

DNMT2 antibody (pAb)

Catalog No: 39205

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.5 µg/µl

Quantity: 100 µg

Molecular Weight: 42 kDa

Background: DNMT2 – DNA methyltransferase 2 (or DNMT 2) 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).

DNMT2 lacks the large N-terminal regulator domain common to other eukaryotic methyltransferases, but it possess a catalytic domain. It is involved in the methylation of aspartic acid transfer RNA at position 38, giving the enzyme the alternative name of TRDMT1. Although referred to as a DNA methyltransferase, DNMT2 does not methylate DNA, but instead is the first RNA cytosine methyltransferase to be identified.

Immunogen: This DNMT2 antibody was raised against a peptide corresponding to amino acid residues 39-53 and 361-376 of mouse DNMT2.

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

Application Notes:

Validated Applications:

WB: 2 - 4 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all Blotto solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Antibodies in solution can be stored at -80°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.

