

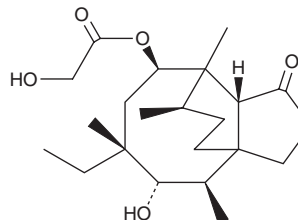
PRODUCT INFORMATION



Dihydropleuromutilin

Item No. 26497

CAS Registry No.: 42302-24-9
Formal Name: 2-hydroxy-acetic acid,
(3aS,4R,5S,6R,8R,9R,9aR,10R)-6-ethyldecahydro-5-
hydroxy-4,6,9,10-tetramethyl-1-oxo-3a,9-propano-
3aH-cyclopentacycloocten-8-yl ester
Synonym: (+)-19,20-Dihydropleuromutilin
MF: C₂₂H₃₆O₅
FW: 380.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Dihydropleuromutilin is supplied as a solid. A stock solution may be made by dissolving the dihydropleuromutilin in the solvent of choice, which should be purged with an inert gas. Dihydropleuromutilin is soluble in the organic solvent DMSO.

Description

Dihydropleuromutilin is a semisynthetic antibiotic and derivative of pleuromutilin (Item No. 19452).¹ It is active against *S. aureus*, *M. hominis*, *M. gallisepticum*, and *M. hyorhinis* bacteria (MICs = 0.5, 0.3, 0.3, and 2 µg/ml, respectively).

Reference

1. Egger, H. and Reinshagen, H. New pleuromutilin derivatives with enhanced antimicrobial activity. II. Structure-activity correlations. *J. Antibiot. (Tokyo)* **29(9)**, 923-927 (1976).

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, 11/07/2022

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