

## HDAC5

(Histone Deacetylase 5, Class IIa)

- **CAT#:** HDAC5
- **Enzyme:** Human HDAC5
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

| Compounds      | IC50 (nM) |
|----------------|-----------|
| Trichostatin A | 15.7      |
| SAHA           | 400       |

- **Scientific Information:** Q9UQL6 (<http://www.expasy.org/uniprot/Q9UQL6>)

### Reference Compound IC50 for HDAC5

