

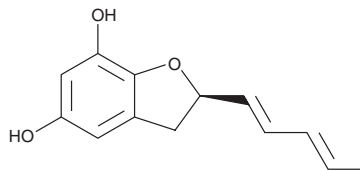
PRODUCT INFORMATION



Asperfuran

Item No. 29989

CAS Registry No.: 129277-10-7
Formal Name: (2R)-2,3-dihydro-2-(1E,3E)-1,3-pentadien-1-yl-5,7-benzofurandiol
Synonym: (-)-Asperfuran
MF: C₁₃H₁₄O₃
FW: 218.3
Purity: ≥95%
UV/Vis.: λ_{max}: 229 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Fungi/*Aspergillus* sp.



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

Laboratory Procedures

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

Description

Asperfuran is a fungal metabolite originally isolated from *A. oryzae* that has antifungal and anticancer activities.¹ It inhibits the growth of 23 fungi in a disc assay when used at a concentration of 50 µg per disc. Asperfuran inhibits chitin synthase with an IC₅₀ value of 300 µM. It inhibits proliferation of HeLa S3 and L1210 cancer cells (IC₅₀s = 25 µg/ml for both).

Reference

1. Pfefferle, W., Anke, H., Bross, M., *et al.* Asperfuran, a novel antifungal metabolite from *Aspergillus oryzae*. *J. Antibiot. (Tokyo)* **43(6)**, 648-654 (1990).

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, 07/20/2021

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