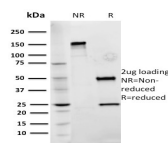


Fibronectin Monoclonal Mouse Antibody (Fn-3)



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; FNZ; 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 July 18, 2025 at 05:34:20 AM. Visit product page to check for updated information before use.

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