

HDAC9

(Histone deacetylase 9, MEF2-interacting transcription repressor MITR, HDAC7B, Class IIa)

- **CAT#:** HDAC9
- **Enzyme:** Human HDAC9
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	79.5
SAHA	165

- **Scientific Information:** Q9UKV0 (<http://www.expasy.org/uniprot/Q9UKV0>)

Reference Compound IC50 for HDAC9

