

TESK1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # T05-34G

Lot # L1894-4

Product Description

Recombinant full-length human TESK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_006285](#).

Gene Aliases

(None)

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.

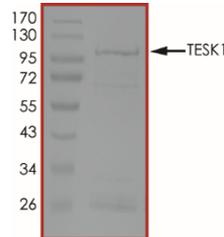
Scientific Background

TESK1 or testis-specific kinase 1 is a serine/threonine protein kinase that contains an N-terminal protein kinase domain and a C-terminal proline-rich domain which is most closely related to those of the LIM motif-containing protein kinases (LIMKs). TESK1 protein can phosphorylate myelin basic protein and histone in vitro and plays an important role at and after the meiotic phase of spermatogenesis. TESK1 is mainly expressed in testicular germ cells (1).

References

1. Toshima, J. et.al: Identification and characterization of a novel protein kinase, TESK1, specifically expressed in testicular germ cells. J. Biol. Chem. 270: 31331-31337, 1995.

Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **106 kDa**.

TESK1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	T05-34G
Lot #	L1894-4
Purity	>70%
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.