

SMARCC2 / BAF170 antibody (pAb)

Catalog Nos: 61471, 61472

RRID: AB_2614974

Isotype: IgG

Application(s): ICC, IF, WB

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 170 kDa

Background: **SMARCC2 / BAF170** (SWI/SNF Related, Matrix Associated, Actin Dependent Regulator Of Chromatin, Subfamily C, Member 2) is involved in transcriptional activation and repression of select genes by chromatin remodeling (alteration of DNA-nucleosome topology). Can stimulate the ATPase activity of the catalytic subunit of these complexes. May be required for CoREST dependent repression of neuronal specific gene promoters in non-neuronal cells. Also involved in vitamin D-coupled transcription regulation via its association with the WINAC complex, a chromatin-remodeling complex recruited by vitamin D receptor (VDR), which is required for the ligand-bound VDR-mediated transrepression of the CYP27B1 gene. Belongs to the neural progenitors-specific chromatin remodeling complex (npBAF complex) and the neuron-specific chromatin remodeling complex (nBAF complex). During neural development a switch from a stem/progenitor to a post-mitotic chromatin remodeling mechanism occurs as neurons exit the cell cycle and become committed to their adult state. The transition from proliferating neural stem/progenitor cells to post-mitotic neurons requires a switch in subunit composition of the npBAF and nBAF complexes. As neural progenitors exit mitosis and differentiate into neurons, npBAF complexes which contain ACTL6A/BAF53A and PHF10/BAF45A, are exchanged for homologous alternative ACTL6B/BAF53B and DPF1/BAF45B or DPF3/BAF45C subunits in neuron-specific complexes (nBAF). The npBAF complex is essential for the self-renewal/proliferative capacity of the multipotent neural stem cells. The nBAF complex along with CREST plays a role regulating the activity of genes essential for dendrite growth (By similarity).

Immunogen: This antibody was raised against a peptide within the C-terminal region of mouse SMARCC2 / BAF170.

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:

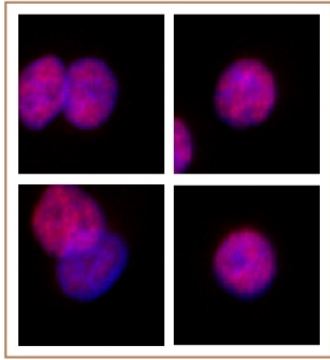
ICC/IF: 1:2,000 dilution

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.

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.



SMARCC2 / BAF170 antibody (pAb) tested by Immunofluorescence.

Formaldehyde fixed HeLa cells stained with SMARCC2 / BAF170 antibody at a 1:2,000 dilution.



SMARCC2 / BAF170 antibody (pAb) tested by Western blot.

Nuclear extract of K-562 cells (30 µg) probed with SMARCC2 / BAF170 antibody (1:500 dilution).