

PRODUCT INFORMATION



Fumiquinazoline D

Item No. 23848

CAS Registry No.: 140715-86-2
Formal Name: (1'R,2S,3'S,5'S,9aS)-1,4',5',9a-tetrahydro-1',2-dimethyl-spiro[9H-imidazo[1,2-a]indole-9,3'-[1,5](iminomethano)[3H][1,4]oxazepino[3,4-b]quinazoline]-3,7',13'(1'H,2H)-trione

MF: C₂₄H₂₁N₅O₄

FW: 443.5

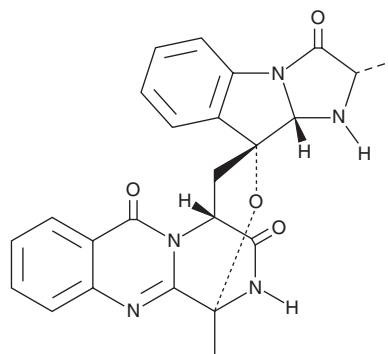
Purity: ≥85%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Aspergillus fumigatus*



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

Laboratory Procedures

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

Description

Fumiquinazoline D is a fungal metabolite originally isolated from *A. fumigatus* that has activity against Gram-positive and Gram-negative bacteria (MICs = 8-16 µg/ml) as well as *F. solani* and *C. albicans* fungi (MICs = 32 and 64 µg/ml, respectively).^{1,2}

References

1. Numata, A., Takahashi, C., Miyamoto, T., *et al.* Structures of cytotoxic substances and new quinazoline derivatives produced by a fungus from a saltwater fish. *Tennen Yuki Kagobutsu Toronkai Koen Yoshishu* **33**, 723-730 (1991).
2. Zhang, H., Liu, R., Yang, J., *et al.* Bioactive alkaloids of *Aspergillus fumigatus*, an endophytic fungus from *Astragalus membranaceus*. *Chem. Nat. Compd.* **53**(4), 802-805 (2017).

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, 08/23/2023

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