

MBD4 antibody (pAb)

Catalog No: 39217

RRID: AB_2793187

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Concentration: 0.5 µg/µl

Molecular Weight: 64 kDa

Background: 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. MBD4 (methyl-CpG binding domain protein 4) is a thymine glycosylase that recognizes the product of deamination at methyl-CpG sites, as a part of DNA repair system. MBD4 is able to bind the hemimethylated DNA or methyl-CpG TpG mismatches.

Immunogen: This MBD4 antibody was raised against a mixture of synthetic peptides corresponding to amino acid residues 268-282 and 337-352 of human MBD4.

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

Application Notes:

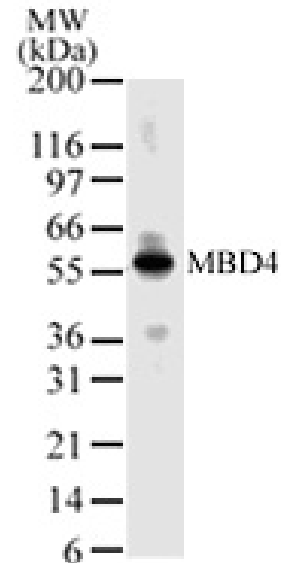
Applications Validated by Active Motif:

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.



MBD4 pAb tested by Western blot.

The analysis was performed using HL-60 whole-cell extract and MBD4 pAb at a 2 µg/ml dilution.