

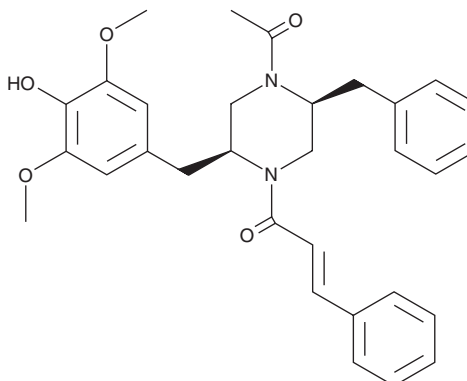
PRODUCT INFORMATION



Hancockiamide B

Item No. 44770

CAS Registry No.: 2758265-35-7
Formal Name: (2E)-1-[(2S,5S)-4-acetyl-2-[(4-hydroxy-3,5-dimethoxyphenyl)methyl]-5-(phenylmethyl)-1-piperazinyl]-3-phenyl-2-propen-1-one
MF: C₃₁H₃₄N₂O₅
FW: 514.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus hancockii*



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

Laboratory Procedures

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

Description

Hancockiamide B is an N-cinnamoylated piperazine fungal metabolite that has been found in *A. hancockii* and has antibacterial and anticancer activities.¹ It is active against *B. subtilis* (MIC = 50 µg/ml). Hancockiamide B is cytotoxic to NS-1 murine myeloma cells (MIC = 25 µg/ml) but not neonatal foreskin fibroblasts at 100 µg/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, 11/24/2025

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