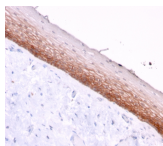


CD44v4/5 Monoclonal Mouse Antibody (3D2)



Product Description

This antibody recognizes an epitope encoded by exon v9 on the variant portion of human CD44. The CD44 molecule belongs to a family of cellular adhesion molecules found on a wide range of normal and malignant cells in epithelial, mesothelial and hemopoiesis tissues. CD44 is a single gene with 20 exons, of which 10 are normally expressed to encode the basic CD44 (H-CAM) molecule. The additional 10 exons (v1 to v10) are only expressed by alternative splicing of the nuclear RNA. The expression of specific cell adhesion molecule CD44 splice variants has been reported to be associated with metastasis in certain human malignancies.

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 3075, Anti-CD44v4/5 (3D2)

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

Product attributes

Antibody number	#3075
Antibody reactivity (target)	CD44v4/5
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3D2
Isotype	IgG1, kappa
Molecular weight	80-95 kDa
Synonyms	LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Epican; Extracellular Matrix Receptor III (ECM III); GP90 Lymphocyte Homing Adhesion Receptor; HCAM; HCELL; hematopoietic cell E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Hermes Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian blood group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDU3; MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)
Human gene symbol	CD44
Entrez gene ID	960
SwissProt	P16070
Unigene	502328
Immunogen	Recombinant human full-length CD44v3-10 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Cervix or Tongue Squamous Cell Carcinoma
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
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	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 January 6, 2026 at 07:28:24 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd44v4-5-monoclonal-mouse-antibody-3d2/>