

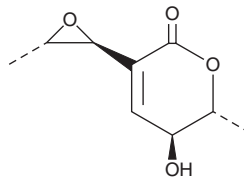
PRODUCT INFORMATION



Aspyrone

Item No. 31389

CAS Registry No.: 17398-00-4
Formal Name: (5S,6R)-5,6-dihydro-5-hydroxy-6-methyl-3-[(2S,3S)-3-methyl-2-oxiranyl]-2H-pyran-2-one
MF: $C_9H_{12}O_4$
FW: 184.2
Purity: $\geq 98\%$
Supplied as: A solid
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

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

Description

Aspyrone is a polyketide fungal metabolite that has been found in *Aspergillus* and has diverse biological activities.^{1,2} It is active against a panel of 13 fungi when used at a concentration of 20 $\mu\text{g/ml}$ and a panel of 21 bacteria in a disc assay when used at a concentration of 100 μg per disc.¹ Aspyrone (10-1,000 mg/L) is nematocidal against *P. penetrans*.²

References

1. Torres, M., Balcells, M., Sala, N., *et al.* Bactericidal and fungicidal activity of *Aspergillus ochraceus* metabolites and some derivatives. *Pestic. Sci.* **53**(1), 9-14 (1999).
2. Kimura, Y., Nakahara, S., and Fujioka, S. Aspyrone, a nematocidal compound isolated from the fungus, *Aspergillus melleus*. *Biosci. Biotech. Biochem.* **60**(8), 1375-1376 (1996).

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