

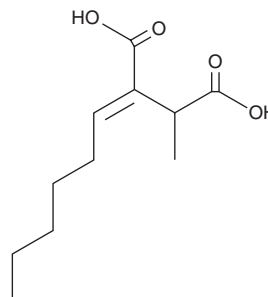
PRODUCT INFORMATION



Piliformic Acid

Item No. 31756

CAS Registry No.: 98985-76-3
Formal Name: 2E-hexylidene-3-methyl-butanedioic acid
MF: C₁₁H₁₈O₄
FW: 214.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Piliformic acid is a fungal metabolite that has been found in *N. pseudotrichia* and has diverse biological activities.^{1,2} It is cytotoxic to BC-1 human breast cancer cells (IC₅₀ = 5 µg/ml).² Piliformic acid is active against *L. braziliensis* amastigotes (IC₅₀ = 78.5 µM). It is also active against the plant pathogenic fungi *C. gloeosporioides* (MIC = 292 µM).¹

References

1. Elias, L.M., Fortkamp, D., Sartori, S.B., *et al.* The potential of compounds isolated from *Xylaria* spp. as antifungal agents against anthracnose. *Braz. J. Microbiol.* **49(4)**, 840-847 (2018).
2. Cota, B.B., Tunes, L.G., Maia, D.N.B., *et al.* Leishmanicidal compounds of *Nectria pseudotrichia*, an endophytic fungus isolated from the plant *Caesalpinia echinata* (Brazilwood). *Mem. Inst. Oswaldo Cruz* **113(2)**, 102-110 (2018).

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, 08/11/2023

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