

## HDAC4

(Histone Deacetylase 4, Class IIa)

- **CAT#:** HDAC4
- **Enzyme:** Human HDAC4
- **Substrate:** Fluorogenic HDAC Class2a Substrate (Boc-Lys(trifluoroacetyl)-AMC), 50  $\mu$ M
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Trichostatin A	7,300
SAHA	76,000

- **Scientific Information:** P56524 (<http://www.expasy.org/uniprot/P56524>)

### Reference Compound IC50 for HDAC4

