

A GloMax[®] 20/20 Luminometer Method for Bright-Glo[™] Luciferase Assay System



INTRODUCTION

The GloMax[®] 20/20 Luminometer in combination with the Bright-Glo[™] Luciferase Assay System 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 (1,2).

The Bright-Glo[™] Luciferase Assay System has been developed specifically to maximize the sensitivity of the assay reagent while providing a luminescent signal half-life of approximately 25 minutes. The light signal can be measured 2 minutes after adding assay reagents. The Bright-Glo[™] Reagent is widely used in the pharmaceutical and biotechnology industries. The Bright-Glo[™] Reagent is compatible with commonly used culture media for mammalian cells containing 0–10% serum (RPMI 1640, MEM α , DMEM and Ham's F12).

The superior performance of the GloMax[®] 20/20 combined with the effectiveness of Bright-Glo[™] Reagent permits detection of very low levels of luciferase activity. The GloMax[®] 20/20 can detect as little as 1×10^{-20} moles luciferase enzyme using the Bright-Glo[™] Assay System. Measurements are linear from 1×10^{-20} to 1×10^{-12} moles luciferase or 8 orders of magnitude (Figure 1). All tests were conducted using the Bright-Glo[™] Luciferase Assay System (Cat.# E2620) and purified recombinant firefly luciferase enzyme (Cat.# E1701).

MATERIALS REQUIRED

- GloMax[®] 20/20 Luminometer
- 1.5 mL microcentrifuge tubes
- Bright-Glo[™] Luciferase Assay System (Cat.# E2610, E2620, E2650)
- p200 pipette and pipette tips

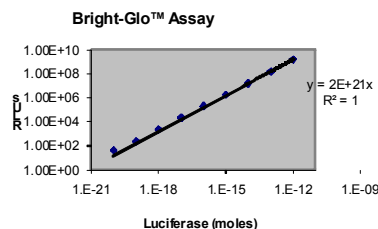


Figure 1. Bright-Glo[™] Assay was performed on the GloMax[®] 20/20 Luminometer using the Bright-Glo[™] Luciferase Assay System and recombinant luciferase.

EXPERIMENT PROTOCOL

1. Reagent Preparation

Bright-Glo[™] Substrate: Use as supplied. Store at -20°C , where it is stable for up to 6 months. The substrate also may be stored at 4°C for up to 1 month.

Bright-Glo[™] Buffer: Use as supplied. Store below 25°C .

Bright-Glo[™] Reagent: Transfer the contents of one bottle of Bright-Glo[™] Buffer to one bottle of Bright-Glo[™] Substrate. Mix by inversion until the substrate is thoroughly dissolved. Use reconstituted reagent on the same day it is prepared or divide into working aliquots and store at -20°C for up to 2 weeks.

Note: The temperature of the Bright-Glo[™] Reagent 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°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.

2. Instrument Setup

- Turn ON the GloMax[®] 20/20. A five-minute warm-up period is recommended but not necessary.
- Touch "Run Promega Protocol" from the "Protocols" menu.
- Select "BrightGlo" from the list of Promega protocols. The "Parameters" screen appears next with pre-programmed settings that are optimized for the Bright-Glo[™] Assay.
- Touch "OK" to go to the "Home" Screen.

3. Sample Analysis

- Remove the cell cultures from the incubator. **Note:** For maximum reproducibility, equilibrate cell cultures to room temperature before adding reagent.
- Add a volume of the Bright-Glo[™] Reagent equal to that of the culture medium.
- Wait a minimum of five minutes to allow for sufficient cell lysis. Then transfer the sample to a 1.5 mL microcentrifuge tube for analysis.
- Insert the tube into the GloMax[®] 20/20, and touch "Measure Luminescence" to begin measurement.

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.

CONTACT INFORMATION

Toll-Free: (800) 356-9526
Fax: (800) 356-1970

www.promega.com

Email: custserv@promega.com

Mailing Address:

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