

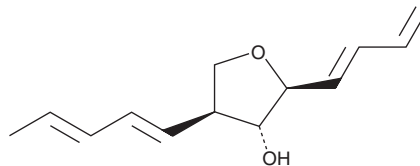
PRODUCT INFORMATION



Aureonitol

Item No. 32925

CAS Registry No.: 71774-51-1
Formal Name: (2S,3R,4S)-2-(1,3-butadien-1-yl)tetrahydro-4-(1,3-pentadien-1-yl)-3-furanol
MF: C₁₃H₁₈O₂
FW: 206.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Chaetomium* sp.



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

Laboratory Procedures

Aureonitol is supplied as a solid. A stock solution may be made by dissolving the aureonitol in the solvent of choice, which should be purged with an inert gas. Aureonitol is soluble in methanol and DMSO.

Description

Aureonitol is a fungal metabolite that has been found in *Chaetomium* and has antiviral activity.¹ It inhibits replication of the influenza A subtype H3N2 with an EC₅₀ value of 30 nM at a multiplicity of infection (MOI) of 0.01. It also inhibits replication of laboratory-adapted and clinical isolates of H3N2 (EC₅₀s = 100 and 312 nM, respectively), as well as clinical isolates of the influenza A subtype H1N1 (EC₅₀ = 417 nM). Aureonitol inhibits hemagglutination activity of H3N2 and H1N1.

Reference

1. Sacramento, C.Q., Marttorelli, A., Fintelman-Rodrigues, N., *et al.* Aureonitol, a fungi-derived tetrahydrofuran, inhibits influenza replication by targeting its surface glycoprotein hemagglutinin. *PLoS One* **10(10)**, e0139236 (2015).

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