

Synthetic peptide: mouse DNMT2

(DNA methyltransferase 2)

Cat. No. C16000095

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 DNMT2 protein. This peptide was used as immunogen to raise the Diagenode crude serum directed against mouse DNMT2 (Cat. No. C15310095). The same peptide can be used to block this antibody in negative control experiments. For more information about DNMT2: see overview in the Technical Data Sheet of the corresponding Diagenode crude serum (Cat. No. C15310095).

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