

Piericidin A

NADH-ubiquinone oxidoreductase inhibitor

Antibiotic. Specific and potent inhibitor of NADH-ubiquinone oxidoreductase (complex I) that binds to ubiquinone binding site(s). Inhibits both mitochondrial and bacterial NADH ubiquinone oxidoreductase. Anticancer agent that prevents up-regulation of GRP78, and exhibits cytotoxic activity for etoposide resistant cancer cells under glucose-deprived conditions.

Citations: 13

[View Online »](#)

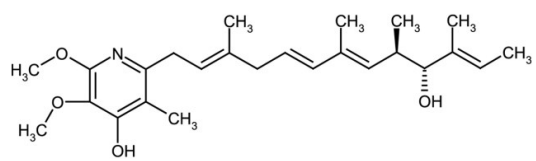
Ordering Information

[Order Online »](#)

ALX-380-235-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Piericidin A1, Shaoguanmycin B, Antibiotic MT 1882-I, Antibiotic SN 198E, Antibiotic IT 143D

Appearance Pale yellow oil.

CAS 2738-64-9

Couple Target NADH-ubiquinone oxidoreductase

Couple Type Inhibitor

Formula $C_{25}H_{37}NO_4$

MW 415.6

Purity ≥95% (HPLC)

RTECS YD4588000

Solubility Soluble in 100% ethanol, methanol, dimethyl formamide or DMSO. Poorly soluble in water.

Source Isolated from *Streptomyces* sp. MST-AS5364.

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