

## SMARCA2 / BRM antibody (mAb)

---

**Catalog Nos:** 39805, 39806

**RRID:** AB\_2615072

**Clone:** 1H7A10

**Isotype:** IgG2a

**Application(s):** ChIP, ICC, IF, WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Rat

**Concentration:** 1 µg/µl

**Molecular Weight:** 190 kDa

---

**Background:** SWI/SNF related, matrix associated, actin dependent regulator of chromatin, subfamily a, member 2 (SMARCA2), also known as BRM, is a member of the SWI/SNF family of ATP-dependant chromatin remodeling proteins. SMARCA2 is the catalytic subunit of a multi-subunit protein complex that utilizes the energy from ATP hydrolysis to alter chromatin structure and activate transcription.

**Immunogen:** This SMARCA2 / BRM antibody was raised against a recombinant protein corresponding to amino acids 48-214 of mouse BRM.

**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:**

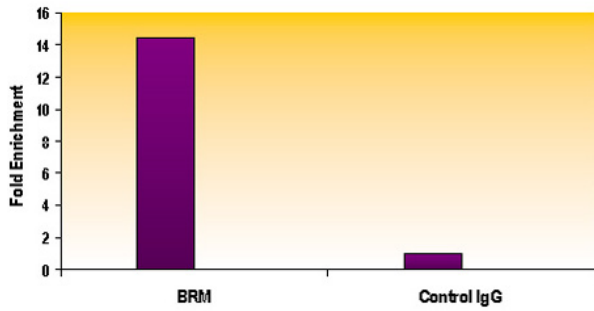
Applications Validated by Active Motif:

ChIP: 10 µg per ChIP

ICC/IF: 0.5 µg/ml dilution

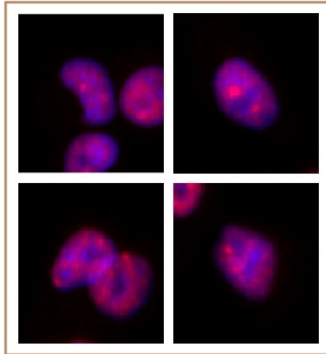
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This 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.



SMARCA2 / BRM antibody (mAb) tested by ChIP.

Chromatin IP performed using the ChIP-IT<sup>®</sup> Express Kit (Catalog No. 53008) and HeLa Chromatin ( $1.5 \times 10^6$  cell equivalents per ChIP) using 10  $\mu$ g of SMARCA2 / BRM antibody or the equivalent amount of rat IgG as a negative control. Real time, quantitative PCR (RT-qPCR) was performed on DNA purified from each of the ChIP reactions using a primer pair specific for the CD44 gene. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control IgG using the ddCT method.



**SMARCA2 / BRM antibody (mAb) tested by Immunofluorescence.**

Formaldehyde fixed HeLa cells stained with SMARCA2 / BRM antibody at a 0.5  $\mu$ g/ml dilution.