

PRODUCT INFORMATION



Filipin II

Item No. 25521

CAS Registry No.: 38620-77-8
Formal Name: 4S,6S,8S,10R,12R,14R,16S,27S-octahydroxy-3R-hexyl-17,28R-dimethyl-oxacyclooctacos-17E,19E,21E,23E,25E-pentaen-2-one

MF: C₃₅H₅₈O₁₀

FW: 638.8

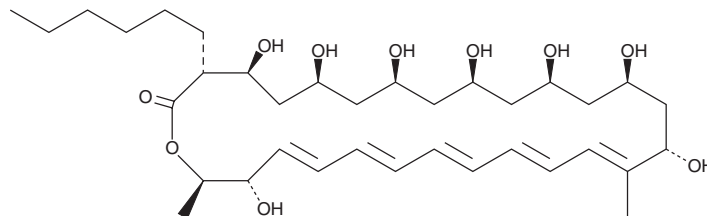
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item origin: Bacteria/*Streptomyces* sp.



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

Laboratory Procedures

Filipin II is supplied as a solid. A stock solution may be made by dissolving the filipin II in the solvent of choice. Filipin II is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Filipin II is a polyene and an active component of the polyene antibiotic filipin complex (Item No. 25073).¹ It has broad-spectrum antifungal activity, reducing growth of various fungi, including *B. dermatitidis*, *C. neoformans*, and *H. capsulatum* (MICs = 1-10 µg/ml). Filipin II inhibits ADP-stimulated respiration in rat liver mitochondria when used at a concentration of 3.5 mg/ml.² It also induces hemolysis of liposomes in a cholesterol-independent manner (IC₅₀ = 0.5 µg/ml).³

References

1. Bergy, M.E. and Eble, T.E. The filipin complex. *Biochemistry* **7**(2), 653-659 (1968).
2. Balcavage, W.X., Beale, M., Chasen, B., et al. Effect of filipin on rat-liver and yeast mitochondria. *Biochim. Biophys. Acta.* **162**(4), 525-532 (1968).
3. Sessa, G. and Weissmann, G. Effects of four components of the polyene antibiotic, filipin, on phospholipid spherules (liposomes) and erythrocytes. *J. Biol. Chem.* **243**(16), 4364-4371 (1968).

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, 12/13/2022

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