

**Purified Mouse Anti-human/mouse/dog  $\beta$ III-Tubulin Antibody \*TU-20, monoclonal\***Catalog number: V103230  
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	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Human, mouse, dog, rat, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\beta$ III-Tubulin
Clone	TU-20

**Biological Properties**

---

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

**Applications**

---

Tubulin  $\beta$ -3 chain (sometimes called Tubulin  $\beta$ -4 chain or Tubulin  $\beta$ -III) is a protein with a molecular weight of 59 kDa, expressed in the microtubule, filopodium and extracellular exosome of cells. In *Homo sapiens*, tubulin  $\beta$ -3 chain binds to GTP and netrin receptor. It aids in organismal processes, in particular, mitotic cell cycle, neuron differentiation and microtubule cytoskeleton organization. Tubulin  $\beta$ -3 chain has been thought to be involved with essential functions such as GTPase activity. Tubulin  $\beta$ -3 chain is clinically significant because abnormalities in its function have been closely linked to diseases like cortical dysplasia, complex with other brain malformations 1 (CDCBM1) and congenital fibrosis of extraocular muscles 3A (CFEOM3A). Congenital fibrosis of extraocular muscles 3A, an autosomal dominant inherited disorder characterized by superior rectus atrophy and specific learning disability, has in particular been of interest to researchers.