

Certificate of Analysis

WHSC1 (NSD2)

Human Wolf-Hirschhorn syndrome candidate 1

(Recombinant enzyme expressed in *E.coli*)

Item # EPI073

Lot # 142157

Product Description: Recombinant human NSD2, Amino acids 825-1208, expressed in *E.coli*. Purified using immobilised metal affinity chromatography.
MW = 44.2kDa.

Tag cleaved by TEV protease.

Formulation: 0.765 mg/ml of enzyme in 25mM Tris/HCl pH8, 100mM NaCl, 2.5mM TCEP, 2.5 mM imidazole, 50% Glycerol. Frozen solution.

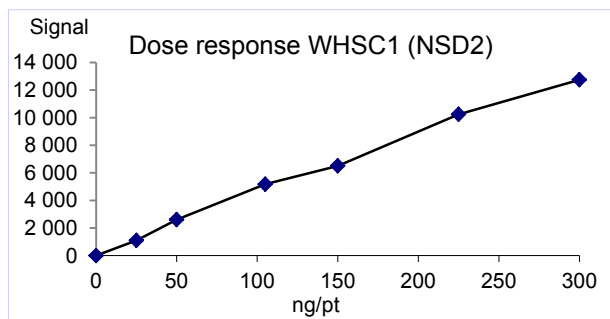
Storage and Stability: Stable for 1 year at -70°C from date of shipment. For maximum recovery of product, centrifuge original vial prior to removing the cap.

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.

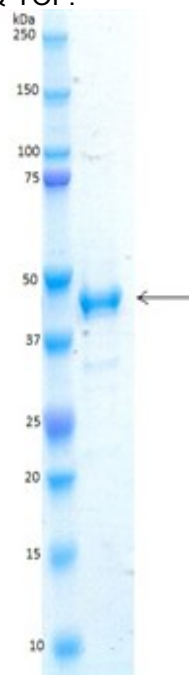
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

HMT Assay: 25-300ng of this lot of enzyme transferred methyl groups from [3H] SAM to core histone 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 WHSC1 (NSD2)

Certificate of Analysis

WHSC1 (NSD2) Assay Protocol

Stock Solutions:

1. **Reaction buffer:** 50mM Tris/HCl pH8, 5mM MgCl₂, 50mM NaCl, 4mM DTT.
2. **WHSC1 (NSD2):** Dilute with reaction buffer. Use 25-300ng per assay point.
3. **Core Histone:** Dilute with reaction buffer to 7500nM.
4. **[3H] SAM:** Dilute with reaction buffer to 500nM.
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 (**25-300ng/pt**) **WHSC1 (NSD2)**.
4. Add 10µl of Core Histone.
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 WHSC1 (NSD2).

Certificate of Analysis

WHSC1 (NSD2) Sequence Information

<u>Protein</u>	Human WHSC1 (NSD2)
<u>Tags</u>	<i>tag cleaved by TEV protease</i>
<u>Accession number</u>	GenBank NP_579877.1

Recombinant WHSC1 (NSD2) amino acid sequence:

1 GHHAHVNSW CFVCSKGGSL LCCESCPAAF HPDCLNIEMP DGSWFCNDCR
51 AGKKLHFQDI IWVKLGNYRW WPAEVCHPKN VPPNIQMKH EIGFPPVFFF
101 GSKDYYWTHQ ARVFPYMEGD RGSRYQGV RG IGRVFKNALQ EAEARFREIK
151 LQREARETQE SERKPPPYKH IKVKNPYGKV QIYTADISEI PKCNCKPTDE
201 NPCGFDSECL NRMLMFECHP QVCPAGEFCQ NQCFTRQYP ETKI IKTDGK
251 GWGLVAKRDI RKGEFVNEYV GELIDEEECM ARIKHAHEND ITHFYMLTID
301 KDRIIDAGPK GNYSRFMNHS CQPNCETLKW TVNGDTRVGL FAVCDIPAGT
351 ELTFNYNLDC LGNEKTVCRC GASNCSGFLG DRPKT

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.