

Recombinant SMYD1 protein

Catalog No: 31405, 31805

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: Full length recombinant SMYD1 protein expressed in Sf9 cells and supplied 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 1 (SMYD1) is a member of the SMYD family of histone methyltransferases that possesses H3K4 methyltransferase activity. It has also been shown to act as a transcriptional repressor. Studies suggest that SMYD1 may play a critical role in cardiomyocyte differentiation and cardiac morphogenesis and that it also may have a function in skeletal muscle development.

Protein Details: Recombinant SMYD1 (accession number NP_938015.1) was expressed in Sf9 cells and contains an N-terminal FLAG-Tag with an observed molecular weight of 58.3 kDa. The recombinant protein is >85% pure by SDS-PAGE.

Application Notes: Recombinant SMYD1 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

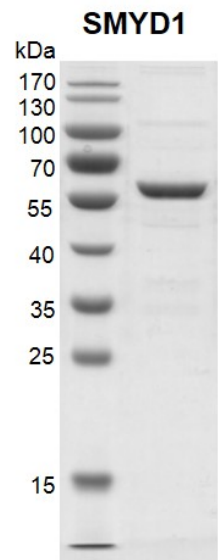
Specific Activity: Not known. May methylate non-histone substrates.

References:

This product was used in the following publications:
Mol. Cell. (2016). 61(2): 247-59. PMID: 26778125.

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 SMYD1 protein gel. SMYD1 protein was run on an SDS-PAGE gel and stained with Coomassie blue.