

NCK Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # N06-30H

Lot # M094-1

Product Description

Recombinant full-length human NCK was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_006153](#).

Gene Aliases

NCK1, MGC12668, NCKalpha

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 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.

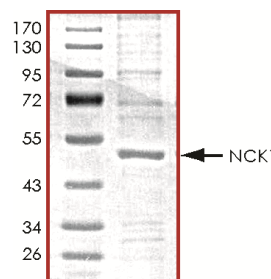
Scientific Background

NCK (or non-catalytic region of tyrosine kinase adaptor protein 1) is an adaptor protein that is located in the cytoplasm and involved in transducing signals from receptor tyrosine kinases to downstream signal recipients such as RAS (1). NCK is a signaling and transforming protein that contains Src homology 2 and 3 (SH2 and SH3) domains. NCK functions by coupling tyrosine phosphorylation via SH2 domains to downstream effectors through SH3 domains. Furthermore, NCK couples tyrosine phosphorylation guidance signals to cytoskeletal events required for the ipsilateral projections of spinal cord neurons and for normal limb movement (2).

References

1. Park, D. et al: Phosphorylation of Nck in response to a variety of receptors, phorbol myristate acetate, and cyclic AMP. Mol. Cell. Biol. 1992; 12 (12): 5816-23.
2. Fawcett, J. P. et al: Nck adaptor proteins control the organization of neuronal circuits important for walking. Proc. Nat. Acad. Sci. 104: 20973-20978, 2007

Purity



The purity was determined to be **>75%** by densitometry. Approx. MW **48kDa**.

NCK Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	N06-30H
Specific Lot Number	M094-1
Purity	>75%
Concentration	0.2 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.