

DNMT1 Protein

Recombinant mouse protein expressed in Sf9 cells

Catalog # D351-31G

Lot # F422-2

Product Description

Recombinant mouse DNMT1 (766-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC053047](#).

Gene Aliases

Dnmt; MCMT; Cxhc9; MTase; Met-1; Dnmt1o; m.Mmul; Met1; MommeD2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

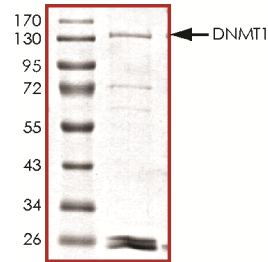
DNMT1 has a role in the establishment and regulation of tissue-specific patterns of methylated cytosine residues. Aberrant methylation patterns are associated with certain human tumors and developmental abnormalities (1). DNMT1 enzyme is crucial for stable expression and function in vivo (2). DNMT1 is essential for maintenance of DNA methylation and the interaction of DNMT1 with the replication machinery is not strictly necessary for maintenance of DNA methylation, but improves its efficiency. DNMT1 is necessary and sufficient to maintain global methylation and aberrant CpG island methylation in human cancer cells.

References

1. Yoder, J. A. et.al: New 5-prime regions of the murine and human genes for DNA (cytosine-5)-methyltransferase. J. Biol. Chem. 271: 31092-31097, 1996.
2. Tucker, K. L. et.al: Complementation of methylation deficiency in embryonic stem cells by a DNA methyltransferase minigene. Proc. Nat. Acad. Sci. 93: 12920-12925, 1996.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity of DNMT1 protein was determined to be **>70%** by densitometry. Approx. MW **132kDa**.

DNMT1 Protein

Recombinant mouse protein expressed in Sf9 cells

Catalog Number	D351-31G
Specific Lot Number	F422-2
Purity	>70%
Concentration	0.05 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.