

Product datasheet

cAMP Assay Kit (Competitive ELISA) ab234585

[2 Images](#)

Overview

Product name	cAMP Assay Kit (Competitive ELISA)
Detection method	Colorimetric
Sample type	Urine, Plasma, Tissue Extracts, Cell Lysate, Cell culture media
Assay type	Quantitative
Sensitivity	0.1 nM
Product overview	cAMP Assay Kit ab234585 uses a sensitive colorimetric competitive ELISA method to quantify cAMP (cyclic AMP) levels in samples from mammals and other species.

The cAMP assay uses a 96-well plate supplied coated with an anti-cAMP antibody.

HRP-labeled cAMP, and free cAMP within a sample, are added to each well of the plate. The cAMP assay is based on the competition between the HRP-labeled cAMP and the free cAMP for the fixed number of cAMP antibody binding sites on the plate.

The more free cAMP is present, the less HRP-labeled cAMP is bound. After a washing step, the amount of bound HRP-labeled cAMP is detected using a colorimetric HRP substrate measured by absorbance at OD 405nm, 650nm, or 740nm. The level of signal is compared to a standard curve for free cAMP, which is run at the same time as the samples.

Unlike other cAMP assays, this kit does not require an acetylation step to achieve the maximum sensitivity.

Notes If you are using fluorometric detection, we recommend [cAMP Assay Kit \(Fluorometric\) \(ab138880\)](#).

Properties

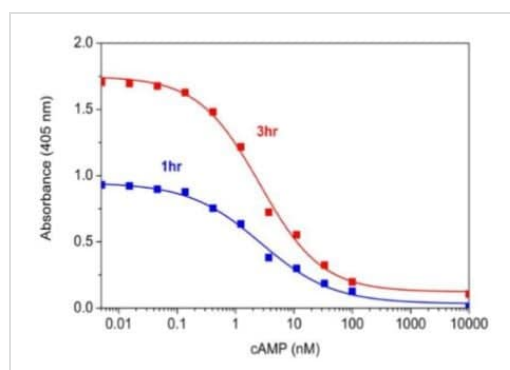
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests	10 x 96 tests
10X Wash Solution	1 x 10ml	1 x 100ml
Anti-cAMP Ab Coated 96-Well Plate	1 unit	10 units
Assay Buffer	1 x 20ml	1 x 100ml
cAMP Standard	1 vial	1 vial
Cell Lysis Buffer	1 x 10ml	1 x 100ml
Green Probe	1 x 10ml	1 x 100ml
HRP-cAMP Conjugate	1 vial	1 vial

Relevance

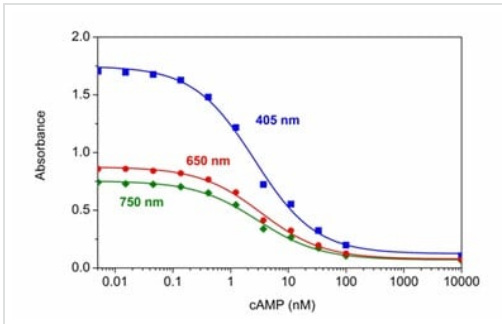
Adenosine 3',5'-cyclic monophosphate (cyclic AMP, cAMP) is an important "second messenger" involved in many physiological processes. It plays a key role as an intracellular second messenger for transduction events that follow a number of extracellular signals.

Images



cAMP dose response was measured with Colorimetric ELISA cAMP Assay Kit (ab234585) in a clear 96-well plate with a microplate reader.

The kit can detect as low as 0.1 nM cAMP in a 100 μ L reaction volume at 405nm after incubation with Green Probe for 1 hour (blue line) and 3 hours (red line).



The Absorbance can be read at 405 nm (blue line), 650 nm (red line) or 740 nm (Green line), the data are from the incubation with Green Probe for 3 hours.

cAMP dose response was measured with Colorimetric ELISA cAMP Assay Kit (ab234585) in a clear 96-well plate with a microplate reader.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors