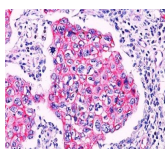


Beta-Catenin Polyclonal Rabbit Antibody (PAb)



Product Description

Beta-catenin associates with the cytoplasmic portion of E-cadherin. The catenin/cadherin complexes play an important role mediating cellular adhesion, including adherens junctions. Beta-catenin is involved in the Wnt signaling pathway as well as other signaling pathways, and is also found in complexes with many different proteins including the tumor suppressor protein APC. Defects in beta-catenin are associated with colorectal cancer, as well as many other cancer types. Beta-catenin is normally localized to the cell membrane, but can translocate to the nucleus in response to certain cell signaling.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0376, Anti-Beta-Catenin (PAb)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#0376
Antibody reactivity (target)	Beta-catenin, Catenin, Beta
Antibody type	Primary
Host species	Rabbit
Clonality	Polyclonal
Clone	PAb
Isotype	IgG, kappa
Molecular weight	92 kDa
Synonyms	Beta-catenin; Catenin beta-1; Catenin (Cadherin associated protein), beta 1; CTNNB
Human gene symbol	CTNNB1
Entrez gene ID	1499
SwissProt	P35222
Unigene	476018
Immunogen	A synthetic peptide from the middle of beta-Catenin (p120) protein
Antibody target cellular localization	Plasma membrane, Nucleus
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified), WB (verified)
Species reactivity	Human, Mouse
Antibody application notes	Immunohistology formalin-fixed 1-2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	HeLa or MCF-7 cells. Breast carcinoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Cell/tissue expression	Epithelial cells
Antibody research areas	Cancer, Cell adhesion, Developmental biology, Signal transduction
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Tumor expression	Breast cancer

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on January 27, 2026 at 09:57:48 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-beta-catenin-p120-rabbit-pab/>