

## HDAC6

(Histone Deacetylase 6, Class IIb)

- **CAT#:** HDAC6
- **Enzyme:** Human HDAC6
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	0.76
SAHA	9.3

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

### Reference Compound IC50 for HDAC6

