

## SIRT1

(Sirtuin 1, protein-lysyl-deacetylase, Class III HDAC)

- **CAT#:** SIRT1
- **Enzyme:** Human SIRT1 (NAD<sup>+</sup>-dependent enzyme)
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), NAD<sup>+</sup>
- **Reference Compound IC50s:**

Compounds	IC50 (μM)
Suramin sodium	2.42
Resveratrol	activator

- **Scientific Information:** Q96EB6 (<http://www.pir.uniprot.org/cgi-bin/upEntry?id=Q96EB6>)

### Reference Compound IC50 for SIRT1

