

# cAMP Assay Kit (Competitive ELISA, Fluorometric) ab138880

**15 References**    **画像数 4**

### 製品の概要

製品名	cAMP Assay Kit (Competitive ELISA, Fluorometric)
検出方法	Fluorescent
サンプルの種類	Urine, Plasma, Tissue Extracts, Cell Lysate, Cell culture media
アッセイタイプ	Competitive
検出感度	1 nM
全工程の試験時間	3h 00m
製品の概要	<p>cAMP Assay Kit ab138880 uses a sensitive fluorometric competitive ELISA method to quantify cAMP (cyclic AMP) levels in samples from mammals and other species.</p> <p>The cAMP assay protocol uses a 96-well plate supplied coated with an anti-cAMP antibody.</p> <p>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.</p> <p>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 fluorometric HRP substrate. The level of signal is compared to a standard curve for free cAMP, which is run at the same time as the samples.</p> <p>Unlike other cAMP assay protocols, this kit does not require an acetylation step to achieve the maximum sensitivity.</p>
特記事項	This product was previously called cAMP Direct Immunoassay Detection Kit (Fluorometric).
試験プラットフォーム	Microplate reader

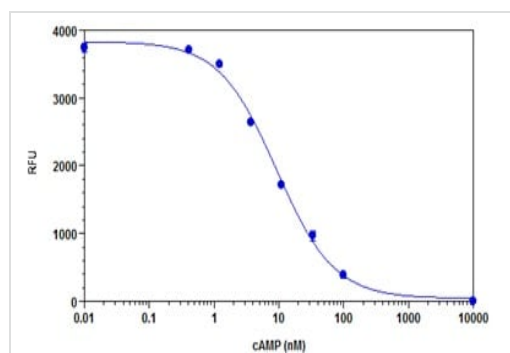
### 製品の特性

**保存方法**      Please refer to protocols.

内容	1 x 96 tests	10 x 96 tests
10X Wash Solution	1 x 10ml	1 x 100ml
AbRed Indicator	1 vial	1 vial
Anti-cAMP-coated 96-well plate	1 unit	1 x 10 units
Assay Buffer	1 x 20ml	1 x 100ml
cAMP Standard	1 x 33μg	1 x 33μg
Cell Lysis Buffer	1 x 10ml	1 x 100ml
HRP-cAMP Conjugate	1 vial	1 vial
Hydrogen Peroxide Solution	1 x 50μl	1 x 250μl
Substrate Buffer	1 x 10ml	1 x 100ml

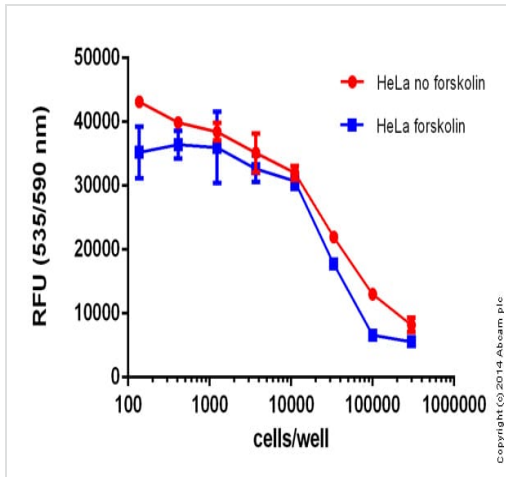
**関連性** Adenosine 3',5'-cyclic monophosphate (cyclic AMP, cAMP) is an important “second messenger” involved in many physiological processes. It plays a key role as an intracellular second messenger for transduction events that follow a number of extracellular signals.

## 画像



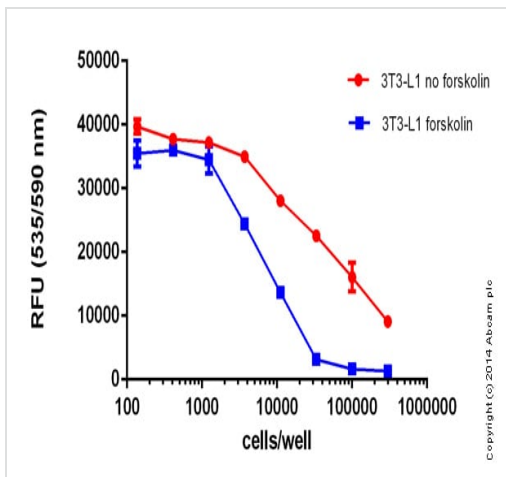
Sample Standard Curve for cAMP

cAMP dose response was measured with ab138880 in a solid black 96-well plate with a fluorescence microplate reader. The kit can detect as low as 1 nM cAMP in a 100 μl reaction volume.



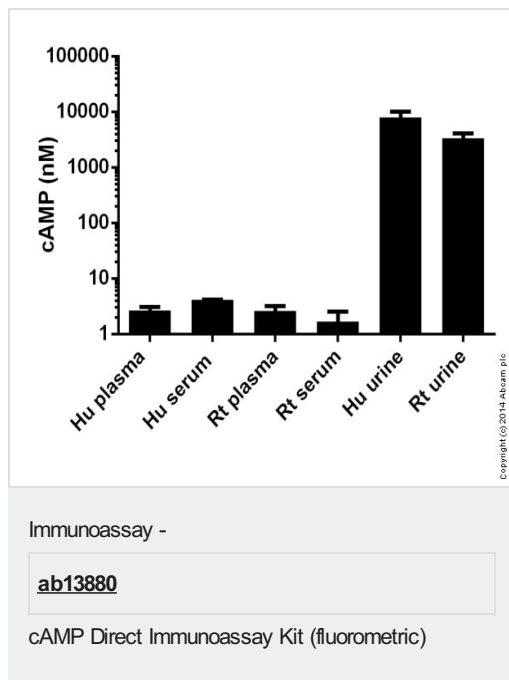
cAMP dose response in various amounts of HeLa cells after and without forskolin treatment (100 micromolar - [ab120058](#))

Immunoassay - 138880 cAMP Direct Immunoassay Kit (Fluorometric)



cAMP dose response in various amounts of 3T3-L1 cells after and without forskolin treatment (100 micromolar - [ab120058](#))

Immunoassay - cAMP Direct Immunoassay (fluorometric) (ab138880)



cAMP concentration (nM) measured in various biological fluids following assay protocol.

**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