

Product Information



β-Catenin Polyclonal Antibody

Item No. 100029 • Lot No. XXXX

- Contents:** This vial contains (100-500 μg of peptide affinity-purified IgG, *lot specific*) in 500 μl TBS, pH 7.4, containing *lot specific* BSA and *lot specific* sodium azide.
- Host:** Rabbit
- Antigen:** Human β-catenin amino acids 43-62. The sequence is 100% identical between human, mouse, and rat. Antigen alignment with other known species sequences:
Human AP S L S G K G N P E E E D V D T S Q V
Pig AP S L S G K G N P E E E D V D T t Q V
Zebrafish AP S L S G K G N P E D D D V D N - Q V
- Cross Reactivity:** (+) Human, mouse, rat, porcine, and bovine β-catenin; other species not tested
- Storage:** ≥6 months at 4°C
- Applications:** Recommended starting dilutions: Western blotting: *lot specific* μg/ml; immunohistochemistry of formalin-fixed paraffin-embedded sections: 1-2 μg/ml; other applications were not attempted and therefore optimal working dilutions should be determined empirically.
- Concentration:** Varies by lot, from 0.2-1.0 mg/ml (100-500 μg/vial). Always 100 ml final working volume for western blotting.

β-Catenin is a multifunctional protein known to be part of the Wnt pathway playing essential roles in development and carcinogenesis.^{1,3} It can act as a regulator of the cell cycle and apoptosis in a variety of different cell systems.^{4,5} The entire primary amino acid sequence is identical among human, rat, and mouse, indicating critical functions of this protein.⁶⁻⁸ Overexpression of β-catenin causes enlarged forebrains and cytoarchitectural distortions in mice.⁹ The reported molecular weight for this protein ranges from 86-97 kDa.^{6,10}

References

1. Yano, H., Hara, A., Takenaka, K., *et al. Neurol. Res.* **22**, 650-656 (2000).
2. Haq, S., Michael, A., Andreucci, M., *et al. Proc. Natl. Acad. Sci. USA* **100(8)**, 4610-4615 (2003).
3. Yasui, K., Li, G., Wang, Y., *et al. Development Growth Differ.* **44**, 467-475 (2002).
4. Olmeda, D., Castel, S., Vilaró, S., *et al. Molecular Biology of the Cell* **14**, 2844-2860 (2003).
5. Jaiswell, A.S., Marlow, B.P., Gupta, N., *et al. Oncogene* **21**, 8414-8427 (2002).
6. Aberle, H., Butz, S., Stappert, J., *et al. J. Cell Sci.* **107**, 3655-3663 (1994).
7. Chung, S.S.W., Lee, W.M., and Cheng, C.Y. *J. Cell. Physiol.* **181**, 258-272 (1999).
8. Nollet, F., Berx, G., Molemans, F., *et al. Genomics* **32**, 413-424 (1996).
9. Chenn, A. and Walsh, C.A. *Cerebral Cortex* **13**, 599-606 (2003).
10. Su, L.-K., Vogelstein, B., and Kinzler, K.W. *Science* **262**, 1734-1737 (1993).



Lane 1: JGC1 cell lysate (30 μg)
Lane 2: RAW 264.7 cell lysate (45 μg)

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/100029

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/20/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com