

Anti-p70S6K

Rabbit Polyclonal Antibody

Catalog # R21-63BR

Lot # R039-6

Cited Applications

- Western blot (1:1000 dilution)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the p70S6K protein

Cross Reactivity

- Western blot p70S6K from human and mouse cells

p70S6K from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 409-413 of human p70 S6K (protein accession number P23443)

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

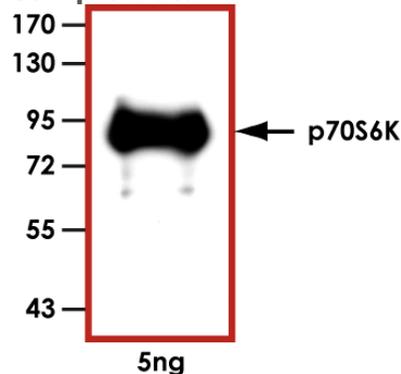
Scientific Background

p70S6K is responsible for the phosphorylation of 40S ribosomal protein S6, and is ubiquitously expressed in human adult tissues (1). p70S6K is activated by serum stimulation and this activation is inhibited by wortmannin and rapamycin. p70S6K activity changes during the cell cycle, and increases 20-fold in G1 cells released from G0 (2). p70S6K activation requires sequential phosphorylation at proline-directed residues in the putative autoinhibitory pseudosubstrate domain, as well as threonine 389, a site phosphorylated by phosphoinositide-dependent kinase 1 (PKD-1).

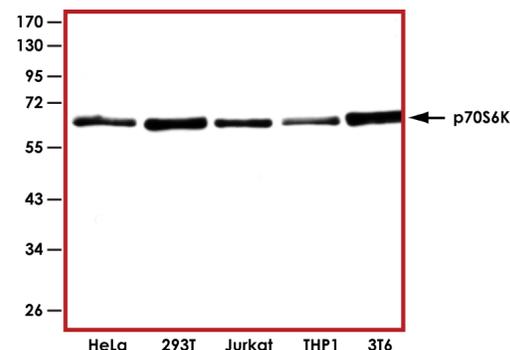
References

1. Ferrari, S. et al: S6 phosphorylation and the p70s6k/p85s6k. Crit Rev Biochem Mol Biol. 1994;29(6):385-413. Review.
2. Edelmann, HM. Et al: Cell cycle regulation of p70 S6 kinase and p42/p44 mitogen-activated protein kinases in Swiss mouse 3T3 fibroblasts. J Biol Chem. 1996 Jan 12;271(2):963-71.

Sample Data



Representative western blot with Anti-p70S6K (1:1000) using 5 ng of human recombinant p70S6K.



Representative western blot with Anti-p70S6K (1:1000) using 30 µg of HeLa, HEK-293T, Jurkat, THP1, and 3T6 cell lysate.

Anti-p70S6K

Rabbit Polyclonal Antibody

Catalog Number

R21-63BR

Specific Lot Number

R039-6

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1 yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.