

DNMT3a (DNA (cytosine-5)-methyltransferase 3a)

CATALOG NO.: DMT-21-125

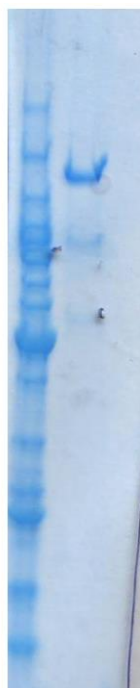
LOT NO.:

DESCRIPTION: Full-length human recombinant DNMT3a (residues 2-912 (C-term.); Genbank Accession # NM_175629; MW = 127 kDa) expressed in *Sf9* insect cells with an N-terminal GST-tag. DNA (cytosine-5)-methyltransferase 3a (DNMT3a) is required for *de novo* DNA methylation and methylates at both CpG and non-CpG sites.

PURITY: >80% by SDS-PAGE.

SUPPLIED AS: ___ µg/µl total protein in 50 mM Tris/HCl pH 8.0, 125 mM NaCl, 3 mM DTT, 10% glycerol (v/v).

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. Freezing and storage of diluted enzyme is not recommended.



Coomassie blue stained SDS-PAGE of 6 µg of RBC DNMT3a. MWs of markers at left are, from top (kDa): 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15 & 10.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.