

## Synthetic peptide: mouse DNMT3B

(DNA methyltransferase 3B)

**Cat. No. C16000044**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide containing a sequence from the N-terminal part of the mouse DNMT3B protein. This peptide was used as immunogen to raise the Diagenode crude serum directed against mouse DNMT3B (Cat. No. C15310044). 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 crude serum (Cat. No. C15310044).

**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:** November 7, 2008

**Last data sheet update:** November 7, 2008