

DNMT3A

(DNA (cytosine-5)-methyltransferase 3A, DNA methyltransferase HsaIII A, DNMT3A)

- **CAT#:** DNMT3A
- **Enzyme:** Human DNMT3A
- **Substrate:** Lambda DNA, 0.0075 mg/ml
SAM, 1 μ M
- **Reaction:**
S-adenosyl-L-[methyl- 3 H]methionine + DNA = S-adenosyl-L-homocysteine + DNA 5-[methyl- 3 H]-cytosine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	271
Sinefungin	430

- **Scientific Information:** Q9Y6K1 (<http://www.expasy.org/uniprot/Q9Y6K1>)

Reference Compound IC50 for DNMT3A

