

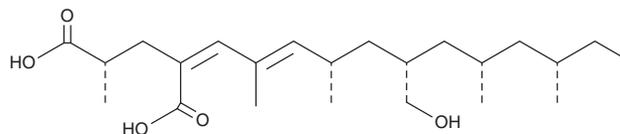
PRODUCT INFORMATION



Radiclonic Acid

Item No. 44992

CAS Registry No.: 49620-14-6
Formal Name: (2Z,4R)-rel-2-[(2E,4R,6R,8R,10R)-6-(hydroxymethyl)-2,4,8,10-tetramethyl-2-dodecen-1-ylidene]-4-methyl-pentanedioic acid
MF: C₂₃H₄₀O₅
FW: 396.6
Purity: ≥95%
Supplied as: An oil
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Radiclonic acid is supplied as an oil. A stock solution may be made by dissolving the radiclonic acid in the solvent of choice, which should be purged with an inert gas. Radiclonic acid is soluble in methanol and DMSO.

Description

Radiclonic acid is a diterpene that has been found in *Talaromyces* species and has diverse biological activities.¹⁻³ It is active against methicillin-resistant *S. aureus* (MRSA; MIC = 3.13 µg/ml).¹ Radiclonic acid is cytotoxic to BIU-87 bladder, Eca-109 esophageal, and BEL-7402 liver cancer cells (IC₅₀s = 12.47, 7.7, and 13.75 µM, respectively).² It stimulates root growth in Chinese cabbage and rice seedlings.³

References

1. Chu, M.-J., Wu, Q.-Q., Liu, B.-T., *et al.* Marine natural products with anti-MRSA activities: Promising candidates for antibacterial drug discovery. *Bioorg. Chem.* **167**, 109242 (2025).
2. Niu, S., Xia, M., Chen, M., *et al.* Cytotoxic polyketides isolated from the deep-sea-derived fungus *Penicillium chrysogenum* MCCC 3A00292. *Mar. Drugs* **17(12)**, 686 (2019).
3. Kinnersley, A.M., Scott, T.C.I., Yopp, J.H., *et al.* Promotion of plant growth by polymers of lactic acid. *Plant Growth Regul.* **9**, 137-146 (1990).

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, 01/29/2026

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