

DOT1L (KMT4), Active

Recombinant human protein expressed in Sf9 cells

Catalog # D344-381G

Lot # L1956-7

Product Description

Recombinant human DOT1L (KMT4) (1-435) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The DOT1L (KMT4) gene accession number is [NM_032482](#).

Gene Aliases

DOT1; KMT4

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.

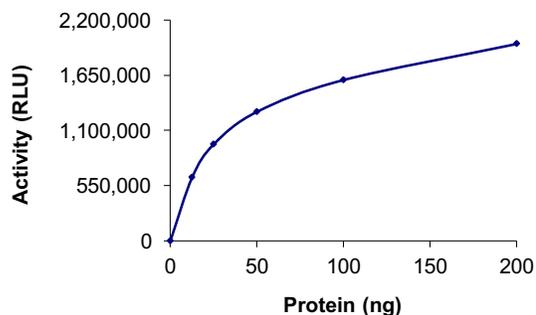
Scientific Background

DOT1-like protein (DOT1L) specifically methylates lysine-79 of histone H3 in the globular domain. DOT1L lacks a SET-domain, which is commonly considered to mediate lysine methylation of histones (1, 2). DOT1 homologs exist in a variety of eukaryotic organisms. The activity of this methyltransferase is involved in cell cycle regulation and leukemogenesis (3).

References

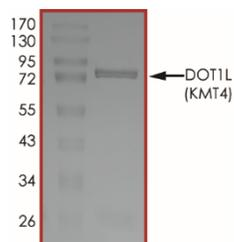
- Feng, Q., et al. Methylation of H3-lysine 79 is mediated by a new family of HMTases without a SET domain. *Curr Biol.* 12(12):1052-8, 2002.
- Min, J., et al. Structure of the catalytic domain of human DOT1L, a non-SET domain nucleosomal histone methyltransferase. *Cell.* 112(5):711-23, 2003.
- Okada, Y., et al. hDOT1L links histone methylation to leukemogenesis. *Cell.* 121(2):167-78, 2005.

Specific Activity



The specific activity of DOT1L (KMT4) was determined to be **14 nmol/min/mg** as per activity assay protocol.

Purity



The purity of DOT1L (KMT4) was determined to be **>90%** by densitometry, approx. MW **78 kDa**.

DOT1L (KMT4), Active

Recombinant human protein expressed in Sf9 cells

Catalog #	D344-381G
Specific Activity	14 nmol/min/mg
Lot #	L1956-7
Purity	>90%
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.

Activity Assay Protocol

Reaction Components

Active Methyltransferase (Catalog #: D344-381G)

Active DOT1L (KMT4) (0.1 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 DOT1L (KMT4) 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 (Reaction Biology Cat. # HMT-35-130)

Nucleosomes (Hela Oligo) diluted in Reaction Buffer to a final concentration of 4 µM.

Assay Protocol

The DOT1L (KMT4) assay is performed using the Methyltransferase-Glo™ Assays kit (Promega, Catalog #: V7601).

- Step 1.** Thaw the active DOT1L (KMT4) 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 DOT1L (KMT4) (Catalog # D344-381G)
 - o 2X Substrate Cocktail: 40 µM of SAM and 4 µM of Nucleosomes (Hela Oligo) 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 DOT1L (KMT4)

Note: A blank control can be set up as outlined in step 3 by replacing the enzyme 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's, 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.