

SET7/9

(Histone-lysine N-methyltransferase SETD7, Histone H3-K4 methyltransferase SETD7, H3-K4-HMTase SETD7, SET domain-containing protein 7, SET7/9, Lysine N-methyltransferase 7, KIAA1717, KMT7, SET7, SET9)

- **CAT#:** SET7/9
- **Enzyme:** Human SET7/9
- **Substrate:** Histone H3, 5 μ M
SAM, 1 μ M
- **Reaction:**
S-adenosyl-L-[methyl- 3 H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N 6 -[methyl- 3 H]-L-lysine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	132,000
Sinefungin	9,100

- **Scientific Information:** Q8WTS6 (<http://www.expasy.org/uniprot/Q8WTS6>)

Reference Compound IC50 for SET7

