

Recombinant Smyd2 Y240F protein

Catalog No: 31324

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 2 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant Smyd2 Y240F 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.

Protein Details: Smyd2 (SET and MYND domain containing protein 2) is a lysine methyltransferase for p53 and histone H3 lysine 36 (H3K36). Recombinant Smyd2 Y240F was expressed as the full-length (accession number NP_064582.2) human protein with an N-terminal poly-histidine tag. The recombinant Smyd2 Y240F protein contains a substitution of tyrosine to phenylalanine at amino acid 240. Smyd2 Y240F is catalytically inactive.

Application Notes: Recombinant Smyd2 Y240F 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. This is a version of the protein with a catalytic site mutant, so it is designed for use as a negative control.

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.