

# β-Catenin monoclonal antibody (12F7)

β-Catenin, a vertebrate homologue of *Drosophila armadillo* encoded by the CTNNB1 gene, performs two functions in the cell: mediating cell-cell adhesion by association between E-cadherin and the cytoskeleton, and acting as a downstream transcriptional activator of Wnt signaling, a key player in such functions as embryonic development, the induction of cell polarity, and the determination of cell fate. In the absence of Wnt signaling, glycogen synthase kinase-3β (GSK-3β) phosphorylates any free β-catenin at the amino terminus. This results in adenomatous polyposis coli tumor suppressor protein (APC)-mediated β-catenin degradation via the ubiquitin-proteasome pathway. Wnt signaling stabilizes β-catenin by inhibiting the activity of GSK-3β and preventing β-catenin degradation. β-catenin forms complexes with DNA-binding proteins such as T-cell factor (Tcf) and lymphocyte-enhancer factor (Lef), which promote expression of a variety of target genes, including c-myc and cyclin D1.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

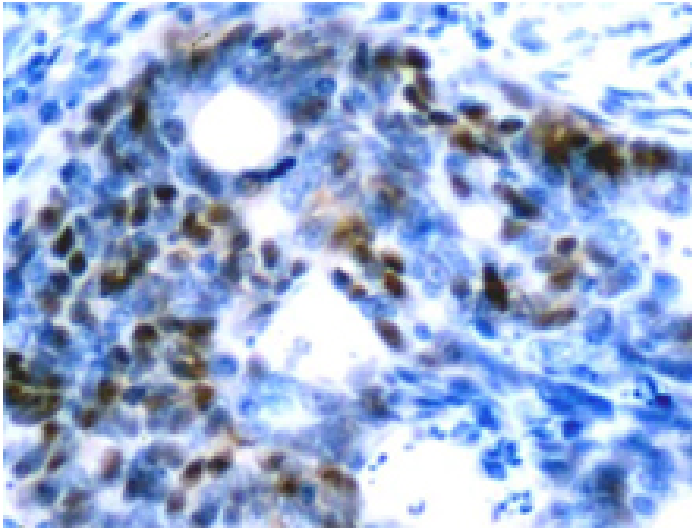
## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ADI-KAM-ST001-E | 100μg |
|-----------------|-------|

Manuals, SDS & CofA

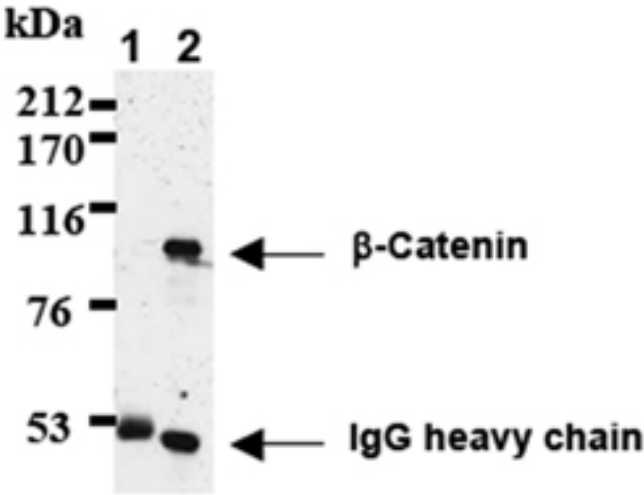
[View Online »](#)



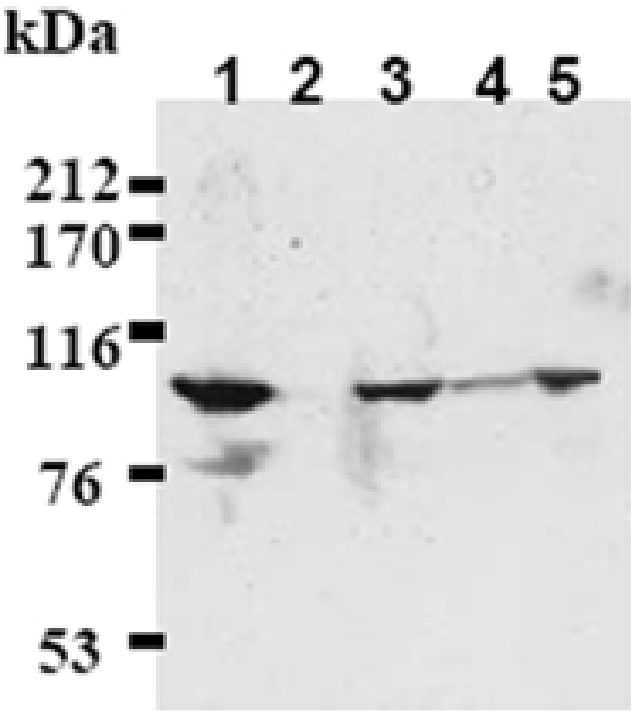
Immunohistochemical detection of colorectal cancer  $\beta$ -catenin on paraffin embedded sections with  $\beta$ -catenin mAb (12F7).



Immunohistochemical detection of gastric epithelium  $\beta$ -catenin on paraffin embedded sections with  $\beta$ -Catenin mAb (12F7).



Immunoprecipitation of Rat-1 cells with normal mouse IgG (1) or  $\beta$ -catenin mAb (12F7) (2). After immunoprecipitated with the antibody, immunocomplex was resolved on SDS-PAGE and immunoblotted with  $\beta$ -catenin mAb (12F7).



Western blot analysis of A431 cells (1), Hela cells (2), NIH/3T3 cells (3), C2C12 cells (4) and Rat-1 cells (5) probed with  $\beta$ -catenin mAb (12F7)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name CTNNB

Application IHC, IP, WB

Application Notes Detects bands of ~97kDa by Western blot.

Clone 12F7

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol.

Host Mouse

Immunogen Recombinant  $\beta$ -catenin (N-terminal fragment).

Isotype IgG1

Purity Detail Protein A affinity purified.

**Recommendation**  
**Dilutions/Conditions** Immunohistochemistry (1-5 $\mu$ g/ml)Immunoprecipitation (2 $\mu$ g/200-300 $\mu$ l cell extract)Western Blot (1 $\mu$ g/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P35222 (human), Q02248 (mouse), Q9WU82 (rat)

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)