

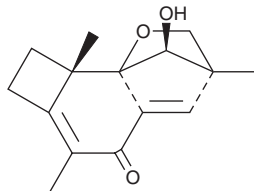
PRODUCT INFORMATION



Radulone A

Item No. 40939

CAS Registry No.: 221374-79-4
Formal Name: (3S,8aS,8bS,9S)-2,3,8a-tetrahydro-9-hydroxy-3,6,8a-trimethyl-3,8b-methano-8bH-cyclobuta[5,6]benzo[1,2-b]pyran-5(7H)-one
MF: $C_{15}H_{18}O_3$
FW: 246.3
Purity: $\geq 70\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Microbial



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

Laboratory Procedures

Radulone A is supplied as a solid. A stock solution may be made by dissolving the radulone A in the solvent of choice, which should be purged with an inert gas. Radulone A is soluble (≥ 10 mg/ml) in ethanol, methanol DMSO, and dichloromethane.

Description

Radulone A is a sesquiterpene fungal metabolite originally isolated from *R. confluens* that has diverse biological activities.¹ It is active against eight bacterial strains (MICs = 41-406 μM). Radulone A inhibits ADP-, collagen-, arachidonic acid-, or phorbol 12-myristate 13-acetate-induced aggregation of isolated human platelets (IC_{50} = 2, 28, 53, and 38 μM , respectively). It induces cytotoxicity in L1210 skin lymphocytic leukemia, HL-60 promyeloblast leukemia, and HeLa cervical cancer cells and COS-7 kidney fibroblasts (IC_{50} s = 1, 2, 20, and 16 μM).

Reference

1. Fabian, K., Lorenzen, K., Anke, T., *et al.* Five new bioactive sesquiterpenes from the fungus *Radulomyces confluens* (Fr.) Christ. *Z. Naturforsch. C. J. Biosci.* **53(11-12)**, 939-945 (1998).

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, 04/09/2025

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