

14-3-3ζ Protein

Full-length recombinant protein expressed in sf9 insect cells

Catalog # Y92-30H

Lot # Q2428-2

Product Description

Recombinant full-length human 14-3-3 zeta was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_003406](#).

Gene Aliases

14-3-3 zeta, YWHAZ, KCIP-1, MGC111427, MGC126532, MGC138156

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

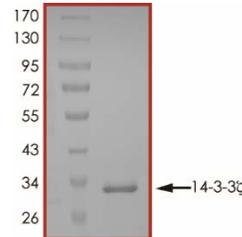
14-3-3ζ (also known as tyrosine 3-monooxygenase/tryptophan 5-monooxygenase activation protein, zeta polypeptide) is a member of the 14-3-3 family of proteins which mediate signal transduction by binding to phosphoserine-containing proteins. 14-3-3ζ protein plays a key role in cancer biology by being an important regulator of major cellular processes such as proliferation, differentiation, senescence and apoptosis (1). 14-3-3ζ protein has been shown to interact with the IRS1 protein, suggesting a role for this protein in regulating insulin sensitivity by interrupting the association between the insulin receptor and IRS1 (2).

References

- Li, Z. et al: Down-regulation of 14-3-3zeta suppresses anchorage-independent growth of lung cancer cells through anoikis activation. *Proc Natl Acad Sci U S A*. 2008; 105(1):162-7.
- Niemantsverdriet, M. et al: Cellular functions of 14-3-3 zeta in apoptosis and cell adhesion emphasize its oncogenic character. *Oncogene*. 2008;27(9):1315-9. 2007.

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 was determined to be >95% by densitometry. Approx. MW 31kDa.

14-3-3ζ Protein

Full-length recombinant protein expressed in sf9 insect cells

Catalog #	Y92-30H
Lot #	Q2428-2
Purity	>95%
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.