

HDAC6

(Histone Deacetylase 6, Class IIb)

- **CAT#:** HDAC6
- **Enzyme:** Human HDAC6
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|----------------|-----------|
| Trichostatin A | 1.11 |
| SAHA | 19.8 |

- **Scientific Information:** Q9UBN7 (<http://www.expasy.org/uniprot/Q9UBN7>)

Reference Compound IC50 for HDAC6

