

Recombinant SMYD3 protein

Catalog No: 31407, 31807

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.2 µg/µl

Source: Human

Buffer Contents: Recombinant SMYD3 protein is supplied at a concentration of 0.2 µg/µl in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

Background: SET and MYND domain-containing protein 3 (SMYD3) is a member of the SMYD family of histone methyltransferases that specifically methylates lysine 4 of histone H3 (H3K4), inducing di- and tri-methylation, but not monomethylation. SMYD3 plays an important role in transcriptional activation as a member of an RNA polymerase complex. It has also been shown to bind DNA containing 5'-CCCTCC-3' or 5'-GAGGGG-3' sequences. SMYD3 expression has been shown to be upregulated in certain cancers.

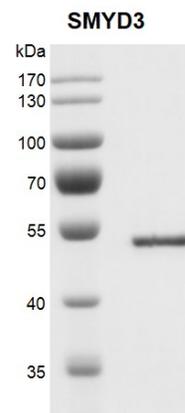
Protein Details: Recombinant SMYD3 (accession number NP_001161212.1) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 50.7 kDa. The recombinant protein is >95% pure by SDS-PAGE.

Application Notes: Recombinant SMYD3 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

Specific Activity: Specifically mediates the methylation of 'Lys-4' and 'Lys-5' of histone H3 and may methylate non-histone substrates.

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.



Recombinant SMYD3 protein gel. SMYD3 protein was run on an SDS-PAGE gel and stained with Coomassie Blue.