

PRODUCT NAME		
MeCP2 polyclonal antibody		
Other names: AUTSX3, MRX16, MRX79, MRXS13, MRXSL, PPMX, RTS, RTT		
<b>Cat. No.</b> C15310106 (CS-106-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A375-001	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse MeCP2 (Methyl-CpG-binding domain Protein 2), using a KLH-conjugated synthetic peptide containing an amino acid sequence from the N-terminal part of the protein.

**Specificity:** Mouse: positive  
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:500 – 1:5,000	Fig 1
Western blotting	1:1,000	Fig 2

**Purity:** Whole antiserum from rabbit containing 0.05% azide.

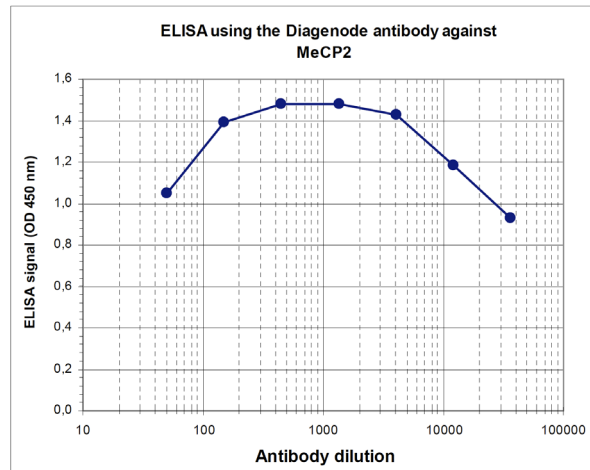
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** April 7, 2010

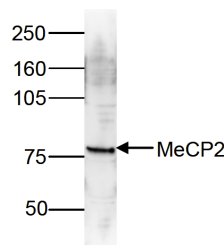
**Target description**

MeCP2 (UniProt/Swiss-Prot entry P51608) is a chromosomal protein with abundant binding sites in chromatin. It belongs to the family of methyl CpG binding proteins which also comprises MBD1, MBD2, MBD3 and MBD4. MeCP2 can bind specifically to methylated promoters, thereby repressing transcription. This transcriptional repression is mediated through interaction with histone deacetylase and the corepressor SIN3A. MeCP2 is essential for development. Mutations in MeCP2 are the cause of several types of mental retardation including Rett syndrome, a progressive neurological disorder that causes mental retardation in females, and mental retardation syndromic X-linked type 13, and may also be involved in Angelman syndrome and susceptibility to some types of autism.



**Figure 1**  
**Determination of the titer**

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse MeCP2 [Cat. No. CS-106-100]. The wells were coated with the peptide used for immunisation of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:136,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against MeCP2**

Nuclear extracts of mouse fibroblasts (NH3T3, 40 µg) were analysed by Western blot using the Diagenode antibody against mouse MeCP2 [Cat. No. CS-106-100] diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.