

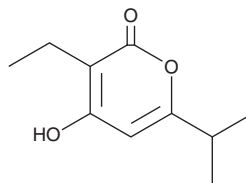
PRODUCT INFORMATION



Germicidin B

Item No. 20591

CAS Registry No.: 150973-78-7
Formal Name: 3-ethyl-4-hydroxy-6-(1-methylethyl)-
2H-pyran-2-one
MF: C₁₀H₁₄O₃
FW: 182.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

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

Description

Germicidin B is a pyranone antibiotic originally derived from *Streptomyces*. It acts as a reversible, autoregulative inhibitor of spore germination and hyphal elongation in the genus *Streptomyces* at concentrations as low as 200 pM.^{1,2} During germination, spores of *Streptomyces* excrete germicidin B along with other germicidin homologs, which inhibit germination of its own arthrospores.^{1,2} At higher concentrations, germicidin can inhibit the porcine brain Na⁺/K⁺-dependent ATPase (ID₅₀ = 100 μM) and prevent the germination of the cress *L. sativum*.¹

References

- Petersen, F. and Zähler, H. Germicidin, an autoregulative germination inhibitor of *Streptomyces viridochromogenes* NRRL B-1551. *J. Antibiot. (Tokyo)* **46(7)**, 1126-1138 (1993).
- Aoki, Y., Matsumoto, D., Kawaide, H., et al. Physiological role of germicidins in spore germination and hyphal elongation in *Streptomyces coelicolor* A3(2). *J. Antibiot. (Tokyo)* **64(9)**, 607-611 (2016).

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, 12/14/2022

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