

Synthetic peptide: human MBD1

(Methyl-CpG-binding domain protein 1)

Cat. No. C16000078

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide corresponding to the N-terminal region of human MBD1.

This peptide was used as immunogen to raise the Diagenode ChIP grade antibody directed against MBD1 (Cat. No. C15410078). The same peptide can be used to block this antibody in negative control experiments.

For more information about MBD1: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15410078).

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