

## HDAC2

(Histone Deacetylase 2, Class I)

- **CAT#:** HDAC2
- **Enzyme:** Human HDAC2
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	21.4
SAHA	389

- **Scientific Information:** Q92769 (<http://www.expasy.org/uniprot/Q92769>)

### Reference Compound IC50 for HDAC2

