

Data Sheet

SET7/9

Human, recombinant, N-terminal His-GST tag
Catalog #: 50019
Lot#: 121012 **Conc.:** 0.26 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% glycerol.

Stability: >6 months at -80°C

References:

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

Description:

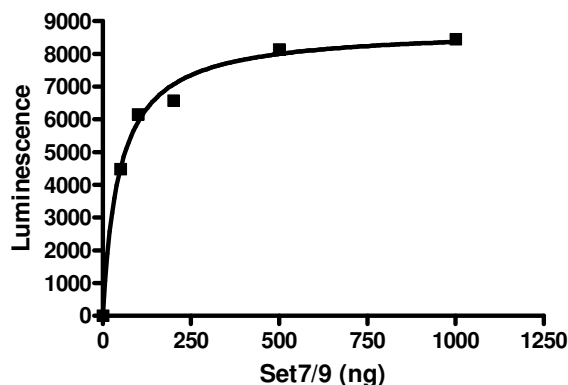
Human recombinant SET domain containing lysine methyltransferase 7 (SETD7), or SET7/9, (GenBank Accession No. NM_030648) a.a. 52-266 (end) with N-terminal His-GST tag, MW = 63 kDa, expressed in a baculovirus infected Sf9 cell expression system.

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.

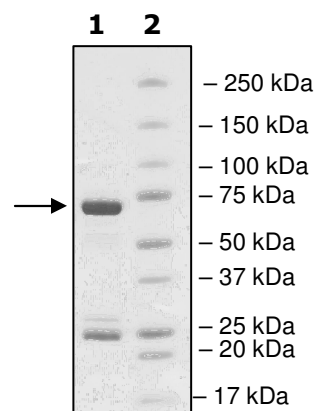
Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:
2.4 µg SET7/9
Lane 2:
Protein Marker

MW: 63 kDa
Purity: ≥62%





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

Amino Acid Sequence:

MHHHHHHGGGSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELG
LEFPNLPYYIDGDVKLTQSMAIIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRIA
YSKDFETLKVDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPM
CLDAFPKLVCFKKRIEAIPQIDKYLKSSKYIAWPLQGQWQATFGGGDHPKSDPENLYFQ
GGGEFFFFDGSTLEGYYVDDALQQGGVYTYEDGGVLQGTYYVDGELNGPAQEYDTDG
RLIFKGQYKDNIRHGVCWIYYPDGGSLVGEVNEDGEMTGEKIAYVYPDERTALYGKFI
DGEMIEGKLATLMSTEEGRPHFELMPGNSVYHFDKSTSSCISTNALLPDPYESERVYV
AESLISSAGEGLFSKVAVGPNTVMSFYNGVRITHQEVDSDRDWALNGNTLSLDEETVID
VPEPYNHVSKYCASLGHKANHSFTPNCIYDMFVHPRFGPIKCIRTLRAVEADEELTVA
YGYDHSPPGKSGPEAPEWYQVELKAFQATQQK

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