

PRODUCT INFORMATION



Rosellichalasin

Item No. 44752

CAS Registry No.: 123452-64-2

Formal Name: (1E,10aS)-13S,13aS,14S,14aR,15aS,15bS-hexahydro-4S,6E,14,14a-tetramethyl-13-(phenylmethyl)-3H-oxacyclododecino[2,3-d]oxireno[f]isoindole-5,9,11(4H,8H,12H)-trione

MF: $C_{28}H_{33}NO_5$

FW: 463.6

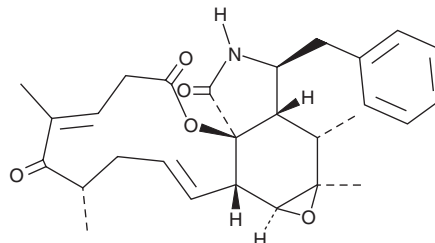
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Rosellichalasin is a fungal metabolite that has been found in *R. necatrix* and has plant growth inhibitory activity.¹ It inhibits lettuce seed germination by 50%, as well as root and hypocotyl growth by 70 and 60%, respectively, when used at a concentration of 100 mg/L.

Reference

1. Kimura, Y., Nakajima, H., and Hamasaki, T. Structure of rosellichalasin, a new metabolite produced by *Rosellinia necatrix*. *Agric. Biol. Chem.* **53**(6), 1699–1701 (1989).

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, 11/18/2025

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