

MeCP2 monoclonal antibody

Other name: AUTSX3, MRX16, MRX79, MRXS13, MRXSL, PPMX, RTS, RTT

Cat. No. C15200225

Type: Monoclonal

Source: Mouse

Lot #: 001

Size: 50 µg/50 µl

Concentration: 1 µg/µl

Specificity: Human, mouse: positive
Other species: not tested

Purity: Affinity purified monoclonal antibody in PBS. Does not contain any preservative.

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

Description: Monoclonal antibody raised in mouse against MeCP2 (Methyl-CpG-binding domain protein 2), using a recombinant protein.

Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1
Immunofluorescence	1:400	Fig 2

Target description

MeCP2 (UniProt/Swiss-Prot entry P51608) is a chromosomal protein with abundant binding sites in the 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.

Results

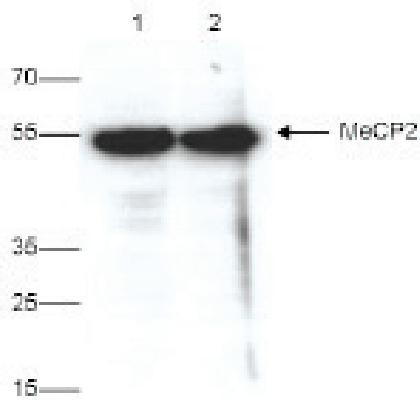


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against MeCP2

Whole cell extracts from HeLa (lane 1) or MEF (lane 2) cells were analysed by Western blot using the Diagenode antibody against MeCP2 (Cat. No. C15200225) diluted 1:1,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

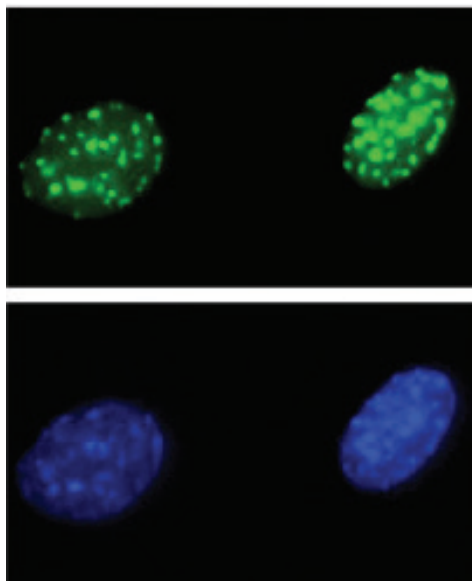


Figure 2. Immunofluorescence using the Diagenode monoclonal antibody directed against MeCP2

HeLa cells were stained with the Diagenode antibody against MeCP2 (Cat. No. C15200225) diluted 1:400 followed by an anti-mouse antibody conjugated to Alexa488 (upper figure). The bottom figure shows staining of the nuclei with DAPI.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: August 21, 2015