

PRODUCT INFORMATION



Aspinolide B

Item No. 40873

CAS Registry No.: 188605-15-4
Formal Name: (2E)-2-butenoic acid, (2R,3S,4E,6R,7S)-3,6,7,8,9,10-hexahydro-6,7-dihydroxy-2-methyl-10-oxo-2H-oxecin-3-yl ester

MF: C₁₄H₂₀O₆

FW: 284.3

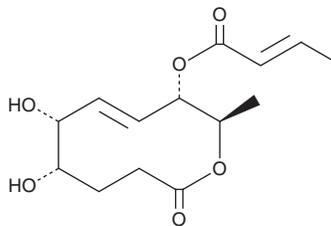
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Nigrospora lacticola*



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

Laboratory Procedures

Aspinolide B is supplied as a solid. A stock solution may be made by dissolving the aspinolide B in the solvent of choice, which should be purged with an inert gas. Aspinolide B is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

Aspinolide B is a polyketide synthase-derived aspinolide that has been found in *T. arundinaceum* and has plant growth regulatory activity.^{1,2} It promotes the growth of *S. lycopersicum* lateral roots when used at a concentration of 50 µg/ml.²

References

1. Cardoza, R.E., McCormick, S.P., Izquierdo-Bueno, I., *et al.* Identification of polyketide synthase genes required for aspinolide biosynthesis in *Trichoderma arundinaceum*. *Appl. Microbiol. Biotechnol.* **106(21)**, 7153-7171 (2022).
2. Malmierca, M.G., Barua, J., McCormick, S.P., *et al.* Novel aspinolide production by *Trichoderma arundinaceum* with a potential role in *Botrytis cinerea* antagonistic activity and plant defence priming. *Environ. Microbiol.* **17(4)**, 1103-1118 (2015).

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, 05/23/2024

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