

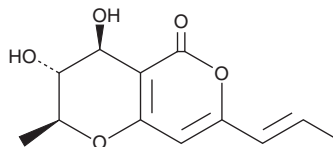
PRODUCT INFORMATION



Radicinol

Item No. 29861

CAS Registry No.: 65647-66-7
Formal Name: (2S,3R,4S)-3,4-dihydro-3,4-dihydroxy-2-methyl-7-(1E)-1-propen-1-yl-2H,5H-pyrano[4,3-b]pyran-5-one
MF: C₁₂H₁₄O₅
FW: 238.2
Purity: ≥95% (may contain up to 5% related metabolites)
UV/Vis.: λ_{max}: 225, 317 nm
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Curvularia lunata*



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

Laboratory Procedures

Radicinol is supplied as an oil. A stock solution may be made by dissolving the radicinol in the solvent of choice, which should be purged with an inert gas. Radicinol is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Radicinol is a phytotoxic fungal metabolite that has been found in *A. chrysanthemi*.¹

Reference

1. Robeson, D.J., Gray, G.R., and Strobel, G.A. Production of the phytotoxins radicinin and radicinol by *Alternaria chrysanthemi*. *Phytochem.* **21(9)**, 2359-2362 (1982).

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