

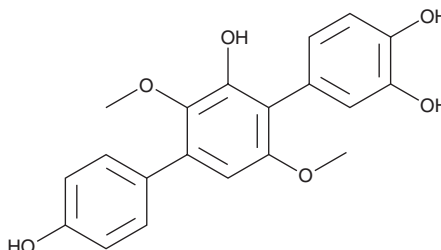
PRODUCT INFORMATION



3-Hydroxyterphenyllin

Item No. 28502

CAS Registry No.: 66163-76-6
Formal Name: 3',6'-dimethoxy-[1,1':4',1''-terphenyl]-2',3,4,4''-tetrol
Synonym: NSC 299113
MF: C₂₀H₁₈O₆
FW: 354.4
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

3-Hydroxyterphenyllin is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxyterphenyllin in the solvent of choice, which should be purged with an inert gas. 3-Hydroxyterphenyllin is soluble in organic solvents such as ethanol and DMSO.

Description

3-Hydroxyterphenyllin is a *p*-terphenyl fungal metabolite originally isolated from *A. candidus* that has diverse biological activities, including antioxidant, antiproliferative, antibacterial, and antiviral properties.¹⁻⁴ It has a 96% scavenging effect on 2,2-diphenyl-1-picrylhydrazyl radicals (DPPH; Item No. 14805) when used at a concentration of 100 µg/ml.² 3-Hydroxyterphenyllin inhibits the growth of HeLa cervical, A549 lung, and HepG2 liver cancer cells (IC₅₀s = 23, 36, and 32 µM, respectively), as well as methicillin-resistant *S. aureus* (MRSA) and *V. vulnificus* bacteria (MIC = 31 µg/ml for both).³ It also inhibits HIV-1 integrase in both coupled and strand transfer assays (IC₅₀s = 2.8 and 12.1 µM, respectively).⁴

References

1. Kurobane, I., Vining, L.C., McInnes, A.G., *et al.* 3-Hydroxyterphenyllin, a new metabolite of *Aspergillus candidus*. Structure elucidation by ¹H and ¹³C nuclear magnetic resonance spectroscopy. *J. Antibiot. (Tokyo)* **32(6)**, 559-564 (1979).
2. Yen, G.-C., Chang, Y.-C., Sheu, F., *et al.* Isolation and characterization of antioxidant compounds from *Aspergillus candidus* broth filtrate. *J. Agric. Food Chem.* **49(3)**, 1426-1431 (2001).
3. 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)**, E348 (2017).
4. 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, 08/16/2023

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