

Recombinant EHMT2 (G9A) H904K protein

Catalog No: 31328

Lot No: 21509001

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 2 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant EHMT2 (G9A) H904K protein supplied at a concentration of 2 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Protein Details: EHMT2 (G9A) H904K is a histone methyltransferase that is specific to lysine 9 of histone H3 (H3K9). Methylation of H3K9 is a signal of transcriptional repression. Recombinant EHMT2 (G9A) H904K was expressed as amino acids 530-1210 of human protein (accession number NP_006700.3) with an N-terminal His-Tag. The recombinant EHMT2 (G9A) H904K protein contains a substitution of histidine for lysine at amino acid 904. EHMT2 (G9A) H904K is catalytically inactive.

Application Notes: Recombinant EHMT2 (G9A) H904K is suitable for use in methyltransferase assays. A recommended starting point for the methyltransferase assay is 50-200 ng of recombinant protein per reaction. The molecular weight of the protein is ~81 kDa. This is a version of the protein with a catalytic site mutant, so it is designed for use as a negative control.

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.