

SUV39H1

(Su(var)3-9 homolog 1; Histone H3-K9 methyltransferase 1)

CATALOG NO.: HMT-11-111

LOT NO.:

DESCRIPTION: Human recombinant SUV39H1 (residues 44-412; Genbank Accession # NM_003173) expressed as an N-terminal fusion protein, with a C-terminal His-tag in *E. coli*. MW = 99.5 kDa. Catalyzes the transfer of methyl groups from S-adenosyl-L-methionine (SAM) to the ε-amino function of protein L-lysine residues, especially to the mono- and dimethylated forms of lysine-9 of histone H3 (H3K9me1, H3K9me2)^{1,2}. A SET-domain lysine-methyltransferase, SUV39H1's activity is the major source of the transcriptionally repressive modification, H3K9me3, in heterochromatin^{1,3}. SUV39H1's role in facultative heterochromatin formation is regulated by SIRT1, both via recruitment and due to an activating deacetylation⁴. DBC1 (deleted in breast cancer 1) disrupts the SUV39H1-SIRT1 complex and also directly inhibits SUV39H1 methyltransferase activity⁵. SIRT1 also regulates global levels of SUV39H1 by inhibiting its polyubiquitination by MDM2⁶. SUV39H1 and SIRT1 are components of the eNoSC complex, which silences rRNA transcription in response to lowered cellular energy status⁷. A multimeric H3K9 methylation complex containing SUV39H1 along with other HMTs (SETDB1, G9a/GLP) has been described⁸. The oncoprotein Evi1, which is inappropriately expressed in AML (acute myeloid leukemia) and MDS (myelodysplastic syndrome), recruits SUV39H1, thus enhancing its transcription repressing activity^{9,10}. Depletion of SUV39H1 by shRNA or inhibition with chaetocin induced reexpression of silenced AML tumour suppressor genes, consistent with its potential as a therapeutic target¹¹.

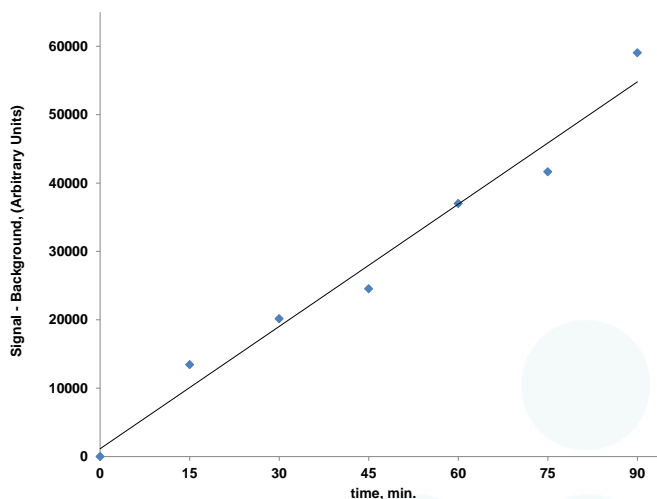
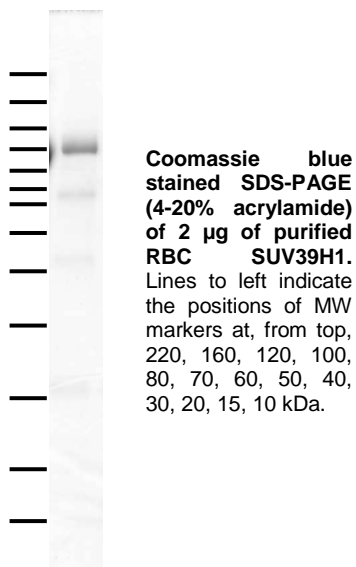
PURITY: >90% by SDS-PAGE.

ASSAY CONDITIONS: RBC's SUV39H1 displays histone methyltransferase activity at enzyme concentrations of 100 nM and above, 30°C, with H3K9-containing substrates, chicken core histones calf thymus histone H3 and H3(1-21) peptide 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), calf thymus histone H3 (5 μM), or H3(1-21) peptide (5 μM) and [³H]-SAM.

SUPPLIED AS: ___ μg/μl in 50 mM Tris/HCl, pH 8.0, 137 mM NaCl, 5 mM MgCl₂, 4 mM DTT, 20% (v/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) S. Rea *et al. Nature* 2000 **406** 593; 2) A.H. Peters *et al. Mol. Cell* 2003 **12** 1577; 3) A.H. Peters *et al. Cell* 2001 **107** 323; 4) A. Vaquero *Nature* 2007 **450** 440; Z. Li *et al. J. Biol. Chem.* 2009 **284** 10361; 6) L. Bosch-Presegue *et al. Mol. Cell* 2011 **42** 210; 7) A. Muruyama *et al. Cell* 2008 **133** 627; 8) L. Fritsch *et al. Mol. Cell* 2010 **37** 46; 9) D. Spensberger & R. Delwel *FEBS Lett.* 2008 **582** 2761; 10) F. Cattaneo & G. Nucifora *J. Cell. Biochem.* 2008 **105** 344; 11) A. Lakshmikuttyamma *et al. Oncogene* 2010 **29** 576



Time course of SUV39H1 methyltransferase reaction in the HotSpotSM assay format. SUV39H1, at 356 nM, was assayed with 5 μM H3(1-21) peptide and 1 μM [³H]-SAM.

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

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com