

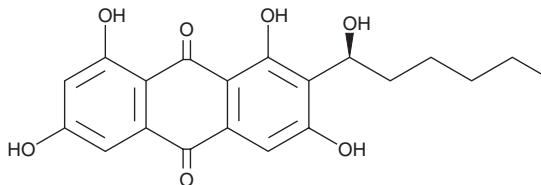
PRODUCT INFORMATION



Averantin

Item No. 28059

CAS Registry No.: 5803-62-3
Formal Name: 1,3,6,8-tetrahydroxy-2-[(1S)-1-hydroxyhexyl]-9,10-anthracenedione
Synonym: (-)-Averantin
MF: C₂₀H₂₀O₇
FW: 372.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 266, 295, 319, 453 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus* sp. WDMH51



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

Laboratory Procedures

Averantin is supplied as a crystalline solid. A stock solution may be made by dissolving the averantin in the solvent of choice, which should be purged with an inert gas. Averantin is soluble in organic solvents such as methanol and DMSO.

Description

Averantin is an anthraquinone fungal metabolite that has been found in *Aspergillus*.^{1,2} It is an intermediate in the biosynthesis of aflatoxin B₁ (Item No. 11293).¹ Averantin is active against *B. subtilis* bacteria and the plant pathogenic fungus *F. solani* *in vitro* (MICs = 16-32 µg/ml).²

References

1. Bennett, J.W., Lee, L.S., Shoss, S.M., *et al.* Identification of averantin as an aflatoxin B₁ precursor: Placement in the biosynthetic pathway. *Appl. Environ. Microbiol.* **39(4)**, 835-839 (1980).
2. Liu, K., Zheng, Y., Miao, C., *et al.* The antifungal metabolites obtained from the rhizospheric *Aspergillus* sp. YIM PH30001 against pathogenic fungi of *Panax notoginseng*. *Nat. Prod. Res.* **28(24)**, 2334-2337 (2014).

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, 08/10/2023

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