

PRODUCT NAME		
MeCP2 polyclonal antibody		
Other names: AUTSX3, MRX16, MRX79, MRXS13, MRXSL, PPMX, RTS, RTT		
Cat. No. C15310088 (CS-088-100)	Type: Polyclonal	Size: 100 µl
Lot #: A131-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human MeCP2 (Methyl-CpG-binding domain protein 2), using 3 different KLH-conjugated synthetic peptides containing an amino acid sequence from the N-terminal, the central and the C-terminal part of the protein, respectively.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:2,000	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 also 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.

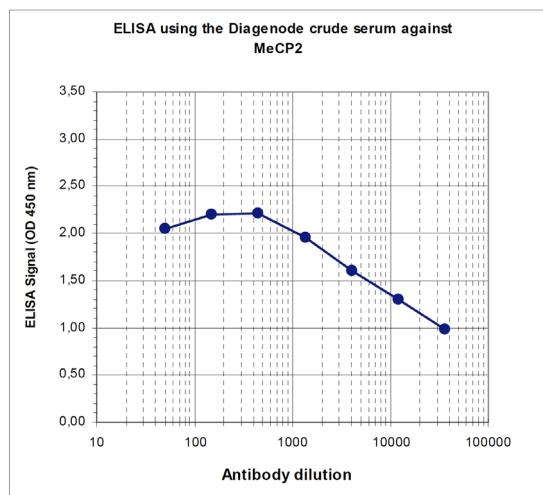


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human MeCP2 (Cat. No. CS-088-100), in antigen coated wells. 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 1:25,000.

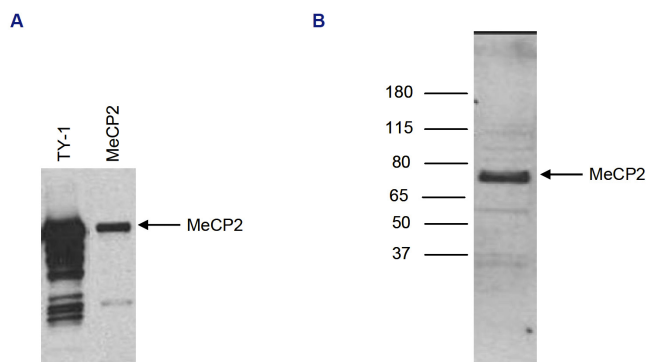


Figure 2
Western blot analysis using the Diagenode antibody directed against MeCP2

Figure 2A: Human osteosarcoma cells (U2OS) were transfected with an expression vector for TY1-tagged MeCP2. The presence of the TY1-MeCP2 in the cell lysates was demonstrated by western blot analysis with the Diagenode antibody against the TY1-tag (Cat. No. MAb-054-050) (lane 1) and with the antibody against MeCP2 (Cat. No. CS-088-100) (lane 2), diluted 1:2,000 in TBST containing 3% milk powder.

Figure 2B: Western blot was performed on nuclear extracts from the human leukemic monocyte lymphoma cell line U937 (60 µg) with the Diagenode antibody against human MeCP2 (Cat. No. CS-088-100), diluted 1:2,000 in TBST containing 3% milk powder. A molecular weight marker (in kDa) is shown on the left. The location of the protein of interest is indicated on the right.