

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 »](#)

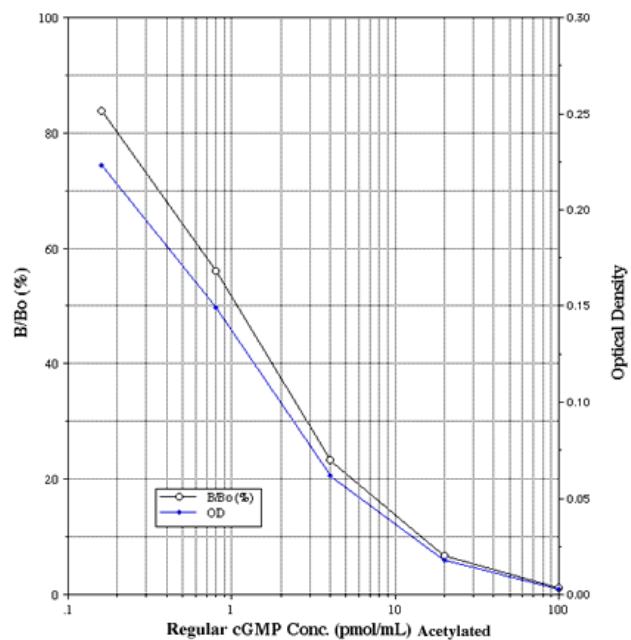
ADI-900-013

96 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



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 IgG Microtiter plate, Conjugate, Antibody, Assay buffer 2, 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)