

Synthetic peptide: mouse DNMT1

(DNA (cytosine-5)-methyltransferase 1)

Cat. No. C16000077

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide corresponding to the N-terminal region of mouse DNMT1.

This peptide was used as immunogen to raise the Diagenode crude serum directed against mouse DNMT1 (cat# CS-077-100). The same peptide can be used to block this antibody in negative control experiments.

For more information about DNMT1: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (cat# CS-077-100).

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 20, 2008

Last data sheet update: August 14, 2008