

SUV39H2 (KMT1B), Active

Full length recombinant protein expressed in Sf9 cells

Catalog # S350-380BG

Lot # P1545-6

Product Description

Recombinant full-length human SUV39H2 (KMT1B) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The SUV39H2 protein accession number is [BC007754](#).

Gene Aliases

KMT1B; FLJ23414

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.

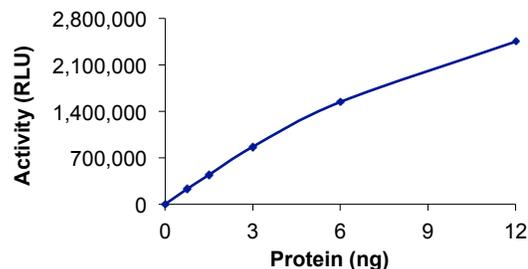
Scientific Background

Homo sapiens suppressor of variegation 3-9 homolog 2 (SUV39H2) is encoded by the SUV39H2 gene and belongs to the histone-lysine methyltransferase family and Suvar3-9 subfamily. It contains an N-terminal chromodomain and a C-terminal SET domain. It trimethylates lysine 9 of histone H3, creating a binding site for the chromo domain of HP1 for epigenetic transcriptional repression. Deregulation of SUV39H2 interferes with mammalian higher-order chromatin organization.

References

1. Rea S, et al: Regulation of chromatin structure by site-specific histone H3 methyltransferases. *Nature*.10;406(6796):593-9, 2000.
2. Aagaard L, et al: Functional mammalian homologues of the Drosophila PEV-modifier Su(var)3-9 encode centromere-associated proteins which complex with the heterochromatin component M31. *EMBO J*. 1;18(7):1923-38, 1999.

Specific Activity



The specific activity of SUV39H2 (KMT1B) was determined to be **92 nmol /min/mg** as per activity assay protocol.

Purity



The purity of SUV39H2 (KMT1B) was determined to be **>70%** by densitometry, approx. MW **67 kDa**.

SUV39H2 (KMT1B), Active

Full length recombinant protein expressed in Sf9 cells

Catalog Number	S350-380BG
Specific Activity	92 nmol/min/mg
Specific Lot Number	P1545-6
Purity	>70%
Concentration	0.05 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.

Activity Assay Protocol

Reaction Components

Active Methyltransferase (Catalog #: S350-380BG)

Active SUV39H2 (KMT1B) (0.05 µg/µl) diluted with Methyltransferase Reaction Buffer and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active SUV39H2 (KMT1B) for optimal results).

Methyltransferase Reaction Buffer

Buffer components: 20mM Tris-HCl, pH 8.0, 50 mM NaCl, 1 mM EDTA, 3 mM MgCl₂, 0.1 mg/ml BSA. Add 1mM DTT prior to use.

MTase-Glo™ Methyltransferase Assay (Promega, Catalog #: V7601)

S-Adenosyl-Methionine (SAM), 1mM
S-Adenosyl-Homocysteine (SAH), 15 µM
Methyltransferase-Glo™ Reagent, 10X
MTase-Glo™ Detection Solution, 1 bottle

Substrate (Catalog #: H12-58)

Histone H3 Peptide (1-21) diluted in distilled H₂O to a final concentration of 1mg/ml.

Assay Protocol

The SUV39H2 (KMT1B) assay is performed using the Methyltransferase-Glo™ Assays kit (Promega, Catalog #: V7601).

- Step 1.** Thaw the active SUV39H2 (KMT1B) and all Methyltransferase-Glo™ Assays kit reagents on ice.
 - Step 2.** Prepare the following working solutions with Methyltransferase Reaction Buffer on ice:
 - o 2X final concentration of Active SUV39H2 (KMT1B) (Catalog # S350-380BG)
 - o 2X Substrate Cocktail: 40 µM of SAM and 100ng/µl of Histone H3 Peptide (1-21) (Catalog # H12-58) in water
 - Step 3.** In a polystyrene 96-well plate, add the following components to bring the initial reaction volume to 20 µl:
 - Component 1.** 10 µl of 2X Substrate Cocktail
 - Component 2.** 10 µl of 2X Active SUV39H2 (KMT1B)
- Note: A blank control can be set up as outlined in step 3 by replacing the substrate working solution with an equal volume of Reaction Buffer.*
- Step 4.** Mix the reaction on an orbital shaker for 2 minutes. Seal the plate with a plate seal and incubate at 37°C for 60 minutes
 - Step 5.** Dilute 10X Methyltransferase-Glo™ Reagent with equal volume of nanopure water, and add 5 µl of the 5X Methyltransferase-Glo™ Reagent to all reaction wells
 - Step 6.** Mix on an orbital shaker for 2 minutes and then incubate at room temperature for 30 minutes.
 - Step 7.** Add 25 µl of MTase-Glo™ Detection Solution to all reaction wells. Mix for 2 minutes and then incubate at room temperature for 30 minutes
 - Step 8.** Read the plate using the KinaseGlo Luminescence Protocol on a GloMax plate reader (Promega; Cat# E7031)
 - Step 9.** Using the SAH standard curve, determine the concentration of SAH produced (nM) and calculate the methyltransferase specific activity as outlined below. For a detailed protocol of how to determine SAH amount from RLU, see MTase-Glo™ Methyltransferase Assay protocol at Promega's website: www.promega.com/protocols

Methyltransferase Specific Activity (SA) (nmol/min/mg)

$$= \frac{[SAH](nM) \times Reaction Volume(\mu l)}{Reaction Time (min) \times Enzyme Amount (mg)} \times 10^{-6}$$

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.