

STK40 (SHIK) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # S54-34G

Lot # T860-4

Product Description

Full-length recombinant human STK40 (SHIK) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_032017](#).

Gene Aliases

RP11-268J15.4; SgK495; SHIK

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

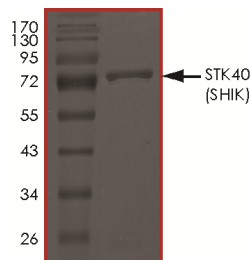
Scientific Background

STK40 is a serine/threonine-protein kinase 40 that functions in cell division and also to promote efficient cytokinesis. STK40 is also called SHIK (sink- homologous inhibitory kinase) which is expressed in all tissues. STK40 also inhibits TNF-induced NF-kappa-B activation and p53-induced transcription (1).

References

1. Huang, J.et.al: Identification of a novel serine/threonine kinase that inhibits TNF-induced NF-kappa-B activation and p53-induced transcription. *Biochem. Biophys. Res. Commun.* 309: 774-778, 2003.

Purity



The purity of STK40 (SHIK) protein was determined to be **>95%** by densitometry. Approx. MW **74 kDa**.

STK40 (SHIK) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	S54-34G
Lot #	T860-4
Purity	>95%
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.