

## DOT1L

(Histone-lysine N-methyltransferase, H3 lysine-79 specific, Histone H3-K79 methyltransferase, H3-K79-HMTase, DOT1-like protein, Lysine N-methyltransferase 4, KIAA1814, KMT4)

- **CAT#:** DOT1L
- **Enzyme:** Human DOT1L
- **Substrate:** Core Histone, 0.05 mg/ml  
SAM, 250 nM
- **Reaction:**  
S-adenosyl-L-[methyl-<sup>3</sup>H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N<sup>6</sup>-[methyl-<sup>3</sup>H]-L-lysine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	540
Sinefungin	104,000

- **Scientific Information:** Q8TEK3 (<http://www.expasy.org/uniprot/Q8TEK3>)

### Reference Compound IC50 for DOT1L

