
Sirtinol

(Catalog No. M45002)

Alternative Names

Naphthol, 2-[(2-Hydroxynaphthalen-1-yl-methylene)amino]-N-(1-phenethyl)benzamide

Description

Sirtinol is a cell permeable specific inhibitor of sirtuin NAD-dependant deacetylases. Sirtinol has been shown to inhibit yeast Sir2 at IC₅₀ = 70 μM, human SIRT2 at IC₅₀ = 40 μM human SIRT1 at IC₅₀ = 131 μM, yet has shown no effect on human HDAC1. In the Arabidopsis plant Sirtinol strongly affected root and vascular tissue development by activating the auxin signaling pathway. Sirtinol has also been shown to induce senescence-like growth arrest in human breast cancer MCF-7 and lung cancer H1299 cells by impaired activation of MAPK.

CAS Number

410536-97-9

Chemical Formula

C₂₆H₂₂N₂O₂

Molecular Weight

394.5

Size

5 mg

Purity

97%

Solubility

Soluble in DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Sirtinol	5 mg	M45002 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.