

PRODUCT INFORMATION

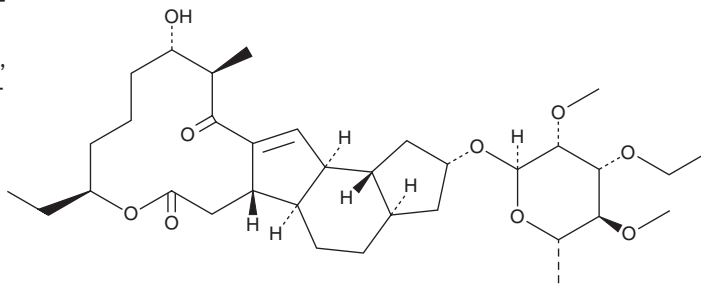


3'-Ethoxy-5,6-dihydro Spinosyn J 17-pseudoaglycone

Item No. 25765

CAS Registry No.: 2055494-09-0
Formal Name: 2-[[6-deoxy-3-O-ethyl-2,4-di-O-methyl- α -L-mannopyranosyl)oxy]-9-ethyl-2R,3,3aR,4,5,5aR,5bS,6,9S,10,11,12,13S,14R,16aS,16bR-hexadecahydro-13-hydroxy-14-methyl-1H-as-indaceno[3,2-d]oxacyclododecin-7,15-dione

MF: C₃₄H₅₄O₉
FW: 606.8
Purity: \geq 95%
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-5,6-dihydro spinosyn J 17-pseudoaglycone is supplied as a solid. A stock solution may be made by dissolving the 3'-ethoxy-5,6-dihydro spinosyn J 17-pseudoaglycone in the solvent of choice, which should be purged with an inert gas. 3'-Ethoxy-5,6-dihydro spinosyn J 17-pseudoaglycone is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

3'-Ethoxy-5,6-dihydro spinosyn J 17-pseudoaglycone is a degradation product of 3'-ethoxy-5,6-dihydro spinosyn J, a component of the insecticide spinosyn J.^{1,2} It has no insecticidal activity against the tobacco budworm up to a concentration of 64 ppm.²

References

1. Kirst, H.A. The spinosyn family of insecticides: Realizing the potential of natural products research. *J. Antibiot. (Tokyo)* **63(3)**, 101-111 (2010).
2. 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, 12/20/2022

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