

Purified Mouse Anti-human/mouse/protozoa γ -Tubulin Antibody *TU-30, monoclonal*Catalog number: V103695
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, protozoa, chicken, cow, rat, pig, plants
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	γ -Tubulin
Clone	TU-30

Biological Properties

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

Applications

Tubulin γ chain is a 34 kDa protein that can be expressed in the γ -tubulin complex and microtubule of cells. Sequencing of tubulin γ chain has exemplified it contains a primary structural unit, the tubulin domain. Tubulin γ chain reacts with GTP. It has been found to be involved in organismal processes, in particular, microtubule nucleation and cytoplasmic microtubule organization. It has been closely linked to key functions like GTPase activity.