

# A Veritas™ Microplate Luminometer Method for Promega Dual-Luciferase® Reporter (DLR) Assay



## 1. INTRODUCTION

The Turner BioSystems Veritas™ Microplate Luminometer in combination with Promega's Dual-Luciferase® Reporter (DLR) Assay kit provides a convenient, rapid, and sensitive procedure for quantifying gene expression. Transcriptional regulation, coupled to the expression of a luciferase reporter gene, is regularly used to study a wide range of biological events in cultured cells. Luciferase is an ideal reporter because of the absence of endogenous luciferase activity in mammalian cells, and the functional enzyme is created immediately upon translation<sup>1,2</sup>.

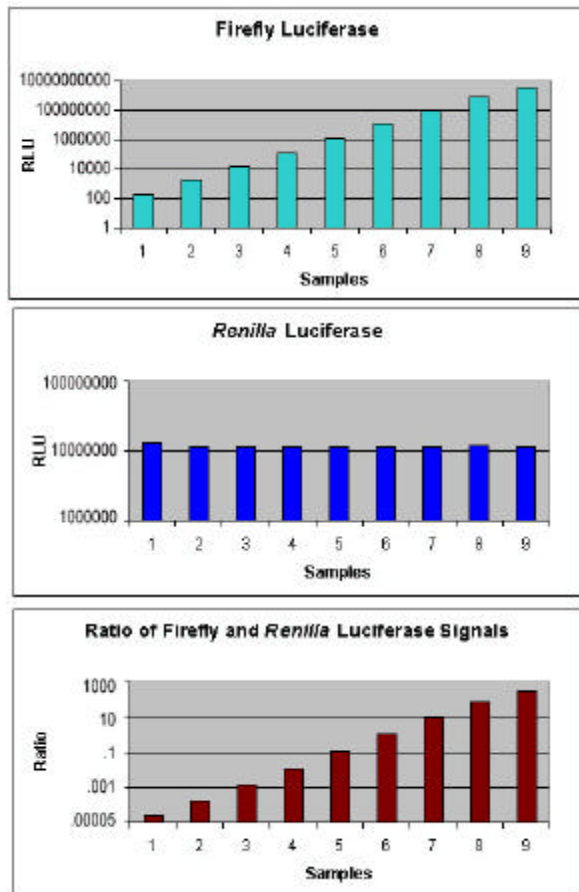
The Dual-Luciferase® Reporter (DLR) Assay System contains two different luciferase reporter enzymes that are expressed simultaneously in each cell. Typically, the experimental reporter is correlated with the effect of specific experimental conditions, while the activity of the co-transfected "control" reporter gene provides an internal control, which serves as the baseline response. Normalizing the experimental reporter gene to the activity of an internal control minimizes the variability caused by differences in cell viability and transfection efficiency. Thus, dual reporter assays allow more reliable interpretation of the experimental data by reducing extraneous influences. The experimental and control luciferase enzymes used in the Dual-Luciferase® Reporter (DLR) Assay have distinct evolutionary origins. The firefly luciferase and the *Renilla* (sea pansy) luciferase can discriminate between their respective bioluminescent substrates and do not cross-activate.

The firefly and *Renilla* substrates have been developed specifically to maximize the sensitivity of the assay reagent. This system is widely used in the life science research because of the superior light generation and high signal to noise ratio. The Dual-Luciferase® Reporter

(DLR) reagents are compatible with commonly used culture media for mammalian cells including RPMI 1640, MEM $\alpha$ , DMEM, and Ham's F12. These reagents can also be used with the Passive Lysis Buffer that comes with the kit and is available separately from Promega (Catalog # E1910).

The sensitivity and wide dynamic range offered by the Veritas™ Microplate Luminometer make it highly suited for the Dual-Luciferase® Reporter (DLR) Assay application. The Veritas™ software facilitates a quick set-up for the Dual-Luciferase® Reporter Assay with a pre-installed template. The internal automatic injectors are easy to use and completely accessible.

The Veritas™ Microplate Luminometer can detect as little as  $1 \times 10^{-19}$  moles firefly luciferase enzyme using Luciferase Assay Reagent II (LAR II) and  $1 \times 10^{-18}$  moles *Renilla* enzyme using Stop & Glo® Reagent. Measurements are linear for more than 8 and 6 orders of magnitude for firefly and *Renilla* substrates, respectively. All tests were conducted using purified recombinant firefly luciferase enzyme (Promega Catalog#E1701) and purified *Renilla* recombinant enzyme (Chemicon Catalog# 4400).



Figures 1—3. DLR assay was performed on the Veritas™ Microplate Luminometer using Promega's Dual-Luciferase® Reporter Assay System with recombinant firefly luciferase ( $1 \times 10^{-19}$  moles to  $1 \times 10^{-11}$  moles) and Renilla luciferase ( $1 \times 10^{-14}$  moles).

## 2. MATERIALS REQUIRED

- ❖ Veritas™ Microplate Luminometer (P/N 9100-002)
- ❖ 96-well plates, white (E&K Scientific EK-25075)
- ❖ Promega Dual-Luciferase Reporter® Assay kit (Promega Catalog #E1980)
- ❖ p200 pipette and pipette tip

## 3. EXPERIMENT PROTOCOL

### 3.1 Reagent Preparation

Luciferase Assay Buffer II and Luciferase Assay Substrate: Use as supplied. Store at  $-20^{\circ}\text{C}$ , where it is stable for up to 6 months. The Luciferase Assay Substrate may also be stored at  $4^{\circ}\text{C}$  for up to one month.

Transfer the contents of one bottle of Luciferase Assay Buffer II into one vial of Luciferase Assay Substrate. Mix by inversion until the substrate is thoroughly dissolved. Use reconstituted Luciferase Assay II Reagent (LAR II) on the same day it is prepared, or aliquot into working volume and store at  $-20^{\circ}\text{C}$  for 1 month or  $70^{\circ}\text{C}$  for up to one year.

Stop & Glo® Substrate and Stop & Glo® Buffer: Use as supplied. Store below  $-20^{\circ}\text{C}$ .

Stop & Glo® Buffer Substrate Solvent: Use as supplied. Store below  $25^{\circ}\text{C}$ .

To make Stop & Glo® Reagent, dilute the 50x Stop & Glo® Substrate to 1x concentration using Stop & Glo® Buffer in a glass or siliconized polypropylene tube. Mix by inversion. Use reconstituted Stop & Glo® Substrate on the same day it is prepared or store at  $-20^{\circ}\text{C}$  for up to two weeks.

Passive Lysis Buffer: To make 1x Passive Lysis Buffer, dilute the 5x Passive Lysis Buffer in DI water. Store below  $25^{\circ}\text{C}$ .

Note: The temperature of the Luciferase Assay Buffer II and Stop & Glo® Buffer should be held constant at room temperature while quantifying luminescence since luciferase activity is temperature dependent. Reagent stored frozen after reconstitution must be thawed below  $25^{\circ}\text{C}$  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 Browse the files in the "DLR" folder and choose the appropriate protocol. For example, if you are using DLR with one injector select "DLR with one injection."

3.2.4 Click on "Options" to select the wells to be read. The default settings on the Promega protocols pre-installed are optimized. However, you may wish to modify the integration time, delay before measurement and injection volume in the

"Other Options" tab. Once you have made your choices, click the "Apply Changes" button to accept changes or the "Save Protocol As" button to save the protocol. You can return to this screen by clicking "Options" from the "Main Dialog Box".

3.2.5 Enter your information into the "Experiment", "Operator", "Plate No.", and "Notes" fields in the "Main Dialog Box".

### 3.3 Sample Analysis

3.3.1 Prepare the 96-well plate containing lysed cell cultures.

**Note:** For maximum reproducibility, equilibrate cell cultures to room temperature before adding reagent.

3.3.2 Prepare the injectors. Place the intake tubing for injector 1 into the bottle of LAR II. Place the intake tubing for injector 2 into the bottle of Stop & Glo<sup>®</sup> Reagent. Prime both injectors using the "Prime" tab located on the "Main Dialog Box".

**Note:** Do not switch injectors. Residual Stop & Glo<sup>®</sup> Reagent will quench the firefly luciferase reporter activity. It is recommended to dedicate a single injector for Stop & Glo<sup>®</sup> Reagent and another injector for LAR II.

3.3.3 Insert the plate into the Veritas<sup>™</sup> Microplate Luminometer and click "Start" to begin assay. RLU values measured by the Veritas<sup>™</sup> Microplate Luminometer will appear in the Excel spreadsheet after all the selected wells in each row have been read. If you encounter an error message, refer to the troubleshooting guide for more information.

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

3.3.5 Remove your plate after the measurements are complete. You can choose the "Reverse Purge" tab to return unused reagent to the bottle. Be sure to flush injectors thoroughly after use.

### 4. REFERENCES

1. Ow, D.W. et al. (1986) Transient and stable expression of the firefly luciferase gene in plant

cells and transgenic plants. *Science* 234, 856—9.

2. De Wet, J.R. et al. (1987) Firefly luciferase gene: structure and expression in mammalian cells, *Mol. Cell. Biol.* 7, 725—37.

### 5. ABOUT PROMEGA CORPORATION

Dual-Luciferase and Stop & Glo are trademarks of Promega Corporation and are registered with the U.S. Patent and Trademark Office. 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

### 6. ABOUT TURNER BIOSYSTEMS

Veritas is a trademark of Turner BioSystems, Inc. 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 Luciferase<sup>®</sup> Assay 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.