

MAD1 antibody (mAb)

Catalog Nos: 39659, 39660

RRID: AB_2793294

Clone: 9B10

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: 85 kDa

Background: MAD1 (MAD1L1, Mitotic arrest deficient 1) is a component of the spindle-assembly checkpoint that prevents the onset of anaphase until all chromosomes are properly aligned at the metaphase plate. MAD1 homodimerizes with MAD2L1 and is present at the kinetochore. Defects in MAD1 are involved in the development and/or progression of various types of cancer.

Immunogen: This MAD1 antibody was raised against a peptide corresponding to amino acids 331-718 of mouse MAD1.

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.