

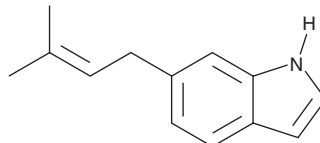
PRODUCT INFORMATION



6-Prenylindole

Item No. 28465

CAS Registry No.: 23158-16-9
Formal Name: 6-(3-methyl-2-buten-1-yl)-1H-indole
MF: C₁₃H₁₅N
FW: 185.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/Streptomyces sp.



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

Laboratory Procedures

6-Prenylindole is supplied as a solid. A stock solution may be made by dissolving the 6-prenylindole in the solvent of choice, which should be purged with an inert gas. 6-Prenylindole is soluble in ethanol, methanol, and DMSO.

Description

6-Prenylindole is a bacterial metabolite that has been found in *Streptomyces* and has antifungal and antimalarial properties.¹ It is active against *A. brassicicola* strain TP-F0423 and *F. oxysporum* f. sp. *tulipae* TU-4-2 (15 and 30 µg/disc in the paper disc assay), and also drug-resistant *P. falciparum* strain K1 (IC₅₀ = 21 µg/ml).²

References

1. Sasaki, T., Igarashi, Y., Ogawa, M., et al. Identification of 6-prenylindole as an antifungal metabolite of *Streptomyces* sp. TP-A0595 and synthesis and bioactivity of 6-substituted indoles. *J. Antibiot. (Tokyo)* **55(11)**, 1009-1012 (2002).
2. Nkunya, M.H., Makangara, J.J., and Jonker, S.A. Prenylindoles from Tanzanian *Monodora* and *Isolona* species. *Nat. Prod. Res.* **18(3)**, 253-258 (2004).

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, 12/05/2022

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