

SUV39H2

(Histone-lysine N-methyltransferase SUV39H2, Suppressor of variegation 3-9 homolog 2, Su(var)3-9 homolog 2, Histone H3-K9 methyltransferase 2, H3-K9-HMTase 2, Lysine N-methyltransferase 1B, KMT1B, SUV92)

- **CAT#:** SUV39H2
- **Enzyme:** Human SUV39H2
- **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	24,000
Sinefungin	103,000
Chaetocin	480

- **Scientific Information:** Q9H511 (<http://www.expasy.org/uniprot/Q9H511>)

Reference Compound IC50 for SUV39H2

