

Okadaic acid . ammonium salt (high purity)

Inhibitor of PP1 and PP2A

Salt form of okadaic acid (Prod. No. ALX-350-003), with slightly greater stability than the free acid after it is put into stock solution (in organic solvents). Potent inhibitor of protein phosphatase 1 (PP1) and 2A (PP2A). Potent tumor promotor and neurotoxin.

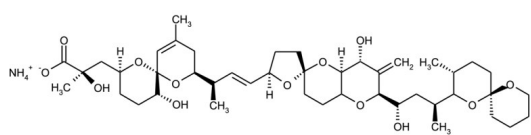
Ordering Information

[Order Online »](#)

ALX-350-010-M001	1mg
ALX-350-010-C100	100µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Halochondrine A . NH ₄ , 9,10-Deepithio-9,10-didehydroacanthifolicin . NH ₄
Appearance	White to off-white solid.
CAS	155716-06-6
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	C ₄₄ H ₆₇ O ₁₃ · NH ₄
MW	804.0 . 18.0
Purity	≥98% (TLC)
Solubility	Soluble in ethanol or methanol.
Source	Isolated from <i>Prorocentrum concavum</i> by preparative flash, medium pressure and high performance liquid chromatography. Salt form generated in aqueous ammonium hydroxide-methanol solution.

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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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