

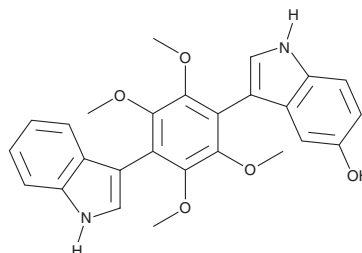
PRODUCT INFORMATION



Petromurin C

Item No. 25089

CAS Registry No.: 194608-29-2
Formal Name: 3-[4-(1H-indol-3-yl)-2,3,5,6-tetramethoxyphenyl]-1H-indol-5-ol
MF: $C_{26}H_{24}N_2O_5$
FW: 444.5
Purity: $\geq 95\%$
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

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

Description

Petromurin C is a bis-indolyl benzenoid fungal metabolite originally isolated from the ascostromata of *P. muricatus*.¹ It is cytotoxic to NS-1 mouse myeloma cells ($IC_{50} = 33 \mu\text{g/ml}$) and has activity against *T. foetus* *in vitro* ($IC_{50} = 100 \mu\text{g/ml}$).²

References

1. Ooike, M., Nozawa, K., Udagawa, S., *et al.* Bisindolylbenzenoids from ascostromata of *Petromyces muricatus*. *Can. J. Chem.* **75**(6), 625-628 (1997).
2. Lacey, H.J., Vuong, D., Pitt, J.I., *et al.* Kumbicins A–D: Bis-Indolyl benzenoids and benzoquinones from an Australian soil fungus, *Aspergillus kumbius*. *Aust. J. Chem.* **69**(2), 152-160 (2015).

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, 10/20/2022

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