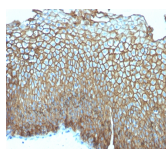


Desmoglein-3 Monoclonal Mouse Antibody (DSG3/2839)



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

Product Description

Recognizes a protein of 130 kDa, identified as Desmoglein-3 (DSG3). This MAbs is highly specific to Desmoglein-3 and does not cross-react with other members of the Desmoglein-family. DSG3 is a calcium-binding transmembrane glycoprotein component of desmosomes in vertebrate epithelial cells. Research has shown that DSG3 has a very high sensitivity (80%) and specificity (100%) in recognizing squamous cell carcinoma (SqCC). Therefore, DSG3 is considered a very important marker for lung SqCC and can be a useful ancillary marker to separate SqCC from other subtypes of lung cancer. Moreover, studies have shown that DSG3 expression in lung SqCC may indicate a poor prognosis.

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 2839, Anti-Desmoglein-3 (DSG3/2839)

Product attributes

Antibody number	#2839
Antibody reactivity (target)	Desmoglein-3
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DSG3/2839
Isotype	IgG1, kappa
Molecular weight of antigen	130 kDa
Synonyms	130kDa pemphigus vulgaris antigen (PVA); Balding (Bal); Cadherin family member 6 (CDHF6); Desmoglein-3 (DSG3)
Human gene symbol	DSG3
Entrez gene ID	1830
SwissProt	P32926
Unigene	1925
Immunogen	Recombinant fragment (around aa 379-491) human DSG3 protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	A431 cells, Skin or Lung Squamous Cell Carcinoma (SCC).
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
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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.
Cell/tissue expression	Epidermal cells
Tumor expression	Squamous cell carcinoma

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 June 23, 2026 at 12:56:48 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/desmoglein-3-monoclonal-mouse-antibody-dsg3-2839/>