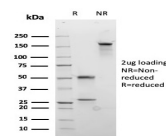


# Desmocollin-2/3 Recombinant Monoclonal Mouse Antibody (rDSC2/3437)



## Product Description

Desmosomes are intercellular adhering junctions that represent cell surface attachment sites for intermediate filament. The desmosome is subdivided into two regions. The plaque region lies adjacent to the plasma, and is believed to contain molecules that attach the intermediate filament cytoskeleton to the desmosome. The core region is composed of transmembrane glycoproteins that are thought to mediate cell-cell adhesion. Desmogleins and desmocollins are the main desmosomal transmembrane proteins. These desmosomal glycoproteins belong to the members of the cadherin family of adhesion molecules. Three different isoforms of both desmogleins and desmocollins have been identified, named as desmoglein 1-3 and desmocollins. Desmosomal cadherins showed differentiation-specific expression in the human epidermis, although the functional significance of this differential expression is not fully understood. Desmocollin-1 can be found in the upper layers. The expression of desmocollin-2 varies in the basal and suprabasal layers. And desmocollin-3 is expressed more evenly throughout the suprabasal layers.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 3437, Anti-Desmocollin-2/3 (rDSC2/3437)**

## Product attributes

Antibody number	#3437
Antibody reactivity (target)	Desmocollin-2/3
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rDSC2/3437
Isotype	IgG1, kappa
Molecular weight	110 kDa (DSC2); 100 kDa (DSC3)
Synonyms	ARVD11; Cadherin family member 2; CDHF2; Desmocollin-2; Desmocollin-3; Desmosomal glycoprotein II and III; Desmosomal glycoprotein II; Desmosomal glycoprotein III; DG2; DGI/III; DSC2; DSC3
Human gene symbol	DSC2
Entrez gene ID	1824
SwissProt	Q02487
Unigene	95612
Immunogen	Recombinant full-length human DSC2 protein
Antibody target cellular localization	Plasma membrane
Verified antibody applications	ELISA
Species reactivity	Human, Mouse, Rat
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	HeLa cells. Skin tissue.
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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	Epidermal cells
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.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 July 1, 2025 at 05:55:29 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/desmocollin-2-3-recombinant-monoclonal-mouse-antibody-rdsc2-3437/>