

MBD2 polyclonal antibody

Other names: DMTase

Cat. No. 15410312

Type: Polyclonal

Source: Rabbit

Lot #: A38-0020

Size: 50 µg/ 42 µl

Concentration: 1.2 µg/µl

Specificity: Human: 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 mouse MBD2 (methyl-CpG binding domain protein 2), using a recombinant protein.

Applications

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

Target description

MBD2 (UniProtKB/Swiss-Prot entry Q9UBB5) belongs to the family of methylated DNA binding proteins. The protein acts as a transcriptional repressor that specifically binds to methylated CpG dinucleotides in promoter sequences. It is part of the MeCP1 complex that also contains the histone deacetylases HDAC1 and HDAC2. On the other hand, MBD2 may be able to demethylate DNA, thus acting as a transcriptional activator. Possibly, MBD2 mediates the effects of DNA methylation in vivo.

Results

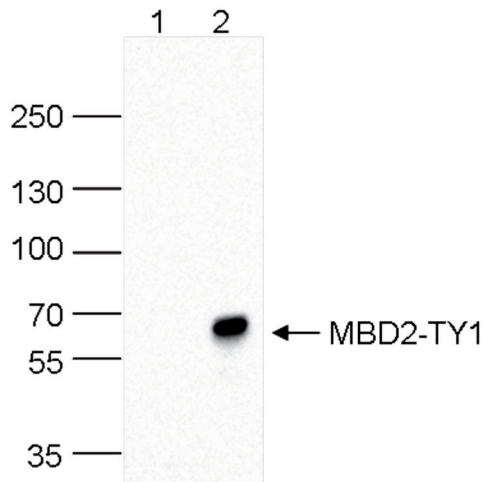


Figure 1. Western blot analysis using the Diagenode antibody directed against MBD2

Whole cell extracts (40 µg) from HeLa cells, transfected with a TY1-tagged MBD2 (lane 2) or an untransfected negative control (lane 1) were analysed by western blot using the Diagenode antibody against MBD2 (Cat. No. 15410312) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the TY1-tagged MBD2 protein 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: November 28, 2014