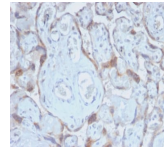


TIMP2 Recombinant Monoclonal Rabbit Antibody (TIMP2/2488R)



Product Description

This antibody recognizes a protein of 21 kDa, identified as tissue inhibitor of metalloproteinases-2 (TIMP-2). It is closely related to TIMP-1 and shows the highest binding affinity to both the latent (pro) and active forms of 72 kDa Type IV collagenase (also known as MMP-2 or gelatinase A). It also has affinity for the active form of 92 kDa Type IV collagenase (also known as MMP-9 or gelatinase B). TIMP-2 inhibits the proteolytic invasiveness of tumor cells and normal placental trophoblast cells.

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 2488, Anti-TIMP2 (TIMP2/2488R)

Product attributes

Antibody number	#2488
Antibody reactivity (target)	TIMP2
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	TIMP2/2488R
Isotype	IgG
Molecular weight	21 kDa
Synonyms	Metalloproteinase inhibitor 2; TIMP metalloproteinase inhibitor 2; TIMP-2; TIMP2; Tissue Inhibitor of Metalloproteinase 2
Human gene symbol	TIMP2
Entrez gene ID	7077
SwissProt	P16035
Unigene	633514
Immunogen	A synthetic peptide from the N-terminus of human TIMP2 protein (aa56-66) (DSGNDIYGNP1)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Cytoplasmic, Secreted (extracellular)
Species reactivity	Cow, Guinea pig, Human, Mouse, Rabbit, Rat, Zebrafish
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user
Positive control	HeLa cells, Normal Placenta, Breast, Colon, Endometrial, Prostate or Ovarian 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
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	Extracellular matrix, Proteases

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, and of LI-COR Bioscience.