

PRODUCT NAME synthetic peptide: mouse DNMT3B			
(DNA (cytosine-5)-methyltransferase 3B)			
Cat. No: sp-076-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide corresponding to the N-terminal region of mouse DNMT3B.

This peptide was used as immunogen to raise the Diagenode ChIP grade antibody directed against mouse DNMT3B (cat# pAb-076-050). The same peptide can be used to block this antibody in negative control experiments.

For more information about DNMT3B: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (cat# pAb-076-100).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: Aliquot upon receipt and store at -20°C. Avoid multiple freeze-thaw cycles.

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

Availability date: August 14, 2008

Last data sheet update: August 20, 2008

Lot #: 001/ Synthesis date: September 15, 2005/ aliquoting date: August 18, 2008