

Data Sheet

SET7/9

Human, recombinant, N-terminal GST tag

Catalog #: 51010

Lot#: 120314 **Conc.:** 1.21 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 16 mM
 Glutathione, 20% Glycerol, and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Ivanov G.S. *et al.*, *Mol. Cell. Biol.* **27 (19)**, 6756-6769 (2007).
2. Couture J.F. *et al.*, *Nat. Struct. Mol. Biol.* **13 (2)**, 140-146 (2006).

Description:

Human recombinant SET7/9 (amino-acids 52-end) with N-terminal GST tag (GenBank Accession No. NM_030648), expressed in an *E. coli* expression system. MW=61 kDa.

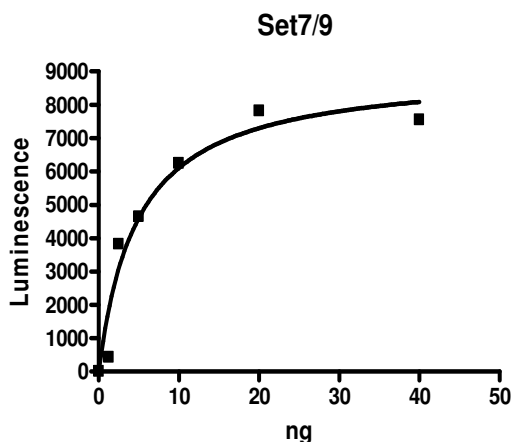
Specific Activity: 0.58 pmol/min/μg.

Assay conditions: 50 μl reaction mix (50 mM TRIS pH8.8, 5 mM MgCl₂, 4 mM DTT, 20 μM S-adenosylmethionine, and 0-10 ng SET7/9) add to the wells coated with the substrate. Incubate at room temperature for 1 hr. Add antibody against methylated K4 residue of histone H3, incubate 1 hr. Then incubate 30 mins with HRP-labeled secondary antibody. Finally, add HRP chemiluminescent substrates and read luminescence.

Application:

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

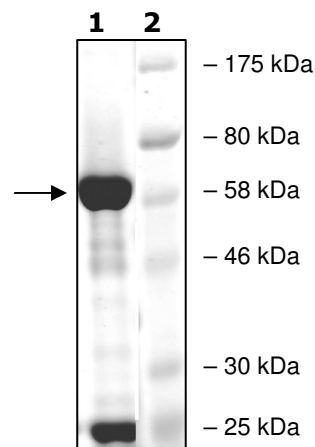
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 5 μg SET7/9
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 61 kDa
Purity:
 ≥55%



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