

PRODUCT NAME DNMT3B polyclonal antibody		
Other names: Dnmt3b, DNA MTase HsallIB, M.HsallIB		
Cat. No. C15310044 (CS-044-100)	Type: Polyclonal	Size: 100 µl
Lot #: A357-001	Source: Rabbit	Concentration: not determined

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 (1).

Specificity: Human and mouse: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:1,000	Fig 2

Purity: Whole antiserum from rabbit containing 0.05% azide.

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.

References:

(1) Peptide design by Andrea Kranz, Dept. of Genomics, Prof. F. Stewart, TU Dresden, Tatzberg 47/49, 01307 Dresden, Germany

Last data sheet update: April 20, 2010

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.

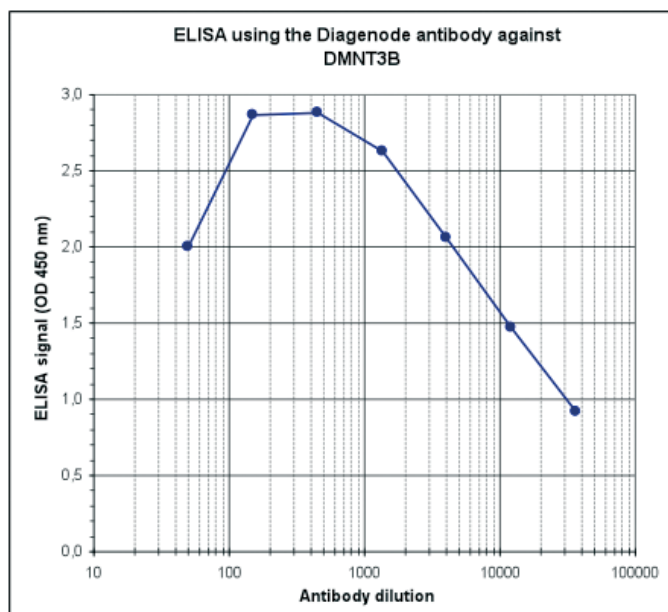


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse DNMT3B (cat# CS-044-100) in antigen coated wells. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:13,500.

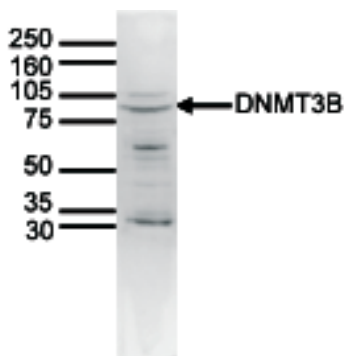


Figure 2
Western blot analysis using the Diagenode antibody directed against DNMT3B

Nuclear extracts from HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against DNMT3B (cat# CS-044-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the MW marker (in kDa) is shown on the left.