

Anabaenopeptin A

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 »](#)

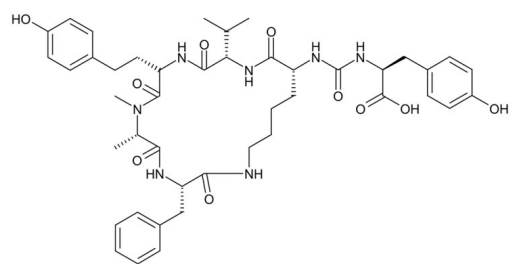
Ordering Information

[Order Online »](#)

ALX-350-183-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.
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{44}H_{57}N_7O_{10}$
Identity	Identity determined by MS.
MW	843.9
Purity	≥95% (HPLC)
RTECS	GU2294250
Solubility	Soluble in methanol.
Source	Isolated from <i>Microcystis aeruginosa</i> .

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)