

## MLL4

(Complex of **MLL4**: Histone-lysine N-methyltransferase MLL4, WW domain-binding protein 7, HRX2, KIAA0304, KMT2D, TRX2, WBP7, **WDR5**: WD repeat-containing protein 5, BMP2-induced 3-kb gene protein, **RBBP5**: Retinoblastoma-binding protein 5, Retinoblastoma-binding protein RBQ-3, **ASH2L**: Set1/Ash2 histone methyltransferase complex subunit ASH2, ASH2-like protein, **DPY30**: Protein dpy-30 homolog, Dpy-30-like protein)

- **CAT#:** MLL4
- **Enzyme:** Human MLL4; 5 component complex
- **Substrate:** Core Histone, 0.05 mg/ml  
SAM, 1  $\mu$ M
- **Reaction:**  
S-adenosyl-L-[methyl-<sup>3</sup>H]methionine + histone L-arginine = S-adenosyl-L-homocysteine + histone [methyl-<sup>3</sup>H]-L-arginine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	8,300
Sinefungin	42,000
BIX01294	>500,000
Chaetocin	310

- **Scientific Information:** **MLL4** Q9UMN6 (<http://www.expasy.org/uniprot/Q9UMN6>)  
**WDR5** P61964 (<http://www.expasy.org/uniprot/P61964>), **RBBP5** Q15291 (<http://www.expasy.org/uniprot/Q15291>), **ASH2L** Q9UBL3 (<http://www.expasy.org/uniprot/Q9UBL3>), **DPY30** Q9C005 (<http://www.expasy.org/uniprot/Q9C005>),

### Reference Compound IC50 for MLL4

