

Data Sheet

NSD1

Human, recombinant, N-terminal GST tag
 Catalog #: 51024
 Lot#: 130524-5B Conc.: 1.02 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 4 mM GSH, 3 mM DTT.

Stability: >12 months at -80°C. Avoid multiple freeze/thaw cycles.

References:

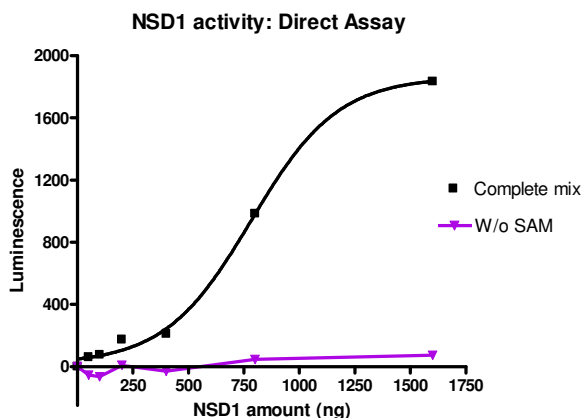
1. Yamada-Okabe, T. and Matsumoto, N.J. *Cell Biochem. Funct.* **26 (2)**, 146-150 (2008).
2. Saugier-veber, P. et al., *Hum. Mutat.* **28 (11)**, 1098-1107 (2007).

Description: Human NSD1 (GenBank Accession No. NM_022455), amino-acids 1849-2094 with N-terminal GST tag, MW = 54 kDa, expressed in an *E. coli* expression system.

Assay Conditions: Assay was performed using the BPS NSD1 Chemiluminescent Assay Kit (Cat# 53010) with 400 μM S-adenosyl-methionine on biotinylated H3 peptide coated plates overnight at room temperature.

Application: Suitable for protein binding studies, as an immunogen, or as a Western blotting control.

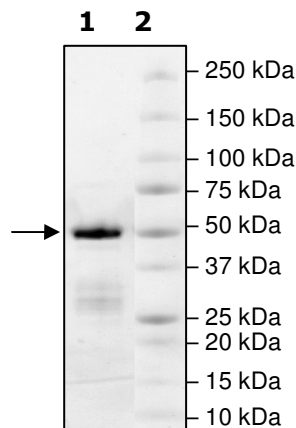
Quality Assurance



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

Lane 1:
 3 μg NSD1
Lane 2:
 Protein Marker

MW: 55 kDa
Purity: ≥76%



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 **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com