

Bub1 antibody (mAb)

Catalog Nos: 39735, 39736

RRID: AB_2793325

Clone: 14H5

Isotype: IgG

Application(s): ICC, IF

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 130 kDa

Background: Bub1 is a serine-threonine protein kinase and an important regulator of cell division. It is essential for the G2/M mitotic spindle assembly/integrity checkpoint as well as critical for chromosome alignment. Bub1 is required for the assembly of a large number of proteins at the kinetochore, including CENP-E, CENP-F, BUB1B, MAD2L1 and PLK1. It plays a role in sister chromatid cohesion and it serves as a substrate for the anaphase-promoting complex. Bub1 mediates cell death in response to chromosome mis-segregation and acts to suppress spontaneous tumorigenesis.

Immunogen: This Bub1 antibody was raised against a recombinant protein corresponding to amino acids 281-419 of human Bub1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 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.