

## Synthetic peptide: human DNMT3A

(DNA methyltransferase 3A)

**Cat. No. C16000086**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide corresponding to amino acids 92 to 106 of human DNMT3A.

This peptide was used as immunogen to raise the Diagenode antibody directed against DNMT3A (Cat. No. C15410086). The same peptide can be used to block this antibody in negative control experiments.

For more information about DNMT3A: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15410086).

**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 20, 2008

**Last data sheet update:** August 14, 2008