

GCN5

(Histone acetyltransferase KAT2A, Lysine acetyltransferase 2A
Histone acetyltransferase GCN5, General control of amino acid synthesis protein 5-like 2
STAF97, GCN5L2, HGCN5, KAT2A)

- **CAT#:** GCN5
- **Enzyme:** Human GCN5
- **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	66,000
Garcinol	64,000
- **Scientific Information:** Q92830 (<http://www.expasy.org/uniprot/Q92830>)

Reference Compound IC50 for GCN5

