

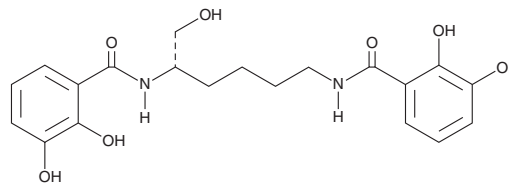
PRODUCT INFORMATION



Myxochelin A

Item No. 26422

CAS Registry No.: 120243-02-9
Formal Name: N,N'-[(1S)-1-(hydroxymethyl)-1,5-pentanediy]bis[2,3-dihydroxy-benzamide]
Synonym: (S)-Myxochelin A
MF: C₂₀H₂₄N₂O₇
FW: 404.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Myxochelin A is supplied as a solid. A stock solution may be made by dissolving the myxochelin A in the solvent of choice, which should be purged with an inert gas. Myxochelin A is soluble in organic solvents such as DMSO, ethanol, and methanol.

Description

Myxochelin A is a microbial metabolite that has been found in *A. disciformis* and has diverse biological activities.¹ It is active against Gram-positive bacteria, including *B. cereus*, *S. aureus*, and *M. luteus*, but not Gram-negative bacteria or fungi in an agar diffusion assay when used at a concentration of 80 µg/disc. Myxochelin A inhibits 5-lipoxygenase (5-LO) activity with an IC₅₀ value of 1.9 µM for the recombinant human enzyme.² It is cytotoxic to 26-L5 colon cancer cells when used at a concentration of 3 µg/ml.³

References

1. Kunze, B., Bedorf, N., Kohl, W., *et al.* Myxochelin A, a new iron-chelating compound from *Angiococcus disciformis* (Myxobacterales). Production, isolation, physico-chemical and biological properties. *J. Antibiot. (Tokyo)* **42**(1), 14-17 (1989).
2. Schieferdecker, S., König, S., Koeberle, A., *et al.* Myxochelins target human 5-lipoxygenase. *J. Nat. Prod.* **78**(2), 335-338 (2015).
3. Miyanaga, S., Obata, T., Onaka, H., *et al.* Absolute configuration and antitumor activity of myxochelin A produced by *Nonomuraea pusilla* TP-A0861. *J. Antibiot. (Tokyo)* **59**(11), 698-703 (2006).

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.

Copyright Cayman Chemical Company, 10/17/2022

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