

PRODUCT INFORMATION



Hancockinone A

Item No. 44772

Formal Name: (R)-3-acetyl-5,6,8,11a-tetrahydroxy-10-(3-methyl-but-2-en-1-yl)-11,11a-dihydro-1H-cyclopenta[b]-anthracene-2,4-dione

MF: C₂₄H₂₂O₇

FW: 422.4

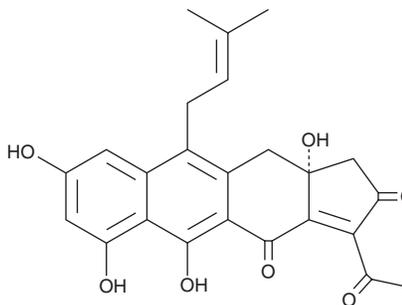
Purity: ≥95%

Supplied as: A 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

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

Description

Hancockinone A is a polycyclic aromatic polyketide that has been found in *A. hancockii* and has antibacterial activity.¹ It contains a unique prenylated 6/6/6/5 tetracyclic skeleton. Hancockinone A is active against *B. subtilis* and *S. aureus* (MIC = 25 µg/ml for both).

Reference

1. Li, H., Shu, S., Kalaitzis, J.A., *et al.* Genome mining of *Aspergillus hancockii* unearths cryptic polyketide hancockinone A featuring a prenylated 6/6/6/5 carbocyclic skeleton. *Org. Lett.* **23(22)**, 8789-8793 (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