

# cGMP ELISA kit

**Most sensitive (0.088 pmol/ml) colorimetric kit for extracellular cGMP quantification.**

This is a colorimetric competitive immunoassay kit for the quantitative determination of extracellular cGMP. Sensitivity is increased by acetylation (reagents included). Absorbance is read at 405 nm. The ready-to-use liquid color-coded reagents provided reduce error.

Cyclic GMP (cGMP), which has been shown to be present at levels typically 10-100 fold lower than cAMP, is formed by the action of the enzyme guanylate cyclase on GTP. Some hormones, such as Acetylcholine, Insulin, and Oxytocin, as well as certain other chemicals like Serotonin and Histamine cause an increase in cGMP levels. Stimulators of guanylate cyclase such as the vasodilators nitroprusside, nitroglycerin, sodium nitrate, and nitric oxide (NO) also stimulate cGMP levels.

**Citations:** 34

[View Online »](#)

## Ordering Information

[Order Online »](#)

**ADI-901-013**

5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive – optional acetylation protocol increases sensitivity 10-fold (0.088 pmol/ml)
- Rapid – reliable results in just 3 hours
- Trusted – regularly cited in peer-reviewed publications
- Reduce errors – ready-to-use color-coded reagents and pre-coated plates
- Simple – efficient and well-established sample handling protocols
- Reliable – consistent lot-to-lot performance
- Save with bulk kit package!

## Handling & Storage

**Use/Stability** Store all components at +4°C, except standard and conjugate at -20°C.

**Shipping** Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Alternative Name** cyclic GMP

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of cGMP in culture supernatants, plasma, serum, saliva, and urine from any species.

**Assay Time** 3 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** GxR Microtiter plate, Conjugate, Antibody, Assay buffer 2 concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution, Acetic anhydride, Triethylamine

**Crossreactivity** cyclic GMP (100%) and <0.001%: GMP, GTP, cyclic AMP, AMP, ATP, cyclic UMP, CTP

**Sensitivity** (non-acetylated) 0.37 pmol/ml; (acetylated) 0.088 pmol/ml (range (non-acetylated) 0.16 – 500 pmol/ml; (acetylated) 0.16 – 100 pmol/ml)

**Species Reactivity** Species independent

**Wavelength** 405 nm

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)