

## Certificate of Analysis

### SETD8

#### Human histone Lysine N-methyltransferase SETD8, active

(Recombinant enzyme expressed in *E.coli*)

Item # EPI057

Lot # 140882

**Product Description:** recombinant human SETD8, amino acids 186-352, expressed in *E.coli*. Purified using immobilised metal affinity chromatography. MW = 19.2kDa.

**Formulation:** 1mg/ml of enzyme in 25mM Tris/HCl pH7.5, 250mM NaCl, 50% Glycerol. 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.

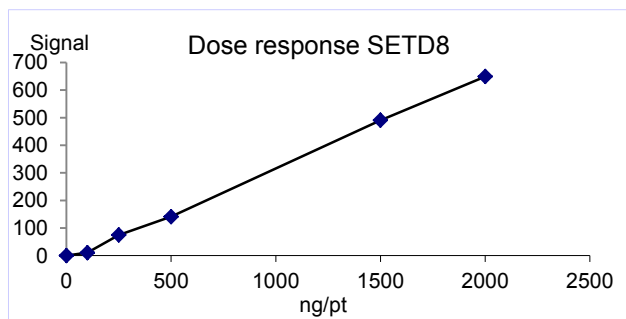
**Tag cleaved by thrombin protease.**

**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.

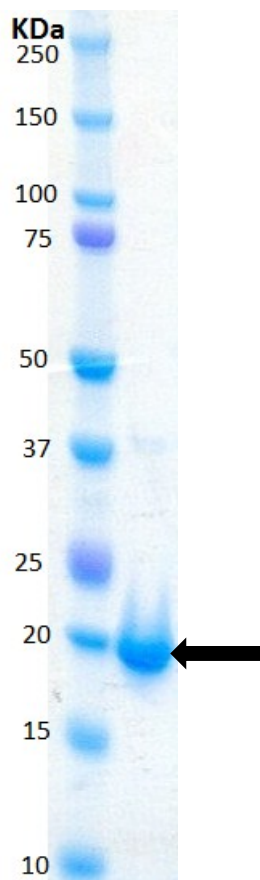
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**HMT Assay:** 100-2000ng of this lot of enzyme transferred methyl groups from [3H] SAM to histone H4 full length 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 SETD8.

## Certificate of Analysis

### SETD8 Assay Protocol

#### Stock Solutions:

1. **Reaction buffer:** 50mM Tris/HCl pH9, 5mM MgCl<sub>2</sub>, 50mM NaCl, 4mM DTT.
2. **SETD8, active:** Dilute with reaction buffer. Use 100-2000ng per assay point.
3. **Histone H4 full length:** Dilute with reaction buffer to 1250mM.
4. **[3H] SAM:** Dilute with SAM solution (1080mM) to 120nM.
5. **Filtration Buffer :** 33mM Citric acid pH2.2

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

1. Add 5µl of 10% DMSO per assay to wells.
2. Add 25µl of [3H] SAM.
3. Add 10µl (**100-2000ng**) **SETD8, active**.
4. Add 10µl of Histone H4 full length.
5. Incubate for 120 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 SETD8.

## Certificate of Analysis

### SETD8 Sequence Information

<b><u>Protein</u></b>	Human SETD8
<b><u>Tags</u></b>	tag cleaved by thrombin protease
<b><u>Accession number</u></b>	GenBank NP_065115.3

#### ***Recombinant SETD8 amino acid sequence:***

1 GSPVRRSSRK SKAELQSEER KRIDELIESG KEEGMKIDLI DGKGRGVIAT  
51 KQFSRGDFVV EYHGDLIEIT DAKKREALYA QDPSTGCYMY YFQYLSKTYC  
101 VDATRETNRL GRLINHSKCG NCQTKLHDID GVPHLILIAS RDIAAGEELL  
151 YDYGDRSKAS IEAHPWLKH

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.