

PRODUCT INFORMATION



Acremine F

Item No. 33228

CAS Registry No.: 863480-61-9
Formal Name: (1S,2R,4S)-5-[(1E)-3-hydroxy-3-methyl-1-buten-1-yl]-2-methyl-5-cyclohexene-1,2,4-triol

Synonym: Acremin F

MF: C₁₂H₂₀O₄

FW: 228.3

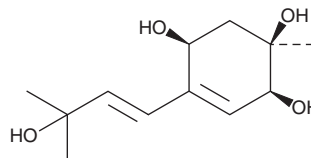
Purity: ≥95%

Supplied as: A residue

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 F is supplied as a residue. A stock solution may be made by dissolving the acremine F in the solvent of choice, which should be purged with an inert gas. Acremine F is soluble in methanol and DMSO.

Description

Acremine F is a fungal metabolite that has been found in *Acremonium*.¹

Reference

1. Assante, G., Dallavalle, S., Malpezzi, L., *et al.* Acremines A–F, novel secondary metabolites produced by a strain of an endophytic *Acremonium*, isolated from sporangiophores of *Plasmopara viticola* in grapevine leaves. *Tetrahedron* **61**(32), 7686-7592 (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, 08/25/2023

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