

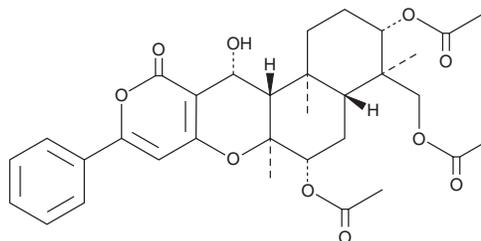
PRODUCT INFORMATION



Phenylpyropene A

Item No. 29594

CAS Registry No.: 189564-20-3
Formal Name: (3S,4R,4aR,6S,6aS,12R,12aS,12bS)-3,6-bis(acetyloxy)-4-[(acetyloxy)methyl]-1,3,4,4a,5,6,6a,12,12a,12b-decahydro-12-hydroxy-4,6a,12b-trimethyl-9-phenyl-2H,11H-naphtho[2,1-b]pyrano[3,4-e]pyran-11-one
MF: C₃₂H₃₈O₁₀
FW: 582.6
Purity: ≥95%
Supplied as: Residue
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

Phenylpyropene A is supplied as a residue. A stock solution may be made by dissolving the phenylpyropene A in the solvent of choice, which should be purged with an inert gas. Phenylpyropene A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Phenylpyropene A is a fungal metabolite originally isolated from *P. griseofulvum* that has enzyme inhibitory and insecticidal activities.¹⁻³ It inhibits acyl-coenzyme A:cholesterol acyltransferase (ACAT; IC₅₀ = 0.8 μM).¹ Phenylpyropene A inhibits diacylglycerol acyltransferase (DGAT) in rat liver microsomes (IC₅₀ = 78.7 μM). It induces 100% mortality in 100% of *M. persicae* when used at a concentration of 5 ppm.³

References

1. Kwon, O.E., Rho, M.C., Song, H.Y., *et al.* Phenylpyropene A and B, new inhibitors of acyl-CoA: Cholesterol acyltransferase produced by *Penicillium griseofulvum* F1959. *J. Antibiot. (Tokyo)* **55(11)**, 1004-1008 (2002).
2. Lee, S.W., Rho, M.C., Choi, J.H., *et al.* Inhibition of diacylglycerol acyltransferase by phenylpyropenes produced by *Penicillium griseofulvum* F1959. *J. Microbiol. Biotechnol.* **18(11)**, 1785-1788 (2008).
3. Horikoshi, R., Goto, K., Mitomi, M., *et al.* Identification of pyripyropene A as a promising insecticidal compound in a microbial metabolite screening. *J. Antibiot. (Tokyo)* **70(3)**, 272-276 (2017).

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, 08/15/2023

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