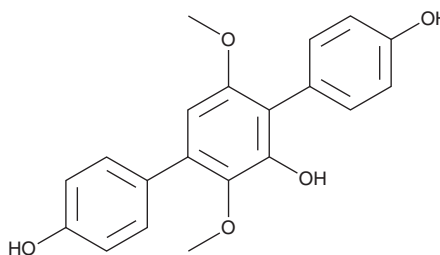


PRODUCT INFORMATION



Terphenyllin Item No. 23480

CAS Registry No.: 52452-60-5
Formal Name: 3',6'-dimethoxy-[1,1':4',1''-terphenyl]-2',4,4''-triol
Synonym: NSC 299114
MF: C₂₀H₁₈O₅
FW: 338.4
Purity: ≥85%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus candidus*



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

Laboratory Procedures

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

Description

Terphenyllin is a flavonoid metabolite originally isolated from *A. candidus* that has antibacterial and cytotoxic properties.^{1,2} It is active against methicillin-resistant *S. aureus* 1 (MRSA-1), MRSA-2, and *V. vulnificus* (MICs = 31.27-31.47 µg/ml) and cytotoxic to HeLa, A549, and HepG2 cells (IC₅₀s = 18.87, 12.33, and 21.2 µM, respectively).² Terphenyllin inhibits HIV-1 integrase in coupled and strand transfer assays with IC₅₀ values of 17.7 µM and 47.7 µM.³

References

1. Marchelli, R. and Vining, L.C. Terphenyllin, a novel *p*-terphenyl metabolite from *aspergillus candidus*. *J. Antibiot. (Tokyo)* **28(4)**, 328-331 (1975).
2. Wang, W., Liao, Y., Tang, C., et al. Cytotoxic and antibacterial compounds from the coral-derived fungus *Aspergillus tritici* SP2-8-1. *Mar. Drugs* **15(11)**, pii: E348 (2017).
3. Singh, S.B., Jayasuriya, H., Dewey, R., et al. Isolation, structure, and HIV-1-integrase inhibitory activity of structurally diverse fungal metabolites. *J. Ind. Microbiol. Biotechnol.* **30(12)**, 721-731 (2003).

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, 08/23/2023

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