

## SIRT3

(NAD-dependent deacetylase sirtuin-3, mitochondrial,  
SIR2-like protein 3, SIR2L3, 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>), 50 μM  
NAD<sup>+</sup>, 500 μM
- **Reference Compound IC50s:**

Compounds	IC50 (μM)
Nicotinamide	35.1

- **Scientific Information:** Q9NTG7 (<http://www.expasy.org/uniprot/Q9NTG7>)

### Reference Compound IC50 for SIRT3

