

## Certificate of Analysis

### NSD1

#### Human Nuclear receptor-binding SET-domain protein 1, active

(Recombinant enzyme expressed in *E.coli*)

Item # EPI041

Lot # 139682

**Product Description:** Recombinant human NSD1, Amino Acids 1810-2120, expressed in *E.coli*. Purified using immobilised metal affinity chromatography.  
MW = 35.6 kDa.

**Formulation:** 1mg/ml of enzyme in 25mM Tris/HCl pH7.5, 125mM NaCl, 1.3mM TCEP, 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.

**Aliases :** ARA267, KMT3B, SOTOS1

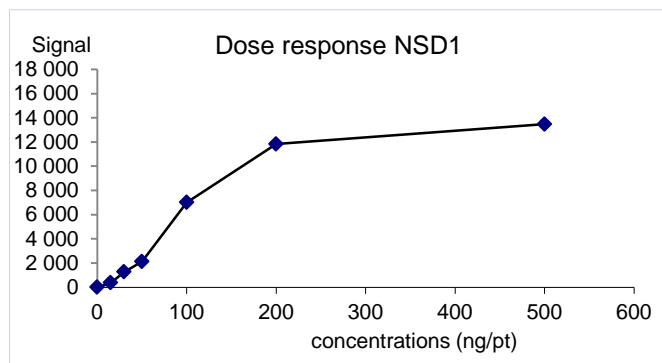
**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and store at -70°C.

**Tag cleaved by TEV protease.**

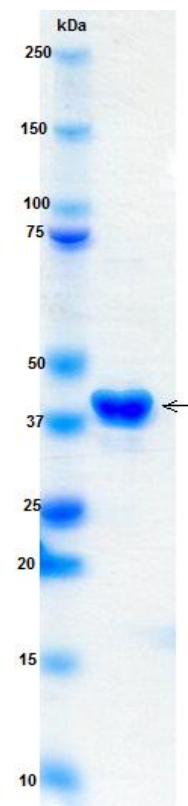
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**HMT Assay:** 15-500ng of this lot of enzyme transferred methyl groups from [3H] SAM to poly nucleosome 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 NSD1.

## Certificate of Analysis

### NSD1 Assay Protocol

#### Stock Solutions:

1. **Reaction buffer:** 50mM Tris/HCl pH9, 0.01% Triton.
2. **NSD1:** Dilute with reaction buffer. Use 15-500ng per assay point.
3. **Polynucleosome:** Dilute with reaction buffer to 15µg/ml.
4. **[3H] SAM:** Dilute with reaction buffer to 100nM.
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 (**15-500ng**) **NSD1**.
4. Add 10µl of 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 NSD1.

## Certificate of Analysis

## NSD1 Sequence Information

<b><u>Protein</u></b>	Human NSD1
<b><u>Tags</u></b>	<i>tag cleaved by TEV protease</i>
<b><u>Accession number</u></b>	GenBank NP_071900.2

**Recombinant NSD1 amino acid sequence:**

1 GARVFPYMEG DVSSKDKMGK GVDGTYKKAL QEAAARFEEL KAQKELRQLQ  
51 EDRKNDKKPP PYKHIKVNRP IGRVQIFTAD LSEIPRCNCK ATDENPCGID  
101 SECINRMLLY ECHPTVCPAG GRCQNQCFSK RQYPEVEIFR TLQRGWGLRT  
151 KTDIKKGEFV NEYVGELIDE EECRARIRYA QEHDITNFYM LTLDKDRIID  
201 AGPKGNYARF MNHCCQPNC E TQKWSVNGDT RVGLFALSDI KAGTELTFFNY  
251 NLECLGNGKT VCKCGAPNCS GFLGVRPKNQ PIATEEKSCK FKKKQQGKRR  
301 TQGEITKERE DE

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.