

Fibronectin Monoclonal Mouse Antibody (TV-1)



Product Description

Fibronectin is a soluble dimeric glycoprotein of 440 kDa, which is present in cells, extracellular matrix, and blood. This MAb reacts with the cellular as well as plasma form of fibronectin. Reportedly, after iv administration, this MAb localizes to tumor vessels where it binds to the underlying basement. The epitope recognized by this antibody is not accessible to the circulating MAb in normal tissues, indicating that it can be used to specifically target tumor vessels in vivo. TV-1 is reportedly useful for delivering vasoactive agents to tumors to induce increased vascular permeability or blood flow prior to treatment with chemotherapeutic drugs or MAbs.

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 0029, Anti-Fibronectin (TV-1)

Duadinat attuibintaa

Call us: 800-304-5357

Product attributes			
Antibody number	#0029		
Antibody reactivity (target)	Fibronectin		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	TV-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	T-cell lymphoma biopsy		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Extracellular matrix		
Species reactivity	Human, Mouse, Pig, Rat		
Expected antibody applications	Flow, surface (published for clone), IHC (frozen) (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, pl +0.0, 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 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. Cancer Res (1995) 55: 2673-2680. (IHC, frozen; WB; in vivo tumor targeting)
- 2. BMC Cancer (2006) 6: 8. (Flow, surface)