

**ASH1L**

(Ash1-Like Protein)

CATALOG NO.: HMT-11-376

LOT NO.:

**DESCRIPTION:** Human recombinant ASH1L (residues 2046-2330; Genbank Accession # NM\_018489; MW = 35.4 kDa) expressed in *E. coli* with an N-terminal His-tag. Catalyzes the transfer of methyl groups from S-adenosyl-L-methionine (SAM) to the ε-amino function of protein L-lysine residues, specifically lysine-36 of histone H3 (H3K36)<sup>1,2</sup>, a mark associated with active transcription. Activity has also been reported at H3K4<sup>3</sup>. A large, multi-domain protein associated with actively transcribed regions of chromatin, ASH1L is the human homolog of *Drosophila* Ash1, a Trithorax group protein. Like its counterpart in *Drosophila*<sup>4</sup>, ASH1L contains a SET histone methyltransferase domain and has been found to play a role in the regulation of Hox gene expression<sup>3,5</sup>. Although the ASH1L SET domain has been shown *in vitro* to methylate histone peptides on lysine-4 of histone H3 (H3K4)<sup>3</sup>, *in vivo* or *in vitro* with nucleosomes as substrate, ASH1L is an H3K36 methyltransferase<sup>1,2,6</sup>. Recruitment of ASH1L by the ncRNA *DBE-T* to the chromosome 4q35 locus associated with FSHD (facioscapulohumeral muscular dystrophy) leads to increased H3K36me2 and inappropriate gene derepression at the FSHD locus<sup>6</sup>. In a possible positive feedback loop, ASH1L increases expression of *DBE-T* itself<sup>6</sup>, suggesting ASH1L's methyltransferase activity or its interaction with *DBE-T* as potential therapeutic targets for FSHD. RBC's ASH1L comprises the catalytic domain (AWS/SET/Post-SET) fused to 6xHis-tag.

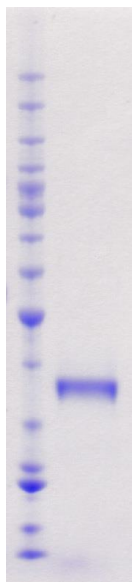
**PURITY:** >80% by SDS-PAGE

**ASSAY CONDITIONS:** RBC's ASH1L displays histone methyltransferase activity at concentrations of 60 nM and 100nM, 60 min. reactions, 30°C, as TCA-precipitated counts in a scintillation/filter plate assay (Multiscreen FB, Topcount), with HeLa oligo or mono/di-nucleosomes, Chicken oligo or mono/di nucleosomes and recombinant tetra or mononucleosomes (0.05 mg/mL as [DNA]). Reaction conditions are: 20 mM Tris, pH 9.0, 1 mM TCEP, 0.01% Triton X-100, 1.25mM MgCl<sub>2</sub>, substrates at concentrations indicated above.

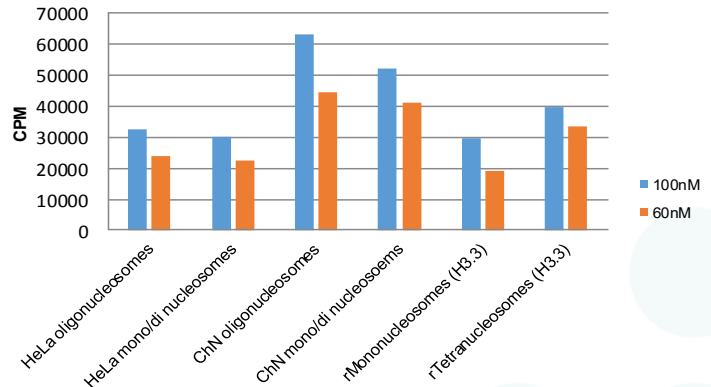
**SUPPLIED AS:** \_\_\_ µg/µL in 100 mM HEPES, pH 7.1, 500 mM NaCl, 0.5 mM TCEP, 5% (v/v) glycerol

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.

**REFERENCES:** 1) Y Tanaka *et al. Gene* 2007 **397** 161; 2) S. An *et al. J. Biol. Chem.* 2011 **286** 8369; 3) G. Gregory & *et al. Mol. Cell. Biol.* 2007 **27** 8466; 4) C. Beisel *et al. Nature* 2002 **419** 857; 5) Y Tanaka *et al. PLOS One* 2011 **6** e28171; 6) D.S. Cagianca *et al. Cell* 2012 **149** 819;



**Coomassie blue stained SDS-PAGE (4-20% acrylamide) of 2 µg of RBC ASH1L.** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Methylation Activity of ASH1L with various types of nucleosomes.** Assays were performed with a scintillation/filter plate assay. 100nM or 60nM ASH1L was incubated 60 min., 30°C with HeLa Oligonucleosomes (Cat. #HMT-35-130), HeLa Mononucleosome (Cat. # HMT-35-123), Chicken Oligonucleosomes (Cat. #HMT-35-177), Chicken Mononucleosomes (Cat. #HMT-35-179), Recombinant Tetranucleosomes (Cat. #HMT-15-367) and Recombinant Mononucleosomes (Cat. #HMT-15-369); 0.05 mg/mL as [DNA]) and 3.25 µM SAM (1:1.6[<sup>3</sup>H]-SAM:unlabeled).

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