

KAT5

(Histone acetyltransferase KAT5, 60 kDa Tat-interactive protein,
Lysine acetyltransferase 5, cPLA(2)-interacting protein, HTATIP, TIP60)

- **CAT#:** KAT5
- **Enzyme:** Human KAT5
- **Substrate:** Histone H2A, 5 μ M
Acetyl CoA, 3 μ M
- **Reaction:**
[Acetyl- 3 H]-Acetyl Coenzyme A + histone L-lysine = Coenzyme A + [Acetyl- 3 H]-L-lysine.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Anacardic Acid	30,000
Garcinol	11,000
Curcumin	25,000

- **Scientific Information:** Q92993 (<http://www.expasy.org/uniprot/Q92993>)

Reference Compound IC50 for KAT5

