

MLL3 Complex (Mixed Lineage Leukemia Protein-3 in complex with WDR5, RbBP5, Ash2L, (DPY-30)₂)

CATALOG NO.: HMT-15-107

LOT NO.:

DESCRIPTION: Human recombinant MLL3 (residues 4689-4911; Genbank Accession # NM_170606; MW = 27 kDa) in complex with human recombinants WDR5 (22-334; NM_017588; 34 kDa), RbBP5 (1-538; NM_005057; 61 kDa), Ash2L (2-534; NM_001105214; 63 kDa) and DPY-30 (1-99; NM_0325742; 13 kDa; two per complex). All proteins were expressed in *E. coli* with N-terminal His-tags. Catalyzes the transfer of methyl groups from S-adenosyl-L-methionine (SAM) to the ε-amino function of protein L-lysine residues, specifically lysine-4 of histone H3 (H3K4). The MLL3 subunit of this preparation comprises the catalytic SET domain¹ as well as the WDR5 interaction motif (*Win*)^{2,3} necessary for assembly of the active complex⁴⁻⁶. Like MLLs 1, 2 and 4, MLL3 complex methylates histone H3 lysine-4 (H3K4), conferring an activating epigenetic mark^{7,8}. MLL3 complex (or MLL4 complex) forms a part of the “activating signal cointegrator-2” (ASC-2) Complex or ASCOM^{9,10}. ASCOM acts as a coactivator of various nuclear receptors¹¹⁻¹⁴ and other transcription factors, including the tumor suppressor p53¹⁵. The chromatin remodeling complex Swi/Snf colocalizes with ASCOM, an interaction mediated by binding of the MLL3 (or MLL4) SET domain to the Swi/Snf subunit INI1¹⁶. ASCOM can be recruited to the *aP2* promoter by PPAR_γ and mice homozygous for MLL3 with a partial SET domain deletion have decreased white fat and a partially impaired potential for adipogenesis¹³. Knockdown of MLLs 3 and 4 diminishes estrogen-induced expression of HOXC10, which is overexpressed in breast cancer¹⁷.

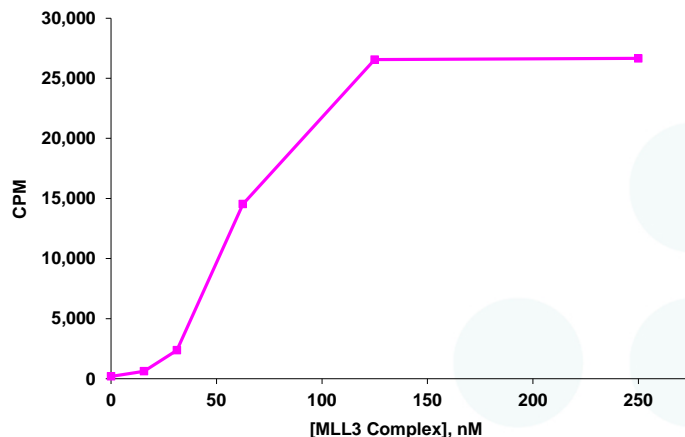
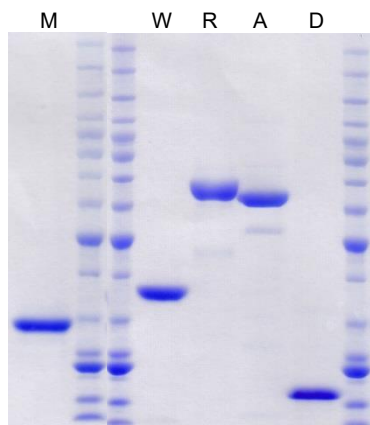
PURITY: >90% by SDS-PAGE.

ASSAY CONDITIONS: RBC's MLL3 Complex displays histone methyltransferase activity at MLL3 concentrations of 10 nM-1 μM, 15-90 min. reactions, 30°C, in the HMT HotSpotSM Assay format, with several H3K4-containing substrates (chicken core histones (0.05 mg/mL), calf-thymus histone H3 (5 μM), H3(1-21) peptide (5 μM)). Reaction conditions are: 50 mM Tris-HCl, pH 8.5, 50 mM NaCl, 5 mM MgCl₂, 1 mM DTT, 1 mM PMSF, protein or peptide substrate at concentrations indicated above, [³H]-SAM and a suitable dilution of “RBC MLL Enhancer” (provided). NOTE: Optimal MLL Enhancer dilution depends on substrate and assay system and must be determined by the user. This typically ranges from 1/10 to 1/100 dilutions, i.e. 10% to 1% of the final reaction volume. MLL3 Complex also displays activity in the absence of Enhancer, in a filter-binding assay with recombinant histone H3.3 as substrate (see Figure below; buffer conditions are the same as for HotSpotSM assays).

SUPPLIED AS: __ μM MLL3 Complex, as defined above, (__ μg/μl total protein) in 20 mM Tris-HCl, pH 7.5, 300 mM NaCl, 1 mM TCEP (tris(2-carboxyethyl)phosphine HCl), 10% (w/v) glycerol, 1 μM ZnCl₂ as determined by OD₂₈₀.

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 are not recommended.

REFERENCES: 1) T.A. Milne *et al. Mol. Cell* 2002 **10** 1107; 2) A. Patel *et al. J. Biol. Chem.* 2008 **283** 32162; 3) J.J. Song & R.E. Kingston *J. Biol. Chem.* 2008 **283** 35258; 4) A. Yokoyama *et al. Mol. Cell. Biol.* 2004 **24** 5639; 5) Y. Dou *et al. Nat. Struct. Mol. Biol.* 2006 **13** 713; 6) A. Patel *et al. J. Biol. Chem.* 2009 **284** 24242; 7) S. Glaser *et al. Development* 2006 **133** 1423; 8) K.I. Ansari & S.S. Mandal *FEBS J.* 2010 **277** 1790; 9) Y.H. Goo *et al. Mol. Cell. Biol.* 2003 **23** 140; 10) S. Lee *et al. Proc.Natl. Acad. Sci. USA* 2006 **103** 15392; 11) M.A. Mahajan & H.H. Samuels *Endocr. Rev.* 2005 **26** 583; 12) Q. Li *et al. Mol. Cell. Biol.* 2007 **27** 8073; 13) J. Lee *et al. Proc.Natl. Acad. Sci. USA* 2008 **105** 19229; 14) D. Kim *et al. Mol. Endocrinol.* 2009 **23** 1556; 15) J. Lee *et al. Proc.Natl. Acad. Sci. USA* 2009 **106** 8513; 16) S. Lee *et al. Mol. Endocrinol.* 2009 **23** 610; 17) K.I. Ansari *et al. J. Mol. Endocrinol.* 2011 doi: 10.1530/JME-11-0078



Methyltransferase Activity of MLL3 Complex. Methylation determined as TCA-precipitable counts in a scintillation/filter plate assay. Reactions were 25 μL, 60 min., 30°C, with 1 μM [³H]-SAM and 1 μM recombinant histone H3.3 (RBC Cat. # HMT-11-134) as substrates.

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