Fibronectin Monoclonal Mouse Antibody (568)

Product Description

The epitope of this antibody is located in the 8th type III repeat in the cell-binding region of fibronectin. Fibronectins are disulfide-linked, dimeric glycoproteins of ~440 kDa. They possess at least four binding sites for collagen, glycosaminoglycans, transglutaminase, and a cell surface receptor. Fibronectins are involved in cell adhesion, tissue organization, and wound healing. Fibronectins are present in basement membranes, interstitial connective tissue matrix, and blood. Cellular fibronectin is widely distributed in the stroma of many malignant tumors. This MAb is excellent for staining of formalin-fixed, paraffin-embedded tissues. Catalog number key for antibody number 0351, Anti-Fibronectin (568)

Call us: 800-304-5357 Email: techsupport@biotium.com

Product attributes

Product attributes				
Antibody number	#0351			
Antibody reactivity (target)	Fibronectin			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	568			
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	High molecular weight proteins secreted by cultivated human fibroblasts			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Extracellular matrix			
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, Immunohistochemistry (formalin-fixed): 0.5-1 ug/mL for 30 minutes at RT, Flow cytometry: 0.5-1 ug/mL for 30 minutes at RT, Flow cytometry: 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris buffer with 1 mM EDTA pH 9.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
Positive control	SW156 cells. 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			
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 research areas	Cell adhesion, Extracellular matrix			

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 October 28, 2025 at 09:02:53 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fibronectin-monoclonal-mouse-antibody-568/