

Acf1 antibody (pAb)

Catalog Nos: 39721, 39722**RRID:** AB_2793321**Isotype:** IgG**Application(s):** ICC, IF, WB**Reactivity:** Drosophila**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 185 kDa

Background: *Drosophila* Acf1 is a component of ACF, the ATP-utilizing chromatin assembly and remodeling factor protein complex. ACF catalyzes the ATP-dependent assembly of periodic nucleosome arrays *in vitro*. The catalytic subunit of ACF is the ISWI ATP-dependant chromatin remodeling protein. ACF is related to CHRAC (Chromatin Accessibility Complex), which also contains Acf1 and ISWI. ACF is involved in the establishment and possibly maintenance of silenced heterochromatic domains in the genome, both at pericentric heterochromatin and at genes repressed by the Polycomb complex.

Immunogen: This Acf1 antibody was raised against recombinant C-terminal fragment of Acf1 from *Drosophila melanogaster*.

Buffer: PBS containing 0.1% sodium azide, 0.2 mg/ml BSA and 35% glycerol. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

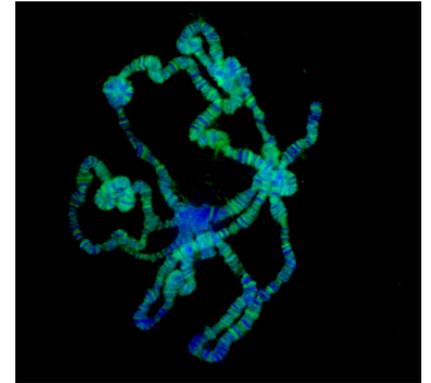
ICC/IF: 1:500 - 1:1,000 dilution

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

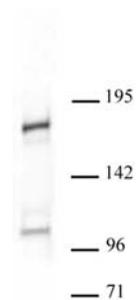
The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

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.



Acf1 pAb tested by immunofluorescence.
Drosophila polytene chromosomes stained with Acf1 pAb (1:1,000 dilution).



Acf1 pAb tested by Western blot.
Drosophila Schneider L2 cells nuclear extract (10 µg) probed with Acf1 pAb (1:500 dilution).