

CBP

(CREB-binding protein, CREBBP)

- **CAT#:** CBP
- **Enzyme:** Human CBP
- **Substrate:** Histone H3, 5 μ M
Acetyl CoA, 3 μ M
- **Reaction:**
[Acetyl-³H]-Acetyl Coenzyme A + histone L-lysine = Coenzyme A + [Acetyl-³H]-L-lysine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Anacardic Acid	20,000
Garcinol	14,000
Curcumin	37,000

- **Scientific Information:** Q92793 (<http://www.expasy.org/uniprot/Q92793>)

Reference Compound IC50 for CBP

