

SCYL3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # S09-30G

Lot # J691-1

Product Description

Recombinant human SCYL3 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC014662](#).

Gene Aliases

PACE-1; PACE1; RP1-97P20.2

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.

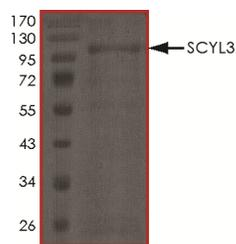
Scientific Background

SCYL3 encodes a protein with a kinase domain and four HEAT repeats which interact with the C-terminal domain of ezrin, an ERM protein which play a role in cell adhesion and migration (1). C-terminal domain of SCYL3 interacts with the C-terminal alpha-helical domain of ezrin. Biochemical assays indicate that SCYL3 had associated rather than intrinsic protein kinase activity. SCYL3 is associated with the cytoplasmic face of the Golgi apparatus. This distribution is dependent upon the presence of the SCYL3 N-terminal myristoylation consensus sequence but is not dependent on an association with ezrin. SCYL3 is highly expressed in breast carcinoma cells and adult colon.

References

- Sullivan, A. et.al: PACE-1, a novel protein that interacts with the C-terminal domain of ezrin. *Exp. Cell Res.* 284: 224-238, 2003.

Purity



The purity of SCYL3 was determined to be **>85%** by densitometry. Approx. MW **110 kDa**.

SCYL3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	S09-30G
Specific Lot Number	J691-1
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.