

Direct cAMP ELISA kit

Most sensitive (0.006 pmol/ml) colorimetric kit for intracellular cAMP quantification.

This is a colorimetric competitive immunoassay kit for the quantitative determination of intracellular cAMP in cells or tissues lysed in 0.1 M HCl, which is used to stop endogenous phosphodiesterase activity and stabilize the released cyclic AMP. The 0.1M HCl treated samples are then analyzed directly in a microtiter plate without extraction, drying and reconstitution. Sensitivity is increased >10-fold by acetylation (reagents included). Absorbance is read at 405 nm. The large signal-to-background ratio offers superior sensitivity compared to competitors. This kit lets you easily assess adenylyl cyclase activation by GPCRs.

Cyclic AMP (cAMP) is one of the most important “second messengers” involved as a modulator of physiological processes. A number of hormones are known to activate cAMP through the action of the enzyme Adenylate cyclase which converts ATP to cAMP. cAMP has been shown to be involved in the cardiovascular and nervous systems, immune mechanisms, cell growth and differentiation, and general metabolism.

Citations: 158

[View Online »](#)

Ordering Information

[Order Online »](#)

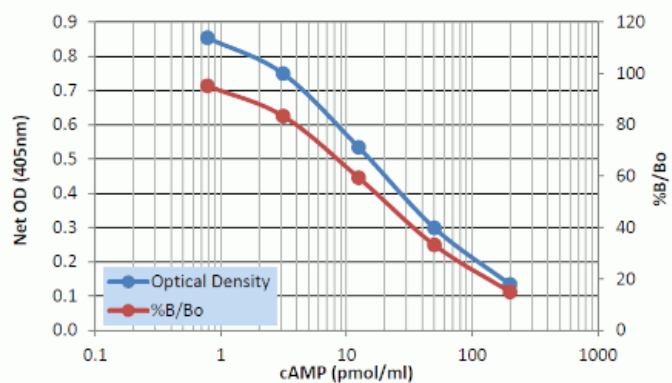
ADI-901-066A

5x96 wells

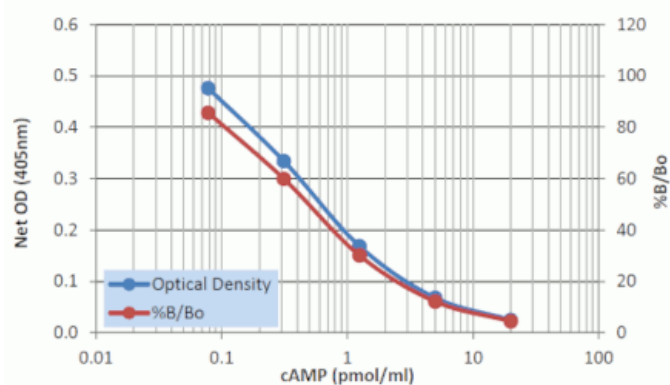
Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive ELISA with an optional acetylation protocol increasing sensitivity >10-fold
- High throughput format with results in <3 hours for up to 39 samples in duplicate
- Widely cited in peer-reviewed publications
- Easy-to-use pre-coated plate and liquid color-coded reagents minimizes errors
- Simple, efficient and well-established sample handling protocols
- Highly reliable with consistent lot-to-lot performance
- Save with bulk kit package!



Typical standard curve used to calculate cAMP concentrations for Non-acetylated assay format.



Typical standard curve used to calculate cAMP concentrations for Acetylated assay format.

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 AMP

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of cAMP in cell lysates and tissue from any species. Cited sample types include tissue culture.

Assay Time 3 hours

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

Contents GxR Microtiter plate, Conjugate, Antibody, 0.1M HCl, Neutralizing reagent, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution, Acetic anhydride, Triethylamine

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

Sensitivity Non-acetylated: 1.18 pmol/ml (range 0.78 – 200 pmol/ml)Acetylated: 0.006 pmol/ml (range 0.078 – 20 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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