

PRODUCT INFORMATION



Elaiophylin

Item No. 15583

CAS Registry No.: 37318-06-2

Synonyms: Azalomycin-B, Gopalamicin,
Salbomycin

MF: $C_{54}H_{88}O_{18}$

FW: 1,025.3

Purity: $\geq 95\%$

UV/Vis.: λ_{\max} : 253 nm

Supplied as: A neat solid

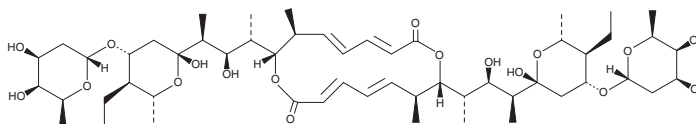
Storage: -20°C

Stability: ≥ 2 years

Special Conditions: Protect from light when in solution.

Item Origin: Bacterium/*Streptomyces hygroscopicus*

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



Laboratory Procedures

Elaiophylin is supplied as a neat solid. A stock solution may be made by dissolving the elaiophylin in the solvent of choice, which should be purged with an inert gas. Elaiophylin is soluble in organic solvents such as ethanol and DMSO. The solubility of elaiophylin in ethanol with sonication is approximately 1 mg/ml and approximately 20 mg/ml in DMSO.

Description

Elaiophylin is a macrodiolide antibiotic that can be isolated from various strains of *Streptomyces*.^{1,2} It displays *in vitro* anti-protozoal activity against both *Plasmodium* and *Trypanosoma* ($\text{IC}_{50}\text{s} = 370$ and 460 ng/ml, respectively) and cytotoxicity against human fetal lung fibroblast MRC-5 cells ($\text{IC}_{50} = 870$ ng/ml).³ Elaiophylin alone has no activity against *Candida*, although it enhances the anti-fungal activity of rapamycin (Item No. 13346).² Elaiophylin also forms stable, long-lasting ion channels in bilayer membranes that are selective for cations.⁴

References

1. Fiedler, H.P., Wörner, W., Zöhner, H., *et al.* Metabolic products of microorganisms. 200 Isolation and characterization of niphithricins A, B, and elaiophylin, antibiotics produced by *Streptomyces violaceoniger*. *J. Antibiot. (Tokyo)* **34**(9), 1107-1118 (1981).
2. Fang, A., Wong, G.K., and Demain, A.L. Enhancement of the antifungal activity of rapamycin by the coproduced elaiophylin and nigericin. *J. Antibiot. (Tokyo)* **53**(2), 158-162 (2000).
3. Otoguro, K., Iwatsuki, M., Ishiyama, A., *et al.* *In vitro* and *in vivo* antiprotozoal activities of bispolides and their derivatives. *J. Antibiot. (Tokyo)* **63**(5), 275-277 (2010).
4. Grigoriev, P.A., Schlegel, R., and Gräfe, U. Cation selective ion channels formed by macrodiolide antibiotic elaiophylin in lipid bilayer membranes. *Bioelectrochemistry* **54**(1), 11-15 (2001).

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, 03/02/2022

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