

MeCP2 antibody (pAb)

Catalog Nos: 39188, 39189

RRID: AB_2793176

Isotype: Serum

Application(s): WB

Reactivity: Human, Mouse

Volumes: 200 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 75 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. MeCP2 (methyl-CpG binding protein 2) can repress transcription by recruiting the Sin3 complex including histone deacetylases (HDACs). MeCP2 is distributed throughout the nucleus in human cells.

Immunogen: This MeCP2 antibody was raised against recombinant protein corresponding to the amino-terminal 320 amino acids of mouse MeCP2.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

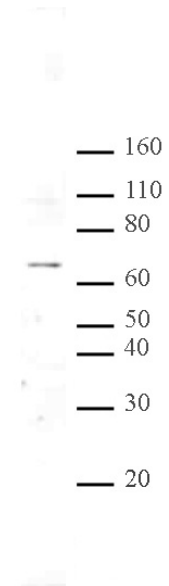
Applications Validated by Active Motif:

WB*: 1:10,000 - 1:50,000 dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



MeCP2 pAb tested by Western blot.

HeLa acid extract (10 µg per lane) was probed with MeCP2 pAb (1:25,000 dilution).