

Recombinant GNMT protein

Catalog No: 81301, 81601

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 3.2 µg/µl

Source: Human

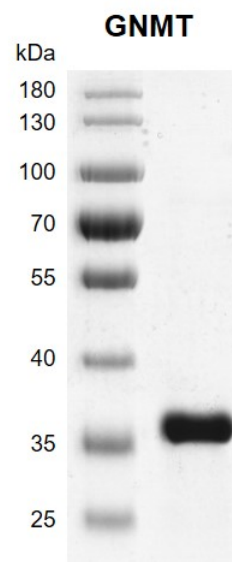
Buffer Contents: Recombinant GNMT protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: GNMT (Glycine N-Methyltransferase) is an enzyme that catalyzes the conversion of S-adenosyl-L-methionine (along with glycine) to S-adenosyl-L-homocysteine and sarcosine. This protein is found in the cytoplasm and acts as a homo-tetramer. Defects in this gene are a cause of GNMT deficiency (hypermethioninemia). Alternative splicing results in multiple transcript variants.

Protein Details: Full length human GNMT protein was expressed in *E. coli* cells (accession number NP_061833.1) with a C-terminal 6×His-Tag. The molecular weight of the protein is 34.9 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

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 arrival.



Recombinant GNMT protein gel
10% SDS-PAGE with Coomassie blue staining

MW: 34.9 kDa

Purity: >95%