

**Purified Mouse Anti-human Tenascin
Antibody *T2H5, monoclonal***Catalog number: V1032320
Unit size: 0.1 mg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Tenascin
Clone	T2H5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IHC(P), IP, WB

Applications

Tenascin (TN) is a 280 kDa protein that can be expressed in the basement membrane, perisynaptic extracellular matrix and extracellular region of cells. It is sometimes called myotendinous antigen, tenascin-C and cytotactin. In humans, TN recognizes syndecan. It upregulates cell population proliferation and gene expression, but it also is a repressor of cell adhesion. TN plays an important role in organismal processes, for instance, mesenchymal-epithelial cell signaling involved in prostate gland development, osteoblast differentiation and extracellular matrix organization. Sequencing of TN has demonstrated it contains 31 conserved structural units: EGF-like 1, EGF-like 2, EGF-like 3, EGF-like 4, EGF-like 5, EGF-like 6, EGF-like 7, EGF-like 8, EGF-like 9, EGF-like 10, EGF-like 11, EGF-like 12, EGF-like 13, EGF-like 14, EGF-like 15, Fibronectin type-III 1, Fibronectin type-III 2, Fibronectin type-III 3, Fibronectin type-III 4, Fibronectin type-III 5, Fibronectin type-III 6, Fibronectin type-III 7, Fibronectin type-III 8, Fibronectin type-III 9, Fibronectin type-III 10, Fibronectin type-III 11, Fibronectin type-III 12, Fibronectin type-III 13, Fibronectin type-III 14, Fibronectin type-III 15, and Fibrinogen C-terminal domain. TN is clinically significant because abnormalities in its function have been thought to be involved with diseases, such as autosomal dominant deafness-56 (DFNA56).