

MCM2 antibody (pAb)

Catalog No: 61717

RRID: AB_2636824

Isotype: IgG

Application(s): WB

Reactivity: Human

Volume: 100 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 115 kDa

Background: MCM2 (Minichromosome Maintenance Complex Component 2) acts as component of the MCM2-7 complex (MCM complex) which is the putative replicative helicase essential for once per cell cycle DNA replication initiation and elongation in eukaryotic cells. The active ATPase sites in the MCM2-7 ring are formed through the interaction surfaces of two neighboring subunits such that a critical structure of a conserved arginine finger motif is provided in trans relative to the ATP-binding site of the Walker A box of the adjacent subunit. The six ATPase active sites, however, are likely to contribute differentially to the complex helicase activity. Required for the entry in S phase and for cell division.

Immunogen: This antibody was raised against a full-length recombinant protein corresponding to human MCM2.

Buffer: Purified IgG in PBS with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

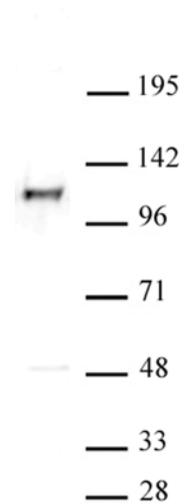
WB*: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*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.



MCM2 antibody (pAb) tested by Western blot.

MCM2 detection by Western blot. The analysis was performed using MCF7 nuclear cell extract (20 µg) and MCM2 antibody at a 1:1,000 dilution.