

# GloSensor™ cAMP Assay

INSTRUCTIONS FOR USE OF PRODUCTS E1171, E1261, E1290, E1291 AND E2301.

Quick  
PROTOCOL

## Live Cell Biosensor for GPCR Studies

The GloSensor™ cAMP Assay provides a simple and powerful new way to measure cAMP levels in cells. cAMP is a key second messenger involved in GPCR signaling, acting through  $G_s$ - and  $G_i$ -coupled proteins. The assay uses a genetically modified form of firefly luciferase containing a cAMP-binding protein moiety. Binding of cAMP causes a conformational change that leads to increased light output. The assay is well suited to HTS and uHTS platforms.

## Simple, Scalable Protocol

The GloSensor™ cAMP Assay is a 'zero step', non-lytic, live-cell assay ideal for kinetic and modulation studies of signaling through cAMP. The simple protocol is easily scalable to your throughput needs.

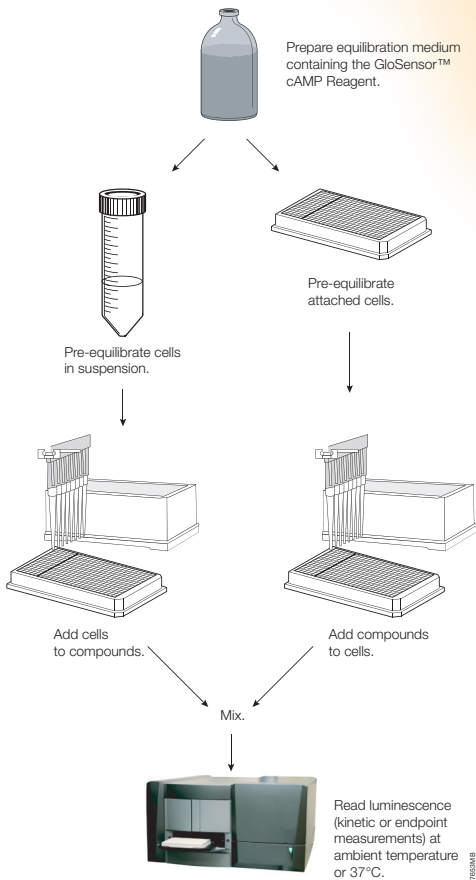
1. **Pre-equilibrate** cells with GloSensor™ cAMP Reagent for ~2 hours.
2. **Treat** cells with specific agonists/antagonists or library compounds.
3. **Measure** luminescence after 15–30 minutes.

## More Information

For the latest information on the GloSensor™ cAMP Assay and the GloSensor™ technology platform, including Frequently Asked Questions, noncommercial materials and more, visit:

[www.promega.com/glosensor](http://www.promega.com/glosensor)

For complete protocol information, see the *GloSensor™ cAMP Assay Technical Manual* #TM076, available at: [www.promega.com/tbs](http://www.promega.com/tbs)



## ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601