

HDAC9

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

- **CAT#:** HDAC9
- **Enzyme:** Human HDAC9
- **Substrate:** Fluorogenic HDAC Class2a Substrate (Boc-Lys(trifluoroacetyl)-AMC), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	7,500
SAHA	141,000

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

Reference Compound IC50 for HDAC9

