

**Purified Mouse Anti-human/mouse/rat  
TPX2 Antibody \*TPX2-01, monoclonal\***

Catalog number: V1032380

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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TPX2
Clone	TPX2-01

**Biological Properties**

---

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

**Applications**

---

Targeting protein for Xklp2 is a 86 kDa protein that can be expressed in the cytosol, intercellular bridge and axon hillock of cells. It is sometimes referred to as Hepatocellular carcinoma-associated antigen 90 and Restricted expression proliferation-associated protein 100. In humans, targeting protein for Xklp2t takes part in processes such as signal transduction by p53 class mediator, G2/M transition of mitotic cell cycle and mitotic spindle organization. It is an integral part of organismal processes, for example, activation of protein kinase activity, mitotic cell cycle and microtubule nucleation. Targeting protein for Xklp2 negatively regulates microtubule depolymerization. It binds with importin- $\alpha$  family protein and protein kinase.