

## SIRT3

(Sirtuin 3, protein-lysyl-deacetylase, Class III HDAC)

- **CAT#:** SIRT3
- **Enzyme:** Human SIRT3 (NAD<sup>+</sup>-dependent enzyme)
- **Substrate:** Fluorogenic peptide from p53 residues 317-320 (QPKK<sub>Ac</sub>), NAD<sup>+</sup>
- **Reference Compound IC50s:**

<b>Compounds</b>	Suramin sodium
<b>IC50 (μM)</b>	273

- **Scientific Information:** Q9NTG7 (<http://www.pir.uniprot.org/cgi-bin/upEntry?id=Q9NTG7>)

### Reference Compound IC50 for SIRT3

