

# PRODUCT INFORMATION



## Alternaric Acid

Item No. 45460

**CAS Registry No.:** 10088-62-7  
**Formal Name:** 4,5-dideoxy-2-C-[(1E)-7-[(6R)-5,6-dihydro-4-hydroxy-6-methyl-2-oxo-2H-pyran-3-yl]-4-methylene-7-oxo-1-hepten-1-yl]-4-ethyl-d-arabinonic acid

**MF:** C<sub>21</sub>H<sub>30</sub>O<sub>8</sub>

**FW:** 410.5

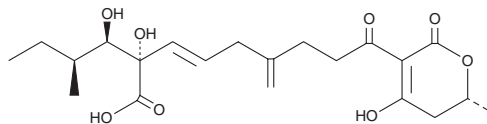
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/*Alternaria* sp.



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

### Laboratory Procedures

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

### Description

Alternaric acid is a phytotoxin that has been found in *A. solani*.<sup>1,2</sup> It inhibits germination and root development of carrot, mustard, radish, and tomato seeds when used at a concentration of 5 µg/ml and induces necrotic lesions in potato shoots.<sup>2</sup> Alternaric acid (0.005-100 µg/ml) is also active against several plant pathogenic fungi.<sup>3</sup>

### References

1. Tabuchi, H., Hamamoto, T., Miki, S., *et al.* Total synthesis and stereochemistry of alternaric acid. *J. Org. Chem.* **59(17)**, 4749-4759 (1994).
2. Brian, P.W., Elson, G.W., Hemming, H.G., *et al.* The phytotoxic properties of alternaric acid in relation to the etiology of plant diseases caused by *Alternaria solani* (Ell. & Mart.) Jones & Grout. *Ann. Appl. Biol.* **39(3)**, 308-321 (1952).
3. Brian, P.W., Curtis, P.J., Hemming, H.G., *et al.* Alternaric acid; a biologically active metabolic product of *Alternaria solani* (Ell. & Mart.) Jones & Grout; its production, isolation and antifungal properties. *J. Gen. Microbiol.* **5(4)**, 619-632 (1951).

#### 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, 05/21/2026

#### 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