

Recombinant SETD7 (SET9) protein

Catalog No: 31319

Lot No: 21509001

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 2 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant SETD7 (SET9) protein supplied in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Protein Details: SETD7 (SET9) is a histone methyltransferase that is specific to lysine 4 of histone H3 (H3K4). Methylation of H3K4 is associated with transcriptional activation. Recombinant SETD7 (SET9) was expressed as the full-length (accession number NP_085151.1) human protein with an N-terminal His-Tag.

Application Notes: Recombinant SETD7 (SET9) 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 ~53 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.