

STK35 (CLIK1) Protein

Recombinant human protein expressed in Sf9 cells

Catalog # S47-35G

Lot # S333-3

Product Description

Recombinant human STK35 (CLIK1) (134-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_080836](#).

Gene Aliases

CLIK1; STK35L1

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.

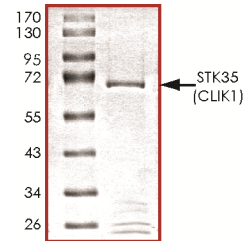
Scientific Background

STK35 is serine/threonine kinase 35 which is a kinase that is predominantly found in the nucleus and interacts with PDLIM1/CLP-36 in the cytoplasm and localize to actin stress fibers. STK35 protein is a regulator of actin stress fibers in nonmuscle cells. STK35 specifically targets STK1 to actin stress fibers and that it may represent a regulator of the actomyosin cytoskeleton in nonmuscle cells (1).

References

1. Vallenius, T. et.al: Clik1: a novel kinase targeted to actin stress fibers by the CLP-36 PDZ-LIM protein. J. Cell Sci. 115: 2067-2073, 2002.

Purity



The purity of STK35 was determined to be **>80%** by densitometry. Approx. MW **70 kDa**.

STK35 (CLIK1) Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	S47-35G
Specific Lot Number	S333-3
Purity	>80%
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.