

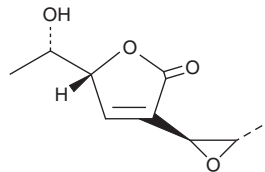
PRODUCT INFORMATION



Asperlactone

Item No. 14435

CAS Registry No.: 76375-62-7
Formal Name: 5R-[(1S)-1-hydroxyethyl]-3-[(2S,3S)-3-methyl-2-oxiranyl]-2(5H)-furanone
MF: C₉H₁₂O₄
FW: 184.2
Purity: ≥98%
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus ochraceus*



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

Laboratory Procedures

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

Description

Asperlactone is a polyketide synthase-derived polyketide that has been found in *A. westerdijkiae* and has anti-inflammatory activity.^{1,2} It inhibits superoxide anion release induced by N-formyl-Met-Leu-Phe (fMLP; Item No. 21495) in isolated human neutrophils when used at a concentration of 10 μM.²

References

1. Bacha, N., Dao, H.P., Atoui, A., *et al.* Cloning and characterization of novel methylsalicylic acid synthase gene involved in the biosynthesis of isoasperlactone and asperlactone in *Aspergillus westerdijkiae*. *Fungal Genet. Biol.* **46(10)**, 742-749 (2009).
2. Hu, H.-C., Li, C.-Y., Tsai, Y.-H., *et al.* Secondary metabolites and bioactivities of *Aspergillus ochraceopetaliformis* isolated from *Anthurium brownii*. *ACS Omega* **5(33)**, 20991-20999 (2020).

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, 08/28/2023

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