

PRODUCT INFORMATION



Fumitremorgin B

Item No. 31173

CAS Registry No.: 12626-17-4

Formal Name: (5aR,6S,12S,14aS)-1,2,3,5a,6,11,12,14a-octahydro-5a,6-dihydroxy-9-methoxy-11-(3-methyl-2-buten-1-yl)-12-(2-methyl-1-propen-1-yl)-5H,14H-pyrrolo[1'',2'':4',5']pyrazino[1',2':1,6]pyrido[3,4-b]indole-5,14-dione

Synonym: Lanosulin

MF: $C_{27}H_{33}N_3O_5$

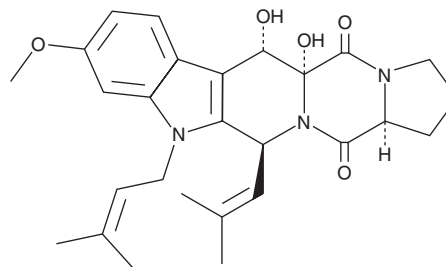
FW: 479.6

Purity: $\geq 95\%$

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

Fumitremorgin B is supplied as a solid. A stock solution may be made by dissolving the fumitremorgin B in the solvent of choice, which should be purged with an inert gas. Fumitremorgin B is soluble in methanol and DMSO.

Description

Fumitremorgin B is a mycotoxin originally isolated from *A. fumigatus*.^{1,2} It induces tremors in and is lethal to mice ($\text{LD}_{50} = 5.4 \text{ mg/kg, i.v.}$). It is also toxic to brine shrimp ($\text{LC}_{50} = 13.6 \text{ }\mu\text{g/ml}$).³ Fumitremorgin B inhibits the proliferation of U937 and PC3 cells ($\text{IC}_{50}\text{s} = 14.12$ and $43.36 \text{ }\mu\text{M}$, respectively).⁴

References

1. Yamazaki, M., Suzuki, S., and Miyaki, K. Tremorgenic toxins from *Aspergillus fumigatus* Fres. *Chem. Pharm. Bull. (Tokyo)* **19**(8), 1739-1740 (1971).
2. Yamazaki, M. and Suzuki, S. Toxicology of tremorgenic mycotoxins, fumitremorgin A and B. *Dev. Toxicol. Environ. Sci.* **12**, 273-282 (1986).
3. Li, X.-J., Zhang, Q., Zhang, A.-L., et al. Metabolites from *Aspergillus fumigatus*, an endophytic fungus associated with *Melia azedarach*, and their antifungal, antifeedant, and toxic activities. *J. Agric. Food Chem.* **60**(13), 3424-3431 (2012).
4. Wang, Y., Li, Z.-L., Bai, J., et al. 2,5-diketopiperazines from the marine-derived fungus *Aspergillus fumigatus* YK-7. *Chem. Biodivers.* **9**(2), 85-93 (2012).

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, 08/25/2023

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