

DNMT3B polyclonal antibody

Other names: Dnmt3b, DNA MTase HsaIIIB, M.HsaIIIB

Cat. No. C15410218

Type: Polyclonal **ChIP-grade**

Source: Rabbit

Lot #: A2343-0040

Size: 50 µg/ 16 µl

Concentration: 3.2 µg/µl

Specificity: Human: positive

Other species: not tested

Purity: Affinity purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Polyclonal antibody raised in rabbit against mouse DNMT3B (DNA methyltransferase 3B), using 3 KLH-conjugated synthetic peptides containing sequences from different parts of the protein.

Applications

	Suggested dilution	Results
ELISA	1:10,000	Fig 1
Western blotting	1:1,000	Fig 2
Immunofluorescence	1:1,000	Fig 3

Target description

DNMT3B (UniProtKB/Swiss-Prot entry Q9UBC3) catalyses the genome wide de novo methylation of CpG residues, which regulates gene expression. DNMT3B is essential for development. DNA methylation on CpG residues is coordinated with methylation of histones. Six different isoforms of DNMT3B, produced by alternative splicing, exist although isoforms 4 and 5 may not be functional due to the absence of two conserved methyltransferase motifs.

Results

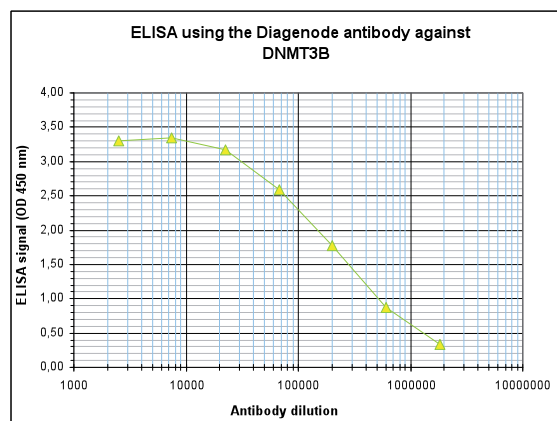


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of Diagenode antibody directed against DNMT3B (Cat. No. C15410218). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:220,000.

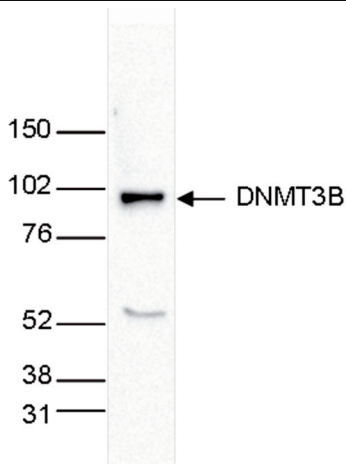


Figure 2. Western blot analysis using the Diagenode antibody directed against DNMT3B

Whole cell extracts (25 µg) from HeLa cells were analysed by Western blot using the Diagenode antibody against DNMT3B (Cat. No. C15410218) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

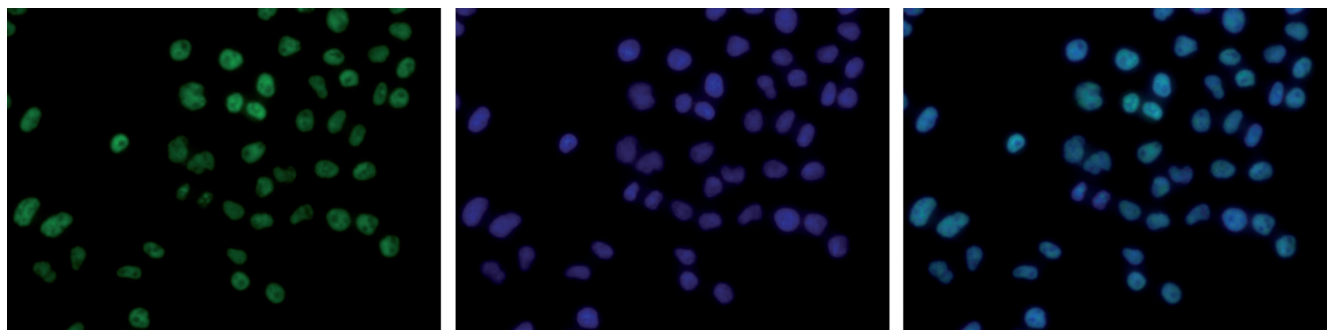


Figure 3. Immunofluorescence using the Diagenode antibody directed against DNMT3B

Human HeLa cells were stained with the Diagenode antibody against DNMT3B (Cat. No. C15410218) and with DAPI. Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labelled with the DNMT3B antibody (left) diluted 1:1,000 in blocking solution followed by an anti-rabbit antibody conjugated to Alexa488. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: April 18, 2014