

## Recombinant Smyd2 protein

---

**Catalog No:** 31323

**Expressed In:** *E. coli*

**Quantity:** 10 µg

**Concentration:** 2 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg recombinant Smyd2 protein supplied at a concentration of 2 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

**Background:** Smyd2 (SET and MYND domain containing protein 2) is a lysine methyltransferase that methylates both histones and non-histone proteins. Specifically, Smyd2 methylates histone H3 Lys4 (H3K4me1) and dimethylates histone H3 Lys36 (H3K36me2). Smyd2 also exhibits methyltransferase activity toward non-histone proteins such as p53 and pRb. Smyd2 monomethylates Lys370 of p53, leading to decreased DNA-binding activity and subsequent inhibition of p53-mediated transcriptional regulation. Smyd2 also monomethylates Lys860 of pRb.

**Protein Details:** Smyd2 (SET and MYND domain containing protein 2) is a lysine methyltransferase for p53 and histone H3 lysine 36 (H3K36). Recombinant Smyd2 was expressed as the full-length (accession number NP\_064582.2) human protein with an N-terminal poly-histidine tag.

**Application Notes:** Recombinant Smyd2 is suitable for use in methyltransferase assays. A recommended starting point for the methyltransferase assay is 50-200 ng of recombinant protein per reaction. The molecular weight of the protein is ~52 kDa.

**Storage and Guarantee:** 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.