

**Purified Mouse Anti-human/mouse/pig  $\alpha/\beta$ -Tubulin dimer Antibody \*TU-08, monoclonal\***

Catalog number: V103080

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, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\alpha/\beta$ -Tubulin dimer
Clone	TU-08

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, ICC

**Applications**

---

Tubulin  $\alpha$ -4A chain (sometimes referred to as Tubulin  $\alpha$ -1 chain, Tubulin H2- $\alpha$ , Testis-specific  $\alpha$ -tubulin or  $\alpha$ -tubulin 1) is a protein with a molecular weight of 58 kDa, located in the cytoplasm and extracellular exosome of cells. In humans, tubulin  $\alpha$ -4A chain has been closely linked to important functions such as GTPase activity. It aids in G2/M transition of mitotic cell cycle. It binds to GTP and protein kinase. It is thought to be essential to organismal processes, for instance, microtubule cytoskeleton organization, ciliary basal body-plasma membrane docking and G2/M transition of mitotic cell cycle. Mutations and abnormalities in tubulin  $\alpha$ -4A chain have been closely linked to a number of diseases, for instance, amyotrophic lateral sclerosis 22, with or without frontotemporal dementia (ALS22). Amyotrophic lateral sclerosis 22, with or without frontotemporal dementia, an autosomal dominant inheritancedisorder characterized by amyotrophic lateral sclerosis and frontotemporal dementia, has in particular been of interest to researchers.