

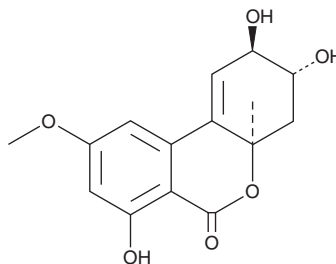
PRODUCT INFORMATION



(-)-Altenuene

Item No. 21130

CAS Registry No.: 889101-41-1
Formal Name: 2R,3R,4aR,4a-tetrahydro-2,3,7-trihydroxy-9-methoxy-4a-methyl-6H-dibenzo[b,d]pyran-6-one
MF: C₁₅H₁₆O₆
FW: 292.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(-)-Altenuene is supplied as a solid. A stock solution may be made by dissolving the (-)-altenuene in the solvent of choice, which should be purged with an inert gas. (-)-Altenuene is soluble in the organic solvent acetone at a concentration of approximately 10 mg/ml.

Description

(-)-Altenuene is a natural mycotoxin that was first isolated from *A. tenuis*. It has antioxidant activity and inhibits cholinesterases.¹ Unlike tenuazonic acid (Item No. 11443), (-)-altenuene does not display toxicity as assessed by the chicken embryo assay.² (-)-Altenuene does not display significant topoisomerase inhibitory activity, unlike alternariol (Item No. 11306).³

References

1. Bhagat, J., Kaur, A., Kaur, R., *et al.* Cholinesterase inhibitor (Altenuene) from an endophytic fungus *Alternaria alternata*: Optimization, purification and characterization. *J. Appl. Microbiol.* **121**(4), 1015-1025 (2016).
2. Griffin, G.F. and Chu, F.S. Toxicity of the *Alternaria* metabolites alternariol, alternariol methyl ether, altenuene, and tenuazonic acid in the chicken embryo assay. *Appl. Environ. Microbiol.* **46**(6), 1420-1422 (1983).
3. Fehr, M., Pahlke, G., Fritz, J.E., *et al.* Alternariol acts as a topoisomerase poison, preferentially affecting the IIa isoform. *Mol. Nutr. Food Res.* **53**(4), 441-451 (2008).

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, 07/01/2025

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