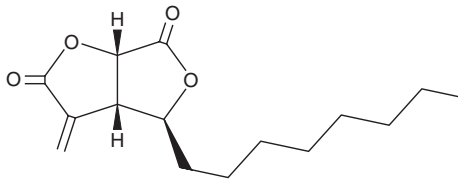


PRODUCT INFORMATION



Avenaciolide Item No. 28128

CAS Registry No.: 26057-70-5
Formal Name: (3aR,4R,6aR)-rel-dihydro-3-methylene-4-octyl-furo[3,4-b]furan-2,6(3H,4H)-dione
Synonym: (±)-Avenaciolide
MF: C₁₅H₂₂O₄
FW: 266.3
Purity: ≥95%
Supplied as: A residue
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungi/*Aspergillus* sp.



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

Laboratory Procedures

Avenaciolide is supplied as a residue. A stock solution may be made by dissolving the avenaciolide in the solvent of choice, which should be purged with an inert gas. Avenaciolide is soluble in ethanol, methanol, dimethyl formamide, and DMSO.

Description

Avenaciolide is a water-insoluble fungal metabolite originally isolated from *A. avenaceus*.¹ It inhibits glutamate transport in isolated rat liver mitochondria ($K_i = 8 \mu\text{M}$).

Reference

1. Johnson, R.N. and Chappell, J.B. The influx and efflux of phosphate in liver mitochondria. *Biochem J.* **116**(4), 37P (1970).

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, 10/11/2022

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