

HDAC1

(Histone deacetylase 1, Class I)

- **CAT#:** HDAC1
- **Enzyme:** Human HDAC1
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	6.8
SAHA	306

- **Scientific Information:** Q13547 (<http://www.expasy.org/uniprot/Q13547>)

Reference Compound IC50 for HDAC1

