

Anabaenopeptin B

Inhibitor of protein phosphatase.

Anabaenopeptins are a highly diverse group of bioactive peptides produced by several genera of cyanobacteria such as *Anabaena*, *Planktothrix*, *Microcystis* and *Nodularia*. These peptides are commonly detected in cyanobacterial blooms along with the well known microcystins. Unique cyclic peptides where lysine is conserved in position 2 and cyclisation is achieved by linkage of the lysine to the C-terminal amino acid. In addition the side chain of a single amino acid is attached via a ureido bond. Bioactive peptides originally isolated from *Anabaena flos-aquae* NRC 525-17 where they were shown to be non-toxic in mice but had dose dependant relaxations in rat aortic preparations with endothelium precontracted with 0.1µM norpinephrine. Anabaenopeptins inhibit protein phosphatase.

Citations: 5

[View Online »](#)

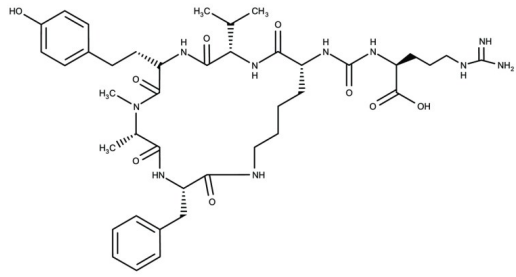
Ordering Information

[Order Online »](#)

ALX-350-184-C100	100µg
------------------	-------

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

Appearance Pale powder or an oil (film), depending on the lot.

CAS 161897-74-1

Couple Target Serine/threonine-protein phosphatase

Couple Type Inhibitor

Formula $C_{41}H_{60}N_{10}O_9$

MW 837.0

Purity ≥95% (HPLC)

RTECS GU2294250

Solubility Soluble in methanol.

Source Isolated from *Microcystis aeruginosa*.

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