Microcystin-RR

PP2A inhibitor

Microcystin-RR (MC-RR) is an Arg-Arg analog of Microcystin-LR (Prod. No. ALX-350-012). Although MC-RR was found to be up to 10-fold less toxic than MC-LR after intra-peritoneal injection in mice, it is both hepatotoxic and carcinogenic. It is a potent inhibitor of protein phosphatase 2A (PP2A).

May require a license for import, please contact us for more information.

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 $\it Microcystis$. Structurally, all microcystins consist of the generalized structure $\rm cyclo(-D-Ala^1-X^2-D-MeAsp^3-Y^4-Adda^5-D-Glu^6-Mdha^7-)$. 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 $_{20}$ 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 mycrocystins to date.

Citations: 42

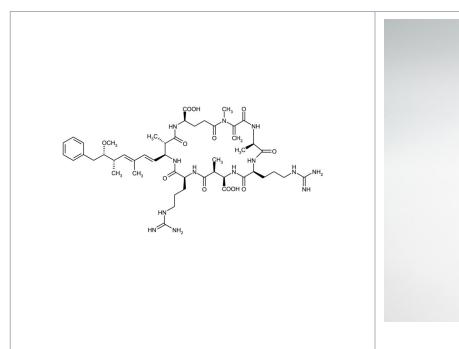
View Online »

Ordering Information

Order Online »

ALX-350-043-M001	1mg
ALX-350-043-C050	50µg
ALX-350-043-C100	100μg
ALX-350-043-C250	250μg
ALX-350-043-C500	500µg

- Hepatotoxin
- Potent inhibitor of PP2A
- Cited in several water qualityrelated research articles





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 MC-RR

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 ≥95% (HPLC)

Solubility Soluble in 80% aqueous methanol.

Source Isolated from *Microcystis aeruginosa*.

Last modified: May 29, 2024

