

SCYL1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # S04-35G

Lot # F528-3

Product Description

Recombinant human SCYL1 (1-556) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The SCYL1 gene accession number is [NM_020680](#).

Gene Aliases

GKLP; HT019; NKTL; NTKL; P105; TAPK; TEIF; TRAP

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

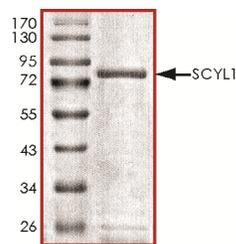
Scientific Background

SCYL1 is a transcriptional regulator belonging to the SCYL1-like family of kinase-like proteins. SCYL1 has a divergent N-terminal kinase domain that is thought to be catalytically inactive, and can bind specific DNA sequences through its C-terminal domain. SCYL1 activates transcription of the telomerase reverse transcriptase and DNA polymerase beta genes. SCYL1 forms multimers following transfection into COS-7 cells (1). A SCYL1 binding protein has been identified that co-localized with SCYL1 in cytoplasm and shows ubiquitous expression (2).

References

1. Kato, M. et.al: Identification and characterization of the human protein kinase-like gene NTKL: mitosis-specific centrosomal localization of an alternatively spliced isoform. *Genomics* 79: 760-767, 2002.
2. Di, Y. et.al: Cloning and characterization of a novel gene which encodes a protein interacting with the mitosis-associated kinase-like protein NTKL. *J. Hum. Genet.* 48: 315-321, 2003.

Purity



The purity of SCYL1 was determined to be **>95%** by densitometry. Approx. MW **85 kDa**.

SCYL1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number S04-35G

Specific Lot Number F528-3

Purity >95%

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.