

HDAC10

(Histone deacetylase 10, Class IIb)

- **CAT#:** HDAC10
- **Enzyme:** Human HDAC10
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	29.5
SAHA	199

- **Scientific Information:** Q969S8 (<http://www.expasy.org/uniprot/Q969S8>)

Reference Compound IC50 for HDAC10

