



Promega

Technical Bulletin

CellTiter-Glo[®] Luminescent Cell Viability Assay

INSTRUCTIONS FOR USE OF PRODUCTS G7570, G7571, G7572 AND
G7573.



www.promega.com

CellTiter-Glo[®] Luminescent Cell Viability Assay

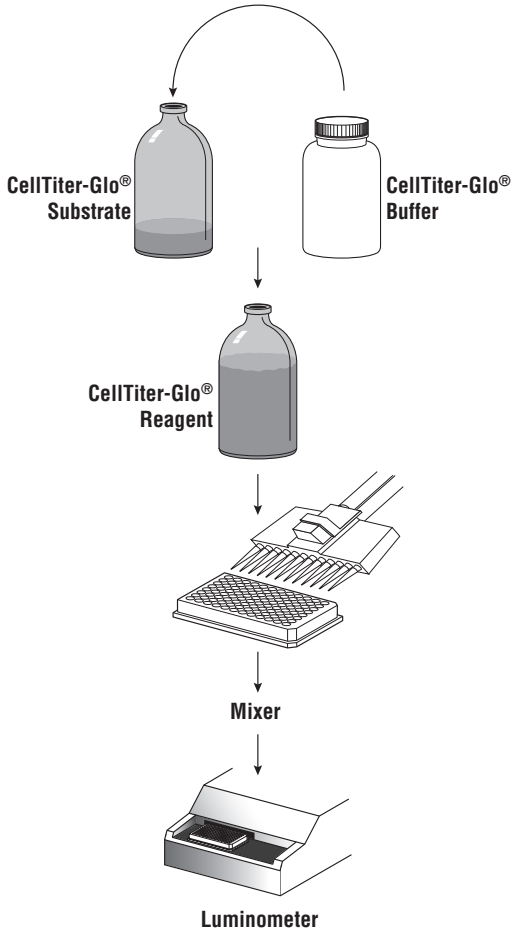
All technical literature is available on the Internet at: www.promega.com/tbs/
Please visit the web site to verify that you are using the most current version of this
Technical Bulletin. Please contact Promega Technical Services if you have questions on use
of this system. E-mail: techserv@promega.com

1. Description.....	1
2. Product Components and Storage Conditions	4
3. Performing the CellTiter-Glo [®] Assay.....	5
A. Reagent Preparation.....	5
B. Protocol for the Cell Viability Assay	6
C. Protocol for Generating an ATP Standard Curve (optional).....	6
4. Appendix	7
A. Overview of the CellTiter-Glo [®] Assay	7
B. Additional Considerations.....	8
C. References.....	10
D. Related Products.....	11

1. Description

The CellTiter-Glo[®] Luminescent Cell Viability Assay^(a-c) is a homogeneous method to determine the number of viable cells in culture based on quantitation of the ATP present, which signals the presence of metabolically active cells. The CellTiter-Glo[®] Assay is designed for use with multiwell-plate formats, making it ideal for automated high-throughput screening (HTS), and cell proliferation and cytotoxicity assays. The homogeneous assay procedure (Figure 1) involves adding a single reagent (CellTiter-Glo[®] Reagent) directly to cells cultured in serum-supplemented medium. Cell washing, removal of medium or multiple pipetting steps are not required.

The homogeneous “add-mix-measure” format results in cell lysis and generation of a luminescent signal proportional to the amount of ATP present (Figure 2). The amount of ATP is directly proportional to the number of cells present in culture in agreement with previous reports (1). The CellTiter-Glo[®] Assay relies on the properties of a proprietary thermostable luciferase (Ultra-Glo[™] Recombinant Luciferase), which generates a stable “glow-type” luminescent signal and improves performance across a wide range of assay conditions. The luciferase reaction for this assay is shown in Figure 3. The half-life of the luminescent signal resulting from this reaction is greater than five hours (Figure 4). This extended half-life eliminates the need for reagent injectors and provides flexibility for continuous or batch-mode processing of multiple plates. The unique homogeneous format reduces pipetting errors that may be introduced during the multiple steps required by other ATP-measurement methods.



3170MA12-0A

Figure 1. Flow diagram showing preparation and use of CellTiter-Glo[®] Assay Reagent.

System Advantages

- **Homogeneous:** “Add-mix-measure” format reduces the number of plate-handling steps to fewer than that required for similar ATP assays.
- **Fast:** Data can be recorded 10 minutes after adding reagent.
- **Sensitive:** Measures cells at numbers below the detection limits of standard colorimetric and fluorometric assays.
- **Flexible:** Can be used with various multiwell formats. Data can be recorded by luminometer or CCD camera imaging device.
- **Robust:** Luminescent signal is very stable, with a half-life >5 hours, depending on cell type and culture medium used.
- **Ability to Multiplex:** Can be used with reporter-gene assays or other cell-based assays from Promega (2,3).

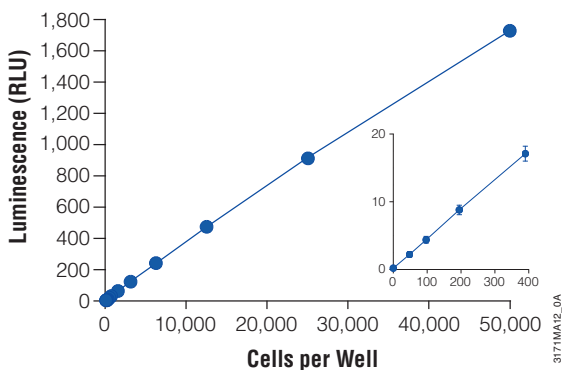


Figure 2. Cell number correlates with luminescent output. A direct relationship exists between luminescence measured with the CellTiter-Glo® Assay and the number of cells in culture over three orders of magnitude. Serial twofold dilutions of Jurkat cells were made in a 96-well plate in RPMI 1640 with 10% FBS according to the methods described in Section 3.B. Luminescence was recorded 10 minutes after reagent addition using a Dynex MLX® Microtiter® plate luminometer. Values represent the mean \pm S.D. of 4 replicates for each cell number. The luminescent signal from 50 Jurkat cells is greater than three standard deviations above the background signal resulting from serum-supplemented medium without cells. There is a linear relationship ($r^2 = 0.99$) between the luminescent signal and the number of cells from 0 to 50,000 cells per well.

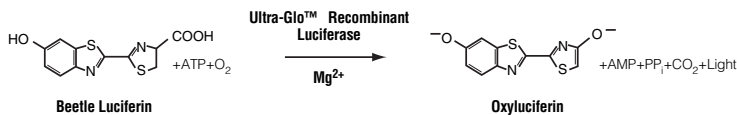


Figure 3. The luciferase reaction. Mono-oxygenation of luciferin is catalyzed by luciferase in the presence of Mg²⁺, ATP and molecular oxygen.

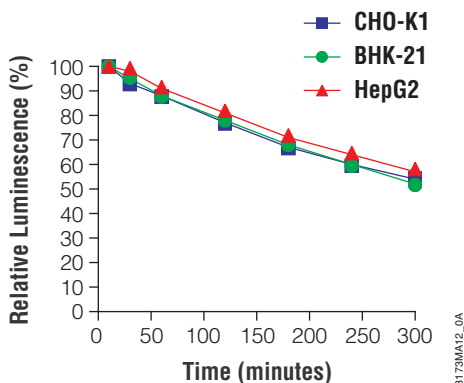


Figure 4. Extended luminescent half-life allows high-throughput batch processing. Signal stability is shown here for three common cell lines. HepG2 and BHK-21 cells were grown and assayed in MEM containing 10% FBS, while CHO-K1 cells were grown and assayed in DME/F-12 containing 10% FBS. CHO-K1, BHK-21 and HepG2 cells, 25,000 cells per well, were added to a 96-well plate. After an equal volume of CellTiter-Glo® Reagent was added, plates were shaken and luminescence monitored over time with the plates held at 22°C. The half-lives of the luminescent signals for the CHO-K1, BHK-21 and HepG2 cells were approximately 5.4, 5.2 and 5.8 hours, respectively.

2. Product Components and Storage Conditions

Product	Size	Cat.#
CellTiter-Glo® Luminescent Cell Viability Assay	10ml	G7570

For Laboratory Use. Substrate is sufficient for 100 assays at 100µl/assay in 96-well plates or 400 assays at 25µl/assay in 384-well plates. Includes:

- 1 × 10ml CellTiter-Glo® Buffer
- 1 vial CellTiter-Glo® Substrate (lyophilized)

Product	Size	Cat.#
CellTiter-Glo® Luminescent Cell Viability Assay	10 × 10ml	G7571

For Laboratory Use. Each vial of substrate is sufficient for 100 assays at 100µl/assay in 96-well plates or 400 assays at 25µl/assay in 384-well plates (1,000 to 4,000 total assays). Includes:

- 10 × 10ml CellTiter-Glo® Buffer
- 10 vials CellTiter-Glo® Substrate (lyophilized)

Product	Size	Cat.#
CellTiter-Glo® Luminescent Cell Viability Assay	100ml	G7572

For Laboratory Use. Substrate is sufficient for 1,000 assays at 100µl/assay in 96-well plates or 4,000 assays at 25µl/assay in 384-well plates. Includes:

- 1 × 100ml CellTiter-Glo® Buffer
- 1 vial CellTiter-Glo® Substrate (lyophilized)

Product	Size	Cat.#
CellTiter-Glo® Luminescent Cell Viability Assay	10 × 100ml	G7573

For Laboratory Use. Each vial of substrate is sufficient for 1,000 assays at 100µl/assay in 96-well plates or 4,000 assays at 25µl/assay in 384-well plates (10,000 to 40,000 total assays). Includes:

- 10 × 100ml CellTiter-Glo® Buffer
- 10 vials CellTiter-Glo® Substrate (lyophilized)

Storage Conditions: For long-term storage, the lyophilized CellTiter-Glo® Substrate and CellTiter-Glo® Buffer should be stored at -20°C. For frequent use, the CellTiter-Glo® Buffer can be stored at 4°C or at room temperature for 48 hours without loss of activity. See product label for expiration date information. Reconstituted CellTiter-Glo® Reagent (Buffer plus Substrate) can be stored at 4°C for 48 hours with ~5% loss of activity, at 4°C for 4 days with ~20% loss of activity or at -20°C for 21 weeks with ~3% loss of activity.

3. Performing the CellTiter-Glo® Assay

Materials to Be Supplied by the User

- opaque-walled multiwell plates adequate for cell culture
- multichannel pipette or automated pipetting station for reagent delivery
- device (plate shaker) for mixing multiwell plates
- luminometer or CCD camera imaging device capable of reading multiwell plates
- **optional:** ATP for use in generating a standard curve (Section 3.C)

3.A. Reagent Preparation

1. Thaw the CellTiter-Glo® Buffer, and equilibrate to room temperature prior to use. For convenience the CellTiter-Glo® Buffer may be **thawed and stored** at room temperature for up to 48 hours prior to use.
2. Equilibrate the lyophilized CellTiter-Glo® Substrate to room temperature prior to use.
3. Transfer the appropriate volume (10ml for the A size or 100ml for the B size) of CellTiter-Glo® Buffer into the amber bottle containing CellTiter-Glo® Substrate to reconstitute the lyophilized enzyme/substrate mixture. This forms the CellTiter-Glo® Reagent.

Note: The entire liquid volume of the CellTiter-Glo® Buffer bottle may be added to the CellTiter-Glo® Substrate vial.

3.A. Reagent Preparation (continued)

4. Mix by gently vortexing, swirling or by inverting the contents to obtain a homogeneous solution. The CellTiter-Glo® Substrate should go into solution easily in less than one minute.

3.B. Protocol for the Cell Viability Assay

1. Prepare opaque-walled multiwell plates with mammalian cells in culture medium, 100µl per well for 96-well plates or 25µl per well for 384-well plates.



Multiwell plates must be compatible with the luminometer used.

2. Prepare control wells containing medium without cells to obtain a value for background luminescence.
3. Add the compound to be tested to experimental wells, and incubate according to culture protocol.
4. Equilibrate the plate and its contents at room temperature for approximately 30 minutes.
5. Add a volume of CellTiter-Glo® Reagent equal to the volume of cell culture medium present in each well (e.g., add 100µl of reagent to 100µl of medium containing cells for a 96-well plate, or add 25µl of reagent to 25µl of medium containing cells for a 384-well plate).
6. Mix contents for 2 minutes on an orbital shaker to induce cell lysis.
7. Allow the plate to incubate at room temperature for 10 minutes to stabilize luminescent signal.

Note: Uneven luminescent signal within standard plates can be caused by temperature gradients, uneven seeding of cells or edge effects in multiwell plates.

8. Record luminescence. (**Note:** Instrument settings depend on the manufacturer. An integration time of 0.25–1 second per well should serve as a guideline.)

3.C. Protocol for Generating an ATP Standard Curve (optional)

It is a good practice to generate a standard curve using the plate on which samples are assayed.

1. Prepare 1µM ATP in culture medium (100µl of 1µM ATP solution contains 10⁻¹⁰ moles ATP).
2. Prepare tenfold serial dilutions of ATP in culture medium (1µM to 10nM; 100µl volumes would contain 10⁻¹⁰ to 10⁻¹² moles of ATP).

3. Prepare a multiwell plate with varying concentrations of standard ATP solution in 100µl medium (25µl for a 384-well plate).
Note: The ATP standard curve (Section 3.C) should be generated immediately prior to addition of the reagent because endogenous ATPase enzymes found in sera may reduce ATP levels.
4. Add a volume of CellTiter-Glo® Reagent equal to the volume of ATP standard present in each well.
5. Mix contents for 2 minutes on an orbital shaker.
6. Allow the plate to incubate at room temperature for 10 minutes to stabilize the luminescent signal.
7. Record luminescence.

4. Appendix

4.A. Overview of the CellTiter-Glo® Assay

The assay system uses the properties of a proprietary thermostable luciferase to enable reaction conditions that generate a stable “glow-type” luminescent signal while simultaneously inhibiting endogenous enzymes released during cell lysis (e.g., ATPases). Release of ATPases will interfere with the accurate measurement of ATP. Historically, firefly luciferase purified from *Photinus pyralis* (LucPpy) has been used in reagents for ATP assays (1,4-7). However, it has only moderate stability in vitro and is sensitive to its chemical environment, including factors such as pH and detergents, limiting its usefulness for developing a robust homogeneous ATP assay. Promega has successfully developed a stable form of luciferase, based on the gene from another firefly, *Photuris pennsylvanica* (LucPpe2), using an approach to select for characteristics that improve performance in ATP assays. The unique characteristics of this mutant (LucPpe2^m) enabled design of a homogeneous single-reagent-addition approach to perform ATP assays on cultured cells. Properties of the CellTiter-Glo® Reagent overcome the problems caused by factors such as ATPases that interfere with the measurement of ATP in cell extracts. The reagent is physically robust and provides a sensitive and stable luminescent output.

Sensitivity and Linearity: The ATP-based detection of cells has been shown to be more sensitive than other methods (8-10). Figure 2 shows that the luminescent signal from 50 Jurkat cells is greater than three standard deviations above the background signal resulting from serum-supplemented medium without cells. There is a linear relationship ($r^2 = 0.99$) between the luminescent signal and the number of cells from zero to 50,000 cells per well in the 96-well format. The luminescence values in Figure 2 were recorded after 10 minutes of incubation at room temperature to stabilize the luminescent signal according to the methods described in Section 3.B. Incubation of the same 96-well plate used in Figure 2 for 300 minutes at room temperature had little

4.A. Overview of the CellTiter-Glo® Assay (continued)

effect on the relationship between the luminescent signal and number of cells ($r^2 = 0.99$). Assay precision may vary up to 10% between 10 minutes and 5 hours for some cell lines. The conditions used to develop the CellTiter-Glo® Assay have resulted in a luminescent signal that is several fold brighter than existing glow-type ATP assay reagents, which require more than one pipetting step.

Speed: The homogeneous procedure to measure ATP using the CellTiter-Glo® Assay is quicker to perform than other ATP assay methods that require multiple steps to extract ATP and measure luminescence. The CellTiter-Glo® Assay also is faster than other commonly used methods to measure the number of viable cells (such as MTT, AlamarBlue or Calcein-AM) that require prolonged incubation steps to enable the cells' metabolic machinery to convert indicator molecules into a detectable signal.

4.B. Additional Considerations

Temperature: The intensity and rate of decay of the luminescent signal from the CellTiter-Glo® Assay depends on the rate of the luciferase reaction. Environmental factors that affect the rate of the luciferase reaction will change the intensity of light output and stability of the luminescent signal. Temperature is one factor that affects the rate of this enzymatic assay and thus the light output. For consistent results, equilibrate assay plates to a constant temperature before performing the assay. Transferring eukaryotic cells from 37°C to room temperature has little effect on the ATP content (5). We have demonstrated that removing cultured cells from a 37°C incubator and allowing them to equilibrate to 22°C for 1-2 hours had little effect on the ATP content. For batch-mode processing of multiple assay plates, take precautions to ensure complete temperature equilibration. Plates removed from a 37°C incubator and placed in tall stacks at room temperature will require longer equilibration than plates arranged in a single layer. Insufficient equilibration may result in a temperature gradient effect between the wells in the center and on the edge of the plates. The temperature gradient pattern also may depend on the position of the plate in the stack.

Chemicals: The chemical environment of the luciferase reaction affects the enzymatic rate and thus luminescence intensity. Differences in luminescence intensity have been observed using different types of culture medium and serum. The presence of phenol red in culture medium should have little impact on luminescence output. Assay of 0.1µM ATP in RPMI medium without phenol red showed ~5% increase in relative light units (RLU) luminescence output compared to RPMI containing the standard concentration of phenol red, whereas RPMI medium containing 2X the normal concentration of phenol red showed a ~2% decrease in RLU.

Solvents for the various chemical compounds tested for cytotoxicity may interfere with the luciferase reaction and thus the light output from the assay. Interference with the luciferase reaction can be determined by assaying a parallel set of control wells containing medium without cells. Dimethylsulfoxide (DMSO), commonly used as a vehicle to solubilize organic chemicals, has been tested at final concentrations up to 2% in the assay and only minimally affects light output.

Plate Recommendations: We recommend using standard opaque-walled multiwell plates suitable for luminescence measurements. Opaque-walled plates with clear bottoms allowing microscopic visualization of cells also may be used; however, these plates will have diminished signal intensity and greater cross talk between wells. Opaque white tape may be used to decrease luminescence loss and cross talk.

Cellular ATP Content: Different cell types have different amounts of ATP, and values reported for the ATP level in cells vary considerably (1,4,11-13). Factors that affect the ATP content of cells may affect the relationship between cell number and luminescence. Anchorage-dependent cells that undergo contact inhibition at high densities may show a change in ATP content per cell at high densities, resulting in a nonlinear relationship between cell number and luminescence. Factors that affect the cytoplasmic volume or physiology of cells also will have an effect on ATP content. For example, oxygen depletion is one factor known to cause a rapid decrease in ATP (1).

Mixing: Optimum assay performance is achieved when the CellTiter-Glo® Reagent is completely mixed with the sample of cultured cells. Suspension cell lines (e.g., Jurkat cells) generally require less mixing to achieve lysis and extract ATP than adherent cells (e.g., L929 cells). Tests were done to evaluate the effect of shaking the plate after adding the CellTiter-Glo® Reagent. Suspension cells cultured in multiwell plates showed only minor differences in light output whether or not the plates were shaken after adding the CellTiter-Glo® Reagent. Adherent cells are more difficult to lyse and show a substantial difference between shaken and nonshaken plates.

Several additional parameters related to reagent mixing include: the force of delivery of CellTiter-Glo® Reagent, sample volume and dimensions of the multiwell. All of these factors may affect assay performance. The degree of reagent mixing required may be affected by the method used to add the CellTiter-Glo® Reagent to the assay plates. Automated pipetting devices using a greater or lesser force of fluid delivery may affect the degree of subsequent mixing required. Complete reagent mixing in 96-well plates should be achieved using orbital plate shaking devices built into many luminometers and the recommended 2-minute shaking time. Special electromagnetic shaking devices using a radius smaller than the well diameter may be required to efficiently mix contents of 384-well plates. The depth of the medium and geometry of the multiwell plates may have an effect on mixing efficiency. We recommend that the end user take these factors into consideration when performing the assay and determine whether a mixing step is necessary for the individual application.

4.C. References

1. Crouch, S.P.M. *et al.* (1993) The use of ATP bioluminescence as a measure of cell proliferation and cytotoxicity. *J. Immunol. Methods* **160**, 81–8.
2. Farfan, A. *et al.* (2004) Multiplexing homogeneous cell-based assays. *Cell Notes* **10**, 2–5.
3. Riss, T., Moravec, R. and Niles, A. (2005) Selecting cell-based assays for drug discovery screening. *Cell Notes* **13**, 16–21.
4. Kangas, L., Grönroos, M. and Nieminen, A.L. (1984) Bioluminescence of cellular ATP: A new method for evaluating cytotoxic agents in vitro. *Med. Biol.* **62**, 338–43.
5. Lundin, A. *et al.* (1986) Estimation of biomass in growing cell lines by adenosine triphosphate assay. *Methods Enzymol.* **133**, 27–42.
6. Sevin, B.U. *et al.* (1988) Application of an ATP-bioluminescence assay in human tumor chemosensitivity testing. *Gynecol. Oncol.* **31**, 191–204.
7. Gerhardt, R.T. *et al.* (1991) Characterization of in vitro chemosensitivity of perioperative human ovarian malignancies by adenosine triphosphate chemosensitivity assay. *Am. J. Obstet. Gynecol.* **165**, 245–55.
8. Petty, R.D. *et al.* (1995) Comparison of MTT and ATP-based assays for the measurement of viable cell number. *J. Biolumin. Chemilumin.* **10**, 29–34.
9. Cree, I.A. *et al.* (1995) Methotrexate chemosensitivity by ATP luminescence in human leukemia cell lines and in breast cancer primary cultures: Comparison of the TCA-100 assay with a clonogenic assay. *AntiCancer Drugs* **6**, 398–404.
10. Maehara, Y. *et al.* (1987) The ATP assay is more sensitive than the succinate dehydrogenase inhibition test for predicting cell viability. *Eur. J. Cancer Clin. Oncol.* **23**, 273–6.
11. Stanley, P.E. (1986) Extraction of adenosine triphosphate from microbial and somatic cells. *Methods Enzymol.* **133**, 14–22.
12. Beckers, B. *et al.* (1986) Application of intracellular ATP determination in lymphocytes for HLA-typing. *J. Biolumin. Chemilumin.* **1**, 47–51.
13. Andreotti, P.E. *et al.* (1995) Chemosensitivity testing of human tumors using a microplate adenosine triphosphate luminescence assay: Clinical correlation for cisplatin resistance of ovarian carcinoma. *Cancer Res.* **55**, 5276–82.

4.D. Related Products

Cell Proliferation Products

Product	Size	Cat.#
CellTiter-Fluor™ Cell Viability Assay* (fluorescent)	10ml	G6080
CellTiter-Blue® Cell Viability Assay (resazurin)	20ml	G8080
CellTiter 96® AQueous One Solution		
Cell Proliferation Assay (MTS, colorimetric)*	200 assays	G3582
CellTiter 96® AQueous Non-Radioactive		
Cell Proliferation Assay (MTS, colorimetric)*	1,000 assays	G5421
CellTiter 96® AQueous MTS Reagent Powder*	1g	G1111
CellTiter 96® Non-Radioactive		
Cell Proliferation Assay (MTT, colorimetric)*	1,000 assays	G4000

*For Laboratory Use. All products available in additional sizes.

Cytotoxicity Assays

Product	Size	Cat.#
CytoTox-Glo™ Cytotoxicity Assay (luminescent)*	10ml	G9290
MultiTox-Glo Multiplex Cytotoxicity Assay (luminescent, fluorescent)*	10 ml	G9270
MultiTox-Fluor Multiplex Cytotoxicity Assay (fluorescent)*	10ml	G9200
CytoTox-Fluor™ Cytotoxicity Assay* (fluorescent)	10ml	G9260
CytoTox-ONE™ Homogeneous Membrane Integrity Assay (LDH, fluorometric)	200–800 assays	G7890
CytoTox-ONE™ Homogeneous Membrane Integrity Assay, HTP	1,000–4,000 assays	G7892
CytoTox 96® Non-Radioactive Cytotoxicity Assay (LDH, colorimetric)*	1,000 assays	G1780

*For Laboratory Use. Most products available in additional sizes.

4.D. Related Products (continued)

Apoptosis Products

Product	Size	Cat.#
Caspase-Glo® 2 Assay*	10ml	G0940
Caspase-Glo® 6 Assay*	10ml	G0970
Caspase-Glo® 3/7 Assay*	2.5ml	G8090
Caspase-Glo® 8 Assay*	2.5ml	G8200
Caspase-Glo® 9 Assay*	2.5ml	G8210
Apo-ONE® Homogeneous Caspase-3/7 Assay	1ml	G7792
DeadEnd™ Fluorometric TUNEL System	60 reactions	G3250
DeadEnd™ Colorimetric TUNEL System	20 reactions	G7360
Anti-ACTIVE® Caspase-3 pAb	50µl	G7481
Anti-Cytochrome C mAb	100µg	G7421
Anti-PARP p85 Fragment pAb	50µl	G7341
Anti-pS ⁴⁷³ Akt pAb	40µl	G7441
Caspase Inhibitor Z-VAD-FMK, 20mM	50µl	G7231
	125µl	G7232

*For Laboratory Use. Most products available in additional sizes.

©U.S. Pat. Nos. 6,602,677 and 7,241,584, Australian Pat. Nos. 754312 and 785294, European Pat. No. 1131441 and other patents pending.

©The method of recombinant expression of *Coleoptera* luciferase is covered by U.S. Pat. Nos. 5,583,024, 5,674,713 and 5,700,673.

©U.S. Pat. No. 7,083,911, Australian Pat. No. 2002255553 and other patents pending.

© 2001, 2004, 2005, 2007, 2008, 2009 Promega Corporation. All Rights Reserved.

Anti-ACTIVE, Apo-ONE, Caspase-Glo, CellTiter 96, CellTiter-Blue, CellTiter-Glo and CytoTox 96 are registered trademarks of Promega Corporation. CellTiter-Fluor, CytoTox-Fluor, CytoTox-Glo, CytoTox-ONE, DeadEnd and Ultra-Glo are trademarks of Promega Corporation.

MLX is a registered trademark of Dynex Technologies, Inc. Microtiter is a registered trademark of Thermo Electron Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.