

## HDAC3

(Histone Deacetylase 3, Class I)

- **CAT#:** HDAC3
- **Enzyme:** Human HDAC3
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	9.7
SAHA	200

- **Scientific Information:** O15379 (<http://www.expasy.org/uniprot/O15379>)

### Reference Compound IC50 for HDAC3

