Beta-Catenin Recombinant Monoclonal Rabbit Antibody (CTNNB1/2030R)



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 2030, Anti-Beta-Catenin (CTNNB1/2030R)

Call us: 800-304-5357

Product attributes				
Antibody number	#2030			
Antibody reactivity (target)	Beta-catenin, Catenin, Beta			
Antibody type	Primary			
Host species	Rabbit			
Clonality	Recombinant Monoclonal			
Clone	CTNNB1/2030R			
Isotype	IgG			
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	Recombinant full-length human β-catenin protein			
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)			
Antibody target cellular localization	Plasma membrane, Nucleus			
Species reactivity	Human			
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissues sections in 10 mM cirate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Western blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user			
	blotting 1-2 ug/mL, Optimal dilution for a specific application			
Positive control	blotting 1-2 ug/mL, Optimal dilution for a specific application			
Positive control Shipping condition	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user			
	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma			
Shipping condition	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when			
Shipping condition Storage Conditions Shelf life	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% 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: 0.2 mg/mL in PBS/0.05% azide, Purified: 0.2 mg/mL in PBS/0.0			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% 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			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Validated in protein array	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) 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 MSA-free: 1 mg/mL in PBS without azide Monospecific Cancer, Cell adhesion, Developmental biology, Signal			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Validated in protein array Antibody research areas	blotting 1-2 ug/mL, Optimal dilution for a specific application should be determined by user HeLa or MCF-7 cells. Breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) 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 Monospecific Cancer, Cell adhesion, Developmental biology, Signal transduction Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in			

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,	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 November 30, 2025 at 04:25:23 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/catenin-beta-recombinant-monoclonal-rabbit-antibody-ctnnb1-2030r/