

CENP-E antibody (mAb)

Catalog Nos: 39619, 39620

RRID: AB_2793278

Clone: 1H12

Isotype: IgG

Application(s): ICC, IF

Reactivity: Human, Mouse

Quantities: 200 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 320 kDa

Background: Centrosome-associated protein E (CENP-E) is a kinesin-like protein and an integral component of kinetochore fibers that link centromeres to spindle microtubules. CENP-E localizes to kinetochores throughout all phases of mitotic chromosome movement. After the onset of anaphase it is released, and by telophase becomes bound to interzonal microtubules of the mitotic spindle. It is likely that CENP-E serves as a chromosome motor during anaphase.

Immunogen: This CENP-E antibody was raised against full-length human CENP-E protein.

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:

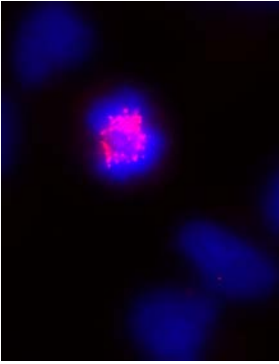
Applications Validated by Active Motif:

ICC/IF: 1 µg/ml dilution

This CENP-E antibody can be used in Western blot to detect CENP-E if the protein has been first enriched by immunoprecipitation.

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.



CENP-E mouse mAb (Clone 1H12) tested by immunofluorescence.

Staining of HeLa cells with CENP-E mouse mAb (Clone 1H12) at a 1 µg/ml dilution (red) and DAPI (blue).

Detection of CENP-E by immunofluorescence.

U2OS cells were stained with CENP-E antibody at a dilution of 1:500. Left panel: DAPI. Middle panel: CENP-E antibody staining. Right panel: merge.

