PRODUCT INFORMATION



Funalenone

Item No. 23491

CAS Registry No.: Formal Name:	259728-61-5 3,4,7,9-tetrahydroxy-2-methoxy- 6-methyl-1H-phenalen-1-one	OH OH
MF:	C ₁₅ H ₁₂ O ₆	
FW:	288.3	
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	но
Stability:	≥4 years	

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

Laboratory Procedures

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

Description

Funalenone is a phenalenone originally isolated from A. niger.¹ It inhibits HIV-1 integrase (IC₅₀ = 10 μ M) and HIV-1 replication in human peripheral blood cells transformed by murine leukemia virus (HPB-M(a); $IC_{50} = 1.7 \mu$ M) but is less cytotoxic to mammalian HPB-M(a) cells ($IC_{50} = 87 \mu$ M).² Funalenone selectively inhibits matrix metalloproteinase 1 (MMP-1; $IC_{50} = 170 \mu$ M) over MMP-2 and MMP-9, which it inhibits by 18.3 and 38.2%, respectively, when used at a concentration of 400 μ M.¹ It also inhibits the bacterial cell wall synthesis enzymes MraY and MurG (IC₅₀ = 25.5 μ M in a membrane plate assay) and inhibits growth of S. aureus with a MIC value of 64 μ g/ml.³

References

- 1. Inokoshi, J., Shiomi, K., Masuma, R., et al. Funalenone, a novel collagenase inhibitor produced by Aspergillus niger. J. Antibiot. (Tokyo) 52(12), 1095-1100 (1999).
- 2. Shiomi, K., Matsui, R., Isozaki, M., et al. Fungal phenalenones inhibit HIV-1 integrase. J. Antibiot. (Tokyo) 58(1), 65-68 (2005).
- 3. Zawadzke, L.E., Wu, P., Cook, L.S., et al. Targeting the MraY and MurG bacterial enzymes for antimicrobial therapeutic intervention. Anal. Biochem. 314(2), 243-252 (2003).

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

uyer 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, 12/16/2022

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