

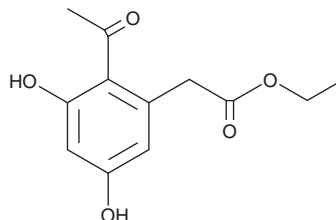
PRODUCT INFORMATION



Curvulin

Item No. 19609

CAS Registry No.: 19054-27-4
Formal Name: 2-acetyl-3,5-dihydroxybenzeneacetic acid, ethyl ester
MF: C₁₂H₁₄O₅
FW: 238.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Protect from light
Item Origin: Fungus/*Curvularia* sp.



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

Laboratory Procedures

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

Description

Curvulin is a phytotoxin first isolated from several species of the mold *Curvularia*. It is reported to inhibit microtubule assembly and has also been shown to inhibit iNOS expression.^{1,2}

References

1. Tadross, P.M., Virgil, S.C., and Stoltz, B.M. Aryne acyl-alkylation in the general and convergent synthesis of benzannulated macrolactone natural products: An enantioselective synthesis of (-)-curvularin. *Org. Lett.* **12(7)**, 1612-1614 (2010).
2. Yao, Y., Hausding, M., Erkel, G., *et al.* Sporogen, S14-95, and S-curvularin, three inhibitors of human inducible nitric-oxide synthase expression isolated from fungi. *Mol. Pharmacol.* **63(2)**, 383-391 (2003).

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, 08/09/2023

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