

HDAC7

(Histone deacetylase 7a, HDAC7A, Class IIa)

- **CAT#:** HDAC7
- **Enzyme:** Human HDAC7
- **Substrate:** Fluorogenic HDAC Class2a Substrate (Boc-Lys(trifluoroacetyl)-AMC), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	4,000
SAHA	105,000

- **Scientific Information:** Q8WUI4 (<http://www.expasy.org/uniprot/Q8WUI4>)

Reference Compound IC50 for HDAC7

