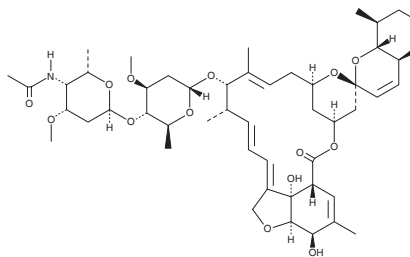


PRODUCT INFORMATION



Eprinomectin B_{1a} Item No. 28130

CAS Registry No.: 133305-88-1
Formal Name: (4''R)-4''-(acetylamino)-5-O-demethyl-4''-deoxy-avermectin A_{1a}
MF: C₅₀H₇₅NO₁₄
FW: 914.1
Purity: ≥99%
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

Eprinomectin B_{1a} is supplied as a solid. A stock solution may be made by dissolving the eprinomectin B_{1a} in the solvent of choice, which should be purged with an inert gas. Eprinomectin B_{1a} is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Eprinomectin B_{1a} is a major component (>90%) of the antiparasitic compound eprinomectin, which also contains eprinomectin B_{1b} (<10%; Item No. 27994) and belongs to the avermectin family of insecticides and anthelmintics.¹

Reference

1. Kvaternick, V., Kellermann, M., Knaus, M., *et al.* Pharmacokinetics and metabolism of eprinomectin in cats when administered in a novel topical combination of fipronil, (S)-methoprene, eprinomectin and praziquantel. *Vet. Parasitol.* **202(1-2)**, 2-9 (2014).

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, 11/23/2022

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