

# PRODUCT INFORMATION

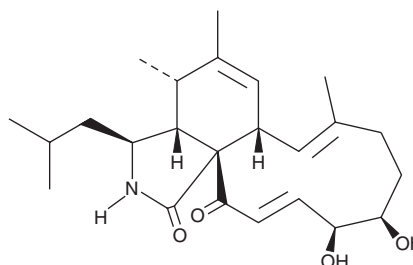


## Aspochalasin D

Item No. 23864

**CAS Registry No.:** 71968-02-0  
**Formal Name:** (3S,3aR,4S,6aS,7E,11R,12S,13E,15aS)-3,3a,4,6a,9,10,11,12-octahydro-11,12-dihydroxy-4,5,8-trimethyl-3-(2-methylpropyl)-1H-cycloundec[d]isoindole-1,15(2H)-dione

**MF:** C<sub>24</sub>H<sub>35</sub>NO<sub>4</sub>  
**FW:** 401.5  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

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

### Description

Aspochalasin D is a co-metabolite originally isolated from *A. microcysticus* with aspochalasin A, B, and C, that was initially thought to be inactive.<sup>1</sup> It has antibacterial activity against Gram-positive and Gram-negative bacteria at a concentration of 1 mg/ml.<sup>2</sup> Aspochalasin D is more cytotoxic, *via* apoptosis, to Ba/F3-V12 cells in an IL-3-free medium than in an IL-3-containing medium (IC<sub>50</sub>s = 0.49 and 1.9 µg/ml, respectively).<sup>3</sup>

### References

1. Keller-Schierlein, W. and Kupfer, E. Metabolites of microorganisms. The Aspochalasin A, B, C, and D. *Helvetica Chimica Acta* **62(5)**, 1501-1524 (1979).
2. Gebhardt, K., Schimana, J., Höltzel, A., *et al.* Aspochalamins A-D and aspochalasin Z produced by the endosymbiotic fungus *Aspergillus niveus* LU 9575 I. Taxonomy, fermentation, isolation and biological activities. *J. Antibiot. (Tokyo)* **57(11)**, 707-714 (2004).
3. Tomikawa, T., Shin-Ya, K., Kinoshita, T., *et al.* Selective cytotoxicity and stereochemistry of aspochalasin D. *J. Antibiot. (Tokyo)* **54(4)**, 379-381 (2001).

#### 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, 10/06/2022

#### 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