

14-3-3θ Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # Y84-30N

Lot # H2631-7

Product Description

Recombinant full-length human tag-free 14-3-3θ was expressed in E. coli cells. The gene accession number is [NM_006826](#).

Gene Aliases

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

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 0.25mM DTT, 0.1mM PMSF, and 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.

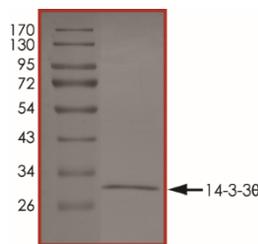
Scientific Background

14-3-3θ (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), 14-3-3θ plays an important role in controlling apoptosis (1). Induced expression of 14-3-3θ protein has been reported in patients with amyotrophic lateral sclerosis. Additionally, 14-3-3θ has been observed to mediate nucleocytoplasmic shuttling of the N protein (coronavirus nucleocapsid protein) which causes severe acute respiratory syndrome (2).

References

- 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.
- 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.

Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **29kDa**.

14-3-3θ Protein

Full-length recombinant protein expressed in E. coli cells

Catalog #	Y84-30N
Lot #	H2631-7
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.