

## EZH2

(**EZH2**: Polycomb protein EZH2, Enhancer of zeste homolog 2, ENX-1, KMT6, **EED**: WD protein associating with integrin cytoplasmic tails 1, **SUZ12**: Suppressor of zeste 12 protein homolog, **RBAP48**: Histone-binding protein RBBP4, **AEBP2**: Zinc finger protein AEBP2)

- **CAT#:** EZH2
- **Enzyme:** Human EZH2; 5 component complex
- **Substrate:** Histone H3, 5  $\mu$ M  
SAM, 250 nM
- **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	46,000
Sinefungin	280,000

- **Scientific Information:** **EZH2** Q15910 (<http://www.expasy.org/uniprot/Q15910>)  
**EED** O75530 (<http://www.expasy.org/uniprot/O75530>), **SUZ12** Q15022 (<http://www.expasy.org/uniprot/Q15022>), **RBBP4** Q09028 (<http://www.expasy.org/uniprot/Q09028>)  
**AEBP2** Q6ZN18 (<http://www.expasy.org/uniprot/Q6ZN18>)

### Reference Compound IC50 for EZH2

