

Nodularin

Inhibitor of PP1 and PP2A

Nodularia spumigena is a cyanobacterium often found in drinking water in developing countries and capable of accumulating in fish and seafood. It produces nodularin, a genotoxic and hepatotoxic monocyclic pentapeptide. Nodularin inhibits protein phosphatase 1 (PP1) ($IC_{50}=1.8nM$), protein phosphatase 2A (PP2A) ($IC_{50}=0.026nM$), and to a lesser extent protein phosphatase 2B (PP2B) ($IC_{50}=8.7\mu M$). Nodularin is similar to microcystin-LR (Prod. No. ALX-350-012) but with increased water solubility. Nodularin is capable of inducing oxidative DNA damage by oxidation of purines as well as apoptosis in HepG2 cells.

May require a license for import, please [contact us](#) for more information.

Citations: 35

[View Online »](#)

Ordering Information

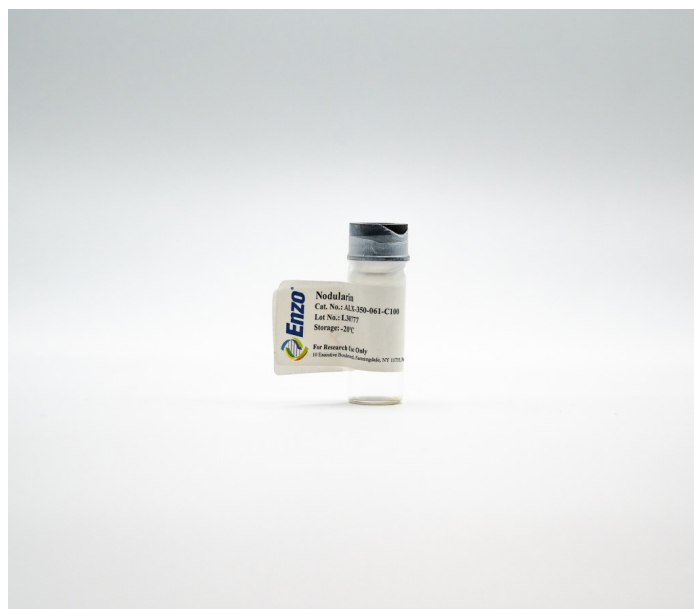
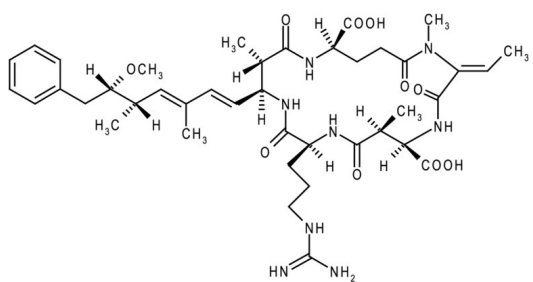
[Order Online »](#)

ALX-350-061-M001	1mg
ALX-350-061-C100	100µg
ALX-350-061-C250	250µg

Manuals, SDS & CofA

[View Online »](#)

- Hepatotoxin
- Potent inhibitor of PP1 and PP2A
- Useful for cytotoxicity and environmental studies



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

Appearance Dry residue containing traces of monobasic potassium phosphate.

CAS 118399-22-7

Couple Target Serine/threonine-protein phosphatase

Couple Type Inhibitor

Formula $C_{41}H_{60}N_8O_{10}$

Identity Identity determined by MS.

MW 825.0

Purity ≥95% (HPLC)

RTECS GU2294250

Solubility Soluble in methanol:water (1:1).

Source Isolated from *Nodularia spumigena*.



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)