

SCYL2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # S03-35G

Lot # A1243-4

Product Description

Recombinant human SCYL2 (1-520) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_017988](#).

Gene Aliases

CVAK104, KIAA1360

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.

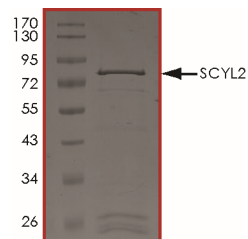
Scientific Background

SCYL2 or SCY1-like 2 (*S. cerevisiae*) is an evolutionarily conserved protein involved in SNARE sorting that is essential for normal embryonic development (1). SCYL2 serves as a regulatory factor for Vpu (viral protein U) which reducing the extent of Vpu phosphorylation and consequently enhancing BST2-mediated viral restriction (2).

References

1. Borner GH. et.al: CVAK104 is a novel regulator of clathrin-mediated SNARE sorting. *Traffic*. 2007 Jul;8(7):893-903.
2. Miyakawa K. et.al: Interferon-induced SCYL2 limits release of HIV-1 by triggering PP2A-mediated dephosphorylation of the viral protein Vpu. *Sci Signal*. 2012 Oct 9;5(245):ra73. doi: 10.1126/scisignal.2003212.

Purity



The purity of SCYL2 protein was determined to be **>85%** by densitometry. Approx. MW **84 kDa**.

SCYL2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	S03-35G
Lot #	A1243-4
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.