

MBD4 recombinant protein

Other names: MED1

Cat. No. C23010108

Source: SF9 cells

Lot #: 001

Size: 10 µg/ 20 µl

Concentration: 0.5 µg/µl

Specificity: Human

Purity: Purified using FPLC, >95% purity as determined by SDS-PAGE.

Storage buffer: 20 mM Tris-Cl pH 7.6, 1 mM EDTA, 0.15 M NaCl, 10% glycerol, 0.5 mM PMSF and 1 mM DTT.

Storage: Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

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

Description: Active full length recombinant human N-terminal His tagged MBD4 (Methyl-CpG Binding Domain Protein 4) protein, produced in SF9 insect cells.

Protein description

MBD4 (UniProt/Swiss-Prot entry O95243) belongs to the family of methyl CpG binding proteins that specifically bind to methylated CpG dinucleotides. MBD4 probably functions to mediate the biological consequences of CpG methylation. Further, MBD4 has thymine glycosylase activity and is involved in DNA mismatch repair. It is specific for G/T mismatches within methylated and unmethylated CpG sites. Mutations of MBD4 are associated with tumors with primary microsatellite-instability, a type of genomic instability that is caused by a defective DNA mismatch repair.

Quality control

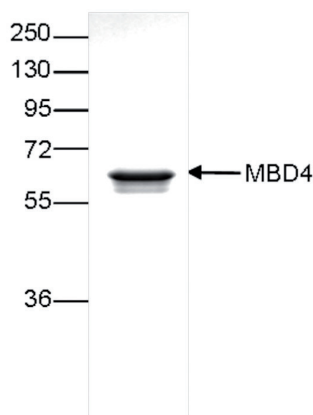


Figure 1.

SDS page of the MBD4 recombinant protein. 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: August 19, 2015