

## Certificate of Analysis

### DOT1L

**Human Disruptor of Telomeric silencing 1-like, active**

**(Recombinant enzyme expressed in *E.coli*)**

**Item # EPI018**

**Lot # 139683**

**Product Description:** Human recombinant DOT1L, amino acids 1-420, expressed in *E.coli*. Purified using immobilised metal affinity chromatography.  
MW = 47.9kDa.

Tag cleaved by TEV.

**Aliases :** KMT4, KIAA1814

**Formulation:** 1mg/ml of enzyme in 25mM Hepes/NaOH pH7.4, 250mM NaCl, 50% Glycerol, 2mM DTT. Frozen solution.

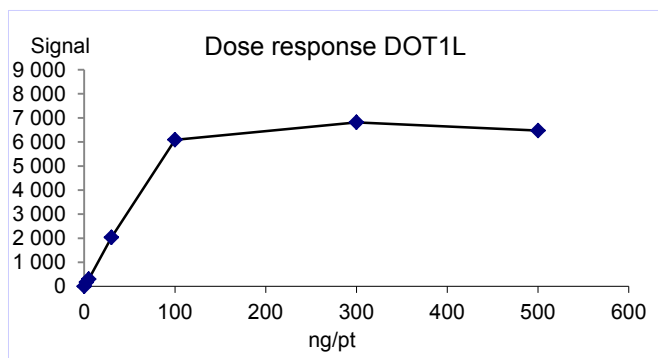
**Storage and Stability:** Stable for 1 year at -70°C from date of shipment. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and store at -70°C.

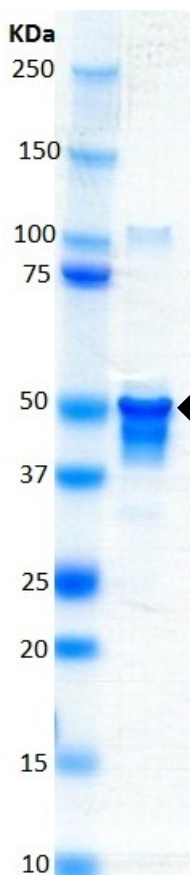
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**HMT Assay:** 2.5-500ng of this lot of enzyme transferred methyl groups from [3H] SAM to HeLa Polynucleosome in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



**MS:** Size was confirmed by mass spectrometry using a Q-TOF.



**SDS-PAGE and Coomassie Stain:** Purity was assessed by SDS-PAGE and Coomassie blue staining using 4µg of DOT1L.

## Certificate of Analysis

### DOT1L Assay Protocol

#### Stock Solutions:

1. **Reaction buffer:** 20 mM Tris/HCl pH8.0, 100mM NaCl, 2mM DTT.
2. **DOT1L, active:** Dilute with reaction buffer. Use 2.5-500ng per assay point.
3. **HeLa Polynucleosome:** Dilute with reaction buffer to 12.5µg/ml.
4. **[3H] SAM:** Dilute with reaction buffer to 200nM.
5. **Filtration Buffer :** 33mM Citric acid pH2.2

#### Assay Procedure (96 well plate format):

1. Add 5µl of 10% DMSO per assay to each well.
2. Add 25µl of [3H] SAM.
3. Add 10µl (**2.5-500ng**) **DOT1L, active**.
4. Add 10µl of HeLa Polynucleosome.
5. Incubate for 15 minutes at 22°C.
6. Stop the reaction by adding 500µl of citric acid, then filter on a GF/B Filter. Wash 3 times with filtration buffer.
7. Dry and add scintillation cocktail.
8. Read in a scintillation counter. Compare the signal of enzyme samples with that of a background sample that contains all assay components except the enzyme DOT1L.

## Certificate of Analysis

### DOT1L Sequence Information

**Protein** Human DOT1L  
**Tags** Tag cleaved by TEV  
**Accession number** GenBank NP\_115871.1

#### ***Recombinant DOT1L amino acid sequence:***

```
1 GMGEKLELRL KSPVGAEP AV YPWPLPVYDK HHDA AHEIIE TIRWVCEEIP
51 DLKLAMENYV LIDYDTKSFE SMQRLCDKYN RAIDSIHQLW KGTTQPMKLN
101 TRPSTGLLRH ILQQVYNHSV TDPEKLNNYE PFSPEVYGET SFDLVAQMID
151 EIKMTDDDLF VDLGSGVGQV VLQVAAATNC KHHYGV EKAD IPAKYAETMD
201 REFRKWMKWY GKKAHEYTL E RGDFLSEWR ERIANTSVIF VNNFAFGPEV
251 DHQLKERFAN MKEGGRI VSS KPFAPLNFRI NSRNLSDIGT IMRVVELSPL
301 KGSVSWTGKP VSYYLHTIDR T ILENYFSSL KNPKLREEQE AARRRQ QRES
351 KSNAATPTKG PEGKVAGPAD APMDSGAEEE KAGAATVKKP SPSKARKKKL
401 NKKGRKMAGR KRGRPKMNT A
```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services.