

SIRT5 Sirtuin Assay Services

SIRT5 [Sirtuin Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

SIRT5

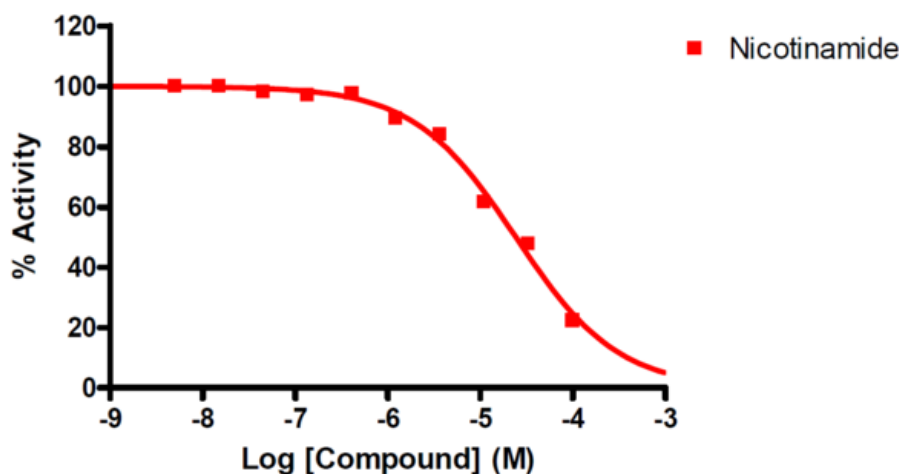
Target Overview

Enzyme	Human SIRT5 (NAD ⁺ -dependent enzyme)
Synonyms	NAD-dependent protein deacylase sirtuin-5, Regulatory protein SIR2 homolog 5, SIR2-like protein 5, SIR2L5, Class III HDAC

Assay Properties

Readout	Enzymatic activity
Substrates	Fluorogenic peptide (Ac-Lys-succ-AMC), 50 μ M NAD ⁺ , 500 μ M
Reference Compound IC50s	Compounds IC50 (μ M) Nicotinamide 24.2
Scientific Information	Q9NXA8 (http://www.uniprot.org/uniprot/Q9NXA8)
Screening Location	Malvern, USA

Reference Compound IC50 for SIRT5



Let's Discover Together.

© Dec 2023 **Reaction Biology**. All rights reserved.

https://reactionbiology.entwicklungsserver.de/datasheet/sirt5_hdac_malvern

