

## Recombinant beta-Catenin protein

**Catalog No:** 82067, 82667

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Source:** Human

**Buffer Contents:** Recombinant beta-Catenin protein is supplied in 25 mM HEPES pH 7.5, 300 mM NaCl, 20% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

**Background:** **beta-Catenin** is a key downstream component of the canonical Wnt signaling pathway. In the absence of Wnt, forms a complex with AXIN1, AXIN2, APC, CSNK1A1 and GSK3B that promotes phosphorylation on N-terminal Ser and Thr residues and ubiquitination of CTNNB1 via BTRC and its subsequent degradation by the proteasome. In the presence of Wnt ligand, CTNNB1 is not ubiquitinated and accumulates in the nucleus, where it acts as a coactivator for transcription factors of the TCF/LEF family, leading to activate Wnt responsive genes.

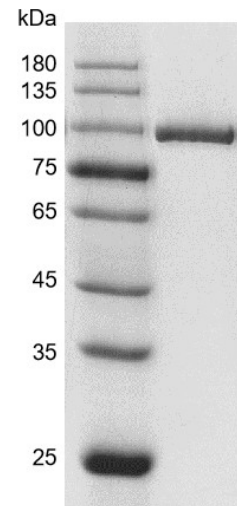
**Protein Details:** Recombinant beta-Catenin protein was expressed in a baculovirus expression system as the full-length protein (accession number NP\_001091679.1) with an N-terminal FLAG tag. The molecular weight of beta-Catenin is 86.79 kDa.

**Application Notes:** This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

### beta-Catenin



### Recombinant beta-Catenin protein

10% SDS-PAGE Coomassie staining

MW: 86.79 kDa,

Purity: ≥ 95%