

Sirtinol

SIRT inhibitor

Specific cell permeable inhibitor of the sirtuin family of NAD-dependent deacetylases (ySir2: IC_{50} =48 μ M; hSIRT1: IC_{50} =60 μ M; hSIRT2: IC_{50} =58 μ M) with no effect on human HDAC1. Reported to inhibit Sir2p transcriptional silencing activity *in vivo* (IC_{50} =25 μ M) and NAD-dependent histone deacetylase activity of purified recombinant yeast Sir2p (IC_{50} =70 μ M) and hSIRT2 (IC_{50} =40 μ M) *in vitro*.

Citations: 10

[View Online »](#)

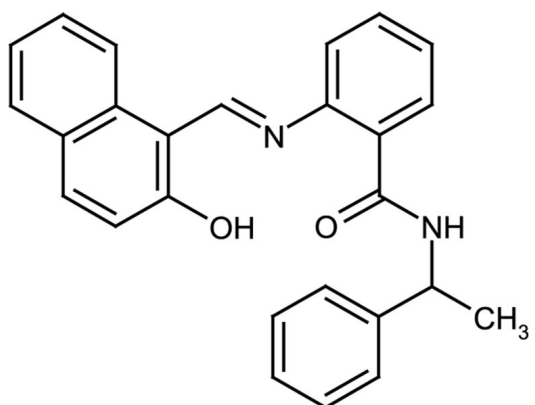
Ordering Information

[Order Online »](#)

ALX-270-308-M001	1mg
ALX-270-308-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Relatively unstable in aqueous solution, especially at acidic pH.
Handling	Protect from moisture. Protect from light. Store under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Sir Two Inhibitor Naphthol, 2-[(2-Hydroxynaphthalen-1-yl-methylene)amino]-N-(1-phenethyl)benzamide
Appearance	Off-white to yellow crystals.
CAS	410536-97-9
Couple Target	SIRT
Couple Type	Inhibitor
Formula	$C_{26}H_{22}N_2O_2$
Identity	NMR
MW	394.5
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol or dimethyl formamide.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)