

Synthetic peptide: mouse MeCP2

[Methyl-CpG-binding domain Protein 2]

Cat. No. C16000106

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 MeCP2 protein. This peptide was used as immunogen to raise the Diagenode crude serum directed against mouse MeCP2 (Cat No. C15310106). The same peptide can be used to block this antibody in negative control experiments. For more information about MeCP2: see overview in the Technical Data Sheet of the corresponding Diagenode crude serum (Cat No. C15310106).

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