

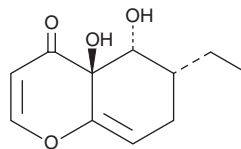
PRODUCT INFORMATION



Agistatin D

Item No. 29169

CAS Registry No.: 144096-47-9
Formal Name: (4aS,5R,6R)-6-ethyl-4a,5,6,7-tetrahydro-4a,5-dihydroxy-4H-1-benzopyran-4-one
MF: C₁₁H₁₄O₄
FW: 210.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Fungus A 6239



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

Laboratory Procedures

Agistatin D is supplied as a solid. A stock solution may be made by dissolving the agistatin D in the solvent of choice, which should be purged with an inert gas. Agistatin D is soluble in acetone and DMSO.

Description

Agistatin D is a pyranacetal fungal metabolite that has been found in the soil fungus FH-A 6239.¹

Reference

1. Göhrt, A., Grabley, S., Thiericke, R., *et al.* Secondary metabolites by chemical screening. Part 29. Agistatines, novel pyranacetals from the fungus FH-A 6239. *Liebigs Ann.* **4**, 627-633 (1996).

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