

cGMP complete ELISA kit

Most sensitive (0.043 pmol/ml) and complete colorimetric kit for intra- and extracellular cGMP quantification.

This is a colorimetric competitive immunoassay kit for the quantitative determination of extracellular cGMP diluted in buffer or intracellular cGMP in cells or tissues lysed in 0.1 M HCl. 0.1M HCl is used to stop endogenous phosphodiesterase activity and stabilize the released cyclic GMP. The 0.1M HCl treated samples are then analyzed directly in a microtiter plate without extraction, drying and reconstitution. 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-900-164

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Versatile and convenient kit to assay both intracellular and extracellular cGMP
- High sensitivity with an optional acetylation protocol which increases sensitivity 10-fold (0.059 pmol/ml)
- Reliable and consistent lot-to-lot performance
- Higher throughput format with results in just 3 hours for up to 39 samples in duplicate
- Simple, efficient and well-established sample handling protocols

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 cell lysates, culture supernatants, plasma, serum, saliva, tissue, 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, 0.1M HCl, Assay buffer 2, Neutralizing reagent, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution, Triethylamine, Acetic anhydride

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

Sensitivity Assay Buffer : 0.42 pmol/ml (non-acetylated) 0.043 pmol/ml (acetylated) HCl: 0.604 pmol/ml (non-acetylated) 0.059 pmol/ml (acetylated) (range 0.8 – 500 pmol/ml (non-acetylated); 0.08 – 50 pmol/ml (acetylated))

Species Reactivity Species independent

Wavelength 405 nm



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-
be@enzolifesciences.com](mailto:info-be@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)