

MBD2 antibody (pAb)

Catalog Nos: 39547, 39548

RRID: AB_2615025

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.79 µg/µl

Molecular Weight: 48 kDa

Background: MBD2 (methyl-CpG binding domain protein 2) 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. MBD2 is a transcriptional repressor which is associated with histone deacetylases (HDAC) in the Sin3 and the MeCP1 complexes in mammalian cells. The HDAC complexes, also containing MBD3 and MeCP2, are involved in packaging the genomic DNA into inactive chromatin.

Immunogen: This MBD2 antibody was raised against a synthetic peptide corresponding to human MBD2.

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:

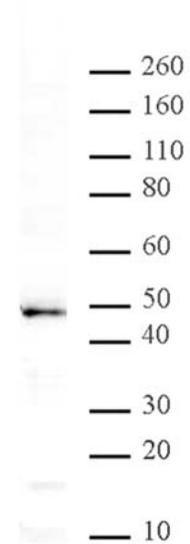
Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 dilution

Individual optimization may be required.

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.



MBD2 pAb tested by Western blot.

The analysis was performed using HeLa nuclear extract and MBD2 pAb at a 1:500 dilution.