

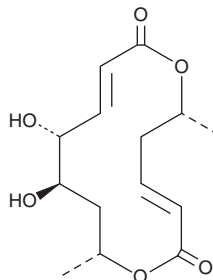
PRODUCT INFORMATION



Colletodiol

Item No. 29945

CAS Registry No.: 21142-67-6
Formal Name: (3E,6R,9E,11R,12R,14R)-11,12-dihydroxy-6,14-dimethyl-1,7-dioxacyclotetradeca-3,9-diene-2,8-dione
Synonym: (+)-Colletodiol
MF: C₁₄H₂₀O₆
FW: 284.3
Purity: ≥70%
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

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

Description

Colletodiol is a fungal metabolite that has been found in *D. grovesii* and has immunosuppressant and antiviral activities.^{1,2} It inhibits concanavalin A- or LPS-induced proliferation of isolated mouse splenocytes (IC₅₀s = 12 and 5 µg/ml, respectively).¹ Colletodiol inhibits influenza A viral replication in HeLa-IAV-Luc cells.²

References

1. Fujimoto, H., Nagano, J., Yamaguchi, K., et al. Immunosuppressive components from an ascomycete, *Diplogelasinospora grovesii*. *Chem. Pharm. Bull.* **46(3)**, 423-429 (1998).
2. Lai, W., Wang, S., and Ye, X. Colletodiol inhibits the replication of influenza A virus WSN/H1N1 by reducing the activity of viral RNA polymerase. *Wei Sheng Wu Xue Bao* **53(12)**, 1334-1339 (2013).

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