

Data Sheet

G9a (EHMT2)

Human, recombinant, N-terminal GST tag

Catalog#: 51001

Lot# 130827 Conc. 0.27 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
 124 mM NaCl, 2.4 Mm KCl, 3 mM DTT and
 10% glycerol.

Stability: >6 months at -70°C . *Avoid
 freeze/thaw cycles.*

References:

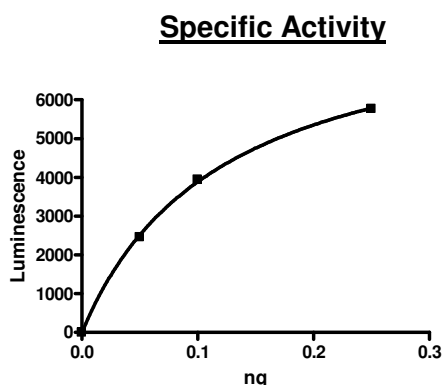
1. Yuan,X. et al., Mol. Cell **27 (4)**, 585-595 (2007).
2. Nishida,M. et al., Genes Cells **12 (7)**, 877-888 (2007).

Description: Human G9a (euchromatic histone-lysine N-methyltransferase 2) (GenBank Accession No. NM_006709) (amino-acids 785-1210), with N-terminal GST tag, MW = 74.6 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 73.7 pmol/min/ μg
 Assay conditions: 50 μl reaction mix (50 mM TRIS pH 8.8, 5mM MgCl₂, 4mM DTT, 0.05% Tween-20, 40 μM S-adenosylmethionine, and G9a (0.05 – 0.5 ng) add to the wells coated with the substrate on a GBioscience Neutravidin black plate. Incubate for 1 hr. Add antibody against methylated K9 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.

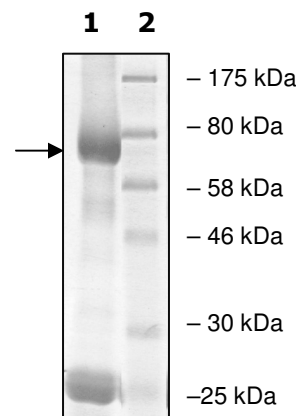
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 5 μg G9a
Lane 2:
 Protein Marker

MW: 74.6 kDa
Purity: $\geq 50\%$



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