

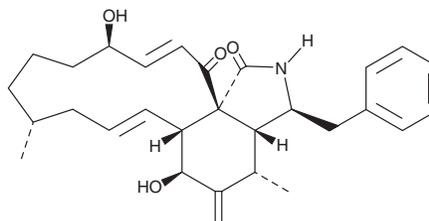
PRODUCT INFORMATION



Deoxaphomin

Item No. 44753

CAS Registry No.: 51053-39-5
Formal Name: (3S,3aR,4S,6S,6aR,7E,10R,14R,15E,17aR)-3,3a,4,5,6,6a,9,10,11,12,13,14-dodecahydro-6,14-dihydroxy-4,10-dimethyl-5-methylene-3-(phenylmethyl)-1H-Ccclotridec[d]isoindole-1,17(2H)-dione
MF: C₂₉H₃₇NO₄
FW: 463.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Drechslera* sp.



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

Laboratory Procedures

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

Description

Deoxaphomin is a polyketide synthase-derived cytochalasin that has been found in *P. semeniperda* and has diverse biological activities.¹⁻⁴ It inhibits anti-mouse IgG-induced capping in isolated mouse splenic lymphocytes by 70% when used at a concentration of 20 μM and inhibits actin filament elongation in C3H fibroblasts by 66% at 2 μM.³ Deoxaphomin (100 μM) inhibits seed germination and radicle growth induced by GR24 (Item No. 13210) in the broomrapes *O. crenata* and *P. ramosa*.² It inhibits the growth of OE21, U373 MG, SK-MEL-28, A549, B16/F10, and Hs 683 cancer cells (IC₅₀s = 0.75, 1.7, 5.9, 2.2, and 0.61 μM, respectively).⁴

References

1. Steyn, P.S. The biosynthesis of polyketide-derived mycotoxins. *J. Environ. Pathol. Toxicol. Oncol.* **11(1)**, 47-59 (1992).
2. Cimmino, A., Fernández-Aparicio, M., Andolfi, A., *et al.* Effect of fungal and plant metabolites on broomrapes (*Orobanche* and *Phelipanche* spp.) seed germination and radicle growth. *J. Agric. Food Chem.* **62(43)**, 10485-10492 (2014).
3. Yahara, I., Harada, F., Sekita, S., *et al.* Correlation between effects of 24 different cytochalasins on cellular structures and cellular events and those on actin in vitro. *J. Cell Biol.* **92(1)**, 69-78 (1982).
4. 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, 12/02/2025

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