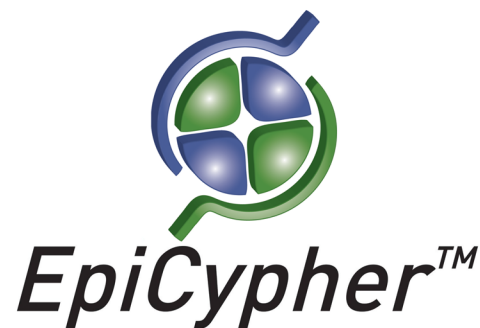


NSD2 / MMSET Catalytic Domain, Recomb. Human

Catalog No. 15-1002
Lot No. 15065001
Pack Size 50 µg



Type HMT
Mol. Wgt. 61 kDa
Expressed In *E. coli*
Epitope Tag GST

Product Description:

NSD2 / MMSET Catalytic Domain, Human Recombinant (WHSC1, accession O96028, amino acids 959-1365), containing an N-terminal GST tag, expressed in *E. coli*. NSD2/MMSET is a SET-domain containing histone methyltransferase, catalyzing the dimethylation of histone H3 at lysine 36. NSD2/MMSET, an oncogene that is overexpressed in several cancers, is thought to drive pathogenesis of t(4;14) positive multiple myeloma.

N.B.-Recombinant NSD2/MMSET requires nucleosomal substrates for activity.

Formulation:

GST-NSD2/MMSET (0.5 µg/µl) in 25 mM Tris pH 8.0, 150 mM NaCl, 1 mM DTT, 1 mM EDTA and 20% glycerol.

Storage and Stability:

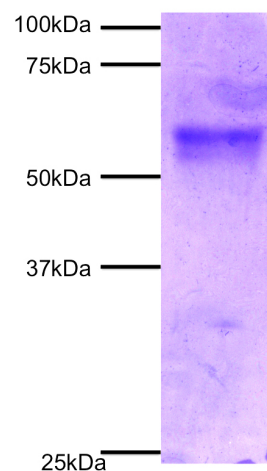
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

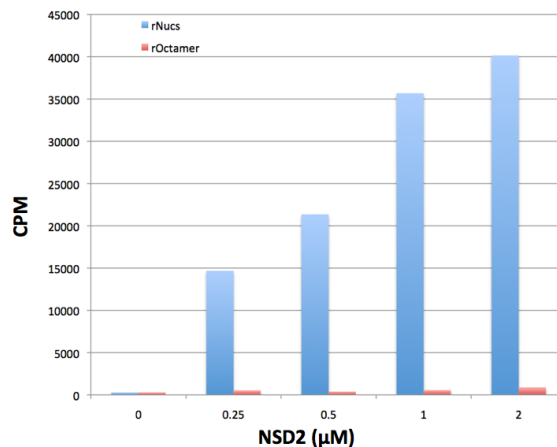
NSD2 / MMSET Catalytic Domain, Human Recombinant is useful for histone H3 methylation experiments, enzyme kinetics and inhibitor screening. Use of 0.5 - 2 µg NSD2 per reaction with HeLa or recombinant nucleosomes as a substrate is recommended.

References:

Kuo AJ et al (2011). *Mol Cell* 44: 609-620.



Protein Gel Data: NSD2 / MMSET Catalytic Domain, Human Recombinant (1 µg) run on a PAGE gel and stained with Coomassie blue. Migration and molecular weights of protein standards are indicated.



Enzyme Activity Data: NSD2 / MMSET Catalytic Domain, Human Recombinant was used in a radioactive methylation assay with either 0.5 µM recombinant nucleosomes or 0.5 µM recombinant histone octamer as the substrate. The reaction was spotted on filter paper and the radioactivity incorporated was counted (CPM). Varying amounts of enzyme were used, as indicated.

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