

HDAC11

(Histone deacetylase 11, HD11, HDA11)

- **CAT#:** HDAC11
- **Enzyme:** Human HDAC11
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	12.1
SAHA	200

- **Scientific Information:** Q96DB2 (<http://www.expasy.org/uniprot/Q96DB2>)

Reference Compound IC50 for HDAC11

