

**FITC Mouse Anti-human/mouse/pig  $\alpha$ -Tubulin Antibody \*TU-01, monoclonal\***

Catalog number: V103120

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, pig, turkey, eisenia, paramecium, nicotiana, yeast
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\alpha$ -Tubulin
Clone	TU-01
Conjugate	FITC

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	FC (QC TESTED), ICC

**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.