

cAMP Assay Kit (Competitive ELISA) ab65355

★★★★★ [2 Abreviews](#) [83 References](#) [画像数 3](#)

医薬用外劇物

製品の概要

製品名	cAMP Assay Kit (Competitive ELISA)
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
アッセイタイプ	Quantitative
検出感度	> 0.002 μ M
検出範囲	0.002 μ M - 0.2 μ M
全工程の試験時間	3h 30m
種交差性	交差種: Mammals, Other species
製品の概要	cAMP Assay Kit ab65355 uses a sensitive colorimetric competitive ELISA method to quantify cAMP (cyclic AMP) levels in samples from mammals and other species. It has a sensitivity of 0.002 μ M.

The cAMP assay protocol uses a 96-well plate supplied coated with Protein G. The Protein G binds an anti-cAMP antibody which is added to the plate.

HRP-labeled cAMP, and free cAMP within a sample, are then 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 450 nm. The level of signal is compared to a standard curve for free cAMP, which is run at the same time as the samples.

This kit uses an acetylation step to achieve the maximum sensitivity.

cAMP assay protocol summary:

- add prepared samples and standards to tubes
- add neutralizing buffer
- add acetylating reagent mix, vortex and incubate for 10 min at room temp
- add assay buffer

- transfer to protein-G 96-well plate
- add cAMP antibody and incubate for 1 hr at room temp
- add cAMP-HRP and incubate for 1 hr at room temp
- wash with assay buffer
- add HRP developer and incubate for 1 hr at room temp
- stop reaction with HCl
- analyze with microplate reader

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K371 cAMP Direct Immunoassay Kit (Colorimetric). K371-100 is the same size as the 100 test size of ab65355.

If you are using fluorometric detection, we recommend **cAMP Assay Kit (Fluorometric)** (**ab138880**).

試験プラットフォーム

Microplate reader

製品の特性

保存方法

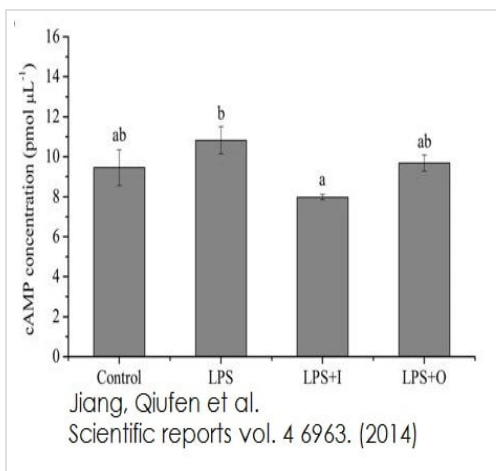
Store at -20°C. Please refer to protocols.

内容	100 tests
Acetylating Reagent A	1 x 750µl
Acetylating Reagent B	1 x 1.5ml
Assay Buffer XXXVI	1 x 25ml
cAMP Standard	1 x 10µl
cAMP-HRP	1 vial
Neutralization Buffer III	1 x 7.5ml
Protein G Coated Plate	1 unit
Rabbit Anti-cAMP Ab	1 vial
TMB Substrate I	1 x 10ml

関連性

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.

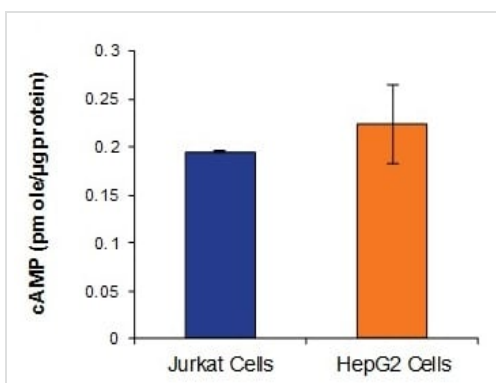
画像



Concentration of cAMP in response to in vitro haemocyte stimulation with LPS and β -AR (B) agonist/antagonist at 24h. Data presented as mean \pm S.D. (N = 6) of cAMP concentration expressed as pmol μL^{-1} . The significant differences among the control and treated groups were subjected to one-way ANOVA, followed by Student-Newman-Keuls multiple range test, which revealed two groups (a to b). The concentration of cAMP with the same letter in common do not differ at the P = 0.05 level of significance.

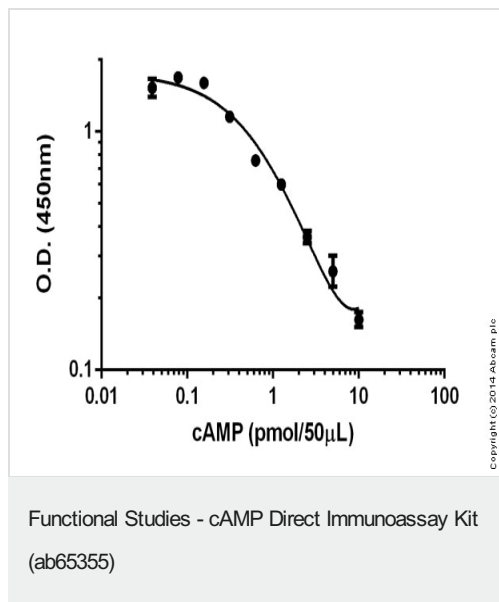
Functional Studies - cAMP Assay Kit (Competitive ELISA) (ab65355)

Jiang, Qiufen et al., Scientific reports vol. 4 6963., Fig 4, doi:10.1038/srep06963



Measurement of cAMP in Jurkat and HepG2 cells with ab65355. Assay performed using the kit protocol.

Functional Studies - cAMP Direct Immunoassay Kit (ab65355)



Standard curve (colorimetric): mean of duplicates (+/-SD)

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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