

---

Product Manual

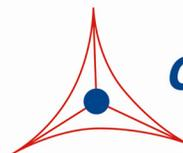
## cGMP ELISA Kit (Chemiluminescent)

### Catalog Numbers

STA-506	96 assays
STA-506-5	5 x 96 assays

**FOR RESEARCH USE ONLY**  
**Not for use in diagnostic procedures**

---



**CELL BIOLABS, INC.**  
*Creating Solutions for Life Science Research*

## **Introduction**

Guanosine 3',5'-cyclic monophosphate (cGMP) is a multi-functional second messenger involved in various cellular activities in many cell and tissue types. It is converted from guanosine triphosphate (GTP) via guanylyl cyclases (GC). It has been shown that the level of cGMP is typically 10-100 fold lower than cAMP in most tissues. cGMP primarily affects cellular activities through four different pathways. These include cGMP-dependent Protein Kinases (PKG/GK), cyclic nucleotide-gated (CNG) channels, cAMP-dependent Protein Kinase (PKA), and Phosphodiesterases.

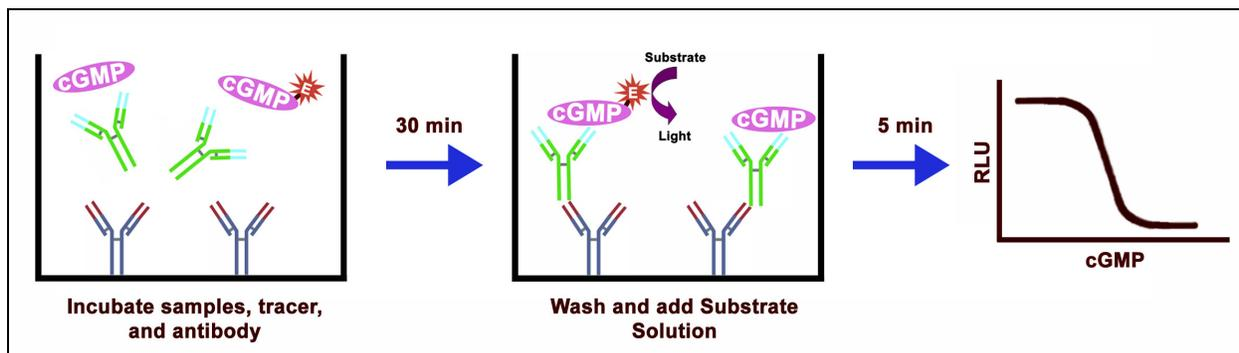
PKGs (PKG I and PKG II) are serine/threonine kinases activated by cGMP. PKG I has several putative targets, many of which are involved in the regulation of smooth muscle cell (SMC) contractility. Other cGMP/PKG I substrates that affect contractility include the myosin-binding subunit (MBS) of MLC Phosphatase and the small GTPase Rho. In addition to its role in smooth muscle relaxation, cGMP/PKG I may also regulate cell survival, proliferation, axon guidance, synaptic plasticity, inflammation, and angiogenesis.

Cell Biolabs' cGMP ELISA Kit is a competitive enzyme immunoassay designed to measure cGMP in cell culture supernatants, plasma, serum, saliva, urine, and cell lysates. The kit selectively measures cGMP levels without any significant cross reactivities to other nucleotides or cyclic nucleotides. The kit has a detection range of 1 pmol/mL to 1000 pmol/mL cGMP.

## **Assay Principle**

An anti-Rabbit IgG polyclonal coating antibody is adsorbed onto a microtiter plate. Cyclic GMP present in the sample or standard competes with Peroxidase cGMP Tracer for plate binding, in the presence of Rabbit Anti-cGMP Polyclonal Antibody.

Following incubation and wash steps, any Peroxidase cGMP Tracer bound to the plate is detected with addition of Chemiluminescent Reagent. The light product formed is inversely proportional to the amount of cGMP present in the sample. This reaction is then measured in a plate luminometer. A standard curve is prepared from cGMP Standard and sample concentration is then determined.



## **Related Products**

1. STA-500: cAMP ELISA Kit (Colorimetric)
2. STA-501: cAMP ELISA Kit (Chemiluminescent)
3. STA-505: cGMP ELISA Kit (Colorimetric)

## **Kit Components**

1. Goat Anti-Rabbit Antibody Coated Plate (Part No. 250101): One strip well 96-well plate.
2. cGMP Standard (Part No. 250501): One 200  $\mu$ L vial provided at 10 mM.
3. Rabbit Anti-cGMP Polyclonal Antibody (Part No. 250502): One 15  $\mu$ L vial.
4. Peroxidase cGMP Tracer Conjugate (Part No. 250503): One 30  $\mu$ L vial.
5. Assay Diluent (Part No. 250005): One 25 mL bottle.
6. Lysis Buffer (Part No. 250006): One 50 mL bottle.
7. 10X Wash Buffer (Part No. 250007): One 50 mL bottle.
8. Chemiluminescent Reagent A (Part No. 250102): One 6 mL amber bottle.
9. Chemiluminescent Reagent B (Part. No. 250103): One 6 mL bottle.

## **Materials Not Supplied**

1. Orbital plate shaker
2. 10  $\mu$ L to 1000  $\mu$ L adjustable single channel micropipettes with disposable tips
3. 50  $\mu$ L to 300  $\mu$ L adjustable multichannel micropipette with disposable tips
4. Multichannel micropipette reservoir
5. Plate Luminometer
6. (optional) Acetylation reagents: acetic anhydride and triethylamine

## **Storage**

Store kit components at 4 °C until their expiration dates. For longer term storage, maintain the Rabbit Anti-cGMP Polyclonal Antibody at -20 °C.

## **Preparation of Reagents**

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- Rabbit Anti-cGMP Polyclonal Antibody: Immediately before use dilute the Rabbit Anti-cGMP Antibody 1:500 with Assay Diluent. Do not store diluted solutions.
- Peroxidase cGMP Tracer Conjugate: Immediately before use dilute the Peroxidase cGMP Tracer Conjugate 1:100 with Assay Diluent. Do not store diluted solutions.

- **Chemiluminescent Reagent:** Immediately before use, mix equal volumes of Chemiluminescent Reagent A with Chemiluminescent Reagent B. Do not store diluted solutions.

### **Preparation of cGMP Standards**

1. Freshly prepare a dilution series of cGMP Standard in the concentration range of 1 mM – 1 nM by diluting the cGMP Standard in Lysis Buffer (Table 1).

<b>Standard Tubes</b>	<b>cGMP Standard (μL)</b>	<b>Lysis Buffer (μL)</b>	<b>cGMP Concentration</b>
1	40	360	1 mM
2	20 of Tube #1	180	100 μM
3	20 of Tube #2	180	10 μM
4	20 of Tube #3	180	1 μM
5	20 of Tube #4	180	100 nM
6	20 of Tube #5	180	10 nM
7	20 of Tube #6	180	1 nM
8	0	180	0

**Table 1. Preparation of cGMP Standard Curve**

### **Preparation of Samples**

1. **Urine, Plasma and Culture Medium Samples:** Urine and plasma may be tested directly or diluted with Lysis Buffer. Culture medium can also be tested with dilutions in Lysis Buffer.
2. **Cell Samples:** Aspirate medium. Add 1 ml of Lysis Buffer for every 35 cm<sup>2</sup> of surface area. Incubate at 4 °C for 20 minutes. Scrape cells off the surface with a cell scraper. Dissociate sample by pipetting up and down until suspension is homogeneous. Transfer to a centrifuge tube and centrifuge at top speed for 10 min. The supernatant can be assayed directly. Protein concentration >1 mg/ml is recommended for reproducible results.
3. **Tissue Samples:** Cyclic nucleotides may be metabolized quickly in tissue, so it is important to rapidly freeze tissues after collection (e.g., using liquid nitrogen). Weigh the frozen tissue and add 5-10 μL of Lysis Buffer per mg of tissue. Homogenize the sample on ice using a Polytron-type homogenizer. Spin at top speed for 5 min and collect the supernatant. The supernatant may be assayed directly.

## **Assay Protocol**

1. Prepare and mix all reagents thoroughly before use.
2. Each cGMP sample, cGMP Standard, and blank should be assayed in duplicate.
3. Add 50  $\mu\text{L}$  of cGMP sample or standard to the Goat Anti-Rabbit Antibody Coated Plate.
4. Add 25  $\mu\text{L}$  of diluted Peroxidase cGMP Tracer Conjugate (see Preparation of Reagents Section) to each tested well.
5. Add 50  $\mu\text{L}$  of diluted Rabbit Anti-cGMP Polyclonal Antibody (see Preparation of Reagents Section) to each tested well.
6. Cover with a Plate Cover and incubate at room temperature for 30 minutes with shaking.
7. Remove Plate Cover and empty wells. Wash microwell strips 5 times with 250  $\mu\text{L}$  1X Wash Buffer per well with thorough aspiration between each wash. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
8. Add 100  $\mu\text{L}$  of Chemiluminescent Reagent (see Preparation of Reagents Section) to each well, including the blank wells. Incubate at room temperature for 5 minutes on an orbital shaker.
9. Read the luminescence of each microwell on a plate luminometer.

## Example of Results

The following figures demonstrate typical cGMP ELISA results. One should use the data below for reference only. This data should not be used to interpret actual results.

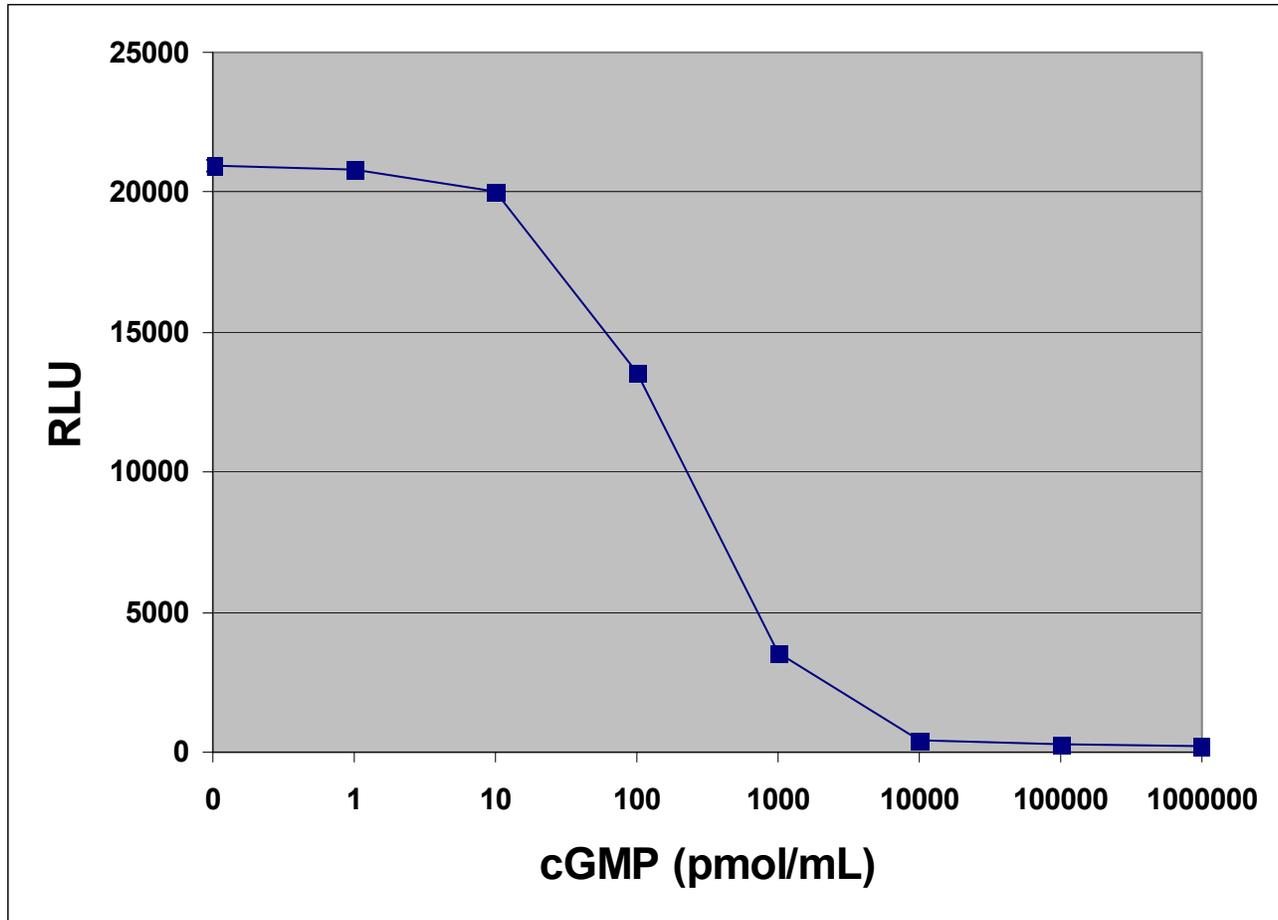


Figure 1: cGMP ELISA Standard Curve

### Cross reactivity of cGMP ELISA Kit

<u>Compounds</u>	<u>Cross Reactivity</u>
cGMP	100%
cAMP	<0.1%
AMP	<0.01%
ADP	<0.01%
ATP	<0.01%
GMP	<0.01%
GTP	<0.01%
CTP	<0.01%

## **References**

1. Hamad, A.M. et al. (2003) *Am. J. Physiol. Lung Cell Mol. Physiol.* **285**:L973.
2. Nakane, M. et al. (1990) *J. Biol. Chem.* **265**:16841.
3. Kuhn, M. (2003) *Circ. Res.* **93**:700.
4. Hood, J. and H.J. Granger. (1998) *J. Biol. Chem.* **273**:23504.
5. Kronemann, N. et al. (1999) *Br. J. Pharmacol.* **126**:349.
6. Suhasini, M. et al. (1998) *Mol. Cell. Biol.* **18**:6983.
7. Arancio, O. et al. (2001) *J. Neurosci.* **21**:143.
8. Arancio, O. et al. (1996) *Cell* **87**:1025.
9. Lev-Ram, V. et al. (1997) *Neuron* **18**:1025.
10. Zhuo, M. et al. (1994) *Nature* **368**:635.
11. Guzik, T.J. et al. (2003) *J. Physiol. Pharmacol.* **54**:469.

## **Warranty**

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

## **Contact Information**

Cell Biolabs, Inc.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2007-2008: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.