

Rsf1 antibody (pAb)

Catalog No: 39579**RRID:** AB_2753149**Isotype:** IgG**Application(s):** WB**Reactivity:** Drosophila**Volume:** 100 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 58 kDa

Background: Rsf-1 (Remodeling and spacing factor 1) is a member of RSF complex. This complex consists of two subunits: Rsf-1 and SNF2H. RSF possesses a chromatin remodeling activity and a chromatin assembly activity. Rsf-1 assembles the nucleosomes randomly with its histone chaperone activity. The resulting nucleosomes are then redistributed by the ATP-utilizing nucleosome mobilization factor SNF2H into a periodic array. Rsf-1 can be involved in transcriptional repression when it is recruited to DNA via the DNA binding of Gal4 or in transcriptional coactivation when associated with hepatitis B virus pX protein (HBX). In response to hepatitis B infection, Rsf-1 can bind to the viral transcription activator pX, which then coactivates the transcription factor NFκB, thus supporting the life cycle of the viral genome. Additionally, Rsf-1 is thought to be a potent oncogene, with elevated expression observed in ovarian carcinomas.

Immunogen: This Rsf1 antibody was raised against a His-Tagged fusion protein corresponding to amino acids 165-350 of *Drosophila* Rsf1.

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

Application Notes:

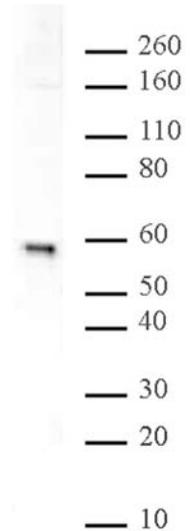
Applications Validated by Active Motif:

WB: 1:1,000 - 1:5,000 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.

**Rsf1 pAb tested by Western blot.**

Schneider's *Drosophila* L2 nuclear extract (10 µg) was probed with Rsf1 pAb (1:1,000 dilution).