

PRODUCT INFORMATION



Aspergillin PZ

Item No. 27722

CAS Registry No.: 483305-08-4
Formal Name: (1R,2S,5R,5aS,5bS,8R,8aS,9R,11aR,13aR)-2,3,4,5,5a,5b,8,8a,9,10,13,13a-dodecahydro-2-hydroxy-5,7,8-trimethyl-9-(2-methylpropyl)-1,5-epoxy-1H-cyclohepta[3,4]benz[1,2-d]isoindole-11,12-dione

Synonym: (+)-Aspergillin PZ

MF: C₂₄H₃₅NO₄

FW: 401.5

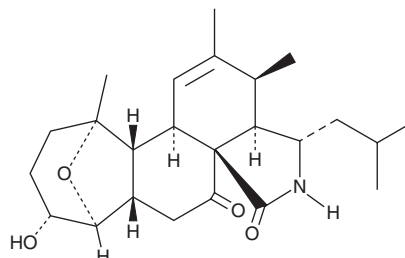
Purity: ≥98%

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

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

Description

Aspergillin PZ is a fungal metabolite originally isolated from *A. awamori*.¹ It induces morphological deformation of *P. oryzae* conidia when used at a concentration of 89 nM. Aspergillin PZ is active against *S. epidermidis* (MIC = 20 μM) but not *S. aureus*, *E. coli*, or *B. cereus* (MICs = >20 μM).² It is cytotoxic to HL-60 promyelocytic leukemia cells (IC₅₀ = 56.61 μM) but not THP-1 acute monocytic leukemia or PC3 prostate cancer cells (IC₅₀s = >80 μM).³

References

1. Zhang, Y., Wang, T., Pei, Y., *et al.* Aspergillin PZ, a novel isoindole-alkaloid from *Aspergillus awamori*. *J. Antibiot. (Tokyo)* **55**(8), 693-695 (2002).
2. Zheng, C.-J., Shao, C.-L., Wu, L.-Y., *et al.* Bioactive phenylalanine derivatives and cytochalasins from the soft coral-derived fungus, *Aspergillus elegans*. *Mar. Drugs* **11**(6), 2054-2068 (2013).
3. Si, Y., Tang, M., Lin, S., *et al.* Cytotoxic cytochalasans from *Aspergillus flavipes* PJ03-11 by OSMAC method. *Tetrahedron Lett.* **59**(18), 1767-1771 (2018).

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, 12/22/2022

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