

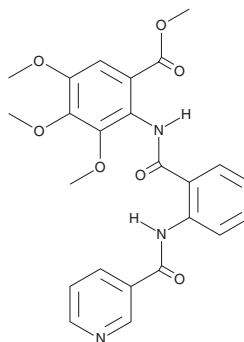
PRODUCT INFORMATION



Preterramide C

Item No. 33097

CAS Registry No.: 81469-77-4
Formal Name: 3,4,5-trimethoxy-2-[[2-[[3-pyridinylcarbonyl)amino]benzoyl]amino]-benzoic acid, methyl ester
MF: C₂₄H₂₃N₃O₇
FW: 465.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus* sp.



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

Laboratory Procedures

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

Description

Preterramide C is a fungal metabolite that has been found in *A. terreus*.¹ It is active against *S. aureus* (MIC = 52.4 μM) but not *E. aerogenes*, *P. aeruginosa*, or *C. albicans* (MICs = >100 μM for all).² Preterramide C inhibits the growth of L5178Y mouse lymphoma cells with an IC₅₀ value of 0.1 μg/ml.³ It also inhibits nitric oxide (NO) production induced by LPS in RAW 264.7 cells (IC₅₀ = 5.48 μM).

References

1. He, F., Bao, J., Zhang, X.-Y., *et al.* Asperterrestide A, a cytotoxic cyclic tetrapeptide from the marine-derived fungus *Aspergillus terreus* SCSGAF0162. *J. Nat. Prod.* **76(6)**, 1182-1186 (2013).
2. Wang, Y., Zheng, J., Liu, P., *et al.* Three new compounds from *Aspergillus terreus* PT06-2 grown in a high salt medium. *Mar. Drugs* **9(8)**, 1368-1378 (2011).
3. Zhou, Y., Mandi, A., Debbab, A., *et al.* New austrialides from the sponge-associated fungus *Aspergillus* sp. *Eur. J. Org. Chem.* **30**, 6009-6019 (2011).

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