

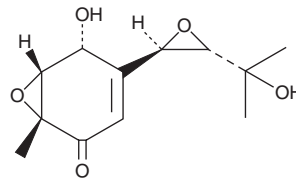
PRODUCT INFORMATION



Acremine I

Item No. 29771

CAS Registry No.: 1110661-29-4
Formal Name: (1R,5R,6R)-5-hydroxy-4-[(2R,3S)-3-(1-hydroxy-1-methylethyl)-2-oxiranyl]-1-methyl-7-oxabicyclo[4.1.0]hept-3-en-2-one
MF: C₁₂H₁₆O₅
FW: 240.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

Acremine I is supplied as a solid. A stock solution may be made by dissolving the acremine I in the solvent of choice, which should be purged with an inert gas. Acremine I is soluble in organic solvents such as ethanol, methanol, dichloromethane, and DMSO.

Description

Acremine I is a fungal metabolite originally isolated from *Acronium byssoides*.¹ It inhibits germination of *P. viticola* sporangia by 23.2 and 32.8% when used at concentrations of 0.5 and 1 mM, respectively.

Reference

1. Arnone, A., Assante, G., Bava, A., et al. Acremine H–N, novel prenylated polyketide metabolites produced by a strain of *Acronium byssoides*. *Tetrahedron* **65**(4), 786-791 (2009).

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