

Product Information



β -Catenin (Phospho-Ser^{33,37}) Polyclonal Antibody

Catalog No. 10009180

Supplied as:	Affinity-purified IgG
Host:	Rabbit
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser ^{33,37} of human β -catenin
Cross Reactivity:	(+) Human β -catenin
Stability:	≥ 1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

β -Catenin is a central component of the cadherin cell adhesion complex and plays an essential role in neural development in the Wntless/Wnt signaling pathway.^{1,2} The role of β -catenin in these processes is thought to be regulated by the sequential phosphorylation of Ser²⁹, Ser³³, Ser³⁷, and Thr⁴¹ by glycogen synthase kinase 3 β (GSK3 β).³ This hyperphosphorylation promotes the ubiquitylation and targeted destruction of β -catenin. Mutations in components of this phosphorylation-regulated process that prevent β -catenin hyperphosphorylation by GSK3 β are strongly associated with cancers.²⁻⁴

References

1. Ding, Y. and Dale, T. Wnt signal transduction: Kinase cogs in a nano-machine? *Trends Biochem. Sci.* **27**, 327-329 (2002).
2. Polakis, P. Wnt signaling and cancer. *Genes Dev.* **14**, 1837-1851 (2000).
3. Liu, C., Li, Y., Semenov, M., *et al.* Control of β -catenin phosphorylation/degradation by a dual-kinase mechanism. *Cell* **108**, 837-847 (2002).
4. Wang, Z.H., Vogelstein, B., and Kinzler, K.W. Phosphorylation of β -catenin at S33, S37, or T41 can occur in the absence of phosphorylation at T45 in colon cancer cells. *Cancer Res.* **63**, 5234-5235 (2003).

Related Products

β -Catenin Polyclonal Antibody - Cat. No. 100029 • 5-Lipoxygenase (Phospho-Ser⁵²³) Polyclonal Antibody - Cat. No. 10007820 • p38 MAPK (Phospho-Thr¹⁸⁰/Tyr¹⁸²) Polyclonal Antibody - Cat. No. 10009177 • MEK 1/2 (Phospho-Ser^{218,222}) Polyclonal Antibody - Cat. No. 10009178 • ERK/MAPK (Phospho-Thr²⁰²/Tyr²⁰⁴) Polyclonal Antibody - Cat. No. 10009179 • CREB (Phospho-Ser¹³³) Polyclonal Antibody - Cat. No. 10009181

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, 06/03/2010

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