

STK32A (YANK1) Protein

Full-length recombinant mouse protein expressed in Sf9 cells

Catalog # Y03-34G

Lot # Z1298-3

Product Description

Full-length recombinant mouse STK32A (YANK1) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_178749](#).

Gene Aliases

YANK1, STK32A, A930015B13Rik

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.

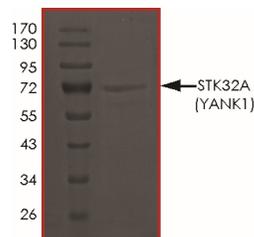
Scientific Background

STK32A or serine/threonine kinase 32A (also known as YANK1) is a highly conserved gene in many species (1). STK32A is a member of the YANK family that also contains STK32B and STK32C. STK32A contains a catalytic domain and metal ion binding domain and is involved in the phosphorylation of serine and threonine residues on various protein substrates. A number of sequence variants have been identified that are shorter with distinct C-terminus (2). The function of this protein is not yet clear.

References

1. Manning G1, et al: The protein kinase complement of the human genome. *Science*. 2002 Dec 6;298(5600): 1912-34.
2. Strausberg RL et al. Generation and initial analysis of more than 15,000 full-length human and mouse cDNA sequences. *Proc. Natl. Acad. Sci. U.S.A.* 2002 Dec;99(26): 16899-16903.

Purity



The purity of STK32A (YANK1) protein was determined to be **>95%** by densitometry. Approx. MW **71 kDa**.

STK32A (YANK1) Protein

Full-length recombinant mouse protein expressed in Sf9 cells

Catalog #	Y03-34G
Lot #	Z1298-3
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.