

Recombinant SMYD5 protein

Catalog No: 31409, 31809

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.3 µg/µl

Source: Human

Buffer Contents: Full length recombinant SMYD5 protein expressed in Sf9 cells at a concentration of 0.3 µg/µl in 25 mM HEPES pH 7.5, 120 mM NaCl, 0.2 mg/ml 3X FLAG peptide and 5% glycerol.

Background: SET and MYND domain-containing protein 5 (SMYD5) is a member of the SMYD family of histone methyltransferases. SMYD5 has been reported to mediate methylation of H4K20 and may methylate non-histone substrates as well.

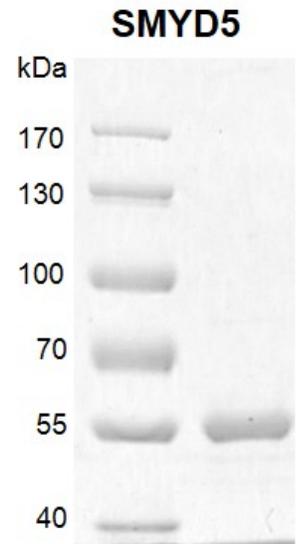
Protein Details: Recombinant Smyd5 (accession number NP_006053.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 49 kDa. The recombinant protein is >93% pure by SDS-PAGE.

Application Notes: Recombinant SMYD5 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

Specific Activity: Specifically mediates the methylation of H4K20 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 SMYD5 protein gel. SMYD5 protein was run on an SDS-PAGE gel and stained with Coomassie Blue.