

DNMT3B antibody (mAb)

Catalog No: 39207

Clone: 52A1018

Isotype: IgG1k

Application(s): ChIP,

Reactivity: Human, Mouse

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Quantity: 100 µg

Molecular Weight: 96 kDa

Background: DNMT3B – DNA methyltransferase (or DNMT 3B) 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.

The DNMT3 family contains two active methyltransferases, DNMT3A & DNMT3B, and one DNMT3-Like protein (DNMT3L). DNMT3B has been shown to be important in the regulation of specific patterns of DNA methylation, specifically *de novo* methylation. 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. Mutations in the DNMT3B gene cause the immunodeficiency-centromeric instability-facial anomalies (ICF) syndrome.

Immunogen: This DNMT3B antibody was raised against a full-length protein corresponding to mouse DNMT3B.

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

Application Notes:

Published Applications:

ChIP

See references for more information. Individual optimization may be required.

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.