

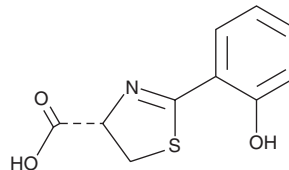
PRODUCT INFORMATION



Dihydroaeruginoic Acid

Item No. 25514

CAS Registry No.: 143209-04-5
Formal Name: (4S)-4,5-dihydro-2-(2-hydroxyphenyl)-4-thiazolecarboxylic acid
Synonym: CGP 52547
MF: C₁₀H₉NO₃S
FW: 223.2
Purity: ≥98%
UV/Vis.: λ_{max}: 221 nm
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

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

Description

Dihydroaeruginoic acid is an antibiotic originally isolated from *P. fluorescens*.¹ It has antimicrobial activity in a disc assay against *R. solani*, *P. ultimum*, *B. cinera*, *S. rolfsii*, *C. gloeosporioides*, *F. oxysporum*, and *S. tritici* fungi as well as *B. subtilis*, *E. herbicola*, and *S. albus* bacteria when used at a concentration of 200 µg/disc.

Reference

1. Carmi, R., Carmeli, S., Levy, E., et al. (+)-(S)-dihydroaeruginoic acid, an inhibitor of *Septoria tritici* and other phytopathogenic fungi and bacteria, produced by *Pseudomonas fluorescens*. *J. Nat. Prod.* **57(9)**, 1200-1205 (1994).

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, 11/30/2022

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