

Recombinant DNMT3L protein

Catalog No: 31414, 31814

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.55 µg/µl

Source: Human

Buffer Contents: Full length recombinant DNMT3L protein is supplied at a concentration of 0.55 µg/µl in 25 mM Tris pH 8.0, 300mM NaCl, 5% Glycerol and 0.04% Triton X-100, 1mM DTT.

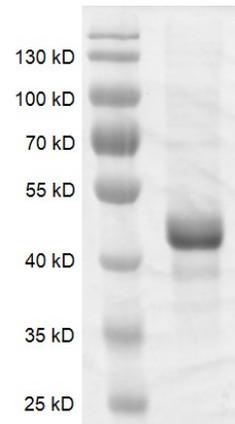
Background: DNMT3L (DNA (Cytosine-5-)-Methyltransferase 3-Like) is a nuclear protein with similarity to the DNMT family of proteins, but does not function as a DNA methyltransferase. It serves as an essential regulator of DNMT3A and DNMT3B, recruiting them to sites for deposition of de novo DNA methylation. DNMT3L is expressed in the presumptive gametes in mammalian embryos and is required for the establishment of genomic imprinting. DNMT3L has a PHD-like domain that allows binding to chromatin through interactions with the amino terminal tail of histone H3 that is devoid of lysine 4 methylation.

Protein Details: Full length recombinant DNMT3L protein (accession number NP_787063.1) was expressed in Sf9 cells and contains an N-terminal 6×His with a molecular weight of 46.6 kDa. The recombinant protein is >80% pure by SDS-PAGE.

Application Notes: Recombinant DNMT3L protein is suitable for use in reconstituting DNMT3 complexes, functional studies.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant DNMT3L protein gel.

Recombinant DNMT3L protein was run on an 8% SDS-PAGE gel and stained with Coomassie blue.