

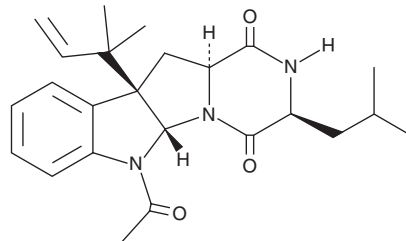
PRODUCT INFORMATION



Verrucofortine

Item No. 28062

CAS Registry No.: 113706-21-1
Formal Name: (3S,5aR,10bR,11aS)-6-acetyl-10b-(1,1-dimethyl-2-propen-1-yl)-6,10b,11,11a-tetrahydro-3-(2-methylpropyl)-2H-pyrazino[1',2':1,5]pyrrolo[2,3-b]indole-1,4(3H,5aH)-dione
Synonyms: Fructigenine B, (-)-Verrucofortine, Verrucozine
MF: C₂₄H₃₁N₃O₃
FW: 409.5
Purity: ≥98%
Supplied as: An oil
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

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

Description

Verrucofortine is a fungal metabolite originally isolated from *P. verrucosum*.¹

Reference

1. Hodge, R.P., Harris, C.M., and Harris, T.M. Verrucofortine, a major metabolite of *Penicillium verrucosum* var. *cyclopium*, the fungus that produces the mycotoxin verrucosidin. *J. Nat. Prod.* **51**(1), 66-73 (1988).

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, 08/10/2023

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