

PRODUCT INFORMATION



(-)-Rugulosin

Item No. 25477

CAS Registry No.: 21884-45-7

Formal Name: [5aR-(5aR*,6S*,13aR*,14S*,17R*,18S*,19S*,20R*)]-1,7,9,15,17,20-hexahydroxy-3,11-dimethyl-5H,6H-6,13a,5a,14-[1,2,3,4]butanetetraylcycloocta[1,2-b:5,6-b']dinaphthalene-5,8,13,16(14H)-tetrone

MF: C₃₀H₂₂O₁₀

FW: 542.5

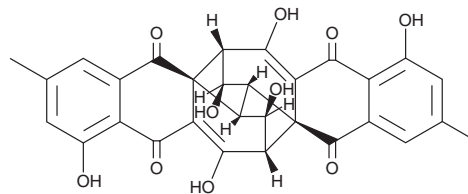
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

(-)-Rugulosin is supplied as a solid. A stock solution may be made by dissolving the (-)-rugulosin in the solvent of choice, which should be purged with an inert gas. (-)-Rugulosin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

(-)-Rugulosin is a fungal metabolite that has been isolated from *Penicillium* species, including *P. islandicum*, which also produces the optical isomer and mycotoxin (+)-rugulosin (Item No. 23558).¹ (-)-Rugulosin is 3-10-fold more cytotoxic to HeLa and L cells compared with (+)-rugulosin.² It is protective against influenza in mice when administered via inhalation but not when administered intraperitoneally with 38.3 and 5% survival rates, respectively.³

References

1. Singh, S.B., Jayasuriya, H., Dewey, R., *et al.* Isolation, structure, and HIV-1-integrase inhibitory activity of structurally diverse fungal metabolites. *J. Ind. Microbiol. Biotechnol.* **30**(12), 721-731 (2003).
2. Umeda, M., Saito, M., and Shibata, S. Comparison of cytotoxic effects of various anthraquinonoids on cultured mammalian cells. *Jpn. J. Exp. Med.* **44**(3), 249-255 (1974).
3. Nakamura, S., Nii, F., Inoue, S., *et al.* Antiinfluenzal effect of rugulosin. *Jpn. J. Microbiol.* **18**(1), 1-7 (1974).

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, 08/23/2023

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