

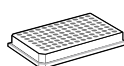
Live-Cell Kinetic Assays

for Time-Course Analysis

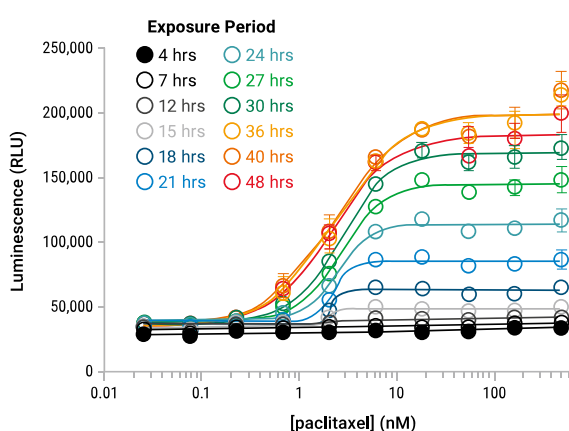
Cell Viability | Apoptosis | Cytotoxicity
Extracellular ATP | Reporter Gene Detection

Transform Your Time-Course Into a One-Plate Experiment

Real-Time Assay

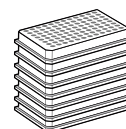


1 Plate



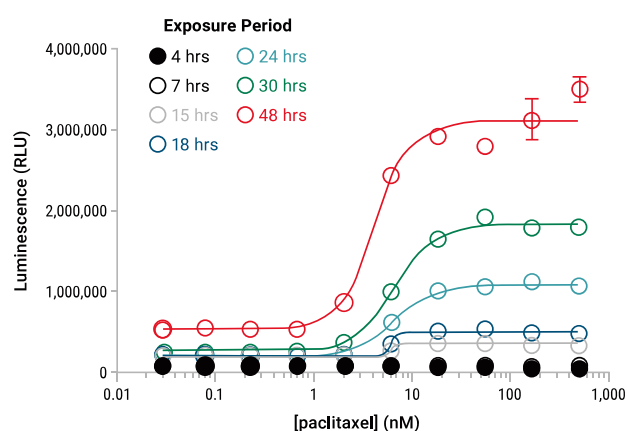
Multiple data points. **One assay plate.**

Endpoint Assay



7 Plates

vs.



Multiple data points. **Multiple assay plates.**

Left panel: Real-time assay. Data for each dose was obtained at multiple time points using a single assay plate with one **RealTime-Glo™ Annexin V Apoptosis Assay** Reagent addition.

Right panel: Endpoint assay. Seven separate plates were used to collect parallel data using an endpoint assay.

Why Use Live-Cell Kinetic Assays?

Live-cell kinetic assays allow the same sample well to be repeatedly measured over multiple time points. **This saves you time and effort, enabling you to collect more informative data in real-time.**

Live-Cell Kinetic Assay Benefits

No more missing time points:

Most endpoint assays lyse the cells to collect data only from a single time point. Live-cell kinetic assays are non-lytic and non-toxic. The cells remain viable, which allows you to detect kinetic changes over a longer time.

Save time, cells, and reagents:

Perform an extended time study with a single plate, instead of the need to have separate plates for each time point. A single-plate assay uses fewer cells and plates, also less culture media and other cell culture tools than multiple plates.

More data from one well with less variability:

Live-cell kinetic assays are non-lytic. This allows additional downstream applications, including multiplexing with other assays to gain more data from a single assay plate. The use of multiple assays from the same cells reduces assay variability in comparison to working with cells in multiple replicate plates.

Live-Cell Kinetic Assay Workflow

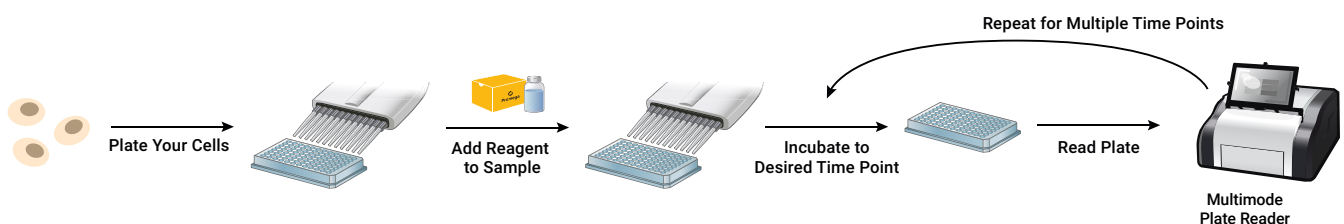
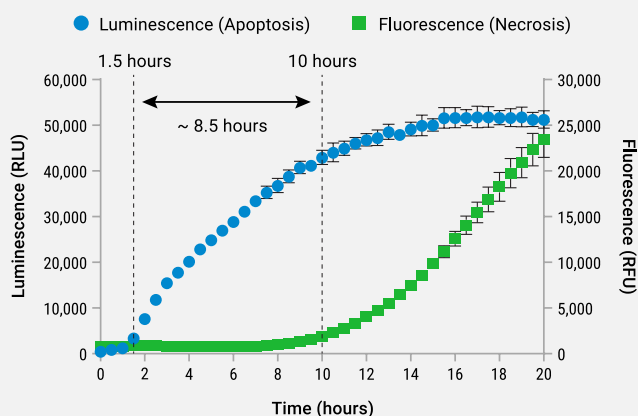


Plate your cells and add the reagent directly to the cells only once. The assay can be performed without further washing or processing steps. Read the plate at each time point, as many times as you need!

RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay

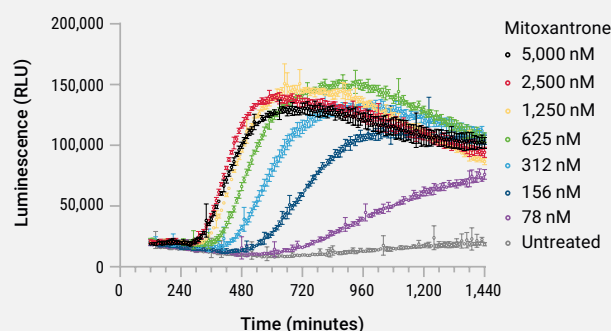
The RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay measures the real-time exposure of phosphatidylserine (PS) on the outer leaflet of cell membranes during the progression of apoptosis. The assay also includes a cell-impermeant, profluorescent DNA dye, which detects necrosis. This simple “add-read-assay” allows to perform real-time kinetics of up to 48 hours.



DLD-1 cells treated with 400 ng/ml rhTRAIL as extrinsic inducer of apoptosis. A time delay between the emergence of PS:Annexin V binding (luminescence) and the loss of membrane integrity (fluorescence) of 8.5 hrs indicate an apoptotic phenotype followed by secondary necrosis.

RealTime-Glo™ Extracellular ATP Assay

The RealTime-Glo™ Extracellular ATP Assay can be used for kinetic monitoring of ATP release from dying, stressed, or activated cells. Extracellular ATP (eATP) functions as a Damage Associated Molecular Pattern (DAMP) signal and is a key biomarker to determine whether a treatment induces Immunogenic Cell Death. The assay is based on an ATP-dependent luciferase reaction and allows real-time kinetics for up to 24 hours.



U937 cells were treated with a dilution of mitoxantrone, an anthracycline shown to induce Immunogenic Cell Death. RealTime-Glo™ Extracellular ATP Assay Reagent was added, and luminescence was collected every 10 minutes for 24 h.

Cell Health Assays Summary

Assay	Marker	Protocol Type	Plate Format	Instrument
RealTime-Glo™ MT Cell Viability Assay	Reducing potential of the cells	Endpoint or continuous readout up to 72 h	96/384/1536	Luminometer
CellTox™ Green Cytotoxicity Assay	DNA	Endpoint or continuous readout up to 72 h	96/384/1536	Fluorometer
RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay	PS, DNA	Continuous readout up to 48 h	96/384/1536	Luminometer (Apoptosis), Fluorometer (Necrosis)
RealTime-Glo™ Extracellular ATP Assay	eATP	Continuous readout up to 24 h	96/384	Luminometer

For more information download our brochure: www.promega.com/CellbasedAssays

Monitoring Bioluminescent Reporter Activity in Real-Time

Live-Cell Detection of NanoLuc® Reporter Activity

Real-time monitoring of reporter activity provides an option to better understand the kinetics of a wide variety of cellular responses. The NanoLuc® luciferase provides a small, sensitive bioluminescent reporter that is compatible with real-time monitoring of gene expression events. Promega offers a collection of NanoLuc® reporter vectors into which you can clone your region of interest, and ready-to-use NanoLuc® reporter vectors for monitoring signaling pathways.

The Nano-Glo® Live-Cell Detection Reagents enable the kinetic measurement of NanoLuc® reporter activity. The reagents differ in their signal brightness and stability and are thus suitable for a variety of experimental goals:

Nano-Glo® Live Cell Assay System

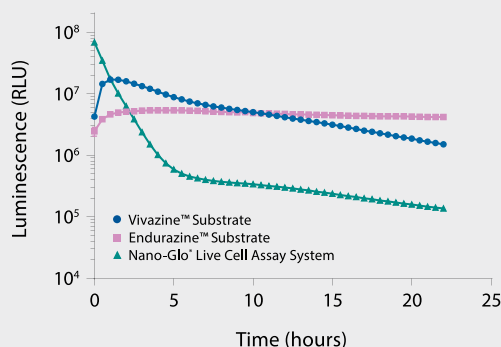
The Nano-Glo® Live Cell Assay System is a single-addition detection reagent, which offers the brightest signal and widest dynamic range for experiments less than 2 hours.

Nano-Glo® Endurazine™ Live Cell Substrate

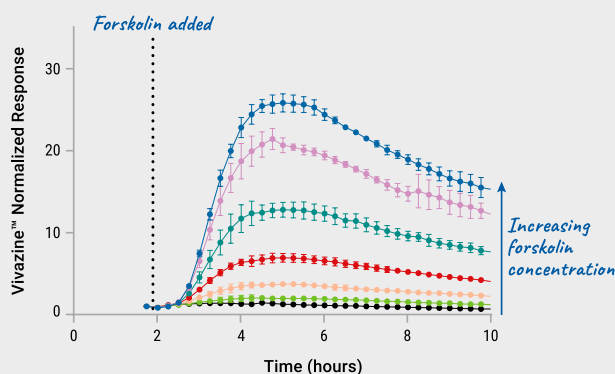
The Nano-Glo® Endurazine™ Live Cell Substrate provides the lowest initial signal intensity but has maximal stability for multi-day experiments.

Nano-Glo® Vivazine™ Live Cell Substrate

The Nano-Glo® Vivazine™ Live Cell Substrate is an intermediate option with increased initial brightness but also increased rate of signal decay compared to the Endurazine™ Substrate.



Comparison of Endurazine™ and Vivazine™ Substrates with the Nano-Glo® Live Cell Assay System for signal intensity and stability. A plasmid encoding NanoLuc® luciferase expressed from the PGK promoter was transiently transfected into HEK293 cells and luminescence signal intensity was measured as indicated.



HEK293 cells were transiently transfected with a construct expressing NlucP from a promoter with cAMP response elements (CRE). Vivazine™ Substrate was added to cells at time zero, and luminescence was measured continuously. After 2 h, varying concentrations of forskolin were added to induce expression of NlucP by increasing the intracellular cAMP concentration. The graph shows a 10 h time-course with per-well normalization of the data.

For more information about Nano-Glo® Live-Cell Reagents, visit: www.promega.com/LiveCellSubstrates

Ordering Information

Live-Cell Cell Health Assays	Quantity	Cat. #
RealTime-Glo™ MT Cell Viability Assay	100 assays	G9711
CellTox™ Green Cytotoxicity Assay	10 ml	G8741
RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay	100 assays	JA1011
RealTime-Glo™ Extracellular ATP Assay	200 assays	GA5010
Live-Cell Detection Reagents	Quantity	Cat. #
Nano-Glo® Live Cell Assay System	100 assays	N2011
Nano-Glo® Endurazine™ Live Cell Substrate	1 ml	N2571
Nano-Glo® Vivazine™ Live Cell Substrate	1 ml	N2581

Number of assays refers to a 96-well plate format.

To learn more about Promega Live-Cell Kinetic Assays, visit: www.promega.com/LiveCellKineticAssays



Cellular Energy Metabolism Assays

- Rapid, selective, and sensitive detection of glucose, lactate, glutamate, glutamine, and glycogen in biological samples
- Highly sensitive, plate-based bioluminescent methods
- Measuring multiple metabolites from a single sample

For more information download our brochure:
www.promega.com/CellularMetabolismAssays



Promega

Promega GmbH

Gutenbergring 10
69190 Walldorf
www.promega.com

Germany/Austria

Phone +49 6227 6906-290
de_techserv@promega.com

Poland

Phone +48 22 531 0668
pl_techserv@promega.com

Central/Eastern Europe

Phone +49 6227 6906-0
de_ceesales@promega.com