

# PRODUCT INFORMATION



## (+)-Anominine

Item No. 32923

**CAS Registry No.:** 120991-21-1  
**Formal Name:** (1S,4R,4aS,8R,8aS)-decahydro-8-(1H-indol-3-ylmethyl)-4,4a-dimethyl-7-methylene-8a-(4-methyl-3-penten-1-yl)-1-naphthalenol

**Synonym:** (+)-Nominine

**MF:** C<sub>28</sub>H<sub>39</sub>NO

**FW:** 405.6

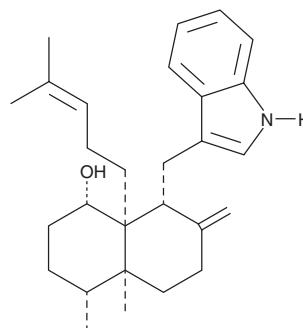
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungi/*Aspergillus* sp.



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

### Laboratory Procedures

(+)-Anominine is supplied as a solid. A stock solution may be made by dissolving the (+)-anominine in the solvent of choice, which should be purged with an inert gas. (+)-Anominine is soluble in methanol and DMSO.

### Description

(+)-Anominine is an indole diterpene fungal metabolite that has been found in *A. nomius* and has insecticidal activity.<sup>1</sup> It induces mortality in 40% of *H. zea* when administered in the diet at 100 ppm and has antifeedant activity in *C. hemipterus* larvae.

### Reference

1. Gloer, J.B. and Rinderknecht, B.L. Nominine: A new insecticidal indole diterpene from the sclerotia of *Aspergillus nomius*. *J. Org. Chem.* **54**(11), 2530-2532 (1989).

#### 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, 11/10/2022

#### 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