

## SMC4 antibody (pAb)

**Catalog Nos:** 61129, 61130

**RRID:** AB\_2793519

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 175 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. SMC4 is the central component of the condensin complex, a complex required for conversion of interphase chromatin into condensed mitotic chromosomes. SMC4 forms a heterodimer with SMC2 and contains a myosin-like ATPase domain that serves as a molecular motor to help organize chromatin. During mitosis, most of the condensin complex is associated with the chromatin.

**Immunogen:** This SMC4 antibody was raised against a peptide from the N-terminus of human SMC4.

**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:

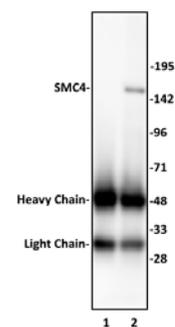
IP: 10 µl per IP

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.

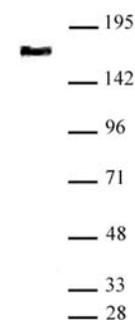
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This 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.



### SMC4 antibody (pAb) tested by Immunoprecipitation.

10 µl of SMC4 antibody was used to immunoprecipitate SMC4 from 500 µg of HeLa nuclear cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the SMC4 antibody at a dilution of 1:1,000.



### SMC4 antibody (pAb) tested by Western blot

Nuclear extract of HeLa cells (20 µg) probed with SMC4 antibody (1:1,000).