

beta Tubulin antibody (mAb)

Catalog Nos: 61157, 61158

RRID: AB_2793531

Clone: 2-28-33

Isotype: IgG1, k

Application(s): IHC, WB

Reactivity: Human, Wide Range Predicted

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 52 kDa

Background: **Beta Tubulin** (β-tubulin) is a structural protein and one of the components of microtubules, cytoskeletal filaments that serve a variety of important roles within the cell, including providing structural stability and performing chromosome segregation.

Immunogen: This beta-tubulin antibody was raised against sarkosyl-resistant filaments from *Strongylocentrotus purpuratus* (sea urchin) sperm axonemes.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.