

Microcystin-RR

(Analytical Standard)

PP2A inhibitor

Microcystin-RR is the Arg-Arg analog of microcystin-LR (Prod. No. ALX-350-431). Hepatotoxic, although found to be up to 10-fold less toxic than microcystin-LR on i.p. injection in mice. Potent inhibitor of protein phosphatase 2A (PP2A). This product can be used as an analytical standard for water testing or for molecular studies when high purity is required.

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

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-Ala¹-X²-D-MeAsp³-Y⁴-Adda⁵-D-Glu⁶-Mdha⁷-). 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 C₂₀ 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: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

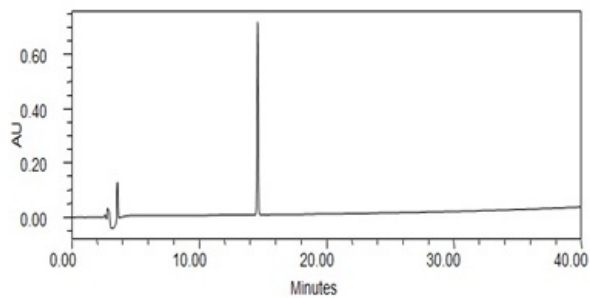
ALX-350-432-C010

10µg

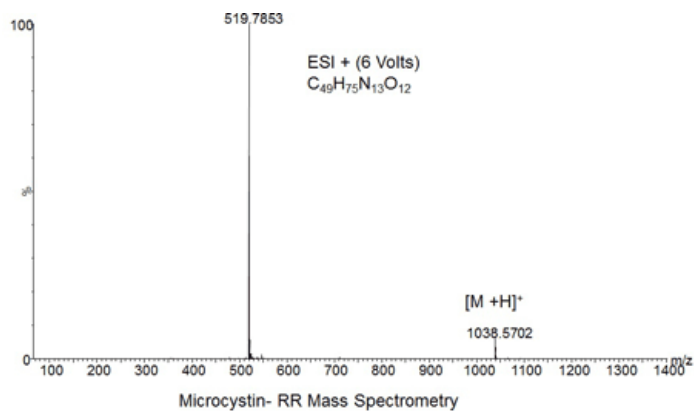
Manuals, SDS & CofA

[View Online »](#)

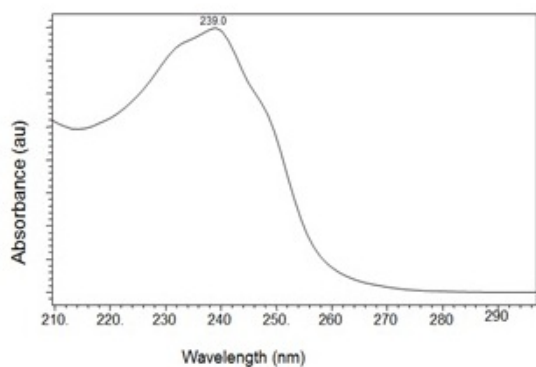
- Highest Purity On the Market!



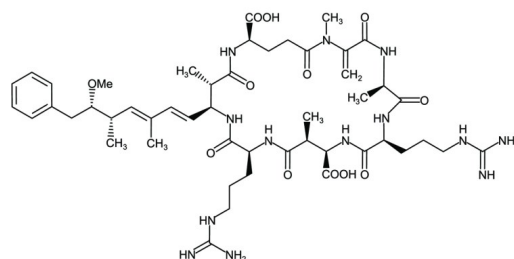
Microcystin- RR Reversed Phase HPLC (238nm)



Microcystin- RR Mass Spectrometry



Microcystin-RR (UV-VIS) at 239nm



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Preparation of stock solution: Thaw vial. Centrifuge vial for maximum product recovery. Use a syringe, draw the desired solvent (suggested solvent is 80% methanol). Make sure that no bubbles are drawn. Inject 1.0ml of desired solvent into the septum to obtain an concentration of 10µg/ml. Make sure that all the solvent has been injected prior to removing the syringe. Gently swirl the vial to get toxin into solution. If required, gently tap the bottom of the vial. Material is ready for analysis. Stock solutions are stable for up to 6 months when stored at -20°C. Unstable at pH >7.7.

Handling Warm to room temperature before opening. Keep sterile. For maximum product recovery after thawing, centrifuge the vial before opening the cap.

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Clear to whitish film adhered to inside of vial.

CAS 111755-37-4

Couple Target Serine/threonine-protein phosphatase

Couple Type Inhibitor

Formula $C_{49}H_{75}N_{13}O_{12}$

Identity Identity determined by MS.

MW 1038.2

Purity ≥99% (HPLC)

Purity Detail



Source

Flash Chromatography followed by HPLC

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Isolated from *Microcystis aeruginosa*.