

PRODUCT INFORMATION



Filipin III

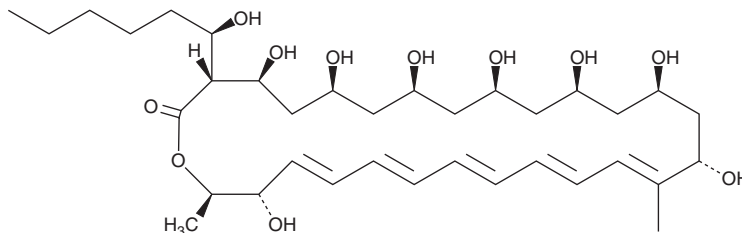
Item No. 70440

CAS Registry No.: 480-49-9
Formal Name: 4S,6S,8S,10R,12R,14R,16S,27S-octahydroxy-3R-(1R-hydroxy-hexyl)17,28R-dimethyl-oxacyclooctacos-17E,19E,21E,21E,23E,25E-pentaen-2-one

MF: $C_{35}H_{58}O_{11}$
FW: 654.8
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 323, 339, 357 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years

Special Conditions: Light sensitive. Do not expose to direct intense light.

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



Laboratory Procedures

Filipin III is supplied as a crystalline solid. A stock solution may be made by dissolving the Filipin III in an organic solvent purged with an inert gas. Filipin III is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Filipin III in these solvents is approximately 2, 5, and 10 mg/ml, respectively. We recommend using the stock solution within 24 hours or it may result in reduced activity.

Filipin III is sparingly soluble in aqueous buffers. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. For maximum solubility in aqueous buffers, Filipin III should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Filipin III has a solubility of approximately 0.4 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Filipin III is a polyene and an active component of the polyene antibiotic filipin complex (Item No. 25073).¹ It is active against various fungi, including *B. dermatitidis*, *C. neoformans*, *H. capsulatum*, *C. albicans*, and *T. mentagrophytes* (MICs = 1-10 μ g/ml). Filipin III induces hemolysis of isolated rabbit erythrocytes (EC_{50} = 0.8 μ g/ml).² It induces apoptosis in HaCaT keratinocytes in a concentration-dependent manner.³ Filipin III binds to cholesterol and has been used to fluorescently label sterols in biological structures.^{4,5} The filipin complex has excitation maxima of 338 and 357 nm and an emission maxima of 480 nm.⁶

References

1. Bergy, M.E. and Eble, T.E. *Biochemistry* **7**(2), 653-659 (1968)
2. Sessa, G. and Weissmann, G. *J. Bio. Chem.* **243**(16), 4364-4371 (1968).
3. Gniadecki, R. *Biochem. Biophys. Res. Commun.* **320**(1), 165-169 (2004).
4. Bittman, R. and Fischkoff, S.A. *Proc. Natl. Acad. Sci. USA* **69**(12), 3795-3799 (1972).
5. Boutté, Y., Men, S., and Grebe, M. *Nat. Protoc.* **6**(4), 446-456 (2011).
6. Drabikowski, W., Lagwinska, E., and Sarzala, M.G. *Biochim. Biophys. Acta* **291**, 61-70 (1973)

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, 04/22/2025

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