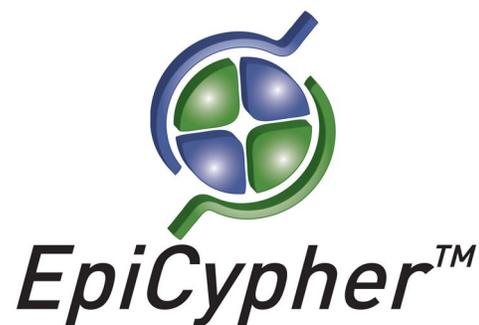
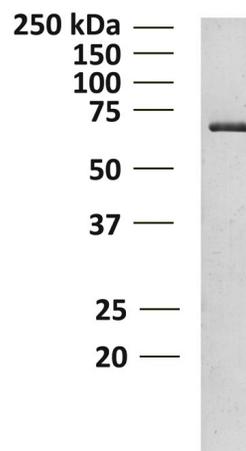


SETD2, Recombinant Human GST-tagged

Catalog No. 15-1011
Lot No. 16244001
Pack Size 50 µg



Type HMT
Mol. Wgt. 68 kDa
Expressed In *E. coli*
Epitope Tag GST



Product Description:

SETD2, Recombinant Human GST-tagged protein (SET2, HYPB, HIF1, KMT3A, accession Q9BYW2, amino acids 1345-1711), containing an N-terminal GST tag, expressed in *E. coli*. SETD2 is a SET-domain containing histone methyltransferase, catalyzing the trimethylation of histone H3 at lysine 36. SETD2 is a candidate tumor suppressor protein and mutated in clear cell renal cell carcinoma.

Protein Gel Data: Recombinant SETD2, Recombinant Human GST-tagged run on a PAGE gel and stained with Coomassie blue. Migration and molecular weight of protein standards are indicated.

Formulation:

SETD2, Recombinant Human GST-tagged (1 µg/µl) in 50mM Tris pH 7.3, 300 mM NaCl, 4mM DTT, 1µM ZnCl₂ and 25% glycerol.

Storage and Stability:

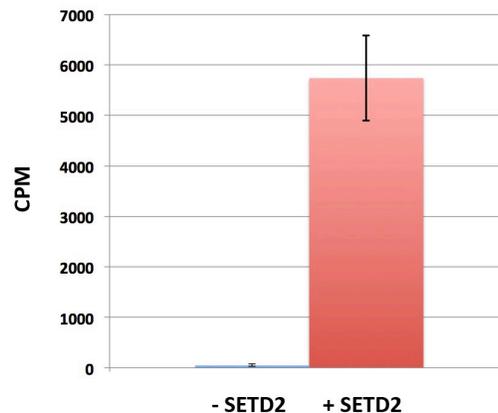
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

SETD2, Recombinant Human GST-tagged is useful for histone H3 methylation experiments, enzyme kinetics and inhibitor screening. Use of 0.5 - 2.5 µg SETD2 per reaction with nucleosomes or recombinant histone H3 as a substrate is recommended.

References:

Hacker KE et al (2016). J Biol Chem DOI: 10.1074/jbc.M116.739375



Enzyme Activity Data: SETD2, Recombinant Human GST-tagged (500 nmol) was used in an HMTase assay with 1 µg Chicken Polynucleosomes (Cat. No. 16-0004) 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.