

## Recombinant KMT5A (SET8) D338A protein

---

**Catalog No:** 31322

**Lot No:** 21509001

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

**Quantity:** 10 µg

**Concentration:** 2 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg recombinant KMT5A (SET8) D338A 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 or SETD8) is a histone methyltransferase specific to lysine 20 of histone H4 (H4K20). Methylation of H4K20 is associated with transcriptional silencing. Recombinant KMT5A (SET8) D338A was expressed as the human protein missing amino acids 13-17 (accession number NP\_065115.3) with an N-terminal His-Tag. The recombinant KMT5A (SET8) D338A protein contains a substitution of aspartic acid to alanine at amino acid 338. KMT5A (SET8) D338A is unable to methylate histone H4.

**Application Notes:** Recombinant KMT5A (SET8) D338A 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. 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.