Direct cGMP ELISA kit

Most sensitive (0.025 pmol/ml) colorimetric kit for intracellular cGMP quantification.

This is a colorimetric competitive immunoassay kit for the quantitation of intracellular cGMP in cells or tissues lysed in 0.1 M HCl, which 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: 61

View Online »

Ordering Information

Order Online »

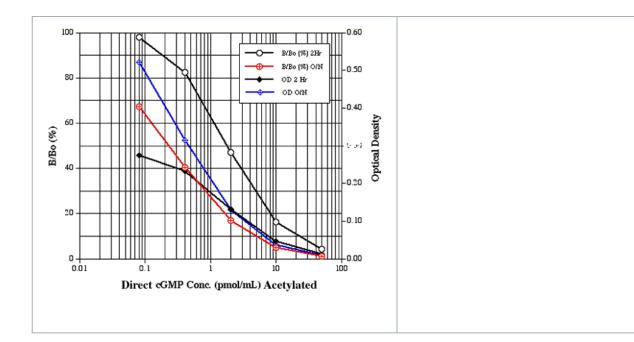
ADI-900-014

96 wells

Manuals, SDS & CofA

View Online »

- Highly sensitive ELISA with an optional acetylation protocol which increases sensitivity 10-fold (0.025 pmol/ml)
- Trusted and widely cited in peerreviewed publications
- Higher throughput format with results for up to 39 samples in duplicate
- Simple, efficient, and wellestablished sample handling protocols
- Reliable and 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 cell lysates

and tissue from any species. Cited sample types include

cerebral spinal fluid and microdialysate.

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,

Neutralizing reagent, 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 format: 0.604 pmol/mlAcetylated 2-hour

format: 0.059 pmol/mlAcetylated overnight format: 0.025

pmol/ml Range: (non-acetylated): 0.8 – 500 pmol/mlRange: (acetylated): 0.08 – 50 pmol/ml)

Species Reactivity Species independent

Wavelength 405 nm



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