

Recombinant KMT5A (SET8) protein

Catalog No: 31321

Lot No: 21509001

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 2 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant KMT5A (SET8) protein supplied in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Protein Details: KMT5A (SET8), also known as PR-Set7, is a histone methyltransferase specific to lysine 20 of histone H4 (H4K20). Methylation of H4K20 is associated with transcriptional silencing. Recombinant KMT5A (SET8) was expressed as the human protein missing amino acids 13-17 (accession number NP_065115.3) with an N-terminal His-Tag.

Application Notes: Recombinant KMT5A (SET8) is suitable for use in methyltransferase assays. A recommended starting point for the methyltransferase assay is 200-500 ng of recombinant protein per reaction. The molecular weight of the protein is ~49 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.