

Stathmin Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # S71-30H

Lot # R111-2

Product Description

Recombinant full-length human Stathmin protein was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [BC014353](#).

Gene Aliases

STMN1; C1orf215; FLJ32206; Lag; LAP18; MGC138869; MGC138870; OP18; PP17; PP19; PR22; SMN

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM 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.

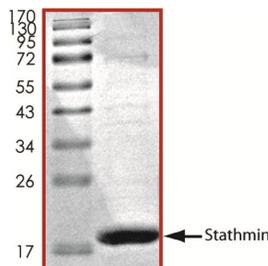
Scientific Background

Stathmin is a ubiquitous, highly conserved protein present in the cytoplasm in a variety of unphosphorylated and phosphorylated forms. Stathmin functions as an intracellular relay integrating regulatory signals of the cellular environment (1). Stathmin is involved in the regulation of the microtubule filament system. Stathmin binds tubulin in a ternary complex, with 2 tubulins for every stathmin and this complex interferes with microtubule dynamics in vitro and in vivo. Overexpression of several stathmin mutants suggest that stathmin induces depolymerization of interphase and mitotic microtubules in its unphosphorylated state, but is inactivated by phosphorylation in mitosis (2).

References

1. Sobel, A. Stathmin: a relay phosphoprotein for multiple signal transduction? Trends Biochem. Sci. 16: 301-305, 1991.
2. Gavet, O. et al: The stathmin phosphoprotein family: intracellular localization and effects on the microtubule network. J. Cell Sci. 111: 3333-3346, 1998.

Purity



The purity of Stathmin was determined to be **>90%** by densitometry. Approx. MW **19kDa**.

Stathmin Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	S71-30H
Specific Lot Number	R111-2
Purity	>90%
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.