

DNMT3B polyclonal antibody

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

Cat. No. C15410076 (pAb-076-005)

Type: Polyclonal **ChIP grade**

Source: Rabbit

Lot #: A16-0042

Size: 50 µg/ 50 µl

Concentration: 1 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 a KLH-conjugated synthetic peptide containing a sequence from the N-terminal part of the protein.

Applications

	Suggested dilution	Results
ChIP	1 µg per ChIP	Fig 1
ELISA	1:1,000	Fig 2

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

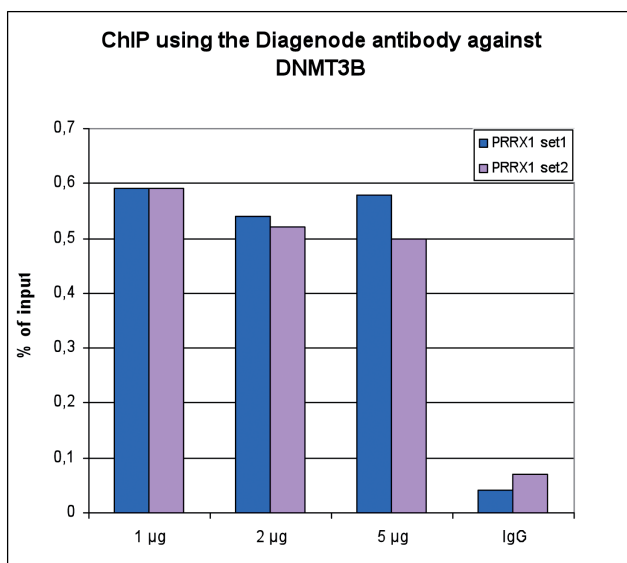


Figure 1. ChIP using the Diagenode antibody against DNMT3B

ChIP assays were performed using GDP control cells, the Diagenode antibody against DNMT3B (Cat. No. pAb-076-005) and optimized PCR primer sets for qPCR. ChIP was performed on sheared chromatin from 1 million cells using the “iDeal ChIP-seq” kit (Cat. No. AB-001-0024). A titration of the antibody consisting of 1, 2, and 5 µg per ChIP experiment was analysed. IgG (5 µg/IP) was used as negative IP control. QPCR was performed with two different primer pairs for the PRRX1 gene, used as a positive control target. Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).

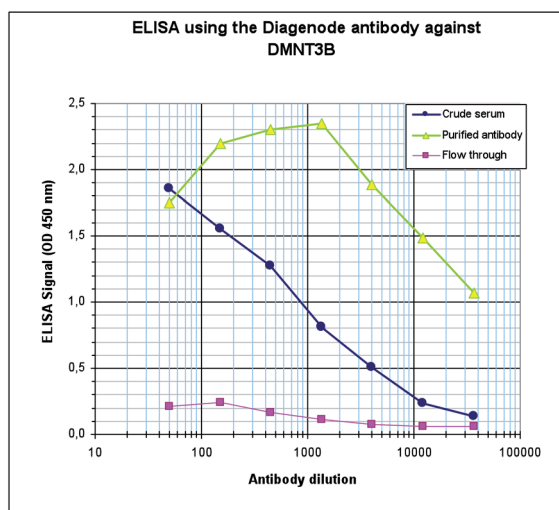


Figure 2. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of Diagenode antibody directed against mouse DNMT3B (Cat. No. pAb-076-005), crude serum and flow through in antigen coated wells. By plotting the absorbance against the antibody dilution (Figure 2), the titer of the antibody was estimated to be 1:32,400.

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: June 1, 2012