

## HDAC1

(Histone deacetylase 1, Class I)

- **CAT#:** HDAC1
- **Enzyme:** Human HDAC1
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	21
SAHA	171

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

### Reference Compound IC50 for HDAC1

