

# cAMP ELISA kit

**Most sensitive (0.027 pmol/ml) colorimetric kit for extracellular cAMP quantification.**

This is a colorimetric competitive enzyme immunoassay kit for the quantitative determination of extracellular cAMP. Sensitivity is increased >10-fold by acetylation (reagents included). Absorbance is read at 405 nm. The ready-to-use liquid color-coded reagents provided reduce error.

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: 48

[View Online »](#)

## Ordering Information

[Order Online »](#)

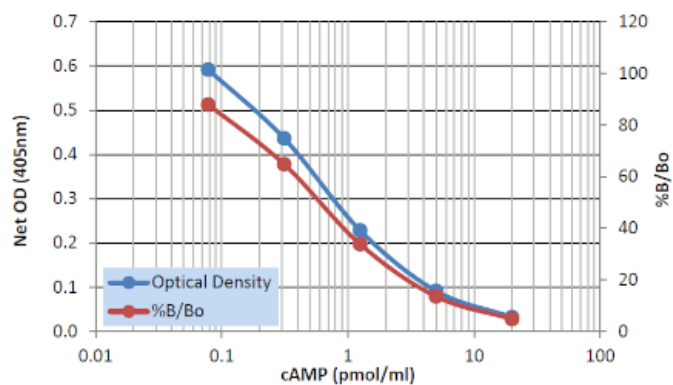
ADI-900-067A

96 wells

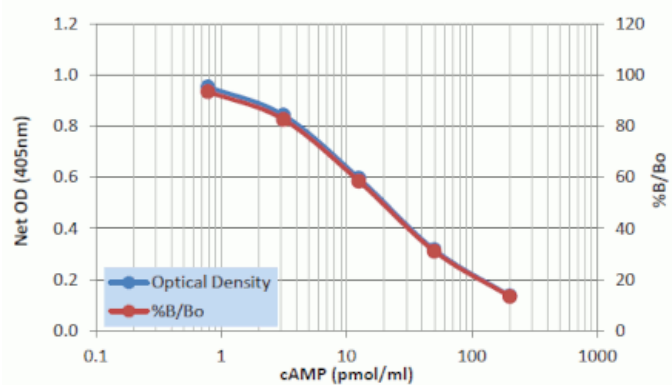
## Manuals, SDS & CofA

[View Online »](#)

- Trusted – regularly cited in peer-reviewed publications
- Sensitive – optional acetylation protocol increases sensitivity >10-fold (0.027 pmol/ml)
- Rapid – results in just 3 hours from up to 39 samples in duplicate
- Easy-to-use – pre-coated plate and liquid color-coded reagents reduce errors
- Simple – efficient and well-established sample handling protocols
- Reliable – consistent lot-to-lot performance



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



Typical standard curve used to calculate cAMP concentrations for Non-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 culture supernatants, serum, and saliva from any species. Cited sample types include cell lysate, tissue, and urine.

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

**Sensitivity** Non-acetylated: 0.49 pmol/ml (range 0.78 – 200 pmol/ml)Acetylated: 0.027 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](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)