

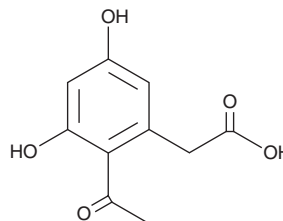
# PRODUCT INFORMATION



## Curvulinic Acid

Item No. 33094

**CAS Registry No.:** 19053-94-2  
**Formal Name:** 2-acetyl-3,5-dihydroxy-benzeneacetic acid  
**MF:** C<sub>10</sub>H<sub>10</sub>O<sub>5</sub>  
**FW:** 210.2  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Curvularia* sp.



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

### Laboratory Procedures

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

### Description

Curvulinic acid is a fungal metabolite originally isolated from *C. siddiqui*.<sup>1</sup> It inhibits *C. bursa-pastoris* seed germination, root growth, and shoot growth by 73.3, 73.5, and 66.7%, respectively, when used at a concentration of 600 µg/ml.<sup>2</sup>

### References

1. Kamal, A., Ahmad, N., Khan, M.A., *et al.* Studies in the biochemistry of microorganisms—I: Curvulin and curvulinic acid, metabolic products of *Curvularia siddiqui*. *Tetrahedron* **18(4)**, 433-436 (1962).
2. Li, E., Yea, Y., Wang, X., *et al.* Herbicidal activity of curvulinic acid isolated from *Nimbya alternantherae*. *Nat. Prod. Commun.* **7(1)**, 51-52 (2012).

#### 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, 08/23/20223

#### 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