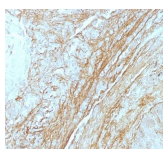


# Tenascin C Monoclonal Mouse Antibody (T2H5)



## 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0392, Anti-Tenascin (T2H5)**

## Product attributes

<b>Antibody number</b>	#0392
<b>Antibody reactivity (target)</b>	Tenascin
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	T2H5
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	210 kDa and 300 kDa
<b>Synonyms</b>	Cytotactin; Glioma-associated-extracellular matrix antigen; GMEG; GP 150-225; Hexabrachion (HXB); J1; Myotendinous antigen; Neuronectin; TNC
<b>Human gene symbol</b>	TNC
<b>Entrez gene ID</b>	3371
<b>SwissProt</b>	P24821
<b>Unigene</b>	143250
<b>Immunogen</b>	Human breast carcinoma
<b>Verified antibody applications</b>	IHC (FPPE) (verified)
<b>Antibody target cellular localization</b>	Extracellular matrix
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL. Does not react with rat; others not known. Immunohistology formalin-fixed 2-4 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	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
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C. Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on December 23, 2024 at 12:04:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-tenascin-t2h5/>