

PRODUCT INFORMATION



Spinosyn A 17-pseudoaglycone

Item No. 25530

CAS Registry No.: 131929-68-5

Formal Name: (2R,3aS,5aR,5bS,9S,13S,14R,16aS,16bR)-2-[(6-deoxy-2,3,4-tri-O-methyl- α -L-mannopyranosyl)oxy]-9-ethyl-2,3,3a,5a,5b,6,9,10,11,12,13,14,16a,16b-tetradecahydro-13-hydroxy-14-methyl-1H-as-indaceno[3,2-d]oxacyclododecin-7,15-dione

MF: $C_{33}H_{50}O_9$

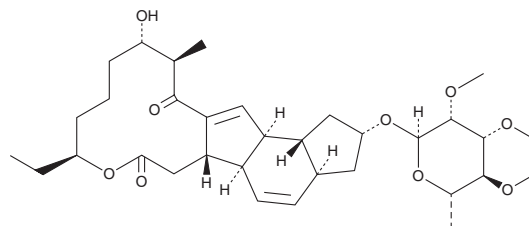
FW: 590.7

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 A 17-pseudoaglycone is supplied as a solid. A stock solution may be made by dissolving the spinosyn A 17-pseudoaglycone in the solvent of choice, which should be purged with an inert gas. Spinosyn A 17-pseudoaglycone is soluble in methanol, ethanol, DMSO, and dimethyl formamide.

Description

Spinosyn A 17-pseudoaglycone is a pseudoaglycone form of the macrocyclic lactone insecticide spinosyn A (Item No. 16528).¹ Unlike spinosyn A, spinosyn A 17-pseudoaglycone is not lethal to tobacco budworm (*H. virescens*) neonate larvae ($LC_{50} = >64$ ppm).

Reference

1. Sparks, T.C., Thompson, G.D., Kirst, H.A., *et al.* Biological activity of the spinosyns, new fermentation derived insect control agents, on tobacco budworm (lepidoptera: noctuidae) larvae. *J. Econ. Entomol.* **91**(6), 1227-1283 (1998).

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, 12/20/2022

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