

DNMT1

(DNA (cytosine-5)-methyltransferase 1, CXXC-type zinc finger protein 9,
DNA methyltransferase HsaI, MCMT, AIM, CXXC9, DNMT)

- **CAT#:** DNMT1
- **Enzyme:** Human DNMT1
- **Substrate:** Poly(dI-dC), 0.001 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	270
Sinefungin	7,400

- **Scientific Information:** P26358 (<http://www.expasy.org/uniprot/P26358>)

Reference Compound IC50 for DNMT1

