Gamma Catenin Recombinant Monoclonal Mouse Antibody (rCTNG/1664)



Product Description

Gamma catenin, also known as JUP or junction plakoglobin, is a member of the catenin family. The catenins (alpha, beta, gamma, and delta) are ubiquitously expressed, cytoplasmic proteins that associate with E-cadherin at cellular junctions. Catenin/cadherin complexes play an important role in mediating cellular adhesion, including adherens junctions.

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 2156, Catenin, Gamma Recombinant Monoclonal Mouse Antibody (rCTNG/1664)

Call us : <u>800-304-5357</u>

Product attributes					
Antibody number	#2156				
Antibody reactivity (target)	Gamma-catenin				
Antibody type	Primary				
Host species	Mouse				
Clonality	Recombinant Monoclonal				
Clone	rCTNG/1664				
Isotype	IgG1, kappa				
Molecular weight	80-87 kDa				
Synonyms	Gamma catenin; JUP, Junction plakoglobin; Catenin gamma				
Human gene symbol	JUP				
Entrez gene ID	3728				
SwissProt	P14923				
Unigene	514174				
Immunogen	Recombinant human gamma catenin protein				
Verified antibody applications	Flow (intracellular) (verified), IF (verified)				
Antibody target cellular localization	Plasma membrane				
Species reactivity	Human				
Species reactivity Antibody application notes	Human ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mil.; Western blot 1-2 ug/mit.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody				
	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/mlilion cells; Immunofluorescence: 1-2 ug/ml.; Western blot 1-2 ug/ml.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect				
Antibody application notes	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody				
Antibody application notes Positive control	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin				
Antibody application notes Positive control Shipping condition	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/mlilion cells; Immunofluorescence: 1-2 ug/ml.; Western blot 1-2 ug/ml.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,				
Antibody application notes Positive control Shipping condition Storage Conditions	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/mlilion cells; Immunofluorescence: 1-2 ug/ml.; Western blot 1-2 ug/ml.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C				
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin Room temperature Store a1 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C 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, Purffied: 0.2 mg/mL in PBS/0.05% BSA/0.05%				
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when				
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life	ELISA: For coating, order Ab without BSA; Flow cytometry: 1-2 ug/million cells; Immunofluorescence: 1-2 ug/mL; Western blot 1-2 ug/ml.; Optimal dilution for a specific application should be determined, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody A431, T47D or MCF-7 cells. Skin Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C 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% BSA/0.05% azide, Purfled, BSA-free: 1 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when stored as recommended Cancer, Cell adhesion, Developmental biology, Signal				

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 November 30, 2025 at 01:39:04 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/catenin-gamma-recombinant-monoclonal-mouse-antibody-rctng-1664/