
DNMT1 Polyclonal Antibody

(Catalog A-1700)

Background

DNA (cytosine-5-)-methyltransferase 1 has a role in the establishment and regulation of tissue-specific patterns of methylated cytosine residues. Aberrant methylation patterns are associated with certain human tumors and developmental abnormalities. Two transcript variants encoding different isoforms have been found for this gene.

Description

DNMT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human

Isotype

IgG

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 760-1109 of DNMT1 (NP_001370.1).

Storage

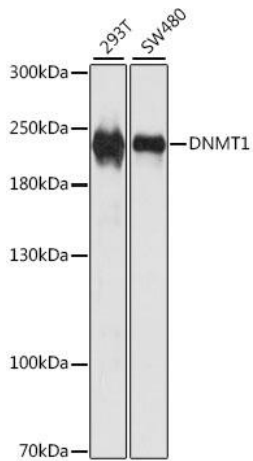
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

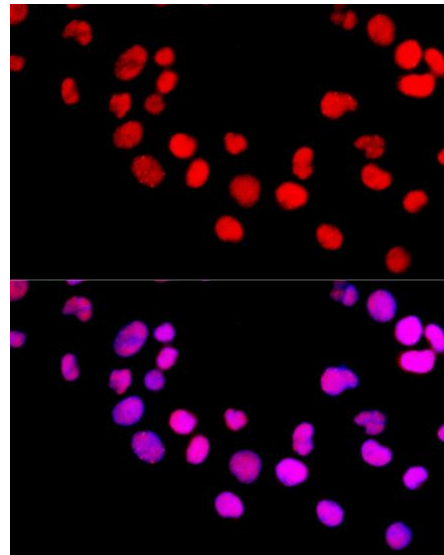
AIM, DNMT, MCMT, CXXC9, HSN1E, ADCADN

Application

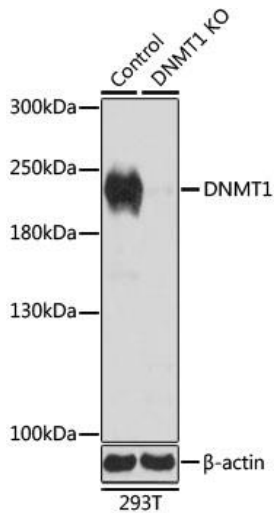
WB, IF, IP; Recommended dilution: WB: 1:500-1:2000, IF:1:50-1:200, IP:1:50-1:100



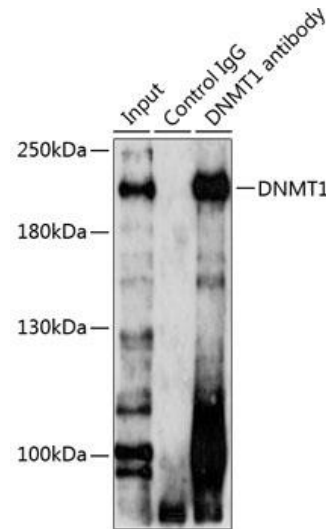
Western blot analysis of extracts of various cell lines, using DNMT1 antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of 293T cells using DNMT1 antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Western blot analysis of extracts from normal (control) and DNMT1 knockout (KO) 293T cells, using DNMT1 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.



Immunoprecipitation analysis of 200ug extracts of 293T cells using 3ug DNMT1 antibody. Western blot was performed from the immunoprecipitate using DNMT1 antibody at a dilution of 1:1000.