

## DNMT1 recombinant protein

**Other names:** DNMT, MCMT, CXXC9, AIM

**Cat. No.** C23010114

**Source:** Sf9 cells

**Lot #:** 001

**Size:** 25 µg/ 50 µl

**Concentration:** 0.5 µg/µl

**Specificity:** Human

**Purity:** Purified using FPLC, >98% purity as determined by SDS-PAGE

**Storage buffer:** 20 mM Tris-Cl pH 7.6, 100 mM EDTA, 0.15 M NaCl, 0.1% NP40 and 10% glycerol.

**Storage:** Store at -80°C; avoid multiple freeze-thaw cycles. Guaranteed stable for 6 months from date of receipt when stored properly.

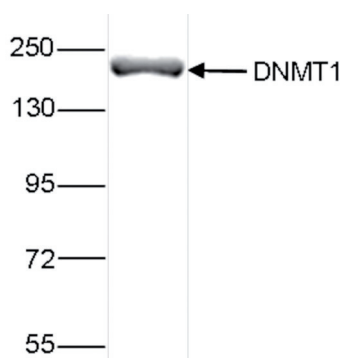
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Full length recombinant human N-terminal His tagged and C-terminal FLAG tagged DNMT1 (DNA (cytosine-5)-methyltransferase 1) protein, produced in Sf9 insect cells.

### Protein description

DNMT1 (UniProt/Swiss-Prot entry P26358) preferentially methylates CpG residues in hemimethylated DNA and is responsible for maintaining methylation patterns established in development. DNA methylation is coordinated with methylation of histones. Inactivation of DNMT1 causes global demethylation and embryonic lethality. Transcriptional repression by DNMT1 is mediated by direct interaction with HDAC2.

### Quality control



**Figure 1.**

SDS page of the DNMT1 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:** February 15, 2016