

MeCP2 polyclonal antibody

Cat. No. C15410217

Type: Polyclonal

Source: Rabbit

Lot #: A34-0020

Size: 50 µg/ 28 µl

Concentration: 1.8 µg/µl

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

Purity: Protein G purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300

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: Polyclonal antibody raised in rabbit against human MeCP2 (Methyl-CpG-binding domain protein 2), using a recombinant protein.

Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1

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

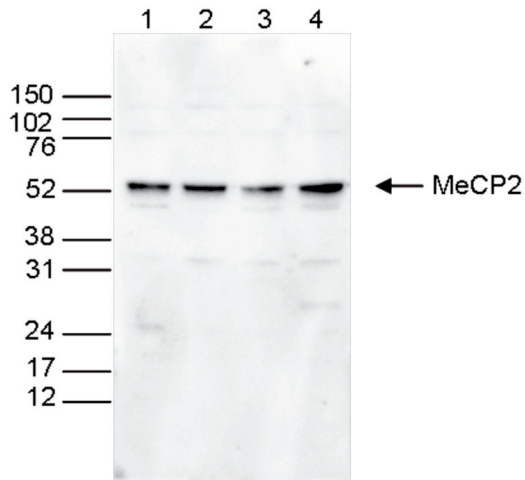


Figure 1. Western blot analysis using the Diagenode antibody directed against MeCP2

Whole cell extracts (25 µg) from HeLa (lane 1), MCF7 (lane 2), K562 (lane 3) and NIH3T3 cells (lane 4) were analysed by Western blot using the Diagenode antibody against MeCP2 (Cat. No. 15410217) 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.

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: January 3, 2014