

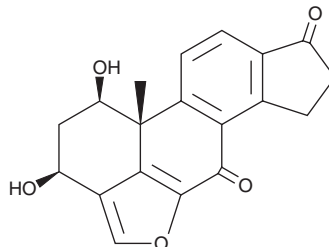
PRODUCT INFORMATION



Demethoxyviridiol

Item No. 27953

CAS Registry No.: 56617-66-4
Formal Name: (1R,3S,11bR)-1,2,3,7,8,11b-hexahydro-1,3-dihydroxy-11b-methyl-cyclopenta[7,8]phenanthro[10,1-bc]furan-6,9-dione
Synonym: Desmethoxyviridiol
MF: C₁₉H₁₆O₅
FW: 324.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Avoid long-term storage in methanol

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

Laboratory Procedures

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

Description

Demethoxyviridiol is a mycotoxin originally isolated from *N. hinnuleum*.¹ It induces lethality in day-old cockerels (LD₅₀ = 4.2 mg/kg). Demethoxyviridiol is also an inhibitor of phosphatidylinositol 3-kinase (PI3K).²

References

1. Cole, R.J., Kirksey, J.W., Springer, J.P., *et al.* Desmethoxyviridiol, a new toxin from *Nodulisporium hinnuleum*. *Phytochemistry* **14**(5-6), 1429-1432 (1975).
2. Dodge, J.A., Sato, M., and Vlahos, C.J. Inhibition of phosphatidylinositol 3-kinase with viridin, demethoxyviridin, viridiol, demethoxyviridiol, virone, wortmannolone, and analogs thereof. *Eli Lilly and Co.* **5,726,167** (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, 12/02/2022

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