

## SIRT2

(Sirtuin 2, protein-lysyl-deacetylase, Class III HDAC)

- **CAT#:** SIRT2
- **Enzyme:** Human SIRT2 (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>	9.8

- **Scientific Information:** Q8IXJ6 (<http://www.pir.uniprot.org/cgi-bin/upEntry?id=Q8IXJ6>)

### Reference Compound IC50 for SIRT2

