

SIRT3 Sirtuin Assay Services

SIRT3 [Sirtuin Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

Target Overview

CAT#	SIRT3
Enzyme	Human SIRT3 (NAD ⁺ -dependent enzyme)
Synonyms	NAD-dependent deacetylase sirtuin-3, mitochondrial, SIR2-like protein 3, SIR2L3, Class III HDAC

Assay Properties

Readout	Enzymatic activity
Substrates	Fluorogenic peptide from p53 residues 379-382 (RHKK _{Ac}), 50 μ M NAD ⁺ , 500 μ M
Reference Compound IC50s	Compounds IC50 (μ M) Nicotinamide 35.1
Scientific Information	Q9NTG7 (http://www.uniprot.org/uniprot/Q9NTG7)
Screening Location	Malvern, USA

Reference Compound IC50 for SIRT3

