

P-Cadherin Monoclonal Mouse Antibody (6A9)

Product Description

This antibody recognizes a protein of 116 kDa, identified as P-Cadherin-1 (CDH3). It is a calcium-dependent cell-cell adhesion glycoprotein comprised of five extracellular cadherin repeats, a transmembrane region and a highly conserved cytoplasmic tail. This gene is located in a six-cadherin cluster in a region on the long arm of chromosome 16 that is involved in loss of heterozygosity events in breast and prostate cancer. In addition, aberrant expression of this protein is observed in cervical adenocarcinomas. Mutations in this gene have been associated with congenital hypotrichosis with juvenile macular dystrophy. 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 1428, Anti-P-Cadherin (6A9)**

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Invest Dermatol (1994) 102:870-877. (WB; IP; adhesion blocking)
2. Cancer (1999) 86(7): 1263-1272. (IHC, FFPE)
3. Cell Comm Adhesion (2002) 9(2): 103-115. (IF; WB; IP)
4. Oncogene (2006) 25: 4595-4604. (IF; WB)

Product attributes

Antibody number	#1428
Antibody reactivity (target)	P-Cadherin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	6A9
Isotype	IgG1, kappa
Molecular weight	116 kDa
Synonyms	Cadherin 3 type 1; Cadherin-3; Cadp; Calcium dependent adhesion protein placental; CDH3; CDHP; HJMD; P-cadherin (Placental); PCAD; Placental cadherin
Human gene symbol	CDH3
Entrez gene ID	1001
SwissProt	P22223
Unigene	191842
Immunogen	A431 trypsinized membranes
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Functional studies (published for clone), IHC (FFPE) (published for clone), IF (published for clone), IP (published for clone), WB (published for clone)
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, Immunofluorescence: 1-2 ug/mL, 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	A431 or PC-3 cells. Pancreas, Placenta or Prostate
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
Cell/tissue expression	Epithelial cells
Tumor expression	Cervical cancer

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
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, LI-COR Bioscience.