

SNFa/BRM Polyclonal Antibody

(Catalog #A-2025)

Background

The chromatin remodeling complex SWI-SNF which is involved in the modulation of gene expression actively relies on two closely related ATPases known as SNF2a/BRM and BRG-1/SNF2b. These proteins are able to assist nuclear receptors for transcriptional activation. They can also regulate cell proliferation by cooperating with p105Rb for repression of E2F-mediated transcription.

Concentration

1 mg/ml

Description

SNFa/BRM Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid in PBS containing 50% glycerol, 0.5% BSA* and 0.02% sodium azide

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P51531

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from the internal region of human Brm

Storage

Store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

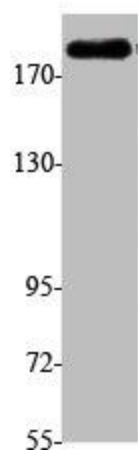
ATP dependent helicase SMARCA2 antibody, BAF190 antibody, BRM antibody, FLJ36757 antibody, Global transcription activator homologous sequence antibody, hBRM antibody, hSNF2a antibody, MGC74511 antibody, SMARCA2 antibody, SNF2 alpha antibody, SNF2/SWI2 like protein 2 antibody, SNF2L2 antibody, SNF2LA antibody, Sth1p antibody, Sucrose nonfermenting 2 like protein 2 antibody

Application

WB, ELISA; Recommended Dilution: WB 1:500-1:2000, ELISA 1:20000

**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*

(kD)



Western Blot analysis of A549 cells using SNFa/BRM
Polyclonal Antibody