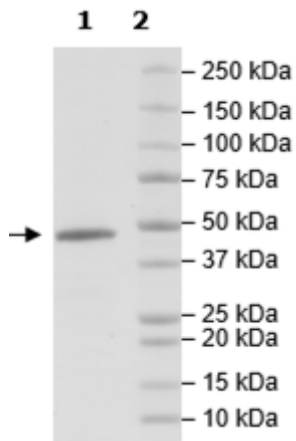


Product Information

Description:	Recombinant human SMYD2 (SET And MYND Domain Containing protein 2), encompassing amino acids 2-433(end). This construct contains an N-terminal FLAG-tag. This protein was affinity purified.
Background:	SMYD2 (SET (su(Var)3-9, Enhancer-of-zeste and Trithorax) And MYND (myeloid, Nervy and DEAF-1) Domain Containing protein 2), also known as KMT3C, is a member of the histone methyltransferase family of proteins, with lysine-specific activity. In addition to its role in histone H3K4 methylation, SMYD2 can also act on non-histone targets, such as p53 and PARP1 (poly-ADP ribose polymerase 1). SMYD2 is found at high levels in many cancer types and has been classified as an oncogene. It regulates cell proliferation and apoptosis, making it a promising cancer therapy target. The use of inhibitors or small interfering RNA can lead to decreased cell cancer viability as SMYD2 is critical for inhibition of TNF (tumor necrosis factor)-induced apoptosis. Targeting SMYD2 alone, or in combination with other therapies, may prove crucial in oncology.
Species:	Human
Construct:	SMYD2 (FLAG-2-433(end))
Concentration:	0.60 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 0.2 mM TCEP, and 80 µg/ml FLAG peptide
MW:	51 kDa
Genbank Accession:	NM_020197
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	0.004 pmole/min/µg
Assay Conditions:	Assay was done according to SMYD2 (KMT3C) Chemiluminescent Assay Kit (#51031) with various concentrations of SMYD2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SMYD2 Activity

