

TSSK3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # S39-34G

Lot # N030-1

Product Description

Recombinant human TSSK3 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_052841](#).

Gene Aliases

SPOGA3; STK22C; STK22D; TSK3

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.

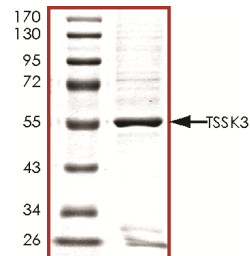
Scientific Background

TSSK3 is a serine/threonine kinase belonging to the subfamily of testis-specific kinases 3 which expressed exclusively in the testis that is thought to play a main role in either germ cell differentiation or mature sperm function (1).

References

1. Visconti, P. E. et.al: Cloning and chromosomal localization of a gene encoding a novel serine/threonine kinase belonging to the subfamily of testis-specific kinases. Genomics 77: 163-170, 2001.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **56 kDa**.

TSSK3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	S39-34G
Specific Lot Number	N030-1
Purity	>85%
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.