

## Data Sheet

### NSD2

Human, recombinant, N-terminal GST tag  
 Catalog #: 51026  
 Lot#: 130516                      Conc.: 0.5 mg/ml

**Formulated in** 50 mM Tris-HCl, pH 8.5,  
 300 mM NaCl, 5 mM DTT, 10 mM  
 Glutathione, and 10% Glycerol.

**Stability:** >6 months at -80°C

**References:**

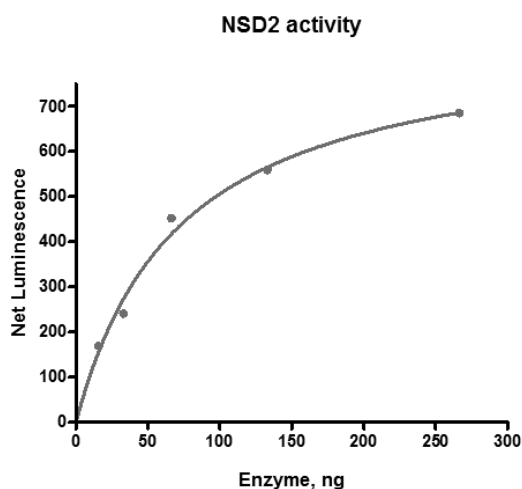
1. Kassambara A. *et al.*, *Biochem. Biophys. Res. Commun.* **379** (4), 840-845 (2009).
2. Brito J.L. *et al.*, *Haematologica* **94** (1), 78-86 (2009).

**Description:** Human NSD2, also known as MMSET and WHSC1, (GenBank Accession No. NM\_133330), amino-acids 941-1240 with N-terminal GST tag, MW = 62 kDa, expressed in *E. coli*.

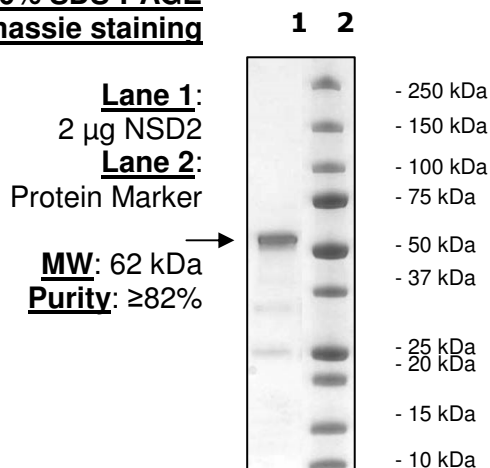
**Specific Activity:** 0.036 pmol/min/μg  
**Assay Conditions:** Add 50 μl reaction mix (100 μM S-adenosylmethionine, and 50-250 ng NSD2 in HMT buffer 7 (BPS #52193)) to wells coated with the substrate. Incubate at 37°C for 2 hr. Add antibody against methylated K36 residue of histone H3, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance



**4-20% SDS-PAGE  
 Coomassie staining**



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Amino acid sequence:**

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYIDGD  
VKLTQSMARIYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLKVDFLSK  
LPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLVCFKKRIEAI  
PQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDLEVLFGPLGSEGDRGSRYYQGVRGIG  
RVFKNALQEAEARFREIKLQREARETQESERKPPPYKHIKVNKPYGKVQIYTADISEIPKC  
NCKPTDENPCGFDSECLNRMLMFECHPQVCPAGEFCQNQCFTKRQYPETKIIKTDGKGW  
GLVAKRDIRKGEFVNEYVGELIDEEECMARIKHAHENDITHFYMLTIDKRIIDAGPKGNYS  
RFMNHSCQPNCE TLKWTVNGDTRVGLFAVCDIPAGTELTFNYNLDCLGNEKTVCRGAS  
NCSGFLGDRPKTSTTLSSSEKGGKTKKKTRRRRAKGEGRQSED

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)