

PRODUCT INFORMATION

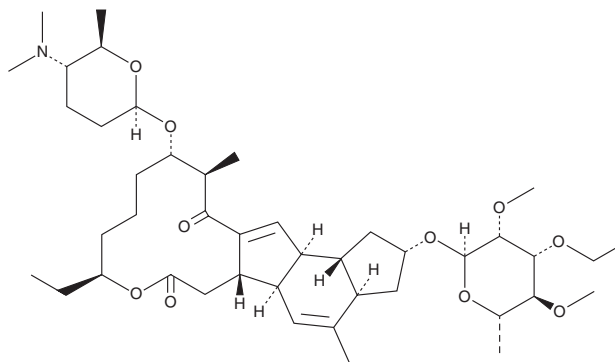


3'-Ethoxy Spinosyn L

Item No. 26664

CAS Registry No.: 187166-15-0
Formal Name: (2S,3aR,5aS,5bS,9S,13S,14R,16aS,16bS)-2-[(6-deoxy-3-O-ethyl-2,4-di-O-methyl- α -L-mannopyranosyl)oxy]-13-[[[(2R,5S,6R)-5-(dimethylamino)tetrahydro-6-methyl-2H-pyran-2-yl]oxy]-9-ethyl-2,3,3a,5a,5b,6,9,10,11,12,13,14,16a,16b-tetradecahydro-4,14-dimethyl-1H-as-indaceno[3,2-d]oxacyclododecin-7,15-dione

Synonym: Spinetoram L
MF: C₄₃H₆₉NO₁₀
FW: 760.0
Purity: \geq 95%
UV/Vis.: λ_{\max} : 245 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

3'-Ethoxy spinosyn L is supplied as a solid. A stock solution may be made by dissolving the 3'-ethoxy spinosyn L in the solvent of choice, which should be purged with an inert gas. 3'-Ethoxy spinosyn L is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

3'-Ethoxy spinosyn L is a minor component of the insecticide spinetoram (Item No. 26623).¹ Formulations containing 3'-ethoxy spinosyn L have been used as insecticides.

Reference

1. Sparks, T.C., Crouse, G.D., Dripps, J.E., *et al.* Neural network-based QSAR and insecticide discovery: Spinetoram. *J. Comput. Aided Mol. Des.* **22**(6-7), 393-401 (2008).

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, 07/26/2023

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