

STK31 Protein

Recombinant protein expressed in Sf9 cells

Catalog # S45-35G

Lot # S203-3

Product Description

Recombinant human STK31 (496-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_031414](#).

Gene Aliases

SGK396; TDRD8

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.

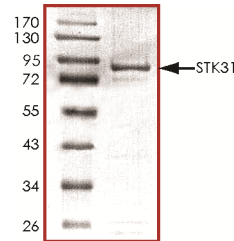
Scientific Background

STK31 is a serine/threonine protein kinase that contains a tudor domain found in RNA-interacting proteins, and a coiled-coil domain. Tudor domains are found in many eukaryotic organisms and have been implicated in protein-protein interactions in which methylated protein substrates bind to these domains. The STK31 gene is conserved across many species and was originally identified as a gene expressed in mouse spermatogonia but not in somatic tissues (1). STK31 has been shown to be specifically expressed in the human testis.

References

1. Wang, P J. et al: An abundance of X-linked genes expressed in spermatogonia. Nature Genet. 27: 422-426, 2001.

Purity



The purity of STK31 was determined to be **>80%** by densitometry. Approx. MW **86 kDa**.

STK31 Protein

Recombinant protein expressed in Sf9 cells

Catalog Number S45-35G

Specific Lot Number S203-3

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.