

DNMT1 antibody (mAb)

Catalog No: 39204**RRID:** AB_2614950**Clone:** 60B1220.1**Application(s):** ChIP, IHC, IP, WB**Reactivity:** Human, Mouse**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Mouse**Isotype:** IgG**Concentration:** 1 µg/µl**Molecular Weight:** 190 kDa

Background: DNMT1 (DNA (cytosine-5)-methyltransferase) proteins are involved in DNA methylation in which a methyl group is added to a cytosine residue on DNA, commonly at the C5 position of a CpG dinucleotide. Three families of DNMTs have been identified: DNMT1, DNMT2, and DNMT3 (comprised of DNMT3A and DNMT3B).

DNMT1 is the most abundant DNMT in somatic cells and shows a preference for methylating hemi-methylated DNA. It is considered to be a maintenance DNA methyltransferase that is important in the maintenance of specific patterns of methylation throughout cellular divisions. Methylation of mammalian DNA has long been recognized to play a major role in a number of cellular functions such as embryonic development, genetic imprinting, X chromosome inactivation and the control of gene expression. DNA methylation is generally associated with transcriptional repression.

Immunogen: This DNMT1 antibody was raised against a peptide corresponding to amino acid residues 637-651 of human DNMT1.

Buffer: PBS containing 0.05% BSA and 0.05% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB: 2 - 4 µg/ml dilution

For DNMT1, we also offer AbFlex® DNMT1 Recombinant Antibody (rAb). For details, see Catalog No. 91175.

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.

DNMT1 antibody (mAb) tested by Western blot.

Detection of DNMT1 by Western blot. The analysis was performed using mouse ES whole-cell extract and DNMT1 antibody (mAb) at a 2 µg/ml dilution.