation:

Donor:

Acceptor:

FRET ▼

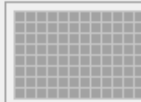
Blue 475 nm ▼

500-550 ▼

580-640 ▼

Coumerin / Fluorescein FRET

▼ Ratiometric



Type:

Excitation:

Donor:

Acceptor:

FRET ▼

AFC 405 nm ▼

415-445 ▼

500-550 ▼

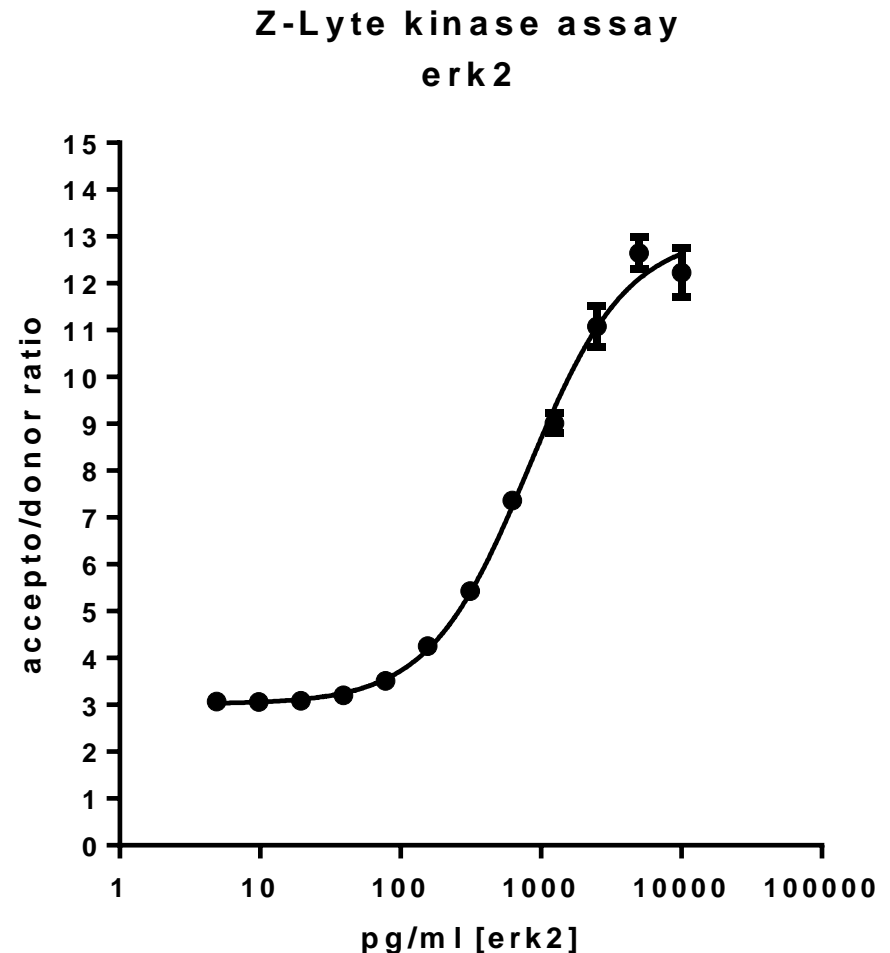
Primary Excitation	Primary Emission	Secondary Excitation	Secondary Emission	Application
475nm	509nm	563nm	582nm	GFP / RFP
447nm	520nm	550nm	573nm	Fluorescein / Rhodamine
409nm	447nm	447nm	520nm	Coumerin / Fluorescein; Life Technology Z-Lyte Assay

FRET: Z-Lyte kinase assay for erk2 activity

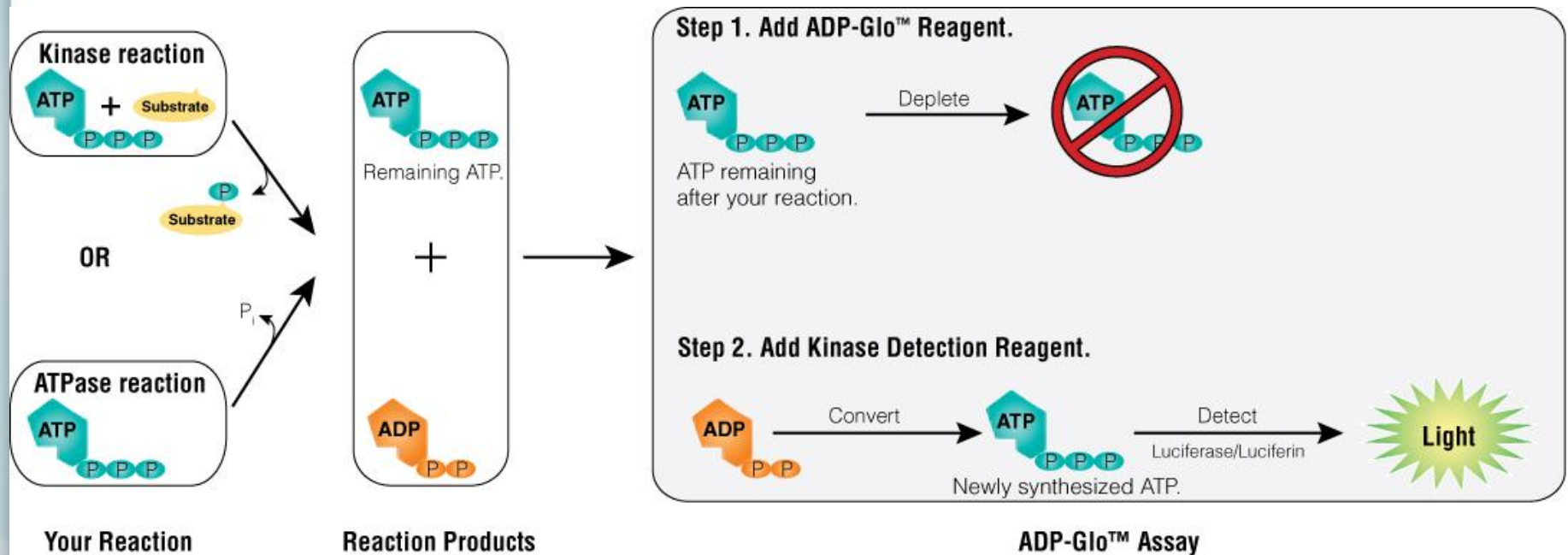
Experiment

- Erk2 kinase activity was determined using the Z-Lyte Kinase assay (Life Technologies) according to manufacturers instructions
- ATP = 0.5 mM
- Activity measurement was taken in a GloMax Discover using the following filter settings

Excitation:	AFC 405 nm
Donor:	415-445
Acceptor:	500-550



ADP-Glo™ Assay and Kinase Profiling Systems



- Measures kinase activity in any purified kinase enzyme
- Direct correlation of kinase activity to luminescence output
- Sensitive and suitable for studying kinases in physiological conditions

ADP-Glo™ Assay and Kinase Profiling Systems

Kinase profiling System – ST-5 Kit components



2 strips



ERK2 GSK3b JNK1 JNK3 p38a p38b p38d p38g

2.5μL Kinase Strip, 50X
(some strips may contain 5μl)



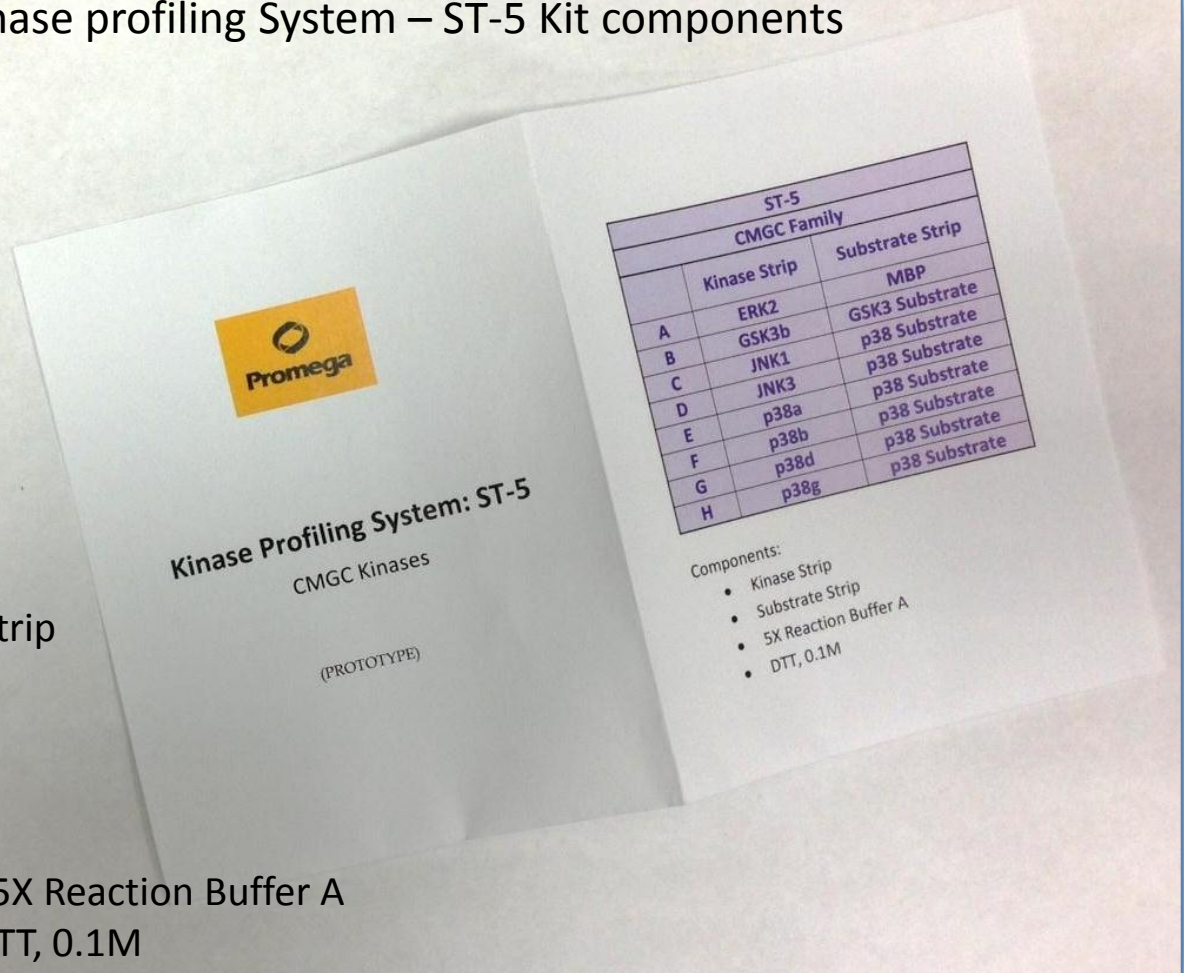
2 strips



37.5μL Substrate/Co-Factor Strip



1.5mL 5X Reaction Buffer A
25μL DTT, 0.1M



Kinase Selectivity Profiling

384-well plate

5 μ l kinase reaction

+

5 μ l ADP-Glo™ Reagent

40 min. Incubation

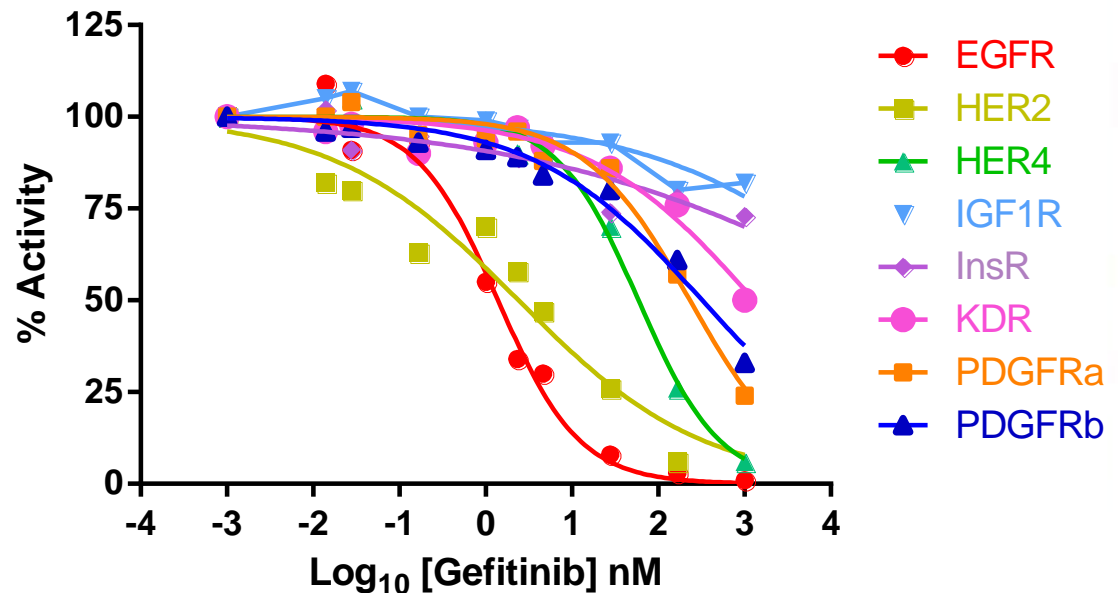
+

10 μ l Kinase Detection Reagent

30-60 min. Incubation

Record Luminescence

Receptor Tyrosine Kinases



Automation Integration

Demonstrated Integration with:

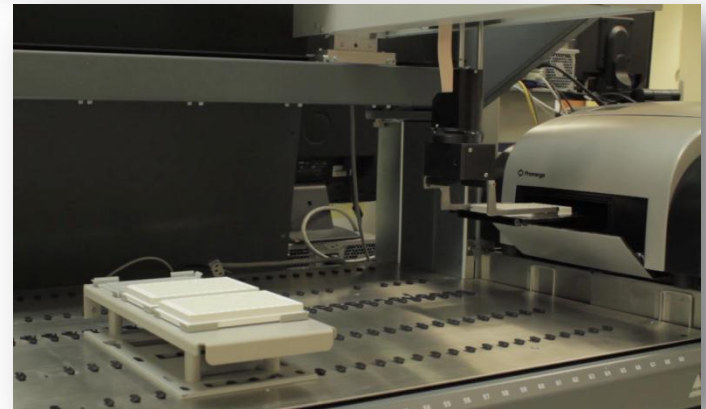
- **Tecan Freedom EVO®** liquid handler
- **Hudson Robotics Solo™** liquid handler and PlateCraneEX™ robotic arm
- Additional platforms planned



3rd party software control of Discover

Data CSV format for **LIMS data** integration

Integrator's Kit (PDF command-line instructions)



GloMax[®] Discover is Designed to be SiLA Compliant

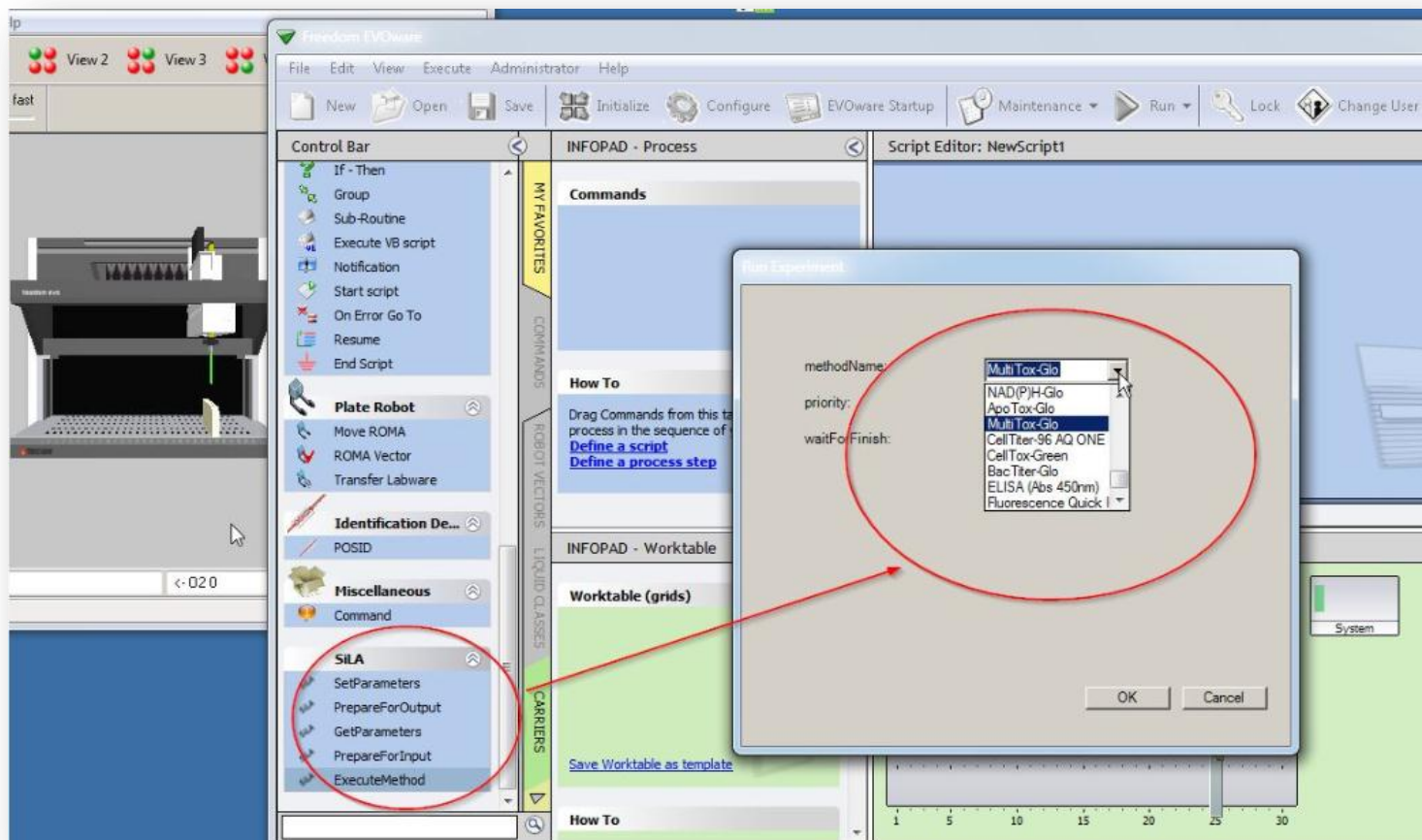
What is SiLA? (SStandards in LLaboratory AAutomation)

- A new industry standard to standardize
 - Device Control and Interfaces
 - Data Capture, Labware, etc.
- Avoids the need for custom software drivers when integrating devices

SiLA Rapid Integration[®]

GloMax[®] Discover SiLA Driver Available in 2014

Demonstrated with Tecan EVOWare[®] 2.6



Promega's SiLA Driver will be available in 2014

GloMax[®] Discover

Integrated	Seamless workflow with Promega Cell and Reporter assays.
Performance	Broad dynamic range, high sensitivity, and low well-to-well cross-talk for more usable data from your experiment.
Easy-to-use	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. Auto-gain adjustment so end-users don't have to worry about it.
Connected to your Workflow	Stand-alone instrument or integrate with automation. Export data to your laboratory network, LIMS, or Cloud. 21CFR Part 11 electronic signature compliant as standard.

Additional information



Request a free demo at
www.promega.com/discover