

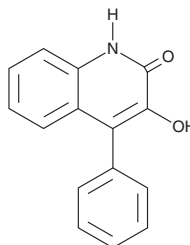
PRODUCT INFORMATION



Viridicatin

Item No. 21763

CAS Registry No.: 129-24-8
Formal Name: 3-hydroxy-4-phenyl-2(1H)-quinolinone
Synonym: NSC 656625
MF: $C_{15}H_{11}NO_2$
FW: 237.3
Purity: $\geq 99\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
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

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

Description

Viridicatin is a fungal metabolite isolated from various *Penicillium* species.^{1,2} This agent exhibits *in vitro* antibiotic activity against *M. tuberculosis* (1:15,000 dilution) but no activity against other tested bacterial strains.²

References

1. Santini, A., Mikusová, P., Sulyok, M., *et al.* *Penicillium* strains isolated from Slovak grape berries taxonomy assessment by secondary metabolite profile. *Mycotoxin Res.* **30**(4), 213-220 (2014).
2. Cunningham, K.G. and Freeman, G.G. The isolation and some chemical properties of viridicatin, a metabolic product of *Penicillium viridicatum* westling. *Biochem J.* **53**(2), 328-332 (1953).

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, 08/08/2023

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