

PRODUCT INFORMATION

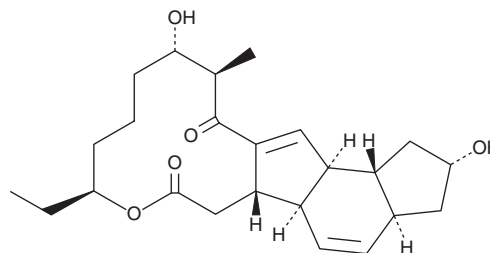


Spinosyn A aglycone

Item No. 27992

CAS Registry No.: 149560-97-4
Formal Name: (2R,3aS,5aR,5bS,9S,13S,14R,16aS,16bR)-9-ethyl-2,3,3a,5a,5b,6,9,10,11,12,13,14,16a,16b-tetradecahydro-2,13-dihydroxy-14-methyl-1H-as-indaceno[3,2-d]oxacyclododecin-7,15-dione

MF: C₂₄H₃₄O₅
FW: 402.5
Purity: ≥95%
UV/Vis.: λ_{max}: 244 nm
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 aglycone is supplied as a solid. A stock solution may be made by dissolving the spinosyn A aglycone in the solvent of choice, which should be purged with an inert gas. Spinosyn A aglycone is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Spinosyn A aglycone is an aglycone form of the insecticide spinosyn A.¹ It is formed by acid hydrolysis of spinosyn A 17-pseudoaglycone (Item No. 25530). Spinosyn A aglycone has no insecticidal activity at concentrations up to 64 ppm.

Reference

1. Creemer, L.C., Kirst, H.A., and Paschal, J.W. Conversion of spinosyn A and spinosyn D to their respective 9- and 17-pseudoaglycones and their aglycones. *J. Antibiot. (Tokyo)* **51**(8), 795-800 (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, 12/19/2022

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