

p70S6K (pT389 + Total) ELISA Kit ab225582

SimpleStep ELISA

6 Images

Overview

Product name p70S6K (pT389 + Total) ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
pT389	6			3.3%
Total	6			4.5%

Inter-assay

Sample	n	Mean	SD	CV%
pT389	3			2.9%
Total	3			4.1%

Sample type Cell Lysate, Tissue Homogenate

Assay type Semi-quantitative

Assay time 1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Mouse, Human

Predicted to work with: Rat 

Product overview Abcam's p70S6K (pT389) and p70S6K (Total) *in vitro* SimpleStep ELISA™ (Enzyme-Linked Immunosorbent Assay) kit is designed for the semi-quantitative measurement of p70S6K (pT389) and Total p70S6K protein in Human and mouse cells.

The SimpleStep ELISA™ employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound

material. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.
It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Microplate

Properties

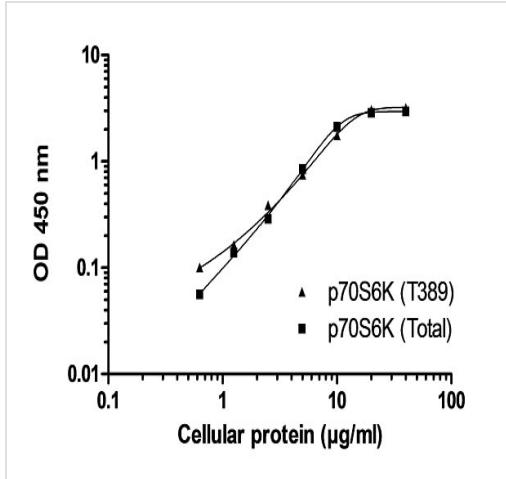
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Wash Buffer PT	1 x 15ml
50X Cell Extraction Enhancer Solution	1 x 1ml
5X Cell Extraction Buffer PTR	1 x 12ml
Lyophilized p38 MAPK alpha Control Lysate	1 vial
p70S6K (pT389) Capture Antibody	1 x 1.5ml
p70S6K (pT389) Detector Antibody	1 x 1.5ml
p70S6K (Total) Capture Antibody	1 x 1.5ml
p70S6K (Total) Detector Antibody	1 x 1.5ml
Plate Seal	1 unit
SimpleStep Pre-Coated 96 Well Microplate (12 x 8 well strips)	1 unit
Stop Solution	1 x 12ml
TMB Substrate	1 x 12ml

Cellular localization

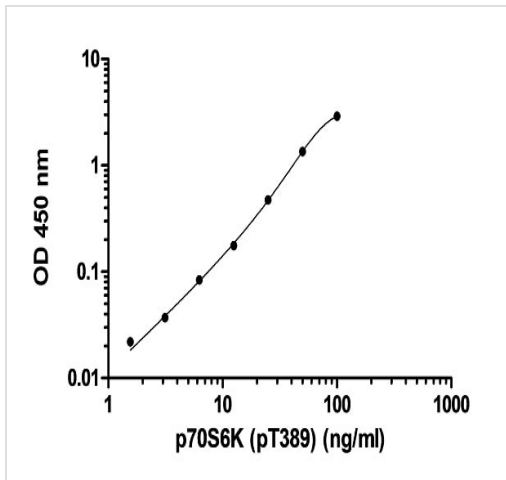
P70 S6 Kinase: Cell junction › synapse › synaptosome By similarity. Mitochondrion outer membrane. Mitochondrion. Note: Colocalizes with URI1 at mitochondrion. Isoform Alpha I: Nucleus. Cytoplasm Isoform Alpha II: Cytoplasm

Images



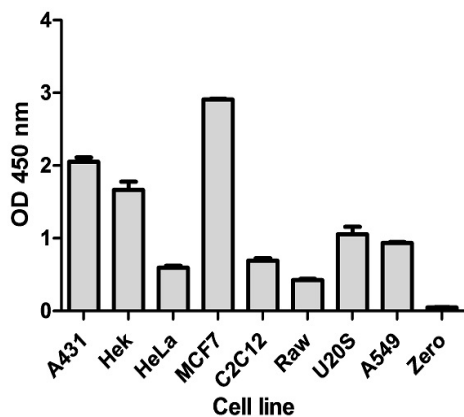
Background-subtracted data values (mean \pm SD) are graphed.

Example of p70S6K (pT389) and p70S6K (Total) cell lysate standard curve



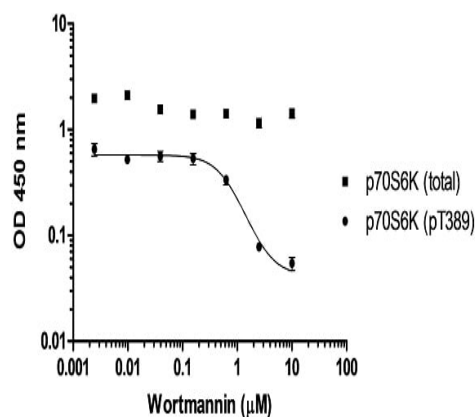
Background-subtracted data values (mean \pm SD) are graphed.

Example of p70S6K (pT389) recombinant protein standard curve



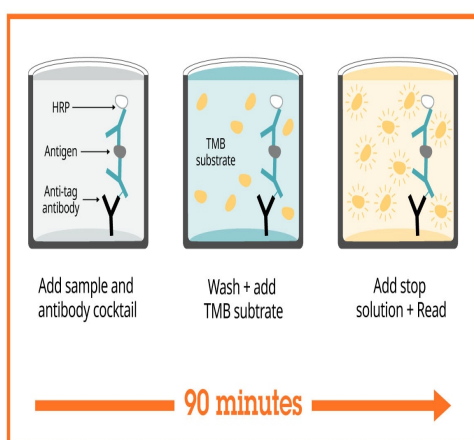
Cell line analysis for Total p70S6K from 100 µg/mL preparations of cell extracts

Data from triplicate measurements (mean \pm SD) are plotted and compared to 1X Cell Extraction Buffer PTR (zero).



Inhibition of p70S6K (pT389) phosphorylation in MCF-7 cells in response to wortmannin treatment

MCF-7 cells were cultured in 96-well tissue culture plates, and treated (2 h) with a dose-range of wortmannin before stimulation with insulin (10 µg/mL for 30 min) and cell lysis. Data from triplicate measurements of p70S6K (pT389) are plotted and compared against total p70S6K protein levels.



Sandwich ELISA - p70S6K (pT389 + Total) ELISA Kit (ab225582)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Get more done with SimpleStep ELISA

**Easy to use**

Single-wash 90-minute protocol

**Flexible**

Matched antibody pairs available

**Precision antibodies**

High sensitivity, specificity and reproducibility

**Scalable**

Now in 10-pack and 384-well formats

To learn more about the advantages of SimpleStep ELISA® kits see [here](#).

Sandwich ELISA - p70S6K (pT389 + Total) ELISA
Kit (ab225582)

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.com/abpromise> or contact our technical team.

Terms and conditions

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