

## HDAC8

(Histone Deacetylase 8, Class I)

- **CAT#:** HDAC8
- **Enzyme:** Human HDAC8
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHK<sub>Ac</sub>K<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	110
SAHA	306

- **Scientific Information:** Q9BY41 (<http://www.expasy.org/uniprot/Q9BY41>)

### Reference Compound IC50 for HDAC8

