

Synthetic peptide: mouse DNMT3B

(DNA (cytosine-5)-methyltransferase 3B)

Cat. No. C16000076

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 No. C15410076). 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 No. C15410076).

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