

PRODUCT INFORMATION



Microcystin-WR

Item No. 43504

CAS Registry No.: 138234-58-9

Formal Name: cyclo[2,3-didehydro-N-methylalanyl-D-alanyl-L-tryptophyl-(3S)-3-methyl-D-β-aspartyl-L-arginyl-(2S,3S,4E,6E,8S,9S)-3-amino-9-methoxy-2,6,8-trimethyl-10-phenyl-4,6-decadienoyl-D-γ-glutamyl]

Synonyms: Cyanoginosin-WR, MC-WR

MF: C₅₄H₇₃N₁₁O₁₂

FW: 1,068.2

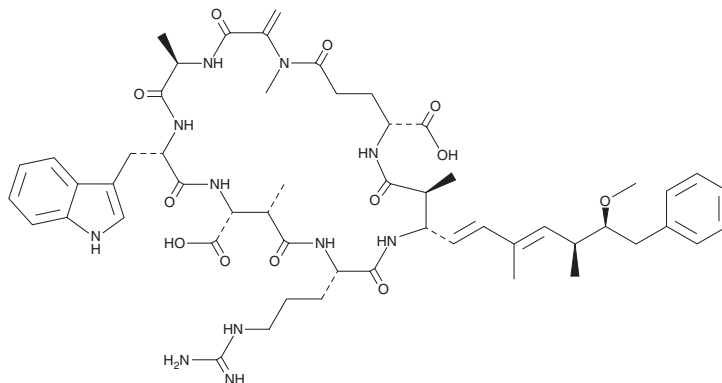
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Microcystis aeruginosa*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Microcystin-WR is supplied as a solid. A stock solution may be made by dissolving the microcystin-WR in the solvent of choice, which should be purged with an inert gas. Microcystin-WR is soluble in methanol:water (1:1).

Description

Microcystin-WR is a cyanotoxin that has been found in *Microcystis* blooms.¹ It is toxic to mice (LD₅₀ = 150-200 µg/kg).

Reference

1. Namikoshi, M., Rinehart, K.L., Sakai, R., *et al.* Identification of 12 hepatotoxins from a Homer Lake bloom of the cyanobacteria *Microcystis aeruginosa*, *Microcystis viridis*, and *Microcystis wesenbergii*: Nine new microcystins. *J. Org. Chem.* **57**(3), 866-872 (1992).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM