

MYST4

(Histone acetyltransferase KAT6B, MOZ, YBF2/SAS3, SAS2 and TIP60 protein 4, Monocytic leukemia zinc finger protein-related factor, KIAA0383, MORF, MOZ2)

- **CAT#:** MYST4
- **Enzyme:** Human MYST4
- **Substrate:** Histone H4, 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	20,000
Garcinol	17,500
Curcumin	14,000

- **Scientific Information:** Q8WYB5 (<http://www.expasy.org/uniprot/Q8WYB5>)

Reference Compound IC50 for MYST4

