

pCAF

(Histone acetyltransferase KAT2B, Lysine acetyltransferase 2B
Histone acetyltransferase PCAF, P300/CBP-associated factor, PCAF, KAT2B)

- **CAT#:** pCAF
- **Enzyme:** Human pCAF
- **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	41,700
Garcinol	69,000

- **Scientific Information:** Q92831 (<http://www.expasy.org/uniprot/Q92831>)

Reference Compound IC50 for pCAF

