

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 μ l, 10 μ 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 α -tubulin and β -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. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

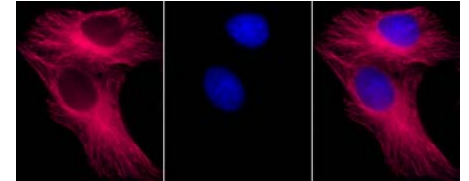
WB*: 1:500 - 1:2,500 dilution

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

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western Blot.

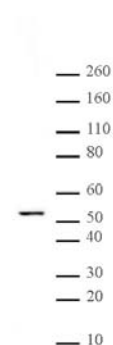
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 μ g/lane) probed with alpha Tubulin mAb (1:5,000 dilution).