

# 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 »](#)

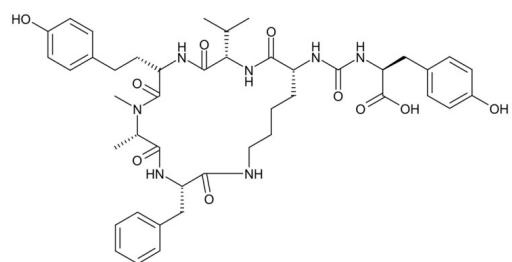
## Ordering Information

[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 *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-  
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)