

PRODUCT INFORMATION



Spinosyn D

Item No. 19158

CAS Registry No.: 131929-63-0

Formal Name: (2S,3aR,5aS,5bS,9S,13S,14R,16aS,16bS)-2-[(6-deoxy-2,3,4-tri-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

Synonyms

A-83543D, Lepicidin D

MF: $C_{42}H_{67}NO_{10}$

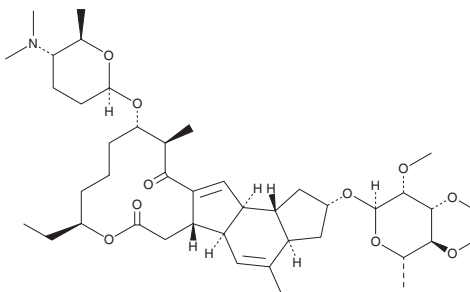
FW: 746.0

Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Spinosyn D is a spinosoid that is a minor component of spinosad, an insecticide based on chemical compounds found in a bacterial species isolated from sugarcane, *S. spinosa*.¹ It acts as an agonist of insect nicotinic acetylcholinesterase receptors.¹ Spinosyn D demonstrates insecticidal activity against *H. virescens* larvae (tobacco budworm) with an LD_{50} value of 0.8 ppm.²

References

1. Duke, S.O., Cantrell, C.L., Meepagala, K.M., *et al.* Natural toxins for use in pest management. *Toxins (Basel)* **2**(8), 1943-1962 (2010).
2. Creemer, L.C., Kirst, H.A., Paschal, J.W., *et al.* Synthesis and insecticidal activity of spinosyn analogs functionally altered at the 2'-,3'- and 4'-positions of the rhamnose moiety. *J. Antibiot. (Tokyo)* **53**(2), 171-178 (2000).

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, 07/13/2023

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