

Purified Mouse Anti-human/mouse/pig α -Tubulin Antibody *TU-01, monoclonal*

Catalog number: V103100

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	α -Tubulin
Clone	TU-01

Biological Properties

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

Applications

Tubulin α chain is a protein with a molecular weight of 42 kDa, located in the microtubule of cells. In humans, tubulin α chain binds with GTP. Tubulin α 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 α chain has shown it contains 2 conserved structural units: tubulin_C and tubulin domain.