

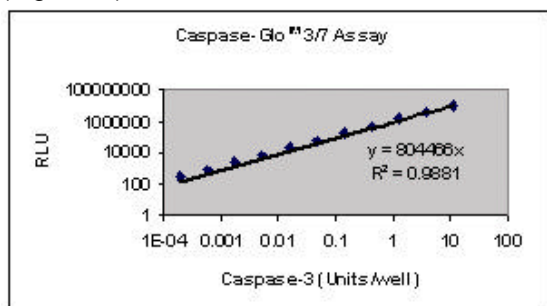
# A Veritas™ Microplate Luminometer Method for Promega's Caspase-Glo™ 3/7 Assay



## 1. INTRODUCTION

Turner BioSystems Veritas™ Microplate Luminometer in combination with Promega's Caspase-Glo™ Assay kit provides a convenient procedure for measuring caspase-3 or -7 activities. A proluminescent caspase-3/7 substrate contains the signature DEVD peptide sequence. Cell lysis occurs following the addition of Caspase-Glo 3/7 Reagent. Lysed cells release Caspase-3 or -7, which cleave the DEVD substrate from the aminoluciferin. Finally, the oxidation of the luciferin by luciferase produces light, which is measured by the Veritas. The amount of light produced is proportional to the amount of caspase-3 or -7 activity.

The ultra-sensitive Veritas enhances the detection limit of the Caspase-Glo 3/7 Assay. The Veritas detects fewer than  $2 \times 10^{-4}$  units/100  $\mu$ L of purified caspase-3 using Caspase-Glo 3/7 Substrate. Measurements are linear from  $1.9 \times 10^{-4}$  units/100  $\mu$ L to 11 units/100  $\mu$ L ( $R^2=0.9881$ ) or for five orders of magnitude (Figure 1.)



*Figure 1. Promega's Caspase-Glo™ Assay on the Veritas™ Microplate Luminometer using BIOMOL's purified caspase-3 diluted in 25 mM HEPES buffer containing 0.1% gelatin. The sample plate incubated for 1 hour at room temperature before measurement.*

## 2. MATERIALS REQUIRED

- Veritas™ Microplate Luminometer (P/N 9100-000)
- 96-well plates, white (E&K Scientific EK-25075)
- Caspase-Glo™ 3/7 Assay kit (Promega Catalog #'s G8090, G8091 and G8092)
- Purified Caspase-3 (Biomol Catalog # SE-169)
- p200 pipette and pipette tips
- Plate shaker

## 3. EXPERIMENT PROTOCOL

### 3.1 Reagent Preparation

**Caspase-Glo™ 3/7 Substrate:** Use as supplied. Store at -20°C.

**Caspase-Glo™ 3/7 Buffer:** Use as supplied. Store at -20°C. Buffer may be thawed and stored at room temperature for 48 hours without loss of activity.

**Caspase-Glo™ 3/7 Reagent:** Transfer the contents of one bottle of Caspase-Glo Buffer to one bottle of Caspase-Glo Substrate. Mix by inversion until the substrate is thoroughly dissolved. Use reconstituted reagent on the same day it is prepared or store at 4°C for 3 days without loss of activity. Reagent stored at -20°C for 4 weeks will give a signal 60% of freshly prepared reagent. Reagent stored at 4°C for 4 weeks will give a signal 75% of freshly prepared reagent.

**Note:** Reagent stored frozen after reconstitution must be thawed below 25°C to ensure reagent performance. Mix well after thawing. The simplest method for thawing is placing the reagent in a water bath at room temperature.

### 3.2 Instrument Setup

3.2.1 Double click on the Veritas icon to start the software.

3.2.2 Click on "Run Promega Protocol" from the "Welcome to Veritas" dialog box.

3.2.3 Open the "Caspase-Glo" template. Enter your information into the "Experiment," "Operator," "Plate No.," and "Notes" fields in the "Main Dialog Box".

3.2.4 Click on "Options" from the "Main Dialog Box" to select the wells to be read and modify the number of runs. Once modified, click the "Apply Changes" button to return to the "Main Dialog Box".

### 3.3 Sample Analysis of Cell Cultures

**Note:** For each set of experimental samples, prepare a negative control and a blank reaction. The blank reaction should consist of Caspase-Glo 3/7 Reagent, vehicle and cell culture medium without cells. The negative control should consist of Caspase-Glo 3/7 Reagent and vehicle-treated cells in cell culture medium. The term "vehicle" refers to the solvent used to dissolve the compound or protein of interest. The experimental samples should consist of Caspase-Glo 3/7 Reagent and treated cells in cell culture medium. Positive and negative controls should be prepared for each plate during multi-batch processing. Use identical cell numbers and volumes for the assay and the negative control samples.

3.3.1 Remove 96-well plates containing cells from the incubator and allow them to equilibrate to room temperature. At the same time, allow the reconstituted Caspase-Glo Reagent to equilibrate to room temperature.

3.3.2 Add a volume of Caspase-Glo Reagent equal to the volume of cell culture medium. For a 96-well plate, typically 100  $\mu$ L of Caspase-Glo Reagent is added to a well containing 100  $\mu$ L of cell culture medium.

3.3.3 Mix gently using a plate shaker at 300–500 rpm for 30 seconds. Incubate at room temperature for 30 minutes to 3 hours. The optimal incubation time will vary according to cell type and should be determined experimentally.

3.3.4 Insert plate into the Veritas and click on "Start" to begin assay. RLU values measured by the Veritas will appear in the Excel spreadsheet

after each selected wells has been read. If you encounter an error message, refer to the *Veritas Operating Manual* - troubleshooting guide for more information.

3.3.5 Once the measurements are complete, you can access Excel to analyze your data.

3.3.6 Remove your plate after measurement.

### 3.4 Sample analysis of purified enzyme

3.4.1 Prepare the blank reactions, positive controls and test samples as described in step 3.3 of this protocol.

3.4.2 Add a volume of Caspase-Glo Reagent equal to the volume of sample material.

3.4.3 Mix gently using a plate shaker at 300–500 rpm for 30 seconds. Incubate at room temperature for 20-60 minutes. Peak luminescence occurs sooner with purified enzymes than cultured cells.

3.4.4 Insert plate into the Veritas and click on "Start" to begin assay. RLU values measured by the Veritas will appear in the Excel spreadsheet after each selected well has been read. If you encounter an error message, refer to the *Veritas Operating Manual* - troubleshooting guide for more information.

3.4.5 Once the measurements are complete, you can access Excel to analyze your data.

3.4.6 Remove your plate after measurement.

## 4. ABOUT PROMEGA CORPORATION

Caspase-Glo is a trademark of Promega Corporation.

Orders for Promega's products may be placed by:

Phone: (800) 356-9526

Fax: (800) 356-1970

E-mail: [custserv@promega.com](mailto:custserv@promega.com)

Mailing Address:

Promega Corporation  
2800 Woods Hollow Rd.  
Madison, WI 53711  
USA

## 5. ABOUT BIOMOL

Orders for Biomol's products may be place by:

Phone: (800) 942-0430  
Fax: (610) 941-9252  
E-mail: [order@biomol.com](mailto:order@biomol.com)

Mailing Address:

Biomol International LP  
5120 Butler Pike  
Plymouth Meeting, PA 19462

## 6. ABOUT TURNER BIOSYSTEMS

Veritas is a trademark of Turner BioSystems.  
Orders for Turner BioSystems' products may be placed by:

Phone: (408) 636-2400 or  
Toll Free: (888) 636-2401 (US and Canada)  
Fax: (408) 737-7919

Web Site: [www.turnerbiosystems.com](http://www.turnerbiosystems.com)

E-Mail: [sales@turnerbiosystems.com](mailto:sales@turnerbiosystems.com)

Mailing Address:

Turner BioSystems, Inc.  
645 N. Mary Avenue  
Sunnyvale, CA 94085 USA

**CAUTION:** The lyophilized Caspase-Glo™ 3/7 Substrate contains dithiothreitol (DTT) and is therefore classified as hazardous. The reconstituted reagent is not known to present any hazards as the concentration of DTT is less than 1%. However, we recommend the use of gloves, lab coats and eye protection when working with these or any chemical reagents. Promega and Turner BioSystems assume no liability for damage resulting from handling or contact with these products.