

SMYD2, Recombinant Human

Catalog No. 15-1006
Lot No. 13141001
Pack Size 25 µg



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

Product Description:

Recombinant full-length human SMYD2 (SET and MYND domain containing protein; also named KMT3C, accession Q9NRG4), containing an N-terminal GST tag, expressed in *E. coli*. SMYD2 is a lysine mono-methyltransferase that has been shown to have activity on histone H3, P53 (Lys370), pRB (Lys380) and HSP90 (Lys616). The SMYD2 locus is amplified in lung carcinoma, serous ovarian carcinoma, and pancreatic ductal adenocarcinoma. SMYD2 expression is linked to a variety of cancers. Elevated levels of SMYD2 in esophageal squamous cell carcinoma patients are correlated with lower rates of

Formulation:

Recombinant GST-SMYD2 (1 µg/µl) in 100mM Tris pH 8.0, 10 mM glutathione and 10% glycerol.

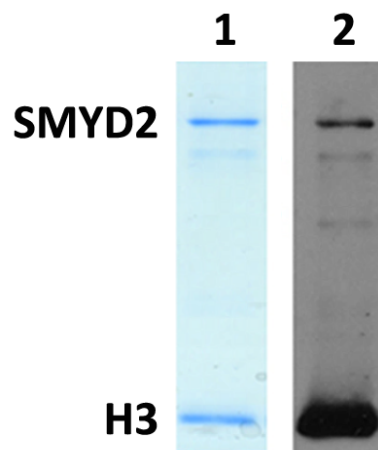
Storage and Stability:

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

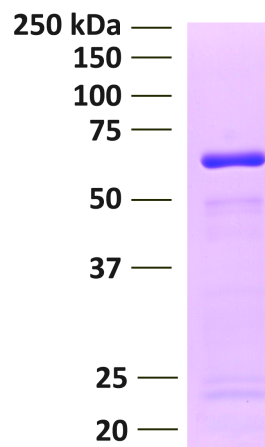
Application Notes:

Recombinant SMYD2, human is useful for histone H3 methylation experiments, enzyme kinetics and inhibitor screening. Use of 1-3 µg SMYD2 per reaction with recombinant histone H3 as a substrate is recommended.

References:



Enzyme Activity Data: Recombinant SMYD2, human (1 µg) was used in a methylation assay with 1 µg recombinant human histone H3 and radioactive SAM and the reaction was run on a PAGE gel. **Lane 1:** Coomassie stain of the gel. **Lane 2:** Autoradiograph. Migration of SMYD2 and H3 is indicated.



Protein Gel Data: Recombinant human SMYD2 (1 µg) run on a PAGE gel and stained with Coomassie blue. Migration and molecular weight of protein standards is indicated.

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