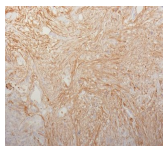


Fibronectin Monoclonal Mouse Antibody (HFN7.1)



Product Description

Fibronectin is a dimeric glycoprotein of 440 kDa, which is present in cells, extracellular matrix, and blood. It possesses at least four binding sites for collagen, glycosaminoglycans, transglutaminase, and a cell surface receptor. Fibronectin is involved in cell adhesion, tissue organization, and wound healing. This MAb is directed against the peptide core and reacts with both the plasma and cellular forms of fibronectin. It blocks the fibronectin-mediated cell attachment not by disrupting the collagen-fibronectin interaction, but by interfering with the attachment of fibronectin to its receptor on the cell surface.

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 0270, Anti-Fibronectin (HFN7.1)

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

Product attributes

Antibody number	#0270
Antibody reactivity (target)	Fibronectin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HFN7.1
Isotype	IgG1, kappa
Molecular weight	220 kDa (monomer); 440 kDa (dimer)
Synonyms	Cold insoluble globulin (CIG); FINC; FN1; FNZ; GFND; GFND2; LETS; Migration stimulating factor (MSF); Ugl-Y3
Human gene symbol	FN1
Entrez gene ID	2335
SwissProt	P02751
Unigene	203717
Immunogen	Human fibronectin purified from serum by affinity chromatography on gelatin-sepharose
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Extracellular matrix
Species reactivity	Human
Expected antibody applications	ELISA (published for clone), Functional studies (published for clone), IF (published for clone), WB (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	SW156 cells or Kidney
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	Cell adhesion, Extracellular matrix
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.

References

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

1. J Biol Chem (1991) 266(14): 23323-23328. (inhibition of integrin binding; epitope mapping)
2. Anat Rec (1996) 245:668-676. (IF, frozen tissue)
3. PNAS USA (2005) 102(17): 5953-5957. (functional studies)
4. Biomaterials (2005) 26: 4523-4531. (IF; functional studies)
5. Hybridoma (2006) 25(4): 202-208. (ELISA; WB)

This datasheet was generated on September 17, 2025 at 02:52:55 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-fibronectin-hfn7-1/>