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 »

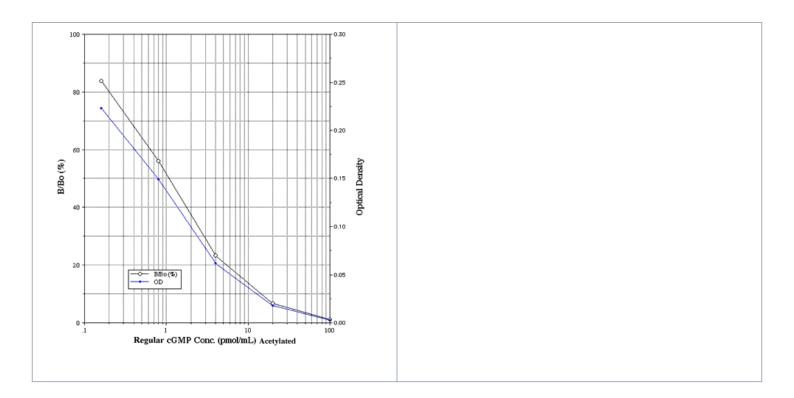
ADI-900-013

96 wells

Manuals, SDS & CofA

View Online »

- Sensitive optional acetylation protocol increases sensitivity 10fold (0.088 pmol/ml)
- Rapid reliable results in just 3 hours
- Trusted regularly cited in peerreviewed publications
- Reduce errors ready-to-use color-coded reagents and precoated plates
- Simple efficient and wellestablished 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 NotesFor 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

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com