

---

## 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 0.09% Sodium azide, 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 (NP\_008823.1).

### Storage

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

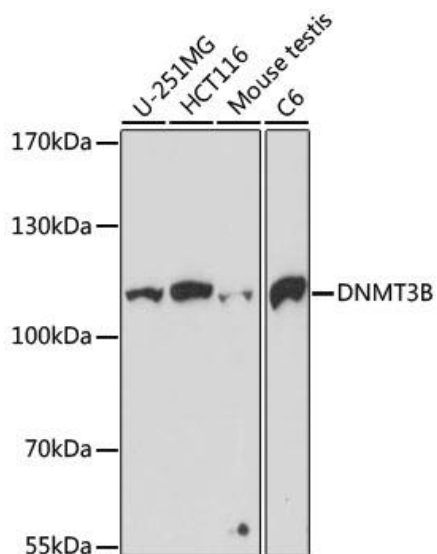
### Alternative Names

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, DNM3B\_HUMAN antibody, Dnmt3b antibody, EC 2.1.1.37 antibody, ICF antibody, ICF1 antibody, M.HsallIB antibody, MGC124407 antibody, RP23-89H14.3 antibody

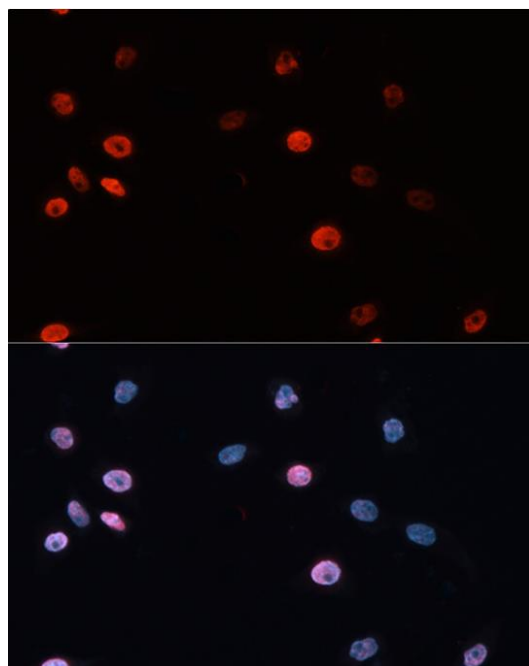
### Application

ELISA, WB, IF/ICC; Recommended dilution: WB: 1:200-1:1000, IF/ICC:1:50-1:200, ELISA - Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.

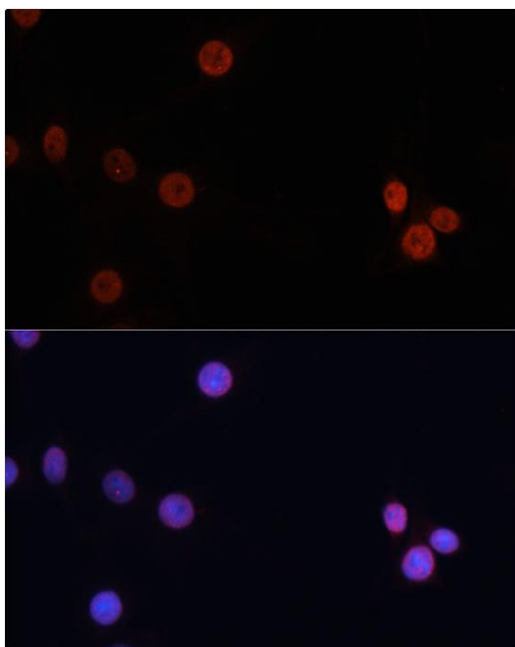
---



Western blot analysis of extracts of various lysates, using DNMT3B Polyclonal Antibody at 1:1000 dilution. 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. Exposure time: 3min.



Immunofluorescence analysis of HeLa cells using DNMT3B Polyclonal Antibody at dilution of 100 (40x lens). Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of C6 cells using DNMT3B Polyclonal Antibody at dilution of 100 (40x lens). Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.