

PRODUCT INFORMATION



Viomellein

Item No. 43200

CAS Registry No.: 55625-78-0

Formal Name: (3R,3'R)-3,3',4,4'-tetrahydro-9',10,10'-trihydroxy-7,7'-dimethoxy-3,3'-dimethyl-[8,8'-bi-1H-naphtho[2,3-c]pyran]-1,1',6,9-tetrone

MF: $C_{30}H_{24}O_{11}$

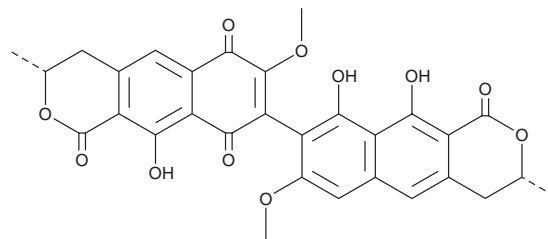
FW: 560.5

Purity: $\geq 70\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Viomellein is a mycotoxin that has been found in *P. viridicatum*.¹ Dietary administration of viomellein (456 mg/kg in feed) induces necrotizing cholangitis in mice. Viomellein is also active against *S. aureus* and *C. albicans* (MICs = 0.78 and 6.25 $\mu\text{g/ml}$).²

References

1. Carlton, W.W., Stack, M.E., and Eppley, R.M. Hepatic alterations produced in mice by xanthomegnin and viomellein, metabolites of *Penicillium viridicatum*. *Toxicol. Appl. Pharmacol.* **38**(2), 455-459 (1976).
2. Ninomiya, A., Masuda, K., Yamada, T., *et al.* Rediscovery of viomellein as an antibacterial compound and identification of its biosynthetic gene cluster in dermatophytes. *Appl. Environ. Microb.* **Online ahead of print**, e0243124 (2025).

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, 04/30/2025

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