

PRODUCT INFORMATION



(-)-Cyclophenol

Item No. 27723

CAS Registry No.: 20007-85-6
Formal Name: (2'S,3'R)-3'-(3-hydroxyphenyl)-
4-methyl-spiro[3H-1,4-
benzodiazepine-3,2'-oxirane]-
2,5(1H,4H)-dione

Synonym: NSC 604990

MF: C₁₇H₁₄N₂O₄

FW: 310.3

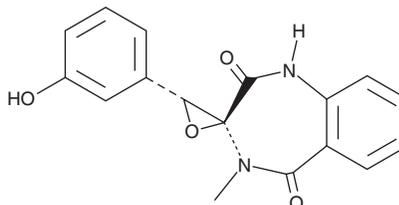
Purity: ≥95%

Supplied as: A solid

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

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

Description

(-)-Cyclophenol is a benzodiazepine alkaloid fungal metabolite originally isolated from *P. viridicatum*.¹ It inhibits protein tyrosine phosphatase 1B (PTP1B; IC₅₀ = 30 μM).²

References

1. Mohammed, Y.S. and Luckner, M. The structure of cyclophenin and cyclophenol, metabolic products from *penicillium cyclopium* westling and *penicillium viridicatum* westling. *Tetrahedron Lett.* **4(28)**, 1953-1958 (1963).
2. Sohn, J.H., Lee, Y.-R., Lee, D.-S., *et al.* PTP1B inhibitory secondary metabolites from marine-derived fungal strains *Penicillium* spp. and *Eurotium* sp. *J. Microbiol. Biotechnol.* **23(9)**, 1206-1211 (2013).

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