

14-3-30 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # Y84-30G

Lot # L326-1

Product Description

Recombinant full-length human 14-3-30 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_006826](#).

Gene Aliases

YWHAQ, 1C5, HS1, 14-3-3

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

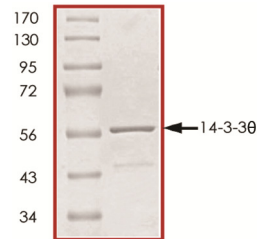
14-3-30 (also known as tyrosine 3-monooxygenase/tryptophan 5-monooxygenase activation protein, theta polypeptide) is a member of the 14-3-3 family of proteins which mediate signal transduction by binding to phosphoserine-containing proteins. Through interaction with ASK1, c-jun NH-terminal kinase, and p38 mitogen-activated protein kinase (MAPK), 14-3-30 plays an important role in controlling apoptosis (1). Induced expression of 14-3-30 protein has been reported in patients with amyotrophic lateral sclerosis. Additionally, 14-3-30 has been observed to mediate nucleocytoplasmic shuttling of the N protein (coronavirus nucleocapsid protein) which causes severe acute respiratory syndrome (2).

References

1. Lau, J.M. et al: The 14-3-3tau phosphoserine-binding protein is required for cardiomyocyte survival. *Mol Cell Biol.* 2007, 27(4):1455-66.
2. The severe acute respiratory syndrome coronavirus nucleocapsid protein is phosphorylated and localizes in the cytoplasm by 14-3-3-mediated translocation. *J Virol.* 2005 Sep;79(17):11476-86.

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

Purity



The purity of 14-3-30 was determined to be **>95%** by densitometry. Approx. MW **56kDa**.

14-3-30 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number Y84-30G

Specific Lot Number L326-1

Purity	>95%
Concentration	0.2 µg/µl
Stability	1 yr 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.