

PKA kinase activity kit

Non-radioactive protein kinase assay
provides rapid and sensitive detection

The PKA Kinase activity kit is an efficient assay that utilizes little sample volume. 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 assay also offers a high signal to background ratio.

PKA (Protein Kinase A), also known as cAMP-dependent protein kinase, phosphorylates serine or threonine residues in target protein in response to elevated levels of cAMP. Inactive PKA exists as a tetrameric protein composed of two regulatory (R) subunits and two catalytic (C) subunits. Activation occurs when two cAMP molecules bind to each R subunit which causes a conformational change that releases the active C subunits.

Citations: 78

[View Online »](#)

Ordering Information

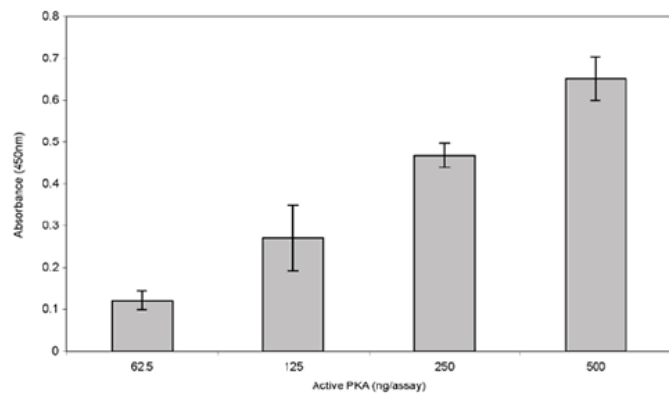
[Order Online »](#)

ADI-EKS-390A	96 wells
--------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Safe and non-radioactive colorimetric measurement of kinase activity
- Flexible format, kinetic and end-point options available
- Fast results in < 4.5 hours
- Efficiency maximized with only 30 µl diluted sample needed per well



Analysis of purified active PKA activity with EKS-390A. Varying quantities of purified active PKA were assayed using the PKA substrate microtiter plate. Assays were incubated for 60 min. at 30°C.

Handling & Storage

Use/Stability Store all components at 4°C, except active kinase at -80°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Protein kinase A, cAMP-dependent protein kinase

Application Activity assay, Colorimetric detection

Application Notes For the quantitative determination of PKA kinase activity in partially purified, purified, or crude enzyme preparations from any species.

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Antibody, Conjugate, Antibody dilution buffer, Kinase assay dilution buffer, ATP, Active Kinase, Wash buffer concentrate, TMB Substrate, Stop solution 2

Species Reactivity Species independent

UniProt ID P17612 (Ca subunit), P22694 (Cβ subunit), P22162 (Cy subunit), P10644 (RIα subunit), P31321 (RIβ subunit), P13861 (RIIα subunit), P31323 (RIIβ subunit)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com