

PRODUCT INFORMATION



Cytochalasin Z2

Item No. 44761

CAS Registry No.: 450398-31-9

Formal Name: (3E,5R,9R,11E,12aS,15S,15aS,16S,18aR)-6,7,8,9,10,12a,15,15a,16,17-decahydro-5-hydroxy-14-(hydroxymethyl)-9,15-dimethyl-16-(phenylmethyl)-2H-oxacyclotetradecino[2,3-d]isoindole-2,18(5H)-dione

MF: $C_{29}H_{37}NO_5$

FW: 479.6

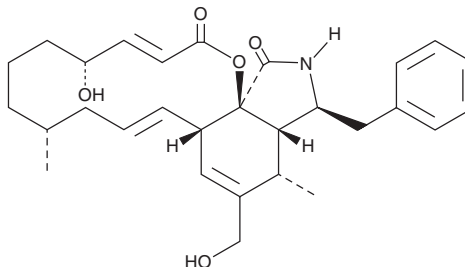
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungi/*Drechslera* sp.



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

Laboratory Procedures

Cytochalasin Z2 is supplied as a solid. A stock solution may be made by dissolving the cytochalasin Z2 in the solvent of choice, which should be purged with an inert gas. Cytochalasin Z2 is soluble (≥ 10 mg/ml) in methanol and DMSO.

Description

Cytochalasin Z2 is a fungal metabolite that has been found in *P. semeniperda* and has anticancer activity.¹ It inhibits the growth of OE21 esophageal squamous cell carcinoma, U373 MG glioblastoma, A549 lung cancer, B16/F10 melanoma, and Hs 683 glioma cancer cells (IC_{50} s = 70, 86, 99, 29, and 94 μM , respectively).

Reference

1. Van Goietsenoven, G., Mathieu, V., Andolfi, A., *et al.* *In vitro* growth inhibitory effects of cytochalasins and derivatives in cancer cells. *Planta Med.* **77**(7), 711-717 (2011).

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, 11/21/2025

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