

SMARCAL1 / HARP antibody (mAb)

Catalog Nos: 61115, 61116

RRID: AB_2614999

Clone: 4B10

Isotype: IgG1

Application(s): IP, WB

Reactivity: Bovine, Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 150 kDa

Background: **SMARCAL1 / HARP** (SWI/SNF-related matrix-associated actin-dependent regulator of chromatin subfamily A-like protein 1, HepA-related protein, DNA-dependent ATPase A) is a member of the SWI / SNF family of ATPases and chromatin remodeling proteins. Unlike other SWI / SNF family members, **SMARCAL1 / HARP** has not been shown to remodel chromatin, but possesses an ATP-dependent annealing helicase activity. SMARCAL1 reanneals single-stranded DNA bubbles in DNA at DNA replication forks that are bound by Replication Protein A (RPA). Defects in SMARCAL1 / HARP are a cause of Schimke immuno-osseous dysplasia (SIOD) SIOD causes spondyloepiphyseal dysplasia (a disorder of bone growth that results in short stature, skeletal abnormalities, and problems with vision and hearing), renal dysfunction and T-cell immunodeficiency.

Immunogen: This SMARCAL1 / HARP antibody was raised against purified bovine SMARCAL1 / HARP protein.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.