

PRODUCT INFORMATION



Cytochalasin C

Item No. 11329

CAS Registry No.: 22144-76-9
Formal Name: (3S,3aR,6S,6aR,7E,10S,12R,13E,15R,15aR)-15-(acetyloxy)-3,3a,6,6a,9,10,12,15-octahydro-6,12-dihydroxy-4,5,10,12-tetramethyl-3-(phenylmethyl)-1H-cycloundec[d]isoindole-1,11(2H)-dione

MF: C₃₀H₃₇NO₆

FW: 507.6

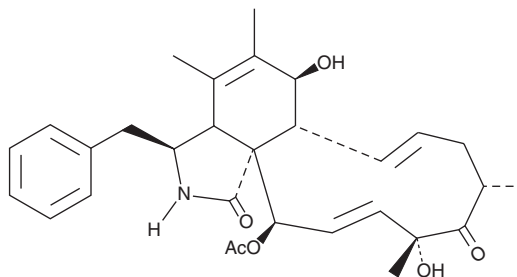
Purity: ≥99%

Supplied as: A powder

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Geniculosporium* sp.



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

Laboratory Procedures

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

Description

Cytochalasin C is an actin polymerization inhibitor and fungal metabolite that has been found in *C. globosum*.^{1,2} It inhibits actin polymerization when used at a concentration of 10 μM.¹ Cytochalasin C is cytotoxic to HeLa cells (IC₅₀ = <0.32 μg/ml).²

References

1. Walling, E.A., Krafft, G.A., and Ware, B.R. Actin assembly activity of cytochalasins and cytochalasin analogs assayed using fluorescence photobleaching recovery. *Arch. Biochem. Biophys.* **264**(1), 321-332 (1988).
2. Umeda, M., Ohtsubo, K., Saito, M., et al. Cytotoxicity of new cytochalasins from *Chaetomium globosum*. *Experientia* **31**(4), 435-438 (1975).

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, 08/11/2023

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