

PRODUCT INFORMATION

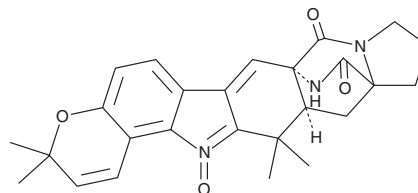


Avrainvillamide

Item No. 31655

CAS Registry No.: 269741-97-1
Formal Name: 3,11,12,13,13aS,14-hexahydro-3,3,14,14-tetramethyl-8H,10H-7aS,12aS-(iminomethano)indolizino[6,7-h]pyrano[3,2-a]carbazole-8,16-dione 15-oxide

Synonym: CJ-17,665
MF: C₂₆H₂₇N₃O₄
FW: 445.5
Purity: ≥90%
UV/Vis.: λ_{max}: 214, 292 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus* sp.



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

Laboratory Procedures

Avrainvillamide is supplied as a solid. A stock solution may be made by dissolving the avrainvillamide in the solvent of choice, which should be purged with an inert gas. Avrainvillamide is soluble in organic solvents such as methanol and DMSO.

Description

Avrainvillamide is a fungal metabolite and monomeric form of stephacidin B (Item No. 31656) that has been found in *A. ochraceus*.^{1,2} It is active against *S. aureus*, *S. pyogenes*, and *E. faecalis* (MICs = 12.5, 12.5, and 25 µg/ml, respectively).² Avrainvillamide inhibits the growth of LNCaP, BT-549, T47D, and MALME-3M cancer cells with GI₅₀ values of 0.24, 0.62, 0.21, and 0.41 µM, respectively.¹

References

1. Wulff, J.E., Herzon, S.B., Siegrist, R., *et al.* Evidence for the rapid conversion of stephacidin B into the electrophilic monomer avrainvillamide in cell culture. *J. Am. Chem. Soc.* **129(16)**, 4898-4899 (2007).
2. Sugie, Y., Hirai, H., Inagaki, T., *et al.* A new antibiotic CJ-17,665 from *Aspergillus ochraceus*. *J. Antibiot.* **54(11)**, 911-916 (2001).

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