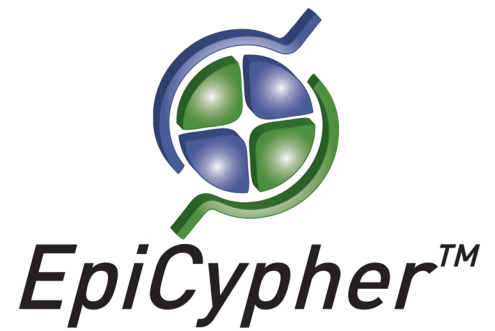


G9a, Recombinant Human, His-Tagged

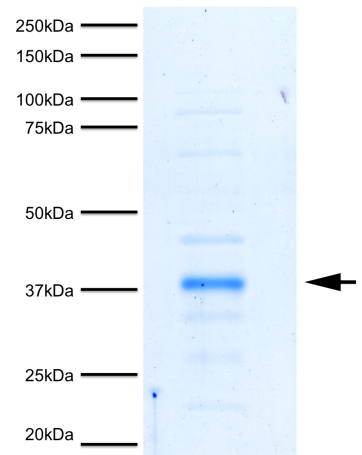
Catalog No. 15-1010
Lot No. 16039001
Pack Size 50 µg



Type HMT
Mol. Wgt. 39 kDa
Expressed In *E. coli*
Epitope Tag GST

Product Description:

Recombinant human G9a protein, (EHMT2 BAT8, KMT1C, accession Q96KQ7, amino acids 913-1193), containing an N-terminal 6xHis tag, expressed in *E. coli*. G9a is a SET-domain containing histone methyltransferase, catalyzing the mono- and di-methylation of histone H3 at lysine 9.



Protein Gel Data: Recombinant G9a, human (1 µg, arrow) run on a PAGE gel and stained with Coomassie blue. Migration and molecular weight of protein standards is indicated.

Formulation:

Recombinant G9a (3.2 mg/ml) in 25 mM Tris-HCl pH 8.0, 150 mM NaCl, 1 mM DTT, 1mM EDTA, and 20% glycerol.

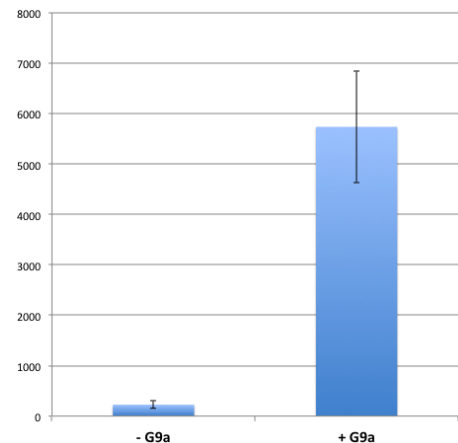
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Recombinant G9a, human is useful for histone H3 methylation experiments, enzyme kinetics and inhibitor screening. Use of 0.1-2 µg G9a per reaction with nucleosomes or recombinant histone H3 as a substrate is recommended.

References:



Enzyme Activity Data: G9a, Recombinant Human, His-Tagged (0.1 µg) was used in an HMTase assay with 0.25 µg HeLa Polynucleosomes (Cat. No. 16-0003) using a standard radiometric filter binding assay protocol.

This product is for *in vitro* research use only and is not intended for use in humans or animals.