

**Purified Mouse Anti-human/mouse/cow  
MAP2 Antibody \*MT-01, monoclonal\***

Catalog number: V1031730

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, cow, cat, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	MAP2
Clone	MT-01

**Biological Properties**

---

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

**Applications**

---

Microtubule-associated protein 2 (MAP-2) is a 200 kDa protein that can be expressed in the dendrite cytoplasm, basal dendrite and proximal neuron projection of cells. In humans, MAP-2 takes part in processes such as cellular protein localization, organelle transport along microtubule and microtubule polymerization. MAP-2 has been thought to be involved with key functions like structural molecule activity. It recognizes  $\tau$  protein, calmodulin and microtubule. MAP-2 plays a role in the downregulation of microtubule polymerization, axon extension and microtubule binding. But it also is a positive regulator of anterograde synaptic vesicle transport and anterograde dense core granule transport. MAP-2 plays an important role in organismal processes, in particular, dendrite morphogenesis, microtubule cytoskeleton organization and dendrite development.