

GloSensor™ cAMP Assay

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

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

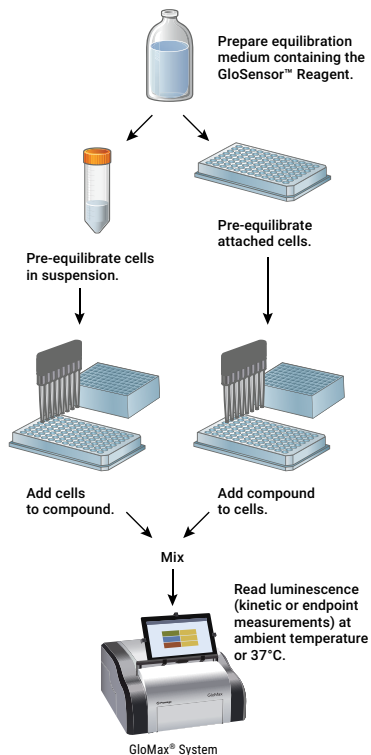
The GloSensor™ cAMP Assay is a 'zero step', nonlytic, live-cell assay ideal for kinetic and modulation studies of signaling through cAMP. The simple protocol is easily scalable to your throughput needs.

- 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

For complete protocol information, see the *GloSensor™ cAMP Assay Technical Manual #TM076*, available at: www.promega.com/protocols



Ordering and Technical Information

www.promega.com • Phone 608-274-4330
Toll free in USA 800-356-9526 • Fax 608-277-2601