

SNF5 antibody (pAb)

Catalog Nos: 39775, 39776

RRID: AB_2614998

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 60 kDa

Background: SNF5 (Sucrose Non Fermenting 5, also called INI1, BAF47 and SMARCB1) is a tumor suppressor protein that is a component of the SWI/SNF chromatin-remodeling complex that regulates gene expression through alterations in chromatin structure. Alterations in chromatin structure facilitated by ATP-dependant chromatin remodeling machines are thought to enhance recognition and binding of trans-acting factors that bind DNA and regulate transcription. SNF5 signals through the p16INK4a pathway to negatively regulate cell cycle progression and mediate chromosome stability. Mutations in SNF5 are a cause of malignant rhabdoid tumors in children.

Immunogen: This SNF5 antibody was raised against a peptide in the C-terminus of SNF5.

Buffer: Purified rabbit IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

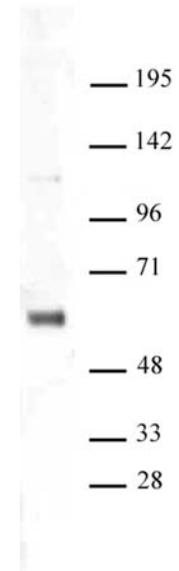
WB*: 1:500 - 1:2,000 dilution

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.



SNF5 antibody (pAb) tested by Western blot. Nuclear extract (20 µg) of Jurkat cells probed with SNF5 antibody (pAb) at a dilution of 1:2,000.