

SETDB1 (KMT1E) Protein

Human recombinant protein expressed in Sf9 cells

Catalog # S346-31G

Lot # B2035-8

Product Description

Recombinant human SETDB1 (KMT1E) (715-end) was expressed by Baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC028671](#).

Gene Aliases

KMT1E, ESET, KG1T, KIAA0067

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

Histone-lysine N-methyltransferase SETDB1 (SETD6) specifically trimethylates 'Lys-9' of histone H3 and plays a central role in the silencing of euchromatic genes. SETDB1 is targeted to histone H3 by TRIM28/TIF1B, a factor recruited by KRAB zinc-finger proteins. Required for activated KRAS-mediated promoter hypermethylation and transcriptional silencing of tumor suppressor genes (TSGs) or other tumor-related genes in colorectal cancer (CRC) cells by forming a corepressor complex. It was also found to be responsible for transcriptionally repressive state of genes in embryonic stem cells (ESCs).

References

- Serra RW, et al: A KRAS-directed transcriptional silencing pathway that mediates the CpG island methylator phenotype. *Elife*. 2014 Mar 12;3:e02313.
- <http://www.uniprot.org/uniprot/Q15047>.

Purity



The purity of SETDB1 (KMT1E) was determined to be **>80%** by densitometry. Approx. MW **~100kDa**.

SETDB1 (KMT1E) Protein

Human recombinant protein expressed in Sf9 cells

| | |
|--------------------|---|
| Catalog # | S346-31G |
| Lot # | B2035-8 |
| Purity | >80% |
| Concentration | 0.1 µg/µl |
| Stability | 1yr at -70°C from date of shipment |
| Storage & Shipping | Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.