

## SETD1A (KMT2F), Active

Recombinant human protein expressed in Sf9 cells

**Catalog # S342-381G**

Lot # L2041-9

### Product Description

Recombinant human SETD1A (1418-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The SETD1A protein accession number is [NM\\_014712](#).

### Gene Aliases

SETD1A; KMT2F; Set1; Set1A

### 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  $-70^{\circ}\text{C}$ . 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.

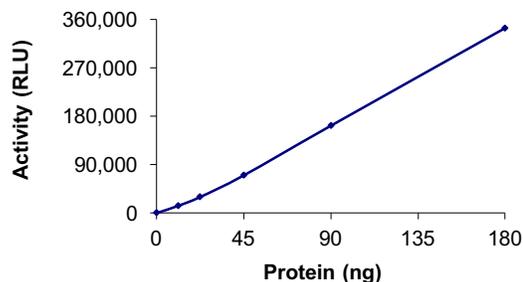
### Scientific Background

As a member of the mixed lineage leukemia (MLL) family of histone methyltransferase, it specifically methylates 'Lys-4' of histone H3, when part of the SET1 histone methyltransferase (HMT) complex, but not if the neighboring 'Lys-9' residue is already methylated. H3 'Lys-4' methylation represents a specific tag for epigenetic transcriptional activation.

### References

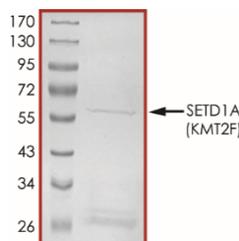
- Li Y, et al: Structural basis for activity regulation of MLL family methyltransferases. *Nature*. 530(7591):447-52. 2016
- <http://www.uniprot.org/uniprot/O15047>

### Specific Activity



The specific activity of SETD1A (KMT2F) was determined to be **500 pmol /min/mg** as per activity assay protocol.

### Purity



The purity of SETD1A (KMT2F) was determined to be **>70%** by densitometry, approx. MW **61 kDa**.

## SETD1A (KMT2F), Active

Recombinant human protein expressed in Sf9 cells

|                    |  |
|--------------------|--|
| Catalog #          | S342-381G  |
| Specific Activity  | 500 pmol/min/mg  |
| Lot #              | L2041-9  |
| Purity             | >70%   |
| Concentration      | 0.05 µg/µl   |
| Stability          | 1yr at $-70^{\circ}\text{C}$ from date of shipment   |
| Storage & Shipping | Store product at $-70^{\circ}\text{C}$ . 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://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 #: S342-381G)

Active SETD1A (KMT2F) (0.05µg/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 SETD1A (KMT2F) for optimal results).

### Methyltransferase Reaction Buffer

Buffer components: 20mM Tris-HCl, pH 8.0, 50 mM NaCl, 1 mM EDTA, 3 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA. Add 1mM DTT prior to use.

### Components of SETD1A (KMT2F) Complex

WDR5 Protein: Catalog #: W325-30H

RbBP7 Protein: Catalog #: R317-30G

ASH2L Protein: Catalog #: A372-30BG

### 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 Reaction Buffer to a final concentration of 20 µM.

## Assay Protocol

The SETD1A (KMT2F) assay is performed using the Methyltransferase-Glo™ Assays kit (Promega, Catalog #: V7601).

- Step 1.** Thaw each active SETD1A (KMT2F) complex component 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 SETD1A (KMT2F) (Catalog # S342-381G) with complex proteins
  - o 2X Substrate Cocktail: 40 µM of SAM and 20 µM of Histone H3 Peptide (1-21) in Reaction Buffer
- 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 SETD1A (KMT2F) complex

*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 RLUs, see MTase-Glo™ Methyltransferase Assay protocol at Promega's website: [www.promega.com/protocols](http://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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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