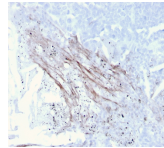


Tenascin C Recombinant Monoclonal Mouse Antibody (TNC/3635)



Product Description

In Western blotting, this antibody reacts with two bands of ~MW of 210 kDa and 300 kDa, identified as two isoforms of Tenascin C. Specificity of this MAb is validated by sequential immunoprecipitation with a PAb against Tenascin C. Tenascin C is a multifunctional, disulfide-linked hexameric extracellular matrix glycoprotein expressed in association with mesenchymal epithelial interactions during development and in the neo-vasculature and stroma of undifferentiated tumors. In adults, it is restricted to certain epithelial-stromal interfaces and increases markedly in hyper-proliferative diseases and in stroma of many neoplasms, including gliomas, breast, squamous and lung carcinomas.

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	#3635
Antibody reactivity (target)	Tenascin C
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	TNC/3635
Isotype	IgG1, kappa
Molecular weight	210 kDa and 300 kDa
Synonyms	Cytotactin; Glioma-associated-extracellular matrix antigen; GMEM; GP 150-225; Hexabrachion (HXB); J1; Myotendinous antigen; Neuronectin; TNC
Human gene symbol	TNC
Entrez gene ID	3371
SwissProt	P24821
Unigene	143250
Immunogen	Recombinant human full-length TNC protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Extracellular matrix
Species reactivity	Human
Positive control	Extracellular matrix in tonsil and blood vessels. Stroma of many tumors such as breast, squamous cell and lung carcinomas. Staining of normal fibroblasts serves as internal positive control (IHC).
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	Breast cancer, Glioma, Lung cancer, Squamous cell carcinoma

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.