

PRODUCT INFORMATION



Lagosin

Item No. 20600

CAS Registry No.: 6834-98-6
Formal Name: (3R,4S,6S,8S,10R,12R,14R,15R,16R,17E,19E,21E,23E,25E,27S,28R)-4,6,8,10,12,14,15,16,27-nonahydroxy-3-[(1R)-1-hydroxyhexyl]-17,28-dimethyl-oxacyclooctacos-17,19,21,23,25-pentaen-2-one
Synonyms: Cogomycin, 14-hydroxy Filipin, Fungichromin, NSC 105388, NSC 276732, Pentamycin

MF: C₃₅H₅₈O₁₂

FW: 670.8

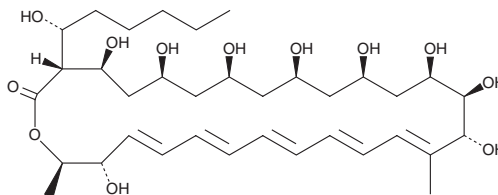
Purity: ≥85%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

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

Description

Lagosin is a polyene macrolide that has been found in *Streptomyces* and has antimicrobial activity.¹⁻³ It is active against *C. albicans*, as well as the plant pathogenic fungus *R. solani* (MICs = 10 and 72 µg/ml, respectively).^{1,2} Lagosin is also active against several strains of *T. vaginalis* that are sensitive or resistant to the antibiotic metronidazole (Item No. 9002409; EC₅₀s = 1.74-2.62 µg/ml after a one-hour incubation).³

References

1. An, D., Wang, X., Li, J., *et al.* The activity of fungichromin against the formation of *Candida albicans* biofilm. *Biol. Pharm. Bull.* **39**(12), 1948-1954 (2016).
2. Shih, H.-D., Liu, Y.-C., Hsu, F.-L., *et al.* Fungichromin: A substance from *Streptomyces padanus* with inhibitory effects on *Rhizoctonia solani*. *J. Agric. Food Chem.* **51**(1), 95-99 (2003).
3. Kranzler, M., Syrowatka, M., Leitsch, D., *et al.* Pentamycin shows high efficacy against *Trichomonas vaginalis*. *Int. J. Antimicrob. Agents* **45**(4), 434-437 (2015).

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, 01/17/2023

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