

PRMT3

(Protein arginine N-methyltransferase 3, Heterogeneous nuclear ribonucleoprotein methyltransferase-like protein 3, HRMT1L3, ANM3)

- **CAT#:** PRMT3
- **Enzyme:** Human PRMT3
- **Substrate:** Histone H3, 5 μ M
SAM, 1 μ M
- **Reaction:**
S-adenosyl-L-[methyl- 3 H]methionine + histone L-arginine = S-adenosyl-L-homocysteine + histone [methyl- 3 H]-L-arginine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	665
Sinefungin	775

- **Scientific Information:** O60678 (<http://www.expasy.org/uniprot/O60678>)

Reference Compound IC50 for PRTMT3

