

**Purified Mouse Anti-human/mouse/rat  $\alpha$ -Tubulin Antibody \*TU-16, monoclonal\***

Catalog number: V103105

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, mouse, rat, pig, nicotiana, dog
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\alpha$ -Tubulin
Clone	TU-16

**Biological Properties**

---

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

**Applications**

---

Tubulin  $\alpha$  chain is a protein with a molecular weight of 42 kDa, located in the microtubule of cells. In humans, tubulin  $\alpha$  chain binds with GTP. Tubulin  $\alpha$  chain has been associated with important functions like GTPase activity. It is an integral part of organismal processes, in particular, microtubule-based process. Sequencing of Tubulin  $\alpha$  chain has shown it contains 2 conserved structural units: tubulin\_C and tubulin domain.