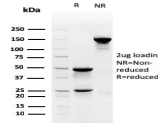


Myofibroblast Marker Monoclonal Mouse Antibody (PR 2D3)



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 pericrypt 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, LI-COR Bioscience.