

DNMT3A, Active

Full length recombinant protein expressed in Sf9 cells

Catalog # D353-380G

Lot # O1059-6

Product Description

Recombinant full-length human DNMT3A was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC043617](#).

Gene Aliases

DNMT3A2; M.Hsa111A

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.

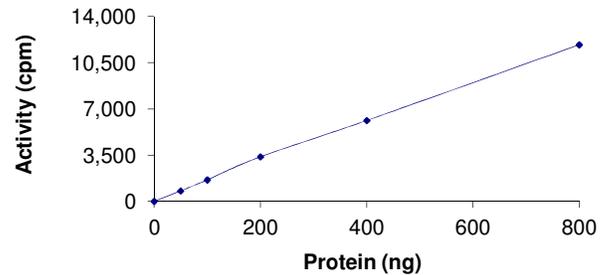
Scientific Background

DNMT3A: DNA (cytosine-5-)-methyltransferase 3 alpha plays important role in genomic imprinting and X chromosome inactivation and is essential for mammalian development. DNMT3A appears to function as a de novo methyltransferase, since it can methylate unmethylated and hemimethylated DNA with equal efficiencies (1). DNMT3A is required for genome wide de novo methylation and is essential for mammalian development (2). DNMT3A is critical for maintaining active chromatin states of genes that are essential for development.

References

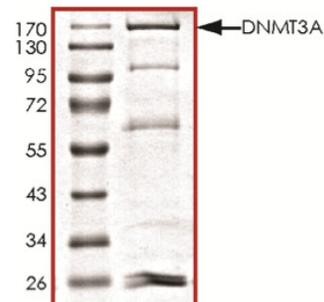
1. Yanagisawa, Y. et al: The human DNA methyltransferases DNMT3A and DNMT3B have two types of promoters with different CpG contents. *Biochim. Biophys. Acta* 1577: 457-465, 2002.
2. Okano, M. et al: DNA methyltransferase Dnmt3a and Dnmt3b are essential for de novo methylation and mammalian development. *Cell* 99: 247-257, 1999.

Specific Activity



The specific activity of DNMT3A was determined to be **250 pmol/min/mg** as per activity assay protocol.

Purity



The purity of DNMT3A was determined to be **>70%** by densitometry, Approx. MW **162 kDa**.

DNMT3A, Active

Recombinant human protein expressed in Sf9 cells

| | |
|---------------------|--|
| Catalog Number | D353-380G |
| Specific Activity | 250 pmol/min/mg |
| Specific Lot Number | O1059-6 |
| Purity | >70% |
| Concentration | 0.1 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. |

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

