

## alpha Tubulin antibody (mAb)

**Catalog Nos:** 39527, 39528

**RRID:** AB\_2793243

**Clone:** 5-B-1-2

**Isotype:** IgG1

**Application(s):** ICC, IF, WB

**Reactivity:** Human, Mouse, Wide Range Predicted

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Ascites

**Host:** Mouse

**Molecular Weight:** 52 kDa

**Background:** Microtubules of the eukaryotic cytoskeleton perform essential and diverse functions and are composed of a heterodimer of  $\alpha$ -tubulin and  $\beta$ -tubulin. Alpha Tubulin expression is reduced in cells deficient in MeCP2, leading to conditions such as Rett syndrome and autistic-spectrum disorders.

**Immunogen:** This alpha Tubulin antibody was raised against total protein purified from sea urchin sperm axonemes.

**Buffer:** Ascites containing 30% glycerol and 0.035% sodium azide.

### Application Notes:

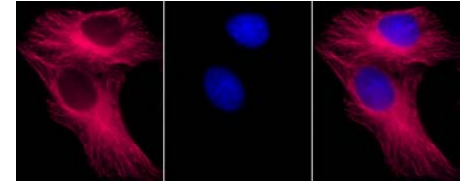
Validated Applications:

WB: 1:500 - 1:2,500 dilution

ICC/IF: 1:1,000 - 1:10,000 dilution

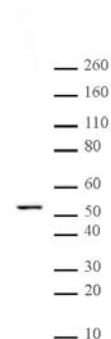
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### alpha Tubulin mAb (Clone 5-B-1-2) tested by immunofluorescence.

Top: HeLa cells stained with alpha Tubulin mAb at 1:500 dilution (red). Middle: Same cells stained with DAPI (blue). Bottom: Both images merged. Staining was carried out using MAX Stain™ Immunofluorescence Tools.



### alpha Tubulin mAb (Clone 5-B-1-2) tested by Western blot.

HeLa whole-cell extract (15  $\mu$ g/lane) probed with alpha Tubulin mAb (1:5,000 dilution).