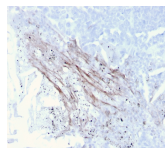


# Tenascin C Recombinant Monoclonal Mouse Antibody (TNC/3635)



## 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](mailto: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; Neurectin; 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 | <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>  |
| 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.