

PRODUCT NAME			
synthetic peptide: C-terminal region of MeCP2			
(Methyl-CpG-binding domain protein 2)			
Cat. No: sp-052-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide derived from the C-terminus of the mouse MeCP2 protein.

This peptide was used as immunogen to raise the Diagenode ChIP-grade antibody directed against MeCP2 (cat# pAb-052-050). 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 our ChIP-grade antibody directed against MeCP2 (cat# pAb-052-050).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: For long storage, aliquot 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 12, 2007

Last data sheet update: November 12, 2007

Lot #: 001/ day of the synthesis: September 8, 2005/ aliquoting day: November 12, 2007