

PRODUCT INFORMATION



Lateropyrone

Item No. 26463

CAS Registry No.: 93752-78-4
Formal Name: 5,9-dihydroxy-2-methyl-4,6-dioxo-4H,6H-pyrano[3',4':4,5]benzo[1,2-b]pyran-8-carboxylic acid, methyl ester

Synonym: Avenacein Y

MF: C₁₅H₁₀O₈

FW: 318.2

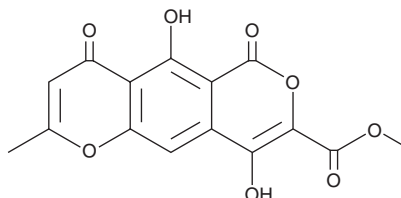
Purity: ≥99%

Supplied as: A solid

Storage: -20°C

Stability: ≥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

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

Description

Lateropyrone is a fungal metabolite produced by *Fusarium* species.¹ It has been found in freshly harvested and stored sugar beet crops in Europe.

Reference

1. Christ, D.S., Märländer, B., and Varrelmann, M. Characterization and mycotoxigenic potential of *Fusarium* species in freshly harvested and stored sugar beet in Europe. *Phytopathology* **101(11)**, 1330-1337 (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, 08/23/2022

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