
DNMT3A Polyclonal Antibody

(Catalog # A-1003)

Background

CpG methylation is an epigenetic modification that is important for embryonic development, imprinting, and X-chromosome inactivation. Studies in mice have demonstrated that DNA methylation is required for mammalian development. This gene encodes a DNA methyltransferase that is thought to function in de novo methylation, rather than maintenance methylation. The protein localizes to the cytoplasm and nucleus and its expression is developmentally regulated. Alternative splicing results in multiple transcript variants encoding different isoforms.

Description

DNMT3A Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

Q9Y6K1

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 400-500 of human DNMT3A (NP_072046.2)

Storage

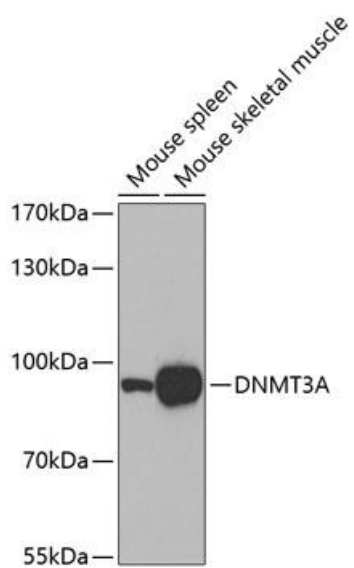
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

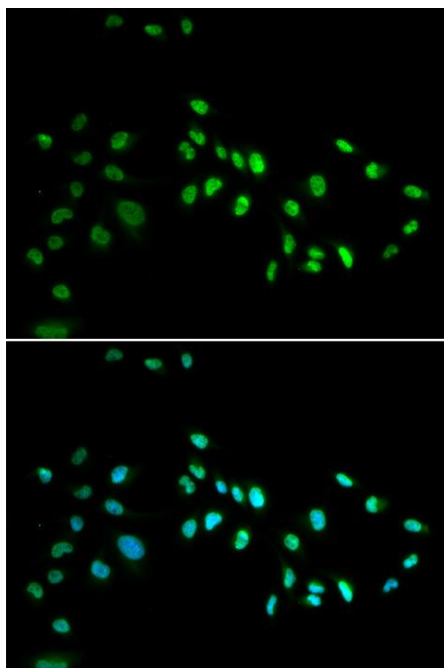
DNA (cytosine 5) methyltransferase 3 alpha antibody, DNA (cytosine-5)-methyltransferase 3A antibody, DNA cytosine methyltransferase 3A2 antibody, DNA methyltransferase 3a antibody, DNA methyltransferase HsaIIIA antibody, DNA MTase HsaI antibody, DNA MTase HsaIIIA antibody, DNMT3A_HUMAN antibody, DNMT 3a antibody, DNMT antibody, DNMT3A2 antibody, M HsaIIIA antibody, MCMT antibody, Methyl CpG binding domain protein 3a antibody, OTTHUMP00000201149 antibody, DNMT3A2, M.HsaIIIA

Application

WB, IF; Recommended dilution, WB 1:500 - 1:2000, IF 1:50 - 1:20



Western blot analysis of extracts of various cell lines, using DNMT3A antibody.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of A549 cells using DNMT3A antibody.