# PRODUCT INFORMATION



# Janthitrem A

Item No. 30166

CAS Registry No.: 73561-89-4

Formal Name: (2R,3S,3aR,4aS,4bS,6aS,9S,9aR,15bS,15cR,

> 17aS)-10,10,12,12,15b,15c-hexamethyl-2-(prop-1-en-2-yl)-3,3a,5,6,6a,7,9,9a,10,12,15, 15b,15c,16,17,17a-hexadecahydro-2H,4bH-

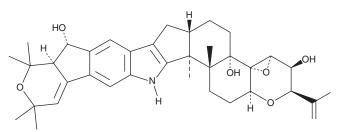
oxireno[2",3":4',4a']chromeno[5',6':6,7] indeno[1,2-b]pyrano[4',3':3,4] cyclopenta[1,2-f]indole-3,4b,9-triol

Synonym: 11,12-Epoxyjanthitrem B

MF:  $C_{37}H_{47}NO_{6}$ FW: 601.8 **Purity:** ≥95% A solid Supplied as: -20°C Storage: Stability: ≥4 years

Item Origin: Fungus/Penicillium sp.

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



## **Laboratory Procedures**

Janthitrem A is supplied as a solid. A stock solution may be made by dissolving the janthitrem A in the solvent of choice, which should be purged with an inert gas. Janthitrem A is soluble in methanol, ethanol, DMSO, and dimethyl formamide.

### Description

Janthitrem A is a mycotoxin that has been found in P. janthinellum. 1 It induces tremors in mice and decreases the time to fall from the rotarod when administered at a dose of 4 mg/kg.<sup>2</sup> Janthitrem A (20 µg/g) also reduces W. cervinata larvae weight gain and food consumption.

### References

- 1. Gallagher, R.T., Latch, G.C., and Keogh, R.G. The janthitrems: Fluorescent tremorgenic toxins produced by Penicillium janthinellum isolates from ryegrass pastures. Appl. Environ. Microbiol. 39(1), 272-273 (1980).
- 2. Babu, J.V., Popay, A.J., Miles, C.O., et al. Identification and structure elucidation of janthitrems A and D from Penicillium janthinellum and determination of the tremorgenic and anti-insect activity of janthitrems A and B. J. Agric. Food Chem. 66(50), 13116-13125 (2018).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the material can be found on our website.

Copyright Cayman Chemical Company, 08/10/2023

### **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