

Data Sheet

G9a (EHMT2)

Human, recombinant, N-terminal GST tag

Catalog#: 51000

Lot#: 140327

Conc.: 0.46 mg/ml

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

Stability: >6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

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 an *E. Coli* expression system.

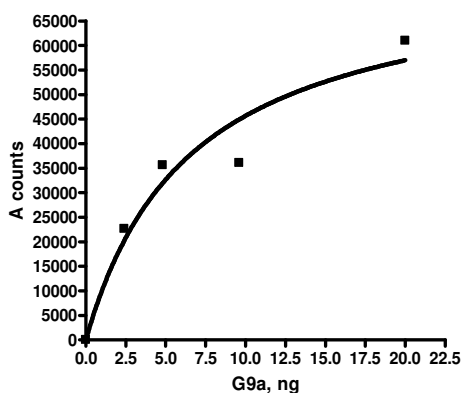
Specific Activity: 22 pmole/min/ μg

Assay Conditions: Enzyme was incubated with 0.5 μM of biotinylated H3 peptide with 40 μM AdoMet in HMT buffer 1A in a 10- μl reaction. 5 μl of Anti-Mouse AlphaLISA acceptor beads (1:250) and 5 μl of Primary Antibody 7 (1:50) were added and incubation continued for 30 min. Streptavidin Donor beads (1:125) were added (10 μl) and mixture was further incubated for 15 min at RT before A-counts were measured.

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

Quality Assurance

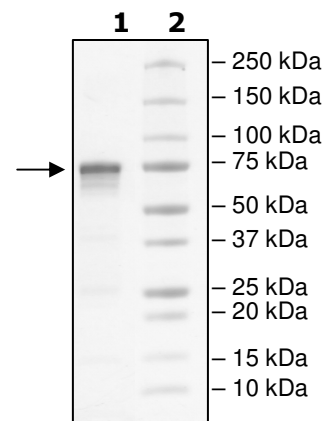
G9a Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg G9a
Lane 2:
Protein Marker

MW: 75.8 kDa
Purity: $\geq 80\%$



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