

Nuclear Extract

(Nuclear Extract of Histone deacetylases)

- **CAT#:** Nuclear Extract
- **Enzyme:** Nuclear Extract from HeLa cells
- **Substrate:** Fluorogenic peptide (K_{Ac}), 50 μ M
- **Reference Compound IC50s:**

Compounds	Trichostatin A
IC50 (nM)	1.3

Reference Compound IC50 for Nuclear Extract

