

Anti-DNMT3L

Rabbit Polyclonal Antibody

Catalog # **D353-63R**

Lot # B3216-50

Cited Applications

ELISA, IP, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the DNMT3L protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 152-164 of human DNMT3L (DNA (cytosine-5-)-methyltransferase-3-like).

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

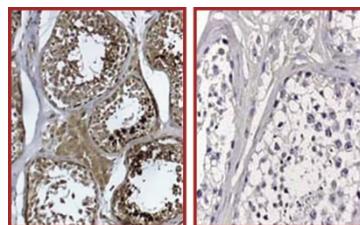
Scientific Background

DNMT3L is an enzymatically inactive regulatory factor that shares sequence homology with the DNA methyltransferase DNMT3A and DNMT3B. DNMT3L is involved in the establishment of DNA methylation patterns during gametogenesis. In mammals, DNA methylation is used for heritable silencing of retrotransposons and imprinted genes and for inactivation of the X chromosome in females (1). DNMT3L has a genomewide role in CpG island methylation beyond genomic imprinting (2). DNMT3L functions as a regulator of methylation at imprinted loci rather than a DNA cytosine methyltransferase because of a lack of catalytic motifs in its sequence.

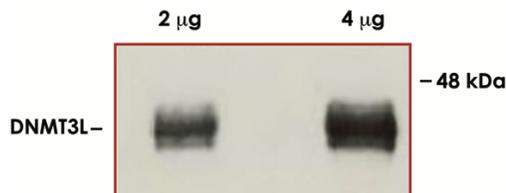
References:

1. Ooi, S.K.T. et al: DNMT3L connects unmethylated lysine 4 of histone H3 to de novo methylation of DNA. *Nature* 448: 714-717, 2007.
2. Smallwood, S. A. et al: Dynamic CpG island methylation landscape in oocytes and preimplantation embryos. *Nature Genet.* 43: 811-814, 2011.

Sample Data



Immunohistochemistry Rabbit Anti-DNMT3L 1:500



Western blot using affinity purified Anti-DNMT3L Antibody (1:1,500) after immunoprecipitation. Transfection of U2OS cells (10cm dish) was accomplished using 3 µg of GAL4-DNMT3L.

Anti-DNMT3L

Rabbit Polyclonal Antibody

Catalog #	D353-63R
Lot #	B3216-50
Purification	Immunoaffinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.