

Purified Mouse Anti-human/mouse/pig β -Tubulin Antibody *TU-13, monoclonal*

Catalog number: V103250

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, plants
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	β -Tubulin
Clone	TU-13

Biological Properties

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

Applications

Tubulin β chain is a 58 kda protein that can be located in the cell body, protein-containing complex and cytoplasmic ribonucleoprotein granule of cells. In humans, Tubulin β chain has been associated with important functions such as GTPase and structural molecule activity, and additionally, plays an important role in synapse organization and G2/M transition of mitotic cell cycle, and additionally, recognizes MHC class I protein, ubiquitin protein ligase and protein-containing complex. It has been found to be involved in organismal processes, in particular, natural killer cell mediated cytotoxicity, spindle assembly and cell division. Tubulin β chain is clinically significant because abnormalities in its function have been thought to be involved with diseases such as congenital symmetric circumferentialSkin creases-1 (CSCSC1) and complex cortical dysplasia with other brain malformations⁶ (CDCBM6). Complex cortical dysplasia with other brain malformations⁶, an autosomal dominant inheritance disorder characterized by congenital cerebellar hypoplasia, cortical dysplasia and retinal dysplasia, has especially been of interest to scientists.