

SMARCA5 antibody (mAb)

Catalog Nos: 39543, 39544

RRID: AB_2615000

Isotype: IgG

Application(s): ChIP, ICC, IF, IP, WB

Reactivity: Human, Mouse

Volumes: 200 µl, 10 µl

Purification: Ascites

Host: Mouse

Molecular Weight: 135 kDa

Background: SMARCA5 is a human homolog of the budding yeast Snf2 protein. SMARCA5 is the catalytic subunit of several chromatin remodeling complexes, such as CHRAC, RSF, ACF, NuRD and NoRC. SMARCA5 is a member of the SWI/SNF superfamily of proteins and members of this family have helicase and ATPase activities and are thought to regulate transcription of certain genes by altering the chromatin structure around those genes. SMARCA5 is a component of the chromatin remodeling and spacing factor RSF (Rsf1), a facilitator of the transcription of class II genes by RNA pol II. It is similar in sequence to the *Drosophila* ISWI chromatin remodeling protein. The protein encoded by this gene is a component of the chromatin remodeling and spacing factor RSF, a facilitator of the transcription of class II genes by RNA polymerase II. The encoded protein is similar in sequence to the *Drosophila* ISWI chromatin remodeling protein.

Immunogen: This antibody was raised against purified human RSF complex which was used to immunize mice.

Buffer: Ascites containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

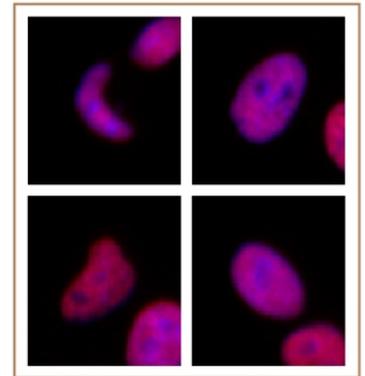
ICC/IF: 1:2,000 dilution

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

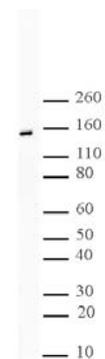
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



SMARCA5 antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with SMARCA5 antibody at a 1:2,000 dilution.



SMARCA5 antibody (mAb) tested by Western blot.

HeLa acid extract (10 µg) was probed with SMARCA5 mAb (1:2,000 dilution).