

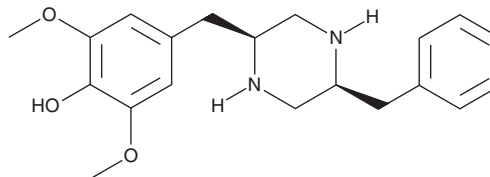
PRODUCT INFORMATION



Hancockiamide D

Item No. 45497

CAS Registry No.: 2758265-37-9
Formal Name: 2,6-dimethoxy-4-[[[(2S,5S)-5-(phenylmethyl)-2-piperazinyl]methyl]-phenol
MF: $C_{20}H_{26}N_2O_3$
FW: 342.4
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungi/*Aspergillus hancockii*



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

Laboratory Procedures

Hancockiamide D is supplied as a crystalline solid. A stock solution may be made by dissolving the hancockiamide D in the solvent of choice, which should be purged with an inert gas. Hancockiamide D is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Hancockiamide D is a fungal metabolite with anticancer and anti-germination activity.¹ It is cytotoxic to NS-1 murine myeloma cells (MIC = 1.6 $\mu\text{g/ml}$). Hancockiamide D inhibits germination in a seed assay using the dicotyledon species *A. thaliana* (MIC = 50 $\mu\text{g/ml}$) but not the monocotyledon species *E. tef* at 100 $\mu\text{g/ml}$.

Reference

1. Li, H., Lacey, A.E., Shu, S., *et al.* Hancockiamides: Phenylpropanoid piperazines from *Aspergillus hancockii* are biosynthesised by a versatile dual single-module NRPS pathway. *Org. Biomol. Chem.* **19**(3), 587-595 (2021).

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, 04/02/2026

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