

p70S6Kb, Unactive

Recombinant full-length human protein expressed in Sf9 cells

Catalog # R22-14G

Lot # Y1093-2

Product Description

Recombinant full-length human p70S6Kb was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC000094](#).

Gene Aliases

RPS6KB2, STK14B, KLS, P70-beta, P70-beta-1, P70-beta-2, S6K2, p70(S6K)-beta, S6K-beta2, SRK

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

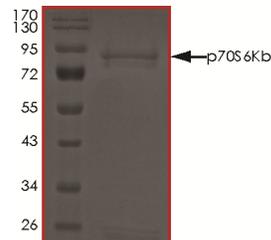
Scientific Background

p70S6Kb is a member of the RSK (ribosomal S6 kinase) family of serine/threonine kinases and is activated by mitogenic stimuli, including growth factors, cytokines, and phorbol esters. p70S6Kb contains 2 nonidentical kinase catalytic domains and phosphorylates the S6 ribosomal protein and eukaryotic translation initiation factor 4B. Phosphorylation of S6 leads to an increase in protein synthesis and cell proliferation. PI3 kinase pathway and mTOR are involved in the activation of p70S6Kb but other pathways can also activate this target protein.

References

- Gout, I. et al: Molecular cloning and characterization of a novel p70 S6 kinase beta containing a proline-rich region. *J. Biol. Chem.* 273: 30061-30064, 1998.
- Saitoh, M. et al: Cloning and characterization of p70(S6K-beta) defines a novel family of p70 S6 kinases. *Biochem. Biophys. Res. Commun.* 253: 470-476, 1998.

Purity



The purity of p70S6Kb was determined to be **>90%** by densitometry. Approx. MW **85kDa**.

p70S6Kb, Unactive

Full length human recombinant protein expressed in Sf9 cells

Catalog #	R22-14G
Lot #	Y1093-2
Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.