

## SETMAR

(Histone-lysine N-methyltransferase SETMAR, SET domain and mariner transposase fusion gene-containing protein, Metnase, Hsmar1, SETMR)

- **CAT#:** SETMAR
- **Enzyme:** Human SETMAR
- **Substrate:** Histone H3, 5  $\mu$ M  
SAM, 1  $\mu$ 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	739
Sinefungin	525

- **Scientific Information:** Q53H47 (<http://www.expasy.org/uniprot/Q53H47>)

### Reference Compound IC50 for SETMAR

