

[D-Asp³]microcystin-LR

[D-Asp³]microcystin-LR is a variant of microcystin-LR isolated from *Microcystis aeruginosa*. In this analog, the D-erythro- β -methylaspartic acid (D-MeAsp) residue at position 3 is replaced with aspartic acid (Asp), hence the designation [D-Asp³]. This microcystin contains leucine (L) at position 2 and arginine (R) at position 4. Like other microcystins, [D-Asp³]microcystin-LR is a potent hepatotoxin.

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

Cyanobacteria are photosynthetic prokaryotes commonly found in freshwater ecosystems. The increasingly frequent occurrence of cyanobacterial blooms in lakes and rivers is closely linked to climate change and human activities, particularly nutrient enrichment. Microcystins are a group of cyclic heptapeptide hepatotoxins produced by several cyanobacterial genera, most notably *Microcystis*. Structurally, all microcystins share a conserved cyclic heptapeptide core with the general structure: cyclo(-D-Ala¹-X²-D-MeAsp³-Y⁴-Adda⁵-D-Glu⁶-Mdha⁷-), where X and Y are variable L-amino acids. D-MeAsp refers to D-erythro- β -methylaspartic acid, and Mdha is N-methyldehydroalanine. Adda (3-amino-9-methoxy-2,6,8-trimethyl-10-phenyl-deca-4,6-dienoic acid) is a unique β -amino acid essential for the toxicity of microcystins. Substitutions at positions 2 and 4 define at least 21 primary microcystin analogs, such as microcystin-LR, while modifications at other positions - including demethylation, oxidation, or amino acid substitutions - have led to the identification of over 90 microcystin variants to date. [D-Asp³]microcystin-LR is one such variant, in which the D-MeAsp at position 3 is replaced by aspartic acid.

- Hepatotoxic

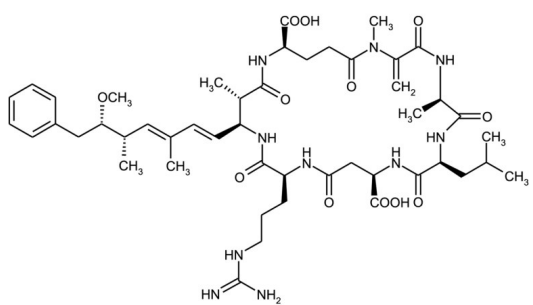
Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-350-173-C025	25 μ g
ALX-350-173-C100	100 μ g



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Microcystin-LR (desmethylated), MC-LR (desmethylated)
Appearance	White to off-white powder or oily film adhered to inside of the vial.
Formula	$C_{48}H_{72}N_{10}O_{12}$
Identity	Identity determined by MS.
MW	981.1
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO or 100% methanol.
Source	Isolated from <i>Microcystis aeruginosa</i> .



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)