

## PRMT6

(Protein arginine N-methyltransferase 6, Heterogeneous nuclear ribonucleoprotein methyltransferase-like protein 6, Histone-arginine N-methyltransferase PRMT6, HRMT1L6, ANM6)

- **CAT#:** PRMT6
- **Enzyme:** Human PRMT6
- **Substrate:** Histone H3, 5  $\mu$ M  
SAM, 1  $\mu$ M
- **Reaction:**  
S-adenosyl-L-[methyl-<sup>3</sup>H]methionine + histone L-arginine = S-adenosyl-L-homocysteine + histone [methyl-<sup>3</sup>H]-L- arginine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	140
Sinefungin	580
BIX01294	>200,000
Chaetocin	>200,000

- **Scientific Information:** Q96LA8 (<http://www.expasy.org/uniprot/Q96LA8>)

### Reference Compound IC50 for PRMT6

