

PKC kinase activity kit

Non-radioactive protein kinase activity assay provides rapid and sensitive detection

The PKC 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.

Protein Kinase C (PKC) is a large superfamily of serine/threonine kinases that mediate essential cellular signals required for activation, proliferation, differentiation and survival. There are at least ten PKC isotypes that are closely related in structure but that have distinct patterns of tissue distribution and function. The PKC isotypes can be subdivided into three classes based on primary structure and biochemical properties. These are: classical or conventional PKC isotypes (cPKC), novel PKC isotypes (nPKC) and atypical PKC isotypes (aPKC).

Citations: 66

[View Online »](#)

Ordering Information

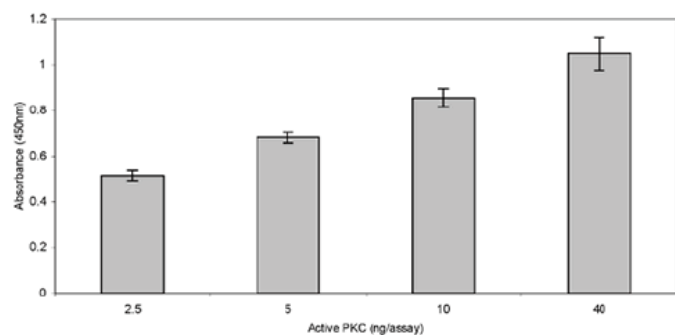
[Order Online »](#)

ADI-EKS-420A	96 wells
--------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Non-radioactive 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 PKC activity with EKS-420A. Varying quantities of purified active PKC were assayed using the PKC 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 C

Application Activity assay, Colorimetric detection

Application Notes For the measurement of PKC kinase activity partially purified, purified, or crude enzyme preparations from any species. Cited sample type includes tissue (ref. 1).

Assay Time 4.5 hours

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

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