

PSKH1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P89-34G

Lot # X574-3

Product Description

Recombinant human PSKH1 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PSKH1 gene accession number is [NM_006742](#).

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.

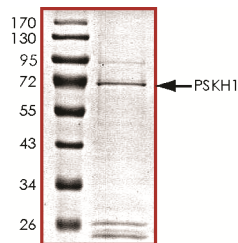
Scientific Background

PSKH1 is a member of the family of protein serine kinase H1 which was originally cloned by homology probing (1). PSKH1 has been localized to the endoplasmic reticulum (ER), the Golgi apparatus, and the plasma membrane (PM). PSKH1 has been shown to play a structural and regulatory role in maintenance of the Golgi apparatus, which is a key organelle within the secretory pathway. PSKH1 is in a cluster of 5 unrelated genes on 16q22.1 chromosome which is transcribed in the opposite direction from that of LCAT which is a member of the same cluster (2).

References

- Hanks, S. K. et.al: Homology probing: identification of cDNA clones encoding members of the protein-serine kinase family. Proc. Nat. Acad. Sci. 84: 388-392, 1987.
- Larsen, F. et.al: A tight cluster of five unrelated human genes on chromosome 16q22.1. Hum. Molec. Genet. 2: 1589-1595, 1993.

Purity



The purity of PSKH1 was determined to be **>70%** by densitometry, approx. MW **72 kDa**.

PSKH1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	P89-34G
Specific Lot Number	X574-3
Purity	>70%
Concentration	0.05 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.