

G9a (EHMT2 (Euchromatic Histone-lysine N-Methyltransferase 2); H3-K9-HMTase 3; KMT1C)

CATALOG NO.: HMT-11-102

LOT NO.:

DESCRIPTION: Human recombinant G9a (residues 786-1210; Genbank Accession # NM_006709) expressed as an N-terminal GST-fusion protein in *E. coli*. MW = 80 kDa. Catalyzes the transfer of methyl groups from S-adenosyl-L-methionine (SAM) to the ε-amino function of protein L-lysine residues (mono-, di- and trimethylation), especially of lysine-9 of histone H3 (H3K9)¹, but with reported activities on H3K27¹, histone H1.4K26², p53 K373³ and other targets (see review⁴). G9a is a SET-domain type histone methyltransferase (HMT), which, in complex with the highly homologous GLP, is the major source of mono- and dimethylated histone H3K9 in euchromatin^{5,6}, marks associated with recruitment of HP1, DNA methylation and gene silencing⁶⁻⁸. A multimeric H3K9 methylation complex containing G9a/GLP along with other HMTs (SETDB1, SUV39H1) has been described⁹. G9a is overexpressed in a variety of cancers and knockdown of G9a/GLP in the MCF7 breast cancer line increases apoptosis³. These results, along with the fact that dimethylation at the G9a/GLP target site, p53 K373, correlates with levels of inactive p53, suggest G9a/GLP inhibition as a potential anti-cancer therapy, especially for tumors expressing wild-type p53³.

PURITY: >80% by SDS-PAGE.

ASSAY CONDITIONS: RBC's G9a displays histone methyltransferase activity at enzyme concentrations of 6.25 nM and above, 30°C, with calf thymus histone H3 or chicken core histones in the HMT HotSpotSM Assay format. Reaction conditions are: 50 mM Tris-HCl, pH 8.5, 50 mM NaCl, 5 mM MgCl₂, 1 mM DTT, 1 mM PMSF, 0.05 mg/mL chicken core histones (0.05 mg/mL) or calf thymus histone H3 (5 μM), [³H]-SAM.

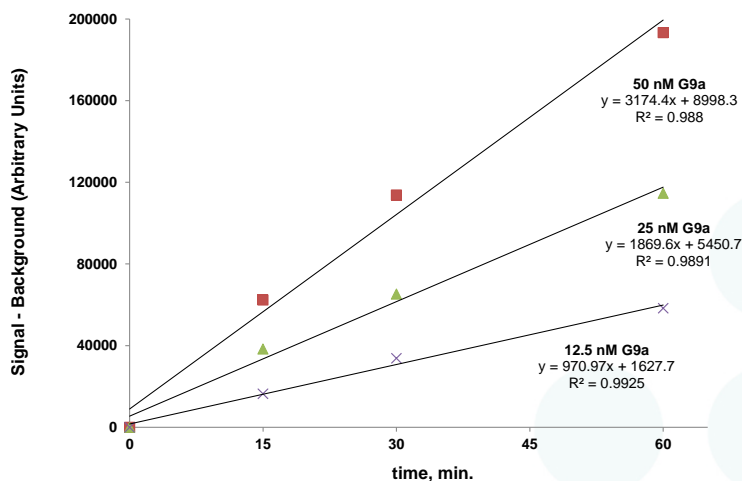
SUPPLIED AS: ___ μg/μl in 25.4 mM Na₂HPO₄, pH 7.4, 4.4 mM KH₂PO₄, 137 mM NaCl, 3 mM DTT, 30% (w/v) glycerol as determined by OD₂₈₀

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. Freezing and storage of diluted enzyme is not recommended.

REFERENCES: 1) M. Tachibana *et al. J. Biol. Chem.* 2001 **276** 25309; 2) P. Trojer *et al. J. Biol. Chem.* 2009 **284** 8395; 3) J. Huang *et al. J. Biol. Chem.* 2010 **285** 9636; 4) Y. Shinkai & M. Tachibana *Genes Dev.* 2011 **25** 781; 5) M. Tachibana *et al. Genes Dev.* 2002 **16** 1779; 6) M. Tachibana *et al. Genes Dev.* 2005 **19** 815; 7) N. Feldman *et al. Nat. Cell Biol.* 2006 **8** 188; 8) M. El Gazzar *et al. J. Biol. Chem.* 2008 **283** 32198; 9) L. Fritsch *et al. Mol. Cell* 2010 **37** 46



Coomassie blue stained SDS-PAGE (4-12% acrylamide) 4 μg of purified RBC G9a. MW marker (left), from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Time courses of G9a methyltransferase reactions in the HotSpotSM assay format. G9a, at the indicated concentrations, was assayed with 0.05 mg/mL chicken core histones and 1 μM [³H]-SAM. Signal/Background ranged from 4.3 (12.5 nM G9a, 15 min.) to 51.2 (50 nM G9a, 60 min.).

This product is not intended for therapeutic or diagnostic use in animals or in humans.