

Recombinant MGMT protein

Catalog No: 81447, 81547

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 0.7 µg/µl

Source: Human

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

Background: MGMT(Methylated-DNA--protein-cysteine methyltransferase)

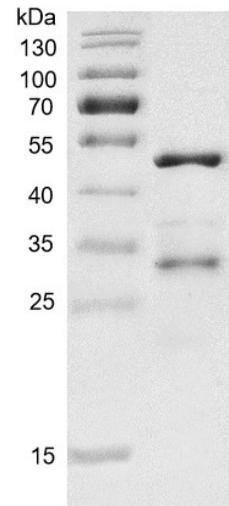
Involved in the cellular defense against the biological effects of O6-methylguanine (O6-MeG) and O4-methylthymine (O4-MeT) in DNA. Repairs the methylated nucleobase in DNA by stoichiometrically transferring the methyl group to a cysteine residue in the enzyme. This is a suicide reaction: the enzyme is irreversibly inactivated.

Protein Details: Recombinant MGMT protein that includes full length of human MGMT protein (accession number NP_002403.3) was expressed in *E. coli* cells with an N-terminal GST tag. The molecular weight of the protein is 48.48 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

MGMT



Recombinant MGMT protein

12.5% SDS-PAGE Coomassie staining

MW: 48.48 kDa

Purity: ≥ 85%