

SMARCA4 / BRG1 antibody (mAb)

Catalog Nos: 39807, 39808

RRID: AB_2614949

Clone: 3G4

Isotype: IgG2b

Application(s): ICC, IF, WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 180 kDa

Background: SWI/SNF related, matrix associated, actin dependent regulator of chromatin, subfamily a, member 4 (SMARCA4), also known as BRG1, is the central catalytic ATPase subunit of numerous chromatin-remodeling complexes, including SWI/SNF. These complexes are able to use energy from ATP hydrolysis to physically disrupt chromatin architecture of target promoters and facilitate the binding of transcription factors to nucleosomal DNA. SMARCA4 is a co-regulator of transcription and has been implicated in the activation and repression of gene expression through the modulation of chromatin in various tissues and physiological conditions. This protein plays a major role in many cellular processes such as DNA replication, repair and recombination. Mammalian SMARCA4 is usually associated with approximately 10-12 BAF subunits or other proteins involved in regulation of gene expression. SMARCA4 is part of the stem cell-specific BAF complex, esBAF, and is involved in regulating stem cell pluripotency.

Immunogen: This SMARCA4 / BRG1 antibody was raised against a recombinant protein corresponding to amino acids 214-279 of mouse BRG1.

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.