

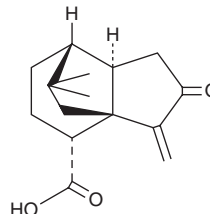
PRODUCT INFORMATION



Terrecyclic Acid

Item No. 28117

CAS Registry No.: 83058-94-0
Formal Name: (3aS,4R,7R,7aR)-octahydro-8,8-dimethyl-3-methylene-2-oxo-3a,7-ethano-3aH-indene-4-carboxylic acid
MF: $C_{15}H_{20}O_3$
FW: 248.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Fungus/*Aspergillus terreus*



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

Laboratory Procedures

Terrecyclic acid is supplied as a solid. A stock solution may be made by dissolving the terrecyclic acid in the solvent of choice. Terrecyclic acid is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Terrecyclic acid is a sesquiterpene originally isolated from *A. terreus* with antibiotic and anticancer activity.¹ It is active against *S. aureus*, *B. subtilis*, and *M. roseus* (MICs = 25, 50, and 25 $\mu g/ml$, respectively). Terrecyclic acid induces a heat shock response, increases levels of reactive oxygen species (ROS), and inhibits NF- κ B activity and cell growth in 3T3-Y9-B12 cells.² *In vivo*, terrecyclic acid (0.1, 1, and 10 mg/ml) reduces the number of ascitic fluid tumor cells in a mouse model of P388 murine leukemia.¹

References

1. Nakagawa, M., Hirota, A., Sakai, H., *et al.* Terrecyclic acid A, a new antibiotic from *Aspergillus terreus*. I. Taxonomy, production, and chemical and biological properties. *J. Antibiot. (Tokyo)* **35**(7), 778-782 (1982).
2. Turbyville, T.J., Wijeratne, E.M.K., Whitesell, L., *et al.* The anticancer activity of the fungal metabolite terrecyclic acid A is associated with modulation of multiple cellular stress response pathways. *Mol. Cancer Ther.* **4**(10), 1569-1576 (2005).

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, 08/11/2023

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