

PRODUCT INFORMATION



Aspterric Acid

Item No. 25919

CAS Registry No.: 67309-95-9
Formal Name: (3R,4R,6aS,9aS)-octahydro-4-hydroxy-7-(1-methylethylidene)-1H-3,9a-methanocyclopent[c]oxocin-4-carboxylic acid

MF: C₁₅H₂₂O₄

FW: 266.3

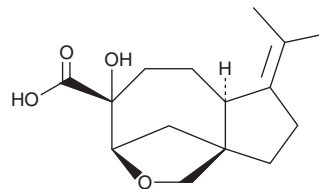
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

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

Description

Aspterric acid is a carotene-type sesquiterpene, plant growth regulator, and inhibitor of pollen development originally isolated from the fungus *A. terreus*.^{1,2} It inhibits pollen development, reduces stem length at first flowering, and increases the time to bolting and first flowering in *A. thaliana* when used at a concentration of 38 μM.¹ Aspterric acid exerts its effects on reproductive growth in *A. thaliana* without inhibiting biosynthesis and transport of the plant growth regulator indole-3-acetic acid (Item No. 16954).²

References

1. Shimada, A., Kusano, M., Takeuchi, S., *et al.* Aspterric acid and 6-hydroxymellein, inhibitors of pollen development in *Arabidopsis thaliana*, produced by *Aspergillus terreus*. *Z. Naturforsch. C.* **57(5-6)**, 459-464 (2002).
2. Shimada, A., Yamane, H., and Kimura, Y. The role of aspterric acid in auxin-regulated reproductive growth of *Arabidopsis thaliana*. *Z. Naturforsch. C.* **63(7-8)**, 554-556 (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, 08/23/2022

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