

PRODUCT NAME			
synthetic peptide: human DNMT3A			
(DNA methyltransferase 3A)			
Cat. No: sp-086-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide corresponding to amino acids 107 to 121 of human DNMT3A.

This peptide was used as immunogen to raise the Diagenode antibody directed against DNMT3A (cat# pAb-086-050). 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# pAb-086-050).

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

Lot #: 001/ Synthesis date: April 13, 2006/ aliquoting date: August 18, 2008