

SMC3 antibody (pAb)

Catalog Nos: 61131, 61132

RRID: AB_2688005

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 150 kDa

Background: The Structural Maintenance of Chromosomes (SMC) family proteins play critical roles in various nuclear events that require structural changes of chromosomes, including mitotic chromosome organization, DNA recombination and repair and global transcriptional repression. SMC3 is a component of the cohesin complex that plays an essential role during chromatid segregation. SMC3 is also involved in DNA recombination and repair, mitotic chromosome organization and microtubule-mediated intracellular transport. SMC3 is acetylated during replication, which is essential for cohesion, and is phosphorylated by ATM subsequent to DNA double strand breakage. SMC3 is frequently elevated in human colon carcinoma and overexpression of the protein transforms fibroblasts.

Immunogen: This SMC3 antibody was raised against a peptide in the C-terminal region of human SMC3.

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

Application Notes:

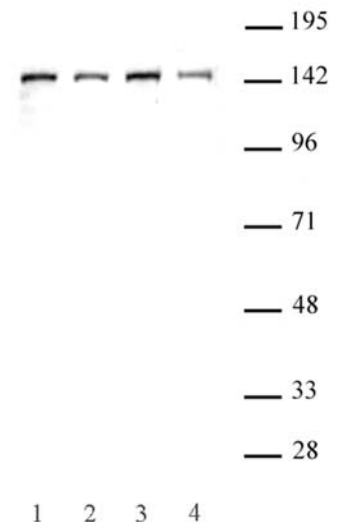
Applications Validated by Active Motif:

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

*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.



SMC3 antibody (pAb) tested by Western blot.
 Detection of SMC3 by Western blot analysis.
 Lane 1: Nuclear extract of HeLa cells (20 μ g).
 Lane 2: Nuclear extract of Hep G2 cells (20 μ g).
 Lane 3: Nuclear extract of HEK293 cells (20 μ g).
 Lane 4: Nuclear extract of U2OS cells (20 μ g).
 All were probed with SMC3 antibody (pAb) at a 1:500 dilution.