

SMARCA1 / SNF2L1 antibody (mAb)

Catalog Nos: 61465, 61466

RRID: AB_2793649

Clone: 2C4

Isotype: IgG2b

Application(s): IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 130 kDa

Background: **SMARCA1 / SNF2L1** (Sucrose Nonfermenting 2-Like Protein 1) is an energy-transducing component of NURF (nucleosome-remodeling factor) and CERF (CECR2-containing-remodeling factor) complexes. Both complexes facilitate the perturbation of chromatin structure in an ATP-dependent manner. SMARCA1 / SNF2L1 is the mammalian ortholog of the *Drosophila* nucleosome-remodeling ATPase ISWI.

Immunogen: This antibody was raised against a recombinant protein corresponding to amino acids 583-784 of human SNF2L1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91225.

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.