

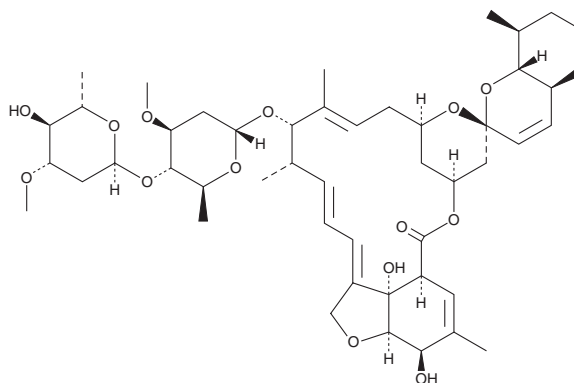
PRODUCT INFORMATION



2-*epi*-Abamectin

Item No. 27993

CAS Registry No.: 106434-14-4
Formal Name: (2S)-5-O-demethyl-avermectin A_{1a}
Synonym: *epi*-Avermectin B_{1a}
MF: C₄₈H₇₂O₁₄
FW: 873.1
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Semi-synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-*epi*-Abamectin is supplied as a solid. A stock solution may be made by dissolving the 2-*epi*-abamectin in the solvent of choice, which should be purged with an inert gas. 2-*epi*-Abamectin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

2-*epi*-Abamectin is a degradation product of abamectin (Item No. 19201).¹ It is toxic to the two-spotted spider mite in a contact assay with an LC₅₀ value of 4 ppm, which is approximately 100-fold less potent than abamectin.²

References

1. Awasthi, A., Razzak, M., Al-Kassas, R., *et al.* Separation and identification of degradation products in abamectin formulation using LC, LTQ FT-MS, H/D exchange and NMR. *Curr. Pharm. Anal.* **8**(4), 415-430 (2012).
2. 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, 12/08/2022

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