

# DNMT2 Polyclonal Antibody

(Catalog # A-1002)

## Background

This gene is specifically involved in the methylation of aspartic acid transfer RNA at the cytosine position 38. Although it is considered part of the family of DNA cytosine methyltransferases, DNMT2 does not methylate CpG in genomic DNA because it lacks the large N-terminal regulatory domain found in other eukaryotic methyltransferases.

## Description

DNMT2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Liquid. PBS with 0.02% sodium azide, pH7.3.

## Specificity

Mouse, Human

## Isotype

IgG

## Uniprot ID

O14717

## Purification

Affinity Purified

## Immunogen

Synthetic Peptide of Human TRDMT1

## Storage

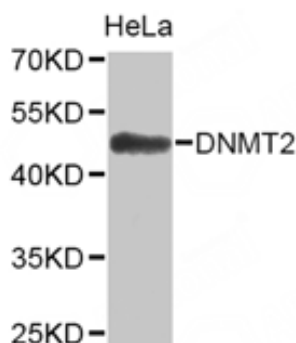
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

## Alternative Names

dDNMT antibody, DmMT 2 antibody, DNA (cytosine-5)-methyltransferase-like protein 2 antibody, DNA 5 cytosine methyltransferase antibody, DNA methyltransferase 2 antibody, DNA methyltransferase homolog HsallP antibody, DNA MTase homolog HsallP antibody, Dnmt 2 antibody, MHsallP antibody, nmt 2 antibody, OTTHUMP00000045198 antibody, PuMet antibody, RNMT 1 antibody, TRDMT 1 antibody, TRDMT\_HUMAN antibody, tRNA (cytosine 5 ) methyltransferase antibody, tRNA (cytosine(38)-C(5))-methyltransferase antibody, tRNA aspartic acid methyltransferase 1 antibody

## Application

IHC, WB; Recommended dilution: WB 1:500 - 1:2000 IHC 1:50 - 1:200



Western blot analysis of extracts of HeLa cellline, using DNMT2 Polyclonal Antibody.