

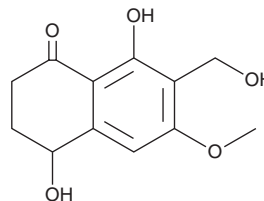
PRODUCT INFORMATION



10-Norparvulenone

Item No. 29452

CAS Registry No.: 618104-32-8
Formal Name: 3,4-dihydro-4,8-dihydroxy-7-(hydroxymethyl)-6-methoxy-1(2H)-naphthalenone
Synonym: (±)-10-Norparvulenone
MF: C₁₂H₁₄O₅
FW: 238.2
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Paraphaeosphaeria* sp.



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

Laboratory Procedures

10-Norparvulenone is supplied as a solid. A stock solution may be made by dissolving the 10-norparvulenone in the solvent of choice, which should be purged with an inert gas. 10-Norparvulenone is soluble in ethanol, methanol, DMSO, and dichloromethane.

Description

10-Norparvulenone is a fungal metabolite originally isolated from *Microsphaeropsis*.¹ It decreases viral sialidase activity in, and increases survival of, MDCK cells infected with the mouse-adapted influenza virus A/PR/8/34 when used at a concentration of 1 µg/ml.

Reference

1. Fukami, A., Nakamura, T., Kim, Y.P., *et al.* A new anti-influenza virus antibiotic, 10-norparvulenone from *Microsphaeropsis* sp. FO-5050. *J. Antibiot. (Tokyo)* **53(10)**, 1215-1218 (2000).

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, 08/16/2023

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