

GLP

(Histone-lysine N-methyltransferase EHMT1, Euchromatic histone-lysine N-methyltransferase 1, G9a-like protein 1, EUHMTASE1, GLP, KIAA1876, KMT1D)

- **CAT#:** GLP
- **Enzyme:** Human GLP
- **Substrate:** Core Histone, 0.05 mg/ml
SAM, 1 μ M
- **Reaction:**
S-adenosyl-L-[methyl- 3 H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N 6 -[methyl- 3 H]-L-lysine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
SAH	460
Sinefungin	13,000
BIX 01294	19,000
Chaetocin	690

- **Scientific Information:** Q9H9B1 (<http://www.expasy.org/uniprot/Q9H9B1>)

Reference Compound IC50 for GLP

