

SMYD2

(SET and MYND domain-containing protein 2, HSKM-B
Lysine N-methyltransferase 3C, KMT3C)

- **CAT#:** SMYD2
- **Enzyme:** Human SMYD2
- **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	526
Sinefungin	15,000

- **Scientific Information:** Q9NRG4 (<http://www.expasy.org/uniprot/Q9NRG4>)

Reference Compound IC50 for SMYD2

