

# PRODUCT INFORMATION



## β-Catenin (human, recombinant)

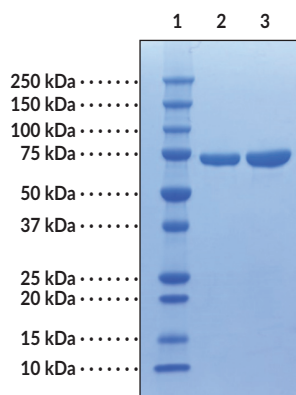
Item No. 28780

### Overview and Properties

|                               |   |
|-------------------------------|---|
| <b>Synonyms:</b>              | Armadillo, Catenin (Cadherin-Associated Protein), β1, 88kDa, Catenin β-1, CTNNB1, EVR7, MRD19, NEDSDV |
| <b>Source:</b>                | Recombinant N-terminal His-TEV-tagged human β-catenin (138-781) purified from <i>E. coli</i>          |
| <b>Amino Acids:</b>           | 138-781   |
| <b>Uniprot No.:</b>           | P35222  |
| <b>Molecular Weight:</b>      | 72.4 kDa  |
| <b>Storage:</b>               | -80°C (as supplied)   |
| <b>Stability:</b>             | ≥1 year   |
| <b>Purity:</b>                | <i>batch specific</i> (≥90% estimated by SDS-PAGE)  |
| <b>Supplied in:</b>           | PBS, pH 7.4, with 1 mM TCEP and 10% glycerol  |
| <b>Protein Concentration:</b> | <i>batch specific</i> mg/ml   |

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: β-Catenin (human, recombinant) (2 μg)  
Lane 3: β-Catenin (human, recombinant) (4 μg)

Representative gel image shown; actual purity may vary between each batch.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/18/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

$\beta$ -Catenin is a transcriptional coactivator that is encoded by the *CTNNB1* gene in humans.<sup>1,2</sup> It is a 781-amino acid protein comprised of an N-terminal domain containing glycogen synthase kinase 3 $\beta$  (GSK3 $\beta$ ) phosphorylation sites, a C-terminal transactivation domain, and a central domain spanning amino acid residues 138-664.<sup>3,4</sup> The central domain consists of 12 armadillo repeats and is required for binding to cadherins, TCF/LEF transcription factors, and adenomatous polyposis coli (APC).  $\beta$ -Catenin has roles in cell adhesion, canonical Wnt signaling, regulation of stem cells, embryonic development, and adult tissue homeostasis, among others.<sup>1,3</sup> In the absence of Wnt, a complex consisting of axin, APC, GSK3 $\beta$ , and casein kinase 1 (CK1), binds to and phosphorylates  $\beta$ -catenin, targeting it for ubiquitination and proteasomal degradation.<sup>1</sup> In the presence of Wnt, phosphorylation of  $\beta$ -catenin is inhibited, allowing  $\beta$ -catenin to translocate into the nucleus, where it interacts with TCF/LEF to activate expression of Wnt target genes. Activating mutations in *CTNNB1* that stabilize  $\beta$ -catenin have been associated with a variety of cancers, including hepatocellular and adrenocortical carcinomas, colorectal cancer, and pilomatricomas.<sup>4-7</sup> Cayman's  $\beta$ -Catenin (human, recombinant) protein can be used for Western blot and ELISA applications.

## References

---

1. MacDonald, B.T., Tamai, K., and He, X. Wnt/ $\beta$ -catenin Signaling: Components, mechanisms, and diseases. *Dev. Cell.* **17(1)**, 9-26 (2009).
2. Nollet, F., Berx, G., Molemans, F., *et al.* Genomic organization of the human  $\beta$ -catenin gene (*CTNNB1*). *Genomics* **32**, 413-424 (1996).
3. Xing, Y., Takemaru, K.-I., Liu, J., *et al.* Crystal structure of a full-length  $\beta$ -catenin. *Structure* **16(3)**, 478-487 (2008).
4. Akiyama, T. Wnt/ $\beta$ -catenin signaling. *Cytokine Growth Factor Rev.* **11**, 273-282 (2000).
5. de La Coste, A., Romagnolo, B., Billuart, P., *et al.* Somatic mutations of the  $\beta$ -catenin gene are frequent in mouse and human hepatocellular carcinomas. *Proc. Natl. Acad. Sci. USA* **95(15)**, 8847-8851 (1998).
6. Clevers, H. Wnt/ $\beta$ -catenin signaling in development and disease. *Cell* **127**, 469-480 (2006).
7. Durand, J., Lampron, A., Mazzuco, T.L., *et al.* Characterization of differential gene expression in adrenocortical tumors harboring  $\beta$ -catenin (*CTNNB1*) mutations. *J. Clin. Endocrinol. Metab.* **96(7)**, E1206-E1211 (2011).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM