

SET8

(Histone-lysine N-methyltransferase SETD8, H4-K20-HMTase SETD8, SET domain-containing protein 8, PR/SET domain-containing protein 07, Lysine N-methyltransferase 5A, KMT5A, PRSET7, SET07)

- **CAT#:** SET8
- **Enzyme:** Human SET8
- **Substrate:** Histone H4, 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	171,000
Sinefungin	203,000

- **Scientific Information:** Q9NQR1 (<http://www.expasy.org/uniprot/Q9NQR1>)

Reference Compound IC50 for SET8

