

CellTiter-Glo® 2.0 Assay

Instructions for Use of Products G9241, G9242 and G9243

Quick Protocol

This Quick Protocol provides instructions for the CellTiter-Glo® 2.0 Assay designed for use with multiwell-plate formats, making the assay ideal for automated high-throughput screening applications.

The CellTiter-Glo® 2.0 Reagent is a ready-to-use reagent based on the original CellTiter-Glo® Luminescent Cell Viability Assay chemistry and eliminates the need to combine buffer with lyophilized substrate when preparing reagent. For detailed instructions, please refer to the *CellTiter-Glo® 2.0 Assay Technical Manual #TM403*, available at:

www.promega.com/protocols/

Reagent Preparation

Store at -30° C to -10° C. For maximum light signal, we recommend long-term storage at less than -65° C. Functional performance (linearity, sensitivity and signal half-life) of the reagent is maintained upon storage at -30° C to -10° C through the expiration date, although light output will decline over time.

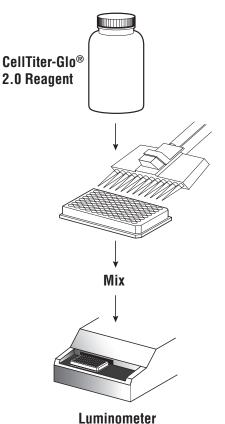
CellTiter-Glo® 2.0 Reagent can be thawed and stored at $+2^{\circ}$ C to $+10^{\circ}$ C for up to 2 months with >85% light output remaining when measuring 1µM ATP in a standard reaction. Do not refreeze the thawed reagent after extended storage above -10° C. Do not dispense CellTiter-Glo® 2.0 Reagent into aliquots due to the risk of ATP contamination.

CellTiter-Glo® 2.0 Reagent is light-sensitive and should be stored in an opaque container.

1. If frozen, thaw CellTiter-Glo® 2.0 Reagent at 4°C overnight.

Notes:

- a. Reagent may also be thawed in a 22°C water bath.
- b. Do not expose the reagent to temperatures above 25°C.
- c. If the CellTiter-Glo 2.0 Reagent is stored at -65°C, do not directly transfer it into a 22°C water bath to thaw. Instead, to avoid potentially cracking the bottle due to the rapid temperature change, leave the material on the bench top for 10-15 minutes and then place it in a 22°C water bath.
- 2. If not at room temperature, equilibrate the CellTiter-Glo® 2.0 Reagent to room temperature by placing the reagent in a 22°C water bath prior to use.
- Note: In a 22°C water bath, 100ml of the thawed reagent (4°C) requires approximately 30 minutes to equilibrate, and 500ml requires approximately 100 minutes to equilibrate.
- 3. Mix gently by inverting the contents to obtain a homogeneous solution. **Note:** Use caution when removing the seal of the CellTiter-Glo® 2.0 Reagent bottle to avoid introducing ATP contamination.





CellTiter-Glo® 2.0 Assay

Instructions for Use of Products G9241, G9242 and G9243

Quick Protocol

Protocol for the Cell Viability Assay

1. Prepare opaque-walled multiwell plates with mammalian cells in culture medium. Volumes and cell number should be optimized for experimental conditions.

Note: Multiwell plates must be compatible with the luminometer used.



- 3. Add test compound to experimental wells and incubate according to your culture protocol.
- 4. Equilibrate the plate and its contents to room temperature for approximately 30 minutes.
- 5. Add a volume of CellTiter-Glo® 2.0 Reagent equal to the volume of cell culture medium present in each well (e.g., for a 96-well plate, add 100µl of CellTiter-Glo® 2.0 Reagent to 100µl of medium containing cells).
- 6. Mix the 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 the luminescent signal.
- 8. Record luminescence.

Notes:

- a. Instrument settings depend on the manufacturer. Use an integration time of 0.25-1 second per well as a guideline.
- b. Uneven luminescent signal within plates can be caused by temperature gradients, uneven seeding of cells or edge effects in multiwell plates.

Sequential Multiplexing of CellTox™ Green Cytotoxicity Assay and CellTiter-Glo® 2.0 Assay

- 1. Completely thaw the CellTox™ Green Dye in a 37°C water bath. Mix the CellTox™ Green Dye using a vortex mixer to ensure homogeneity. A brief centrifugation may be necessary for complete recovery of the CellTox™ Green Dye. Although performance of CellTox™ Green Dye is optimal in black plates, white plates are optimal for sequential multiplexing formats that include a luminescent second measure.
- 2. Add the CellTox™ Green Dye at seeding or dosing so the final CellTox™ Green Dye concentration is 1X (stock = 1,000X) and incubate under desired conditions.
- 3. Equilibrate plate to room temperature.
- 4. Measure fluorescence intensity at 485–500nm_{Ex}/520–530nm_{Em}. Depending upon instrument manufacturer and optical detection configuration, replicate wells may benefit from 1 minute of shaking on an orbital shaker before measuring fluorescence.
- 5. Equilibrate the CellTiter-Glo® 2.0 Reagent to room temperature. After the final fluorescence measurement, add an equal volume of CellTiter-Glo® 2.0 Reagent to each well.
- 6. Place on an orbital shaker for 2 minutes and measure luminescence after 10 minutes.

CellTiter-Glo is a registered trademark of Promega Corporation. CellTox is a trademark of Promega Corporation.

