

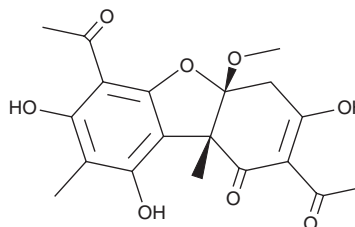
PRODUCT INFORMATION



(-)-Mycousnine

Item No. 31784

CAS Registry No.: 77480-55-8
Formal Name: (4aR,9bS)-2,6-diacetyl-4a,9b-dihydro-3,7,9-trihydroxy-4a-methoxy-8,9b-dimethyl-1(4H)-dibenzofuranone
Synonym: Mycousnin
MF: C₁₉H₂₀O₈
FW: 376.4
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 285 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

(-)-Mycousnine is supplied as a solid. A stock solution may be made by dissolving the (-)-mycousnine in the solvent of choice, which should be purged with an inert gas. (-)-Mycousnine is soluble in the organic solvent DMSO.

Description

(-)-Mycousnine is a microbial metabolite and derivative of usnic acid (Item No. 22585) originally isolated from *M. nawae* that has antibacterial and antifungal activities.^{1,2} It is active against the Gram-positive bacteria *B. subtilis*, *K. rhizophila*, and *S. aureus* (MICs = 4, 8, and 4 g/ml, respectively) but not the Gram-negative bacteria *E. coli*, *S. typhimurium*, and *K. pneumoniae* (MICs = >128 g/ml for all).² (-)-Mycousnine is also active against the fungi *T. mentagrophytes*, *T. rubrum*, and *C. albicans* (MICs = 25, 25, and 100 µg/ml, respectively).¹

References

1. Sassa, T. and Igarashi, M. Structures of (-)-mycousnine, (+)-isomycousnine and (+)-oxymycousnine, new usnic acid derivatives from phytopathogenic *Mycosphaerella nawae*. *Agric. Biol. Chem.* **54(9)**, 2231-2237 (1990).
2. Lee, J., Lee, J., Kim, G.J., et al. Mycousfurans A and B, antibacterial usnic acid congeners from the fungus *Mycosphaerella* sp., isolated from a marine sediment. *Mar. Drugs* **17(7)**, 422 (2019).

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, 08/17/2023

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