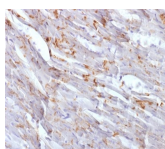


# N-Cadherin / Cadherin-2 / CD325

## Monoclonal Mouse Antibody

### (CDH2/1573)



## Product Description

This antibody recognizes a protein of ~140 kDa, identified as N-Cadherin (NCAD), also known as CD325. N-cadherin is a 140 kDa protein belonging to a family of transmembrane molecules that mediate calcium-dependent intercellular adhesion. Cadherins are involved in controlling morphogenetic movements during development and regulate cell surface adhesion through homotypic adhesion with the same cadherin species. Expression of N-cadherin has been reported on a variety of normal tissues including neuronal, endothelial and muscle cells, and a subpopulation of early hematopoietic progenitor cells. Results aid in the classification of malignant non-carcinomatous neoplasms including mesotheliomas, chordomas, synovial sarcomas, malignant melanomas, epithelioid sarcomas, epithelioid angiosarcomas, clear cell sarcomas as well as serous and endometrioid tumors of the ovary have been demonstrated to be N-cadherin positive, whereas mucinous tumors are negative. Other N-cadherin-positive neoplasms include renal cell carcinomas and some variant breast tumors, including medullary breast carcinomas and sarcomatoid metaplastic breast carcinomas.

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 1573, Anti-Cadherin-2|CD325|N-Cadherin (CDH2/1573)**

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

## Product attributes

Antibody number	#1573
Antibody reactivity (target)	Cadherin-2, CD325, N-Cadherin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CDH2/1573
Isotype	IgG1, kappa
Molecular weight	130-140 kDa
Synonyms	Cadherin-2 N cadherin neuronal; Cadherin-2 type 1; Cadherin-2; Calcium dependent adhesion protein neuronal; CD325; CDH2; CDHN; CDw325; N-Cadherin; NCAD
Human gene symbol	CDH2
Entrez gene ID	1000
SwissProt	P19022
Unigene	464829
Immunogen	Recombinant human CDH2 intracellular domain (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human, Mouse
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Immunohistology (formalin): 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Western blotting 0.5-1 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	HeLa or HUVEC cells. Heart, Pancreas or Cerebral Cortex
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
Antibody research areas	Cancer, Cell adhesion, Developmental biology
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	Metastatic cells

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 September 9, 2025 at 08:50:28 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-n-cadherin-cadherin-2-cd325-cdh21573/>