

PKA Kinase Activity Assay Kit ab139435

53 References 3 图像

概述

产品名称

检测方法

精确度

样品类型

检测类型

检测时间

产品概述

平台

性能

PKA Kinase Activity Assay试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Sample				< 10%

批次间

样品	n	Mean	SD	CV%
Sample				<=12%

Adherent cells, Suspension cells, Tissue Extracts, Purified protein

Enzyme activity

4h 30m

PKA Kinase Activity Kit (ab139435) is a non-radioactive assay providing a safe, rapid and reliable method for the screening of inhibitors or activators of PKA and for quantitating the activity of PKA in purified or partially purified enzyme preparations. This kit is based on a solid phase enzyme-linked immuno-absorbent assay (ELISA) that utilizes a specific synthetic peptide as a substrate for PKA and a polyclonal antibody that recognizes the phosphorylated form of the substrate. The assay is designed for the analysis of PKA activity in the solution phase. This kit gives you the ability to an end-point or kinetic assay read-out in a convenient 96-well plate based assay and offers easy sample handling protocols.

The kit offers the following advantages:

1. Safe - non-radioactive measurement of kinase activity

2. Flexible - kinetic and end-point options available

3. Fast - results in < 4.5 hours

4. Efficient - only 30 µl diluted sample needed per well

Microplate reader

存放说明

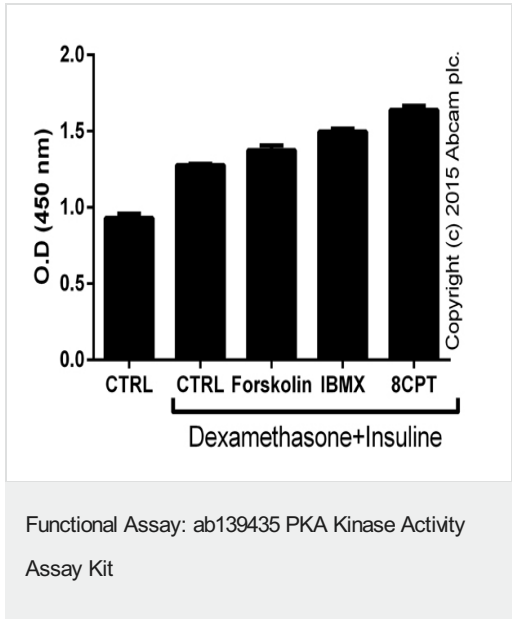
Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 30ml
Active PKA 20 ng/μl	1 x 30μl
Antibody Dilution Buffer	1 x 10ml
Anti-Rabbit IgG: HRP Conjugate	1 x 20μl
ATP	1 x 2mg
Kinase Assay Dilution Buffer	1 x 10ml
PKA Phosphospecific Substrate Antibody	1 x 5ml
PKA Substrate Microtiter Plate	1 unit
Stop Solution 2	1 x 10ml
TMB Substrate	1 x 10ml

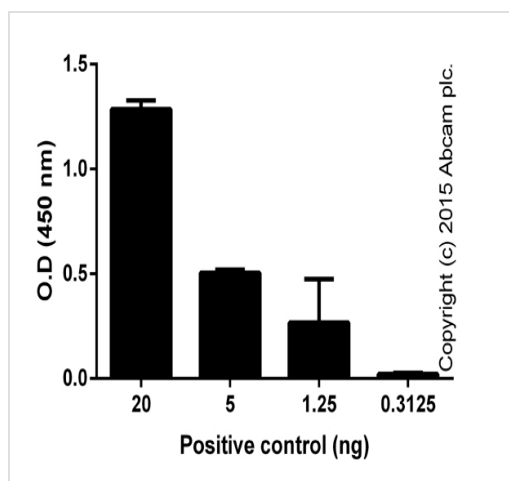
相关性

In cell biology, Protein kinase A (PKA) is a family of enzymes whose activity is dependent on cellular levels of cyclic AMP (cAMP). PKA is also known as cAMP-dependent protein kinase. Protein kinase A has several functions in the cell, including regulation of glycogen, sugar, and lipid metabolism.

图片

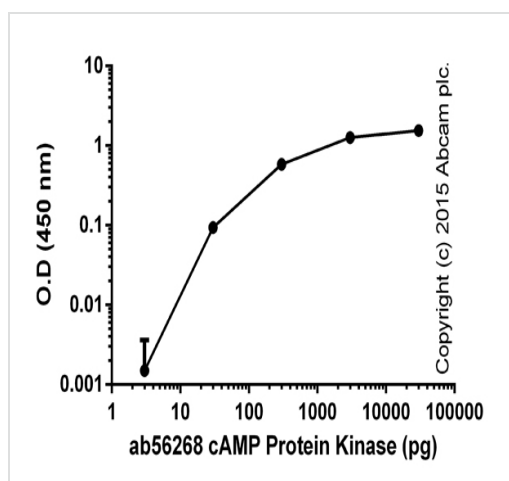


1e7 3T3-L1 cells were treated for 15 minutes with 1 μM dexamethasone ([ab120743](#)) and 1 μg/mL insulin ([ab123768](#)) with the addition of 50 μM forskolin ([ab120058](#)), 0.5 mM IBMX ([ab120840](#)) or 0.1 mM 8-CPT-cAMP ([ab120424](#)). Cell extracts were prepared according to the protocol. Lysates were diluted 50-fold, and tested in duplicates (+/-SD).



Titration of positive control (active PKA), background signal subtracted (duplicates; +/- SD).

Functional Assay: ab139435 PKA Kinase Activity
Assay Kit



Titration of cAMP Protein Kinase (**ab56268**), background signal subtracted (duplicates; +/- SD).

Functional Assay: ab139435 PKA Kinase Activity
Assay Kit

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.cn/abpromise> or contact our technical team.

Terms and conditions

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