

WHSC1L1 (NSD3) Protein

Human recombinant protein expressed in Sf9 cells

Catalog # N378-31G

Lot # B2058-4

Product Description

Recombinant human WHSC1L1 (NSD3) (1038-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC113469](#).

Gene Aliases

NSD3, DC28, FLJ20353, pp14328

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.

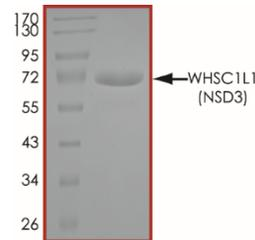
Scientific Background

Histone-lysine N-methyltransferase NSD3 or Wolf-Hirschhorn syndrome candidate 1-like protein 1 (WHSC1L1) is a nuclear receptor binding SET domain (NSD) protein that belongs to a family of three histone methyltransferases including NSD1, NSD2/MMSET/WHSC1, and NSD3/WHSC1L1. Preferentially methylates 'Lys-4' and 'Lys-27' of histone H3. H3 'Lys-4' methylation represents a specific tag for epigenetic transcriptional activation, while 'Lys-27' is a mark for transcriptional repression.

References

- Kim SM, et al: Characterization of a novel WHSC1-associated SET domain protein with H3K4 and H3K27 methyltransferase activity. *Biochem Biophys Res Commun.* 2006 Jun 23;345(1):318-23.
- Morishita M, et al: Cancers and the NSD family of histone lysine methyltransferases. *Biochim Biophys Acta.* 2011 Dec;1816(2):158-63.

Purity



The purity of WHSC1L1 (NSD3) was determined to be **>95%** by densitometry. Approx. MW **~72kDa**.

WHSC1L1 (NSD3) Protein

Human recombinant protein expressed in Sf9 cells

Catalog #	N378-31G
Lot #	B2058-4
Purity	>95%
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.