

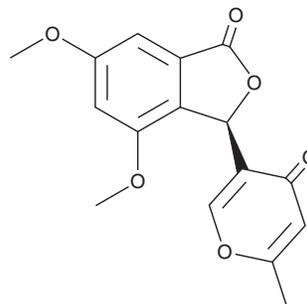
# PRODUCT INFORMATION



## Hancockiamide C

Item No. 44771

**CAS Registry No.:** 2758265-36-8  
**Formal Name:** 1-[(2S,5S)-4-acetyl-2-[(7-methoxy-1,3-benzodioxol-5-yl)methyl]-5-(phenylmethyl)-1-piperazinyl]-3-phenyl-2E-propen-1-one  
**MF:** C<sub>31</sub>H<sub>32</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 512.6  
**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 C is supplied as a solid. A stock solution may be made by dissolving the hancockiamide C in the solvent of choice, which should be purged with an inert gas. Hancockiamide C is soluble in methanol and DMSO.

### Description

Hancockiamide C is an N-cinnamoylated piperazine fungal metabolite that has been found in *A. hancockii* and has anti-germination activity.<sup>1</sup> It inhibits germination in a seed assay using the dicotyledon species *A. thaliana* (MIC = 6.3 µg/ml) but not the monocotyledon species *E. tef* 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, 11/25/2025

### 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