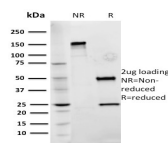


Fibronectin Monoclonal Mouse Antibody (Fn-3)



Product Description

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. Epitope of this MAb is located in the 2nd-3rd type-III repeats of fibronectin. Fibronectins are extracellular matrix glycoproteins that are essential for embryonic development. Fibronectins are also 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 reacts with human cellular fibronectin, but not plasma fibronectin.

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 2848, Anti-Fibronectin (Fn-3)

Product attributes

Antibody number	#2848
Antibody reactivity (target)	Fibronectin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	Fn-3
Isotype	IgG1, kappa
Molecular weight	220 kDa (monomer); 440 kDa (dimer)
Synonyms	Cold insoluble globulin (CIG); FINC; FN1; FN2; GFND; GFND2; LETS; Migration stimulating factor (MSF); Ugi-Y3
Human gene symbol	FN1
Entrez gene ID	2335
SwissProt	P02751
Unigene	203717
Immunogen	FR5 cells, derived by SV40 transformation of human mammary epithelial cells.
Antibody target cellular localization	Extracellular matrix
Species reactivity	Human
Expected antibody applications	IF (published for clone), WB (published for clone)
Positive control	SW156 cells. Kidney.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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 (<i>Bos taurus</i>), 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.

References

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

- Am J Physiol Cell Physiol (2002) 282: C654-C664. (IF; WB)
- Meth Mol Biol (2009) https://doi.org/10.1007/978-1-59745-413-1_18 (IF)
- J Tissue Eng Regen Med. (2020) 14:761-773. (IF, frozen tissue sections)

This datasheet was generated on January 8, 2026 at 03:26:49 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/fibronectin-monoclonal-mouse-antibody-fn-3/>