

## HDAC7

(Histone deacetylase 7a, HDAC7A, Class IIa)

- **CAT#:** HDAC7
- **Enzyme:** Human HDAC7
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	99
SAHA	510

- **Scientific Information:** Q8WUI4 (<http://www.expasy.org/uniprot/Q8WUI4>)

### Reference Compound IC50 for HDAC7

