

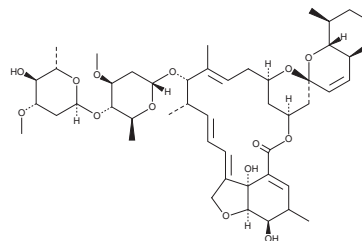
PRODUCT INFORMATION



Δ^2 -Avermectin B_{1a}

Item No. 25119

MF: C₄₈H₇₂O₁₄
FW: 873.1
Purity: ≥95%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A crystalline 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

Δ^2 -Avermectin B_{1a} is supplied as a solid. A stock solution may be made by dissolving the Δ^2 -Avermectin B_{1a} in the solvent of choice, which should be purged with an inert gas. Δ^2 -Avermectin B_{1a} is soluble in organic solvent DMSO.

Description

Δ^2 -Avermectin B_{1a} is a degradation product formed by the reversible base-catalyzed isomerization of ivermectin B_{1a} (Item No. 18768).¹ It is less active than ivermectin against *T. urticae* (LC₉₀s = 0.23 and 0.038 ppm, respectively).

Reference

1. Pivnichny, J.V., Arison, B.H., Preiser, F.A., et al. Base-catalyzed isomerization of avermectins. *J. Agric. Food Chem.* **36**(4), 826-828 (1988).

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, 11/30/2022

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