

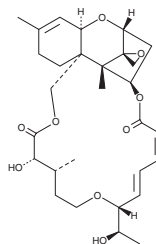
PRODUCT INFORMATION



Roridin A

Item No. 43401

CAS Registry No.: 14729-29-4
Formal Name: 7'R-deoxo-7'-[(1R)-1-hydroxyethyl]-
verrucarin A
MF: $C_{29}H_{40}O_9$
FW: 532.6
Purity: $\geq 70\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Roridin A is a mycotoxin that has been found in *M. verrucaria*.¹⁻⁴ It reduces the growth of *P. falciparum* ($EC_{50} = 0.31$ ng/ml).¹ Roridin A (2.3, 3.9, or 4.1 mg/kg) is toxic to rabbits.⁴ It has been found in agricultural soil samples.

References

1. Isaka, M., Punya, J., Lertwerawat, Y., et al. Antimalarial activity of macrocyclic trichothecenes isolated from the fungus *Myrothecium verrucaria*. *J. Nat. Prod.* **62**(2), 329-331 (1999).
2. Nguyen, M.V., Han, J.W., Kim, H., et al. Curvicollide D, a new modified γ -lactone from the culture broth of *Albifimbria verrucaria* and its antifungal activity against plant pathogenic fungi. *J. Antibiot. (Tokyo)* **75**(9), 514-518 (2022).
3. Nguyen, L.T.T., Jang, J.Y., Kim, T.Y., et al. Nematicidal activity of verrucarin A and roridin A isolated from *Myrothecium verrucaria* against *Meloidogyne incognita*. *Pestic. Biochem. Physiol.* **148**, 133-143 (2018).
4. Habermehl, G.G., Busam, L., Heydel, P., et al. Macrocyclic trichothecenes: Cause of livestock poisoning by the Brazilian plant *Baccharis coridifolia*. *Toxicon.* **23**(5), 731-745 (1985).

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, 05/01/2025

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