

HDAC5

(Histone Deacetylase 5, Class IIa)

- **CAT#:** HDAC5
- **Enzyme:** Human HDAC5
- **Substrate:** Fluorogenic HDAC Class2a Substrate (Boc-Lys(trifluoroacetyl)-AMC), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	2,800
SAHA	27,200

- **Scientific Information:** Q9UQL6 (<http://www.expasy.org/uniprot/Q9UQL6>)

Reference Compound IC50 for HDAC5

