

SIRT2 Sirtuin Assay Services

SIRT2 [Sirtuin Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

Target Overview

CAT#	SIRT2
Enzyme	Human SIRT2 (NAD ⁺ -dependent enzyme)
Synonyms	Sirtuin 2, protein-lysyl-deacetylase, Class III HDAC

Assay Properties

Readout	Enzymatic activity	
Substrates	Fluorogenic peptide from p53 residues 379-382 (RHKK _{Ac}), 50μM NAD ⁺ , 500 μM	
Reference Compound IC50s	Compounds	IC50 (μM)
	Suramin Sodium	9.8
Scientific Information	Q8IXJ6 (http://www.uniprot.org/uniprot/Q8IXJ6)	
Screening Location	Malvern, USA	

Reference Compound IC50 for SIRT2

