

## DNMT3B

(DNA (cytosine-5)-methyltransferase 3B, DNA methyltransferase HsaIIIB, DNMT3B)

- **CAT#:** DNMT3B
- **Enzyme:** Human DNMT3B-full length
- **Substrate:** Lambda DNA, 0.0075 mg/ml  
SAM, 1  $\mu$ 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	122
Sinefungin	140

- **Scientific Information:** Q9UBC3 (<http://www.expasy.org/uniprot/Q9UBC3>)

### Reference Compound IC50 for DNMT3B

