

HDAC4

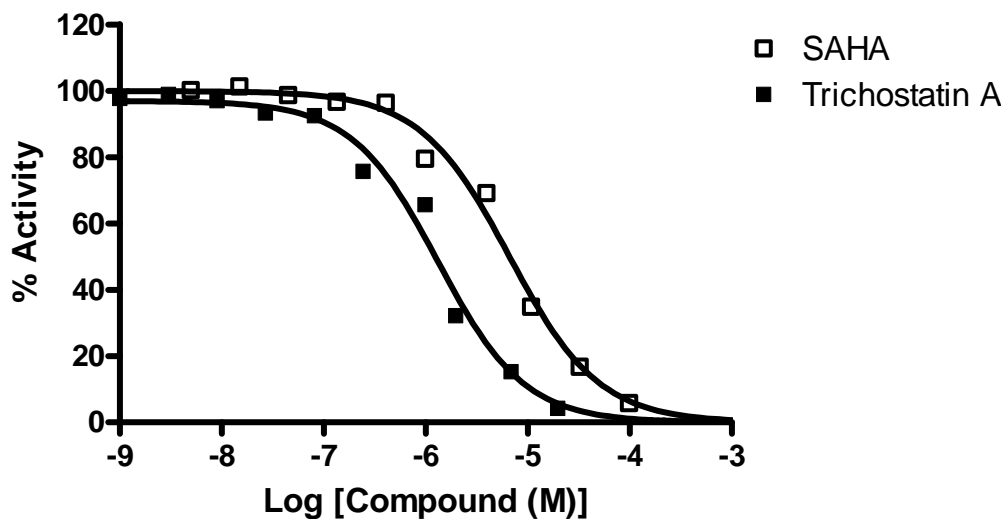
(Histone Deacetylase 4, Class IIa)

- **CAT#:** HDAC4
- **Enzyme:** Human HDAC4
- **Substrate:** Fluorogenic HDAC Substrate Class2a (Acetyl-Lys(trifluoroacetyl)-AMC), 50 μ M
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
Trichostatin A	1.321
SAHA	6.638

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

Control Compound IC50 for HDAC-4



HDAC5

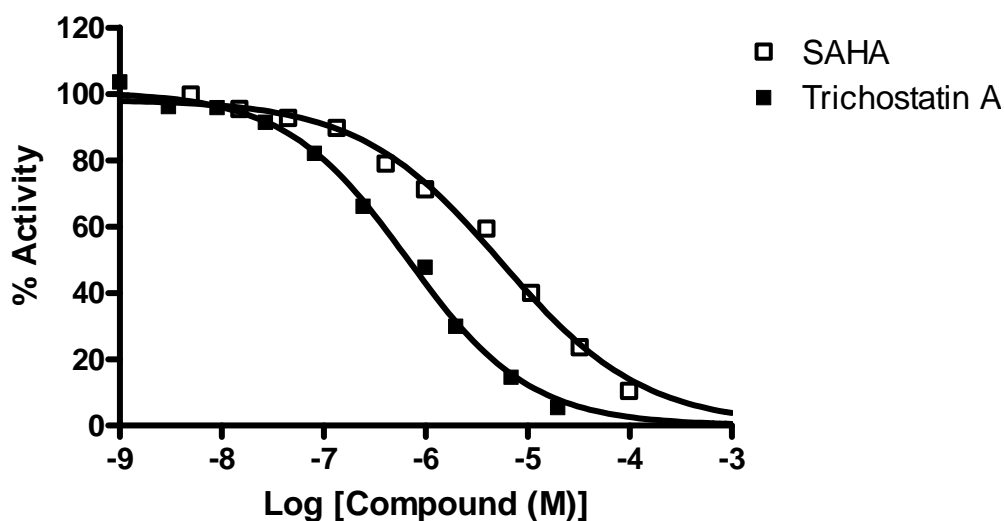
(Histone Deacetylase 5, Class IIa)

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

Compounds	IC50 (μ M)
Trichostatin A	0.662
SAHA	5.511

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

Control Compound IC50 for HDAC-5



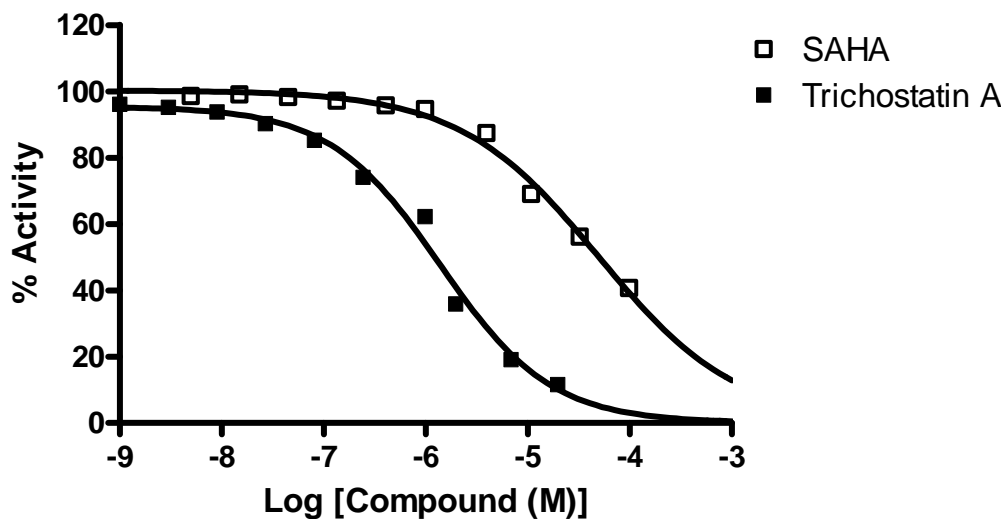
HDAC7

(Histone Deacetylase 7, Class IIa)

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

Compounds	IC50 (μ M)
Trichostatin A	1.386
SAHA	49.48
- **Scientific Information:** Q8WUI4 (<http://www.expasy.org/uniprot/Q8WUI4>)

Control Compound IC50 for HDAC-7



HDAC9

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

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

Compounds	IC50 (μ M)
Trichostatin A	3.282
SAHA	56.28

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

Control Compound IC50 for HDAC-9

