

Recombinant DNMT3B (212-358) protein

Catalog No: 31542, 31942

Lot No: 09115001

Expressed In: *E. coli*

Quantity: 100 µg

Concentration: 0.8 µg/µl

Source: Human

Buffer Contents: Recombinant DNMT3B (212-358) was expressed in *E. coli* cells at a concentration of 0.8 µg/µl in 25 mM Tris pH 7.4, 150 mM NaCl, 5% glycerol.

Background: DNMT3B (DNA (Cytosine-5-)-Methyltransferase 3 Beta) protein is a *de novo* DNA methyltransferase. It is involved in DNA methylation in which a methyl group is added to a cytosine residue on DNA, commonly at the C5 position of a CpG dinucleotide. Methylation of mammalian DNA has long been recognized to play a major role in a number of cellular functions such as embryonic development, genetic imprinting, X chromosome inactivation and control of gene expression. DNA methylation is generally associated with transcriptional repression.

Trimethylation of Lys36 on histone H3 tail (H3K36me3) is associated with DNA methylation and elongation phase of transcription. PWWP domain of DNMT3B reads this epigenetic mark to guide DNA methylation.

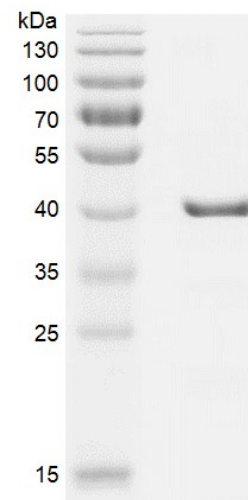
Protein Details: The peptide corresponding to amino acids 212 - 358 that contains the PWWP sequences of DNMT3B (accession number NP_715640.2) was expressed in *E. coli* and contains an N-terminal GST tag with an observed molecular weight of 42.4 kDa. The recombinant protein is >90% pure by SDS-PAGE.

Application Notes: Recombinant DNMT3B (212-358) is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

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.

DNMT3B (212-358)



Recombinant DNMT3B (212-358) gel. DNMT3B (212-358) was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.