

# Cylindrospermopsin

## Protein synthesis inhibitor

Cylindrospermopsin is a tricyclic alkaloid hepatotoxin produced by cyanobacteria. Cylindrospermopsin exhibits a completely different mechanism of toxicity than microcystins by targeting/inhibiting the synthesis of glutathione, proteins, and pyrimidine nucleotides. Cylindrospermopsin is considered to be genotoxic. It was shown to break double stranded DNA and reduce cell viability in HepG2 cells.

Cyanobacteria are photosynthetic prokaryotes mostly present in freshwater ecosystems. The increasingly frequent appearance of cyanobacteria blooms in lakes and rivers is linked to climate changes and human activities. Microcystins are a group of cyclic heptapeptide hepatotoxins produced by a number of cyanobacterial genera. The most notable of which, and namesake, is the widespread genus *Microcystis*. Structurally, all microcystins consist of the generalized structure cyclo(-D-Ala<sup>1</sup>-X<sup>2</sup>-D-MeAsp<sup>3</sup>-Y<sup>4</sup>-Adda<sup>5</sup>-D-Glu<sup>6</sup>-Mdha<sup>7</sup>-). X and Y are variable L-amino acids, D-MeAsp is D-erythro-β-methylaspartic acid and Mdha is N-methyldehydroalanine. Adda is the cyanobacteria unique C<sub>20</sub> β-amino acid 3-amino-9-methoxy-2,6,8-trimethyl-10-phenyl-deca-4,6-dienoic acid. Substitutions of the variable L-amino acids at positions 2 and 4 give rise to at least 21 known primary microcystin analogs and alterations in the other constituent amino acids result in more than 90 reported microcystins to date.

Citations: 38

[View Online »](#)

## Ordering Information

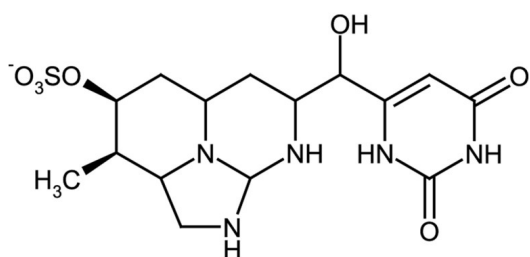
[Order Online »](#)

ALX-350-149-M001	1mg
ALX-350-149-C100	100µg
ALX-350-149-C500	500µg

## Manuals, SDS & CofA

[View Online »](#)

- Hepatotoxic
- Protein synthesis inhibitor
- Cited in several environment-related research articles



## 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** EN-Cylindrospermopsin

**Appearance** Glassy solid.

**CAS** 143545-90-8

**Formula**  $C_{15}H_{21}N_5O_7S$

**Identity** Identity determined by MS.

**MW** 415.4

**Purity** ≥95% (HPLC)

**RTECS** UV9104310

**Solubility** Soluble in DMSO, 100% methanol, or water.

**Source** Isolated from *Cylindrospermopsis raciborskii*.



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)