

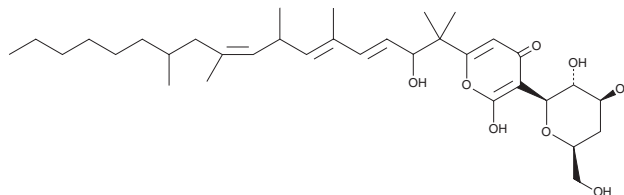
# PRODUCT INFORMATION



## Deoxyfusapyrone

Item No. 27963

**CAS Registry No.:** 156856-32-5  
**Formal Name:** 3-(4-deoxy- $\beta$ -xylo-hexopyranosyl)-2-hydroxy-6-  
[(3E,5E,8Z)-2-hydroxy-1,1,5,7,9,11-hexamethyl-  
3,5,8-heptadecatrien-1-yl]-4H-pyran-4-one  
**MF:** C<sub>34</sub>H<sub>54</sub>O<sub>8</sub>  
**FW:** 590.8  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Fungus/*Fusarium* sp.



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

### Laboratory Procedures

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

### Description

Deoxyfusapyrone is an  $\alpha$ -pyrone fungal metabolite originally isolated from *F. semitectum* and has antifungal activity.<sup>1</sup> It is active against *C. neoformans*, *A. fumigatus*, *A. niger*, and *A. flavus* human mycoses (MICs = 1.56-6.25  $\mu$ g/ml). Deoxyfusapyrone is also active against a variety of filamentous fungi, but not yeast or the bacterium *B. megaterium*, in a disc assay.

### Reference

1. Altomare, C., Perrone, G., Zonno, M.C., *et al.* Biological characterization of fusapyrone and deoxyfusapyrone, two bioactive secondary metabolites of *Fusarium semitectum*. *J. Nat. Prod.* **63(8)**, 1131-1135 (2000).

#### 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