

MBD1 antibody (mAb)

Catalog No: 39215
RRID: AB_2615026

Clone: 100B272.1

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG

Concentration: 1 µg/µl

Molecular Weight: 70 kDa

Background: MBD1 (methyl-CpG binding domain protein 1) is thought to function as a mediator of the biological consequences of DNA methylation. Methylation of mammalian DNA has long been recognized to play a major role in a number of cellular functions such as development and control of gene expression. It is generally associated with the repressive chromatin state. The complex series of events leading to this repressive state involve the coordinated regulation of DNA methyltransferases and two other groups of proteins called the Methyl-CpG binding proteins (MBD proteins) and the Kaiso family of proteins. The MBD family of proteins include MeCP2, MBD1, MBD2, MBD3 and MBD4. MBD1 is recruited to both methylated and non-methylated CpGs via separate domains: the CXXC domain targets the protein to non-methylated CpGs whereas the MBD domain targets the protein to methylated CpG. During S phase of the cell cycle, MBD1 can interact with SETDB1 (H3K9 methyltransferase).

Immunogen: This MBD1 antibody was raised against a synthetic peptide corresponding to human MBD1.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

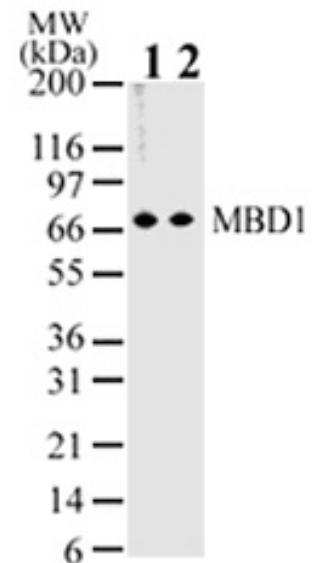
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. 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.



MBD1 antibody tested by Western blot. MBD1 detection by Western blot. The analysis was performed using HeLa whole-cell extract and MBD1 antibody at 2 µg/ml (lane 1) and a 1 µg/ml (lane 2) dilutions.