
DNMT3B Polyclonal Antibody

(Catalog A-1004)

Background

DNMT3B has been demonstrated to methylate both unmethylated and hemimethylated DNA equally and is supposed to mediate de novo methylation together with Dnmt1. Increased activation or amount of Dnmt3b is believed to be involved in carcinogenesis, as well as other genetic and epigenetic diseases.

Description

DNMT3B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9UBC3

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 345-475 of human DNMT3B (Q9UBC3).

Storage

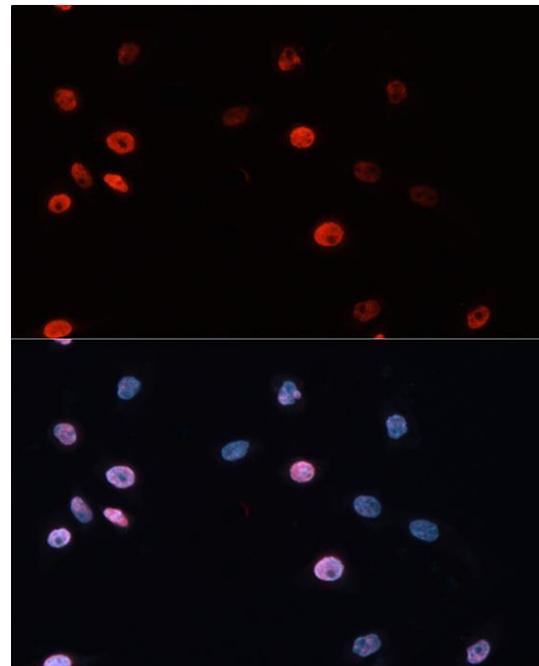
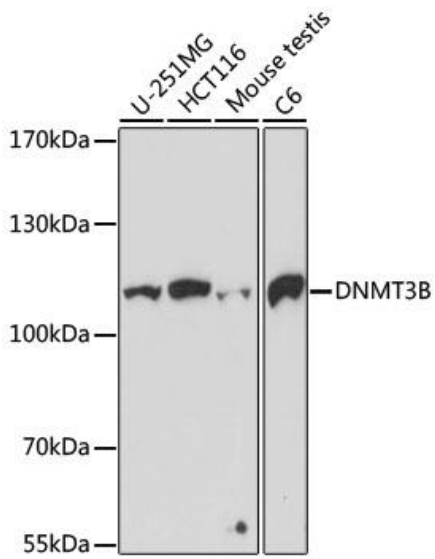
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

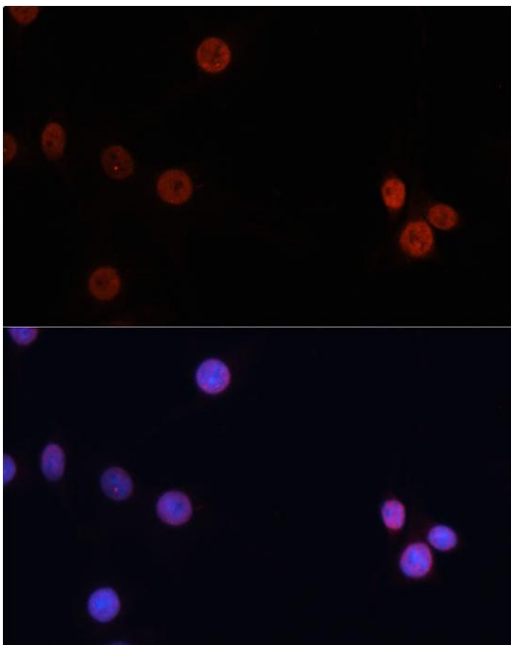
Cytosine 5methyltransferase 3B antibody, DNA (cytosine 5) methyltransferase 3 beta antibody, DNA (cytosine-5)-methyltransferase 3B antibody, DNA antibody, DNA methyltransferase HsallIB antibody, DNA MTase HsallIB antibody, DNMT3B_HUMAN antibody, Dnmt3b antibody, EC 2.1.1.37 antibody, ICF antibody, ICF1 antibody, M.HsallIB antibody, MGC124407 antibody, RP23-89H14.3 antibody

Application

WB, IF; Recommended dilution: WB: 1:200-1:1000, IF:1:50-1:200



Immunofluorescence analysis of HeLa cells using DNMT3B antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of C6 cells using DNMT3B antibody at dilution of 1:100. Blue: DAPI for nuclear staining.