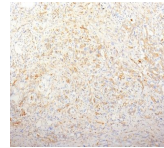


Fibronectin Monoclonal Mouse Antibody (TV-1)



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)**

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)

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, 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 |

| 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 |
| 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, LI-COR Bioscience.