

**Purified Mouse Anti-human/mouse/rat  $\gamma$ -Tubulin GCP2 Antibody \*GCP2-01, monoclonal\***

Catalog number: V103705

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, rat, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\gamma$ -Tubulin GCP2
Clone	GCP2-01

**Biological Properties**

---

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

**Applications**

---

$\gamma$ -Tubulin complex component 2 (also known as  $\gamma$ -ring complex protein 103 kDa, Spindle pole body protein Spc97 homolog or hSpc97) is a protein with a molecular weight of 119 kDa, expressed in the cytosol,  $\gamma$ -tubulin complex and nucleoplasm of cells. In humans,  $\gamma$ -Tubulin complex component 2 is thought to be essential to organismal processes, in particular, microtubule nucleation by interphase microtubule organizing center, cytoplasmic microtubule organization and microtubule nucleation. It binds to  $\gamma$ -tubulin.