

PRODUCT INFORMATION



Harzianum A

Item No. 27740

CAS Registry No.: 156250-74-7

Formal Name: (4 β)-4-[hydrogen (2Z,4E,6E)-2,4,6-octatrienedioate]12,13-epoxy-trichothec-9-en-4-ol

MF: C₂₃H₂₈O₆

FW: 400.5

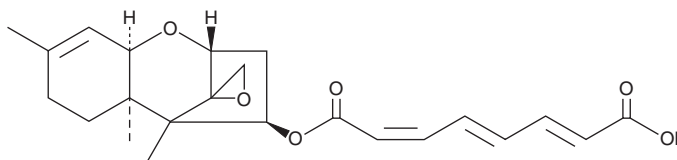
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/*Trichoderma* sp.



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

Laboratory Procedures

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

Description

Harzianum A is a trichothecene fungal metabolite originally isolated from *T. harzianum*.¹ It is cytotoxic to HeLa, MCF-7, and HT-1080 cells (IC₅₀s = 5.07, 10.13, and 0.65 $\mu\text{g}/\text{ml}$, respectively) and exhibits antifungal activity against *C. albicans* and *S. cerevisiae* when used at a concentration of 100 $\mu\text{g}/\text{ml}$.^{1,2}

References

1. Corley, D.G., Miller-Wideman, M., and Durley, R.C. Isolation and structure of harzianum A: A new trichothecene from *Trichoderma harzianum*. *J. Nat. Prod.* **57**(3), 422-425 (1994).
2. Lee, H.B., Kim, Y., Jin, H.Z., et al. A new Hypocrea strain producing harzianum A cytotoxic to tumour cell lines. *Lett. Appl. Microbiol.* **40**(6), 497-503 (2005).

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, 08/10/2023

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