

# PRODUCT INFORMATION



## Fusapyrone

Item No. 20751

CAS Registry No.: 156856-31-4

Formal Name: 3-(4-deoxy- $\beta$ -xylo-hexopyranosyl)-2-hydroxy-6-[(3E,5E,8Z)-2-hydroxy-7-(hydroxymethyl)-1,1,5,9,11-pentamethyl-3,5,8-heptadecatrien-1-yl]-2H-pyran-2-one

MF: C<sub>34</sub>H<sub>54</sub>O<sub>9</sub>

FW: 606.8

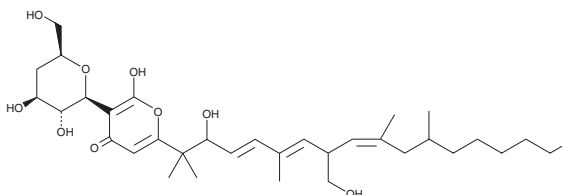
Purity:  $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability:  $\geq 2$  years

Item Origin: Fungi/*Fusarium* sp.



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

### Laboratory Procedures

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

### Description

Fusapyrone is a broad-spectrum antifungal metabolite first isolated from *Fusarium* species. It has been investigated for use in the control of postharvest crop diseases such as inhibiting the growth of ochratoxin-producing strains of *Aspergillus* section *Nigri* in wine grapes.<sup>1</sup>

### Reference

1. Favilla, M., *et al.* Inhibition of species of the *Aspergillus* section *Nigri* and ochratoxin a production in grapes by fusapyrone. *Appl. Environ. Microbiol.* **74**(7), 2248-25253 (2008).

#### 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, 02/13/2019

#### 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