

HDAC4

(Histone Deacetylase 4, Class IIa)

- **CAT#:** HDAC4
- **Enzyme:** Human HDAC4
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	36.1
SAHA	282

- **Scientific Information:** P56524 (<http://www.expasy.org/uniprot/P56524>)

Reference Compound IC50 for HDAC4

