

p70 S6K activity kit

The p70 S6K 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.

p70 S6K is a mitogen stimulated serine/threonine protein kinase that regulates the cell growth and G1 phase cell cycle progression. p70 S6K is activated by serum stimulation and can be inhibited by both wortmannin and rapamycin. Activation of p70 S6K by multi-site phosphorylation requires phosphoinositide 3-kinase (PI3-K)-dependent signal(s) leading to sequential phosphorylation at two key residues: Thr389 and Thr229. Thr229 phosphorylation is mediated by phosphoinositide-dependent kinase 1 (PDK-1), and is dependent upon prior phosphorylation of Thr389. Multiple residues in the p70 S6K pseudosubstrate domain (e.g. Ser411, Thr421, and Ser424) are also subject to activating phosphorylation by upstream kinases.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-EKS-470

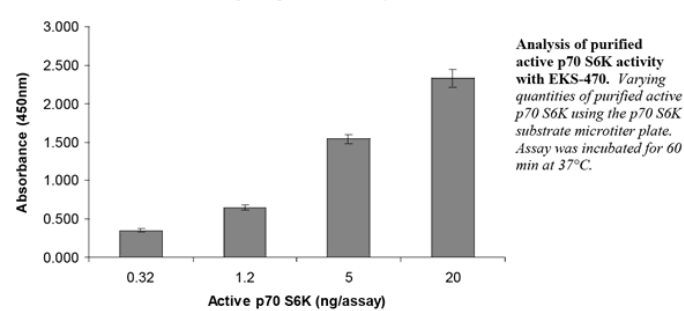
96 wells

Manuals, SDS & CofA

[View Online »](#)

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

Example of p70 S6K Activity Curve



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

Application Activity assay, Colorimetric detection

Application Notes For the measurement of p70 S6K 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 P23443 (S6K1), Q9UBS0 (S6K2)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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