

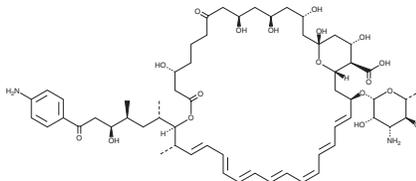
PRODUCT INFORMATION



Levorin Complex

Item No. 26708

CAS Registry No.: 1403-17-4
MF: $C_{59}H_{86}N_2O_{18}$ (for Levorin A₃)
FW: 1,111.3
UV/Vis.: λ_{max} : 360, 380, 403 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Bacterium/*Streptomyces griseus*



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

Laboratory Procedures

Levorin complex is supplied as a solid. A stock solution may be made by dissolving the levorin complex in the solvent of choice, which should be purged with an inert gas. Levorin complex is soluble in organic solvent DMSO (warmed).

Description

Levorin complex is a macrocyclic heptaene antifungal complex originally isolated from *S. griseus*.¹ It is active against *C. albicans*, *T. glabrata*, *A. niger*, *T. mentagrophytes*, and *T. soudanense in vitro* (MICs = 0.03-25 $\mu\text{g/ml}$).² Oral administration of levorin complex reduces mortality in a mouse model of *C. albicans* infection with a protective dose (PD₅₀) value of 25 mg/kg. This product is a mixture of levorin A₀, levorin A₂, and levorin A₃.

References

1. Lechevalier, H., Acker, R.F., Corke, C.T., *et al.* Candicidin, a new antifungal antibiotic. *Mycologia* **45(2)**, 155-171 (1953).
2. Wagman, G.H., Testa, R.T., Patel, M., *et al.* New polyene antifungal antibiotic produced by a species of *Actinoplanes*. *Antimicrob. Agents Chemother.* **7(4)**, 457-461 (1975).

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, 03/23/2023

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