

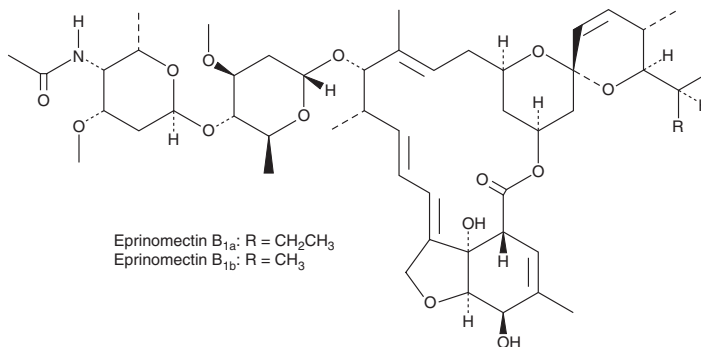
PRODUCT INFORMATION



Eprinomectin

Item No. 28182

CAS Registry No.: 123997-26-2
Formal Name: (4''R)-4''-(acetylamino)-4''-deoxy-avermectin B₁
Synonym: MK-397
MF: C₅₀H₇₅NO₁₄ (for B_{1a})
FW: 914.1 (for B_{1a})
Purity: ≥95% (mixture of B_{1a} and B_{1b})
UV/Vis.: λ_{max}: 246 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

Eprinomectin is supplied as a crystalline solid. A stock solution may be made by dissolving the eprinomectin in the solvent of choice, which should be purged with an inert gas. Eprinomectin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of eprinomectin in these solvents is approximately 30 mg/ml.

Description

Eprinomectin is an avermectin with anthelmintic and insecticidal activities.¹ It is composed of eprinomectin B_{1a} (Item No. 28130) and eprinomectin B_{1b} (Item No. 27994).² Eprinomectin (≥0.08 mg/kg, applied topically) reduces worm burden by 99% or more in bovine models of adult *H. placei*, *O. ostertagi*, *T. axei*, *Cooperia punctata*, *N. helvetianus*, *O. radiatum*, or *D. viviparus* infection.¹ It also reduces parasite burden in bovine models of horn fly (*H. irritans*) or mange mite (*C. bovis*) infestation, as well as in cows naturally infested with lice (*L. vituli*), when applied topically at doses ranging from 0.16 to 0.5 mg/kg.

References

- Shoop, W.L., Egerton, J.R., Eary, C.H., et al. Eprinomectin: A novel avermectin for use as a topical endectocide for cattle. *Int. J. Parasitol.* **26(11)**, 1237-1242 (1996).
- Jiang, X., Hao, H.X., Growney, J.D., et al. Inactivating mutations of *RNF43* confer Wnt dependency in pancreatic ductal adenocarcinoma. *Proc. Natl. Acad. Sci. U.S.A.* **110(31)**, 12649-12654 (2013).

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, 7/25/2023

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