

## WRAD<sub>2</sub> Complex

(WDR5, RbBP5, Ash2L, (DPY-30)<sub>2</sub>)

CATALOG NO.: HMT-14-109

LOT NO.:

**DESCRIPTION:** 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). Members of the WRAD<sub>2</sub> complex are found in the active, assembled histone H3K4 methyltransferase complexes of all human SET1 family members, MLL1<sup>1-3</sup>, MLL2<sup>4</sup>, MLL3<sup>5,6</sup>, MLL4<sup>5,6</sup>, SET1A<sup>6</sup> and SET1B<sup>6</sup>. Binding interactions occur between WDR5 and a “WDR5 interaction motif” (*Win*)<sup>7,8</sup> in SET1 family histone methyltransferases, and between the subunit pairs WDR5-RbBP5, RbBP5-Ash2L and Ash2L-(DPY-30)<sub>2</sub><sup>9</sup>. Addition of the WRAD<sub>2</sub> subunits *in vitro* to recombinant MLL1 (3745-3969) stimulates methyltransferase activity about 600-fold<sup>9</sup>. Thus, WRAD<sub>2</sub> is considered an essential sub-complex of SET1 family H3K4 methyltransferases, necessary for complex assembly and full activity<sup>2,6,9</sup>. Moreover, recent work has shown that the WRAD<sub>2</sub> complex by itself has H3K4 methyltransferase activity, making it the first known non-SET domain/non-Dot1L lysine methyltransferase<sup>9,10</sup>. Interaction between MLL1 and WRAD<sub>2</sub> is required for sequential mono- and dimethylation of H3K4<sup>9</sup> and for methylation of nucleosomal histone H3<sup>10</sup>. The *Win* motif interacts with WDR5 at a site also implicated in binding the N-terminus of histone H3 and peptides based on both *Win* and H3 sequence have been found to disrupt the MLL1-WDR5 interaction and inhibit methyltransferase activity<sup>7,11</sup>. Unlike SET-domain inhibition, targeting of WRAD<sub>2</sub>'s MLL regulatory interactions or its intrinsic methyltransferase activity would represent routes to MLL inhibition that would be specific to the SET1 family.

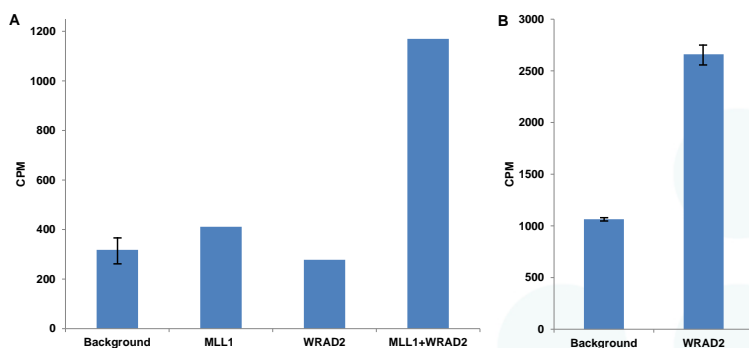
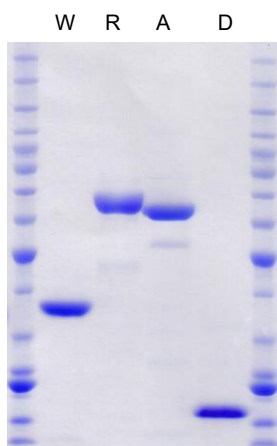
**PURITY:** >90% by SDS-PAGE.

**ASSAY CONDITIONS:** RBC's WRAD<sub>2</sub> Complex displays methyltransferase activity with an H3(1-21) peptide in a reverse phase resin binding assay. Reaction conditions are: 100 mM CHES/KOH, pH 8.5, 5 mM MgCl<sub>2</sub>, 0.8 mM DTT, protein or peptide substrate, [<sup>3</sup>H]-SAM.

**SUPPLIED AS:** \_\_\_ μM WRAD<sub>2</sub> 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 ZnCl<sub>2</sub> as determined by OD<sub>280</sub>.

**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) A. Yokoyama *et al. Mol. Cell. Biol.* 2004 **24** 5639; 2) Y. Dou *et al. Nat. Struct. Mol. Biol.* 2006 **13** 713; 3) M.M. Steward *et al. Nat. Struct. Mol. Biol.* 2006 **13** 852; 4) R. Mo *et al. J. Biol. Chem.* 2006 **281** 15714; 5) S. Lee *et al. Proc. Natl. Acad. Sci USA* 2006 **103** 15392; 6) Y. Cho *et al. J. Biol. Chem.* 2007 **282** 20395; 7) A. Patel *et al. J. Biol. Chem.* 2008 **283** 32162; 8) J.J. Song & R.E. Kingston *J. Biol. Chem.* 2008 **283** 35258; 9) A. Patel *et al. J. Biol. Chem.* 2009 **284** 24242; 10) A. Patel *et al. J. Biol. Chem.* 2011 **286** 3359; 11) H. Karatas *et al. J. Med. Chem.* 2010 **53** 5179



**WRAD<sub>2</sub> Complex Stimulates MLL1 Activity (A) and Displays Intrinsic Methyltransferase Activity (B).** Reactions (10 μl) were incubated 30 min., 30°C and stopped by denaturation with guanidinium HCl. Proteins/peptides were separated from unreacted SAM by binding to a reverse phase resin (C18) and eluted for scintillation counting. Background samples were denatured prior to SAM addition. Both MLL1 and WRAD<sub>2</sub> were 1 μM. **A.** Substrates were 0.05 mg/mL chicken core histones, 1 μM [<sup>3</sup>H]-SAM. **B.** Substrates were 50 μM H3(1-21) peptide and 10 μM [<sup>3</sup>H]-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