

PRODUCT INFORMATION

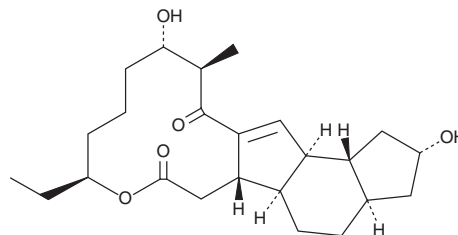


Dihydrospinosyn A aglycone

Item No. 23641

CAS Registry No.: 727695-12-7
Formal Name: (2R,3aR,5aR,5bS,9S,13S,14R,16aS,16bR)-9-ethyl-2,3,3a,4,5,5a,5b,6,9,10,11,12,13,14,16a,16b-hexadecahydro-2,13-dihydroxy-14-methyl-1H-as-indaceno[3,2-d]oxacyclododecin-7,15-dione

MF: C₂₄H₃₆O₅
FW: 404.5
Purity: ≥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

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

Description

Dihydrospinosyn A aglycone is a derivative of spinosyn A aglycone with a reduced double bond at positions 5 and 6.

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, 12/17/2022

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