

TSSK6 (SSTK) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # S23-34G

Lot # Z1372-4

Product Description

Full-length recombinant human TSSK6 (SSTK) protein was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_032037](#).

Gene Aliases

CT72; FKSG82; SSTK; TSSK4; TSSK6

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.

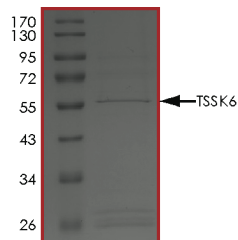
Scientific Background

TSSK6 or testis-specific serine kinase 6 is a member of the CAMK (calcium/calmodulin-dependent) serine/threonine protein kinase family which is required for postmeiotic chromatin remodeling and male fertility (1). Tssk6-null spermatozoa have impaired DNA condensation. TSSK6 is highly expressed in testis. TSSK6 is present in the equatorial segment of human sperm (2).

References

1. Spiridonov, N. et.al: Identification and characterization of SSTK, a serine/threonine protein kinase essential for male fertility. *Molec. Cell. Biol.* 25: 4250-4261, 2005.
2. Hao, Z. et.al: Expression analysis of the human testis-specific serine/threonine kinase (TSSK) homologues: a TSSK member is present in the equatorial segment of human sperm. *Molec. Hum. Reprod.* 10: 433-444, 2004.

Purity



The purity of TSSK6 (SSTK) protein was determined to be **>75%** by densitometry. Approx. MW **57 kDa**.

TSSK6 (SSTK) Protein

Full-length recombinant human protein expressed in Sf9 cells

| | |
|--------------------|--|
| Catalog # | S23-34G |
| Lot # | Z1372-4 |
| Purity | >75% |
| 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.