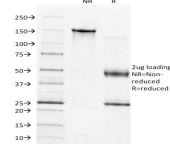


TIMP1 Monoclonal Mouse Antibody (2A5)



Product Description

TIMP-1, TIMP-2, TIMP-3 and TIMP-4 (for tissue inhibitor of metalloproteinases -1, -2, -3 and -4) complex with metalloproteinases such as collagenases, gelatinases and stromelysins, resulting in irreversible inactivation of the metalloproteinase. TIMP-1 is identical to EPA (erythroid-potential activity). PTH has been shown to be a regulator of TIMP-2 in osteoblastic cells. TIMP-3 may be involved in regulating trophoblastic invasion of the uterus as well as in regulating remodeling of the extracellular matrix during the folding of epithelia, and in the formation, branching and expansion of epithelial tubes. TIMP-4 is most highly expressed in heart tissues. Studies have demonstrated that TIMP1 is useful as a biomarker for early detection of colorectal cancer, outperforming CEA. Additionally, TIMP1 studies have demonstrated its role in CRC tumorigenesis, as well as observing its overexpression in metastatic lymph nodes.

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.

References

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

Gut (1998) 43:791-797. (WB; IHC, FFPE)

Product attributes

Antibody number	#1648
Antibody reactivity (target)	TIMP1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	2A5
Isotype	IgG1, kappa
Molecular weight	23 kDa
Synonyms	Collagenase inhibitor; EPO; Erythroid-Potentiating Activity (EPA); Fibroblast collagenase inhibitor; Human Collagenase Inhibitor (HCI); Metalloproteinase inhibitor 1; Tissue Inhibitor of Metalloproteinase 1 (TIMP1)
Human gene symbol	TIMP1
Entrez gene ID	7076
SwissProt	P01033
Unigene	522632
Immunogen	A synthetic peptide (aa 72-82) of human TIMP1 protein (FQALGDAADIR)
Antibody target cellular	Secreted (extracellular)
Species reactivity	Human
Expected antibody applications	IHC (FFPE) (published for clone), WB (published for clone)
Antibody application notes	ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	293 cells. Colon Carcinoma.
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
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
Tumor expression	Colon cancer

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, of LI-COR Bioscience.