

PRODUCT NAME			
synthetic peptide: mouse DNMT2			
(DNA methyltransferase 2)			
Cat. No: sp-095-050	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# CS-095-100). 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# CS-095-100).

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: November 7, 2008

Last data sheet update: November 7, 2008

Lot #: 001/ Synthesis date: August 17, 2007/ aliquoting date: November 5, 2008