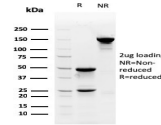


Myofibroblast Marker Monoclonal Mouse Antibody (PR 2D3)



Product Description

Myofibroblasts are a unique group of smooth muscle-like fibroblasts that play an important role in oncogenesis, inflammation, repair, wound contraction and fibrosis. Like smooth muscle (SM) cells, myofibroblasts contain microfilament bundles and express -SM Actin, the Actin isoform that is present in myoepithelial cells and SM cells and especially abundant in vascular SM cells. Myofibroblasts secrete inflammatory and anti-inflammatory cytokines, chemokines, growth factors and lipid and gaseous inflammatory mediators, as well as extracellular matrix proteins and proteases in most organs and tissues. Besides being temporarily present following tissue injuries and fibrocontractive diseases, myofibroblasts are also present under normal conditions in regions such as the skin, pulmonary septa and periodontal ligaments. Stem cell factor and platelet-derived growth factor (PDGF) are two secreted proteins responsible for differentiating myofibroblasts from embryological stem cells. PR 2D3 reacts with a cell membrane component of cells in the pericypt sheath; with smooth muscle cells and myofibroblasts. This antibody is considered to be the gold standard for the identification of myofibroblasts.

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.

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #3069 |
| Antibody reactivity (target) | Myofibroblast Marker |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | PR 2D3 |
| Isotype | IgG1, kappa |
| Molecular weight | 28 kDa |
| Human gene symbol | Not Known |
| Entrez gene ID | Not Known |
| SwissProt | Not Known |
| Unigene | Not Known |
| Immunogen | Crude homogenate of normal human colorectal mucosa |
| Verified antibody applications | WB (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | Myofibroblasts. |
| 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 |
| Cell/tissue expression | Myofibroblasts |

| 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, or LI-COR Bioscience.