

Catenin β Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog # C06-30G

Lot # L2003-4

Product Description

Recombinant full-length human Catenin β was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001904](#).

Gene Aliases

CTNNB, FLJ25606

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.

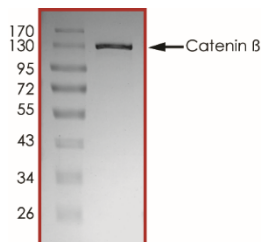
Scientific Background

β-catenins are cytoplasmic proteins that are ubiquitously expressed. These proteins associate with E-cadherin at cellular junctions (1). β-catenin interacts with TCF and LEF transcription factors and is an essential member of the Wntless-Wnt signal transduction pathway. The adenomatous polyposis coli (APC) tumor-suppressor protein, together with Axin and GSK3β, form a Wnt-regulated signaling complex that mediates phosphorylation-dependent degradation of β-catenin by the proteasome. APC and Siah-1 mediate a novel β-catenin degradation pathway linking p53 activation to cell cycle control. Activating mutations in the human β-catenin gene have been found in human colon cancer and melanomas (2).

References

1. Morin, P.J. et al: Activation of beta-catenin-Tcf signaling in colon cancer by mutations in beta-catenin or APC. *Science* 1997; 275(5307):1787-1790.
2. Liu, J. et al: Siah-1 mediates a novel beta-catenin degradation pathway linking p53 to the adenomatous polyposis coli protein. *Mol. Cell.* 2001;7(5):927-36.

Purity



The purity of Catenin β was determined to be **>95%** by densitometry.
Approx. MW **115kDa**.

Catenin β

Full-length recombinant protein expressed in Sf9 cells

Catalog #	C06-30G
Lot #	L2003-4
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.

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.