

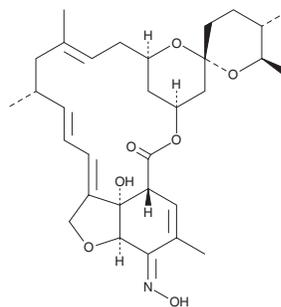
PRODUCT INFORMATION



Milbemycin A₃ oxime

Item No. 27967

CAS Registry No.: 114177-14-9
Formal Name: 5-demethoxy-28-deoxy-6R,28-epoxy-5-(hydroxyimino)-25R-methyl-milbemycin B
MF: C₃₁H₄₃NO₇
FW: 541.7
Purity: ≥95%
UV/Vis.: λ_{max}: 245 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

Milbemycin A₃ oxime is supplied as a crystalline solid. A stock solution may be made by dissolving the milbemycin A₃ oxime in the solvent of choice, which should be purged with an inert gas. Milbemycin A₃ oxime is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of milbemycin A₃ oxime in these solvents is approximately 30 mg/ml.

Description

Milbemycin A₃ oxime is an anthelmintic, a derivative of milbemycin A₃ (Item No. 18360), and a component of milbemycin oxide (Item No. 17164).¹ It reduces the number of microfilariae of the heartworm *D. immitis* in the blood of naturally infested dogs by 85.2% when administered at a dose of 0.05 mg/kg. Formulations containing milbemycin A₃ oxime, in combination with milbemycin A₄ oxime, have been used in the treatment of parasitic infections in dogs.

Reference

1. Tsukamoto, Y., Sato, K., Mio, S., *et al.* Synthesis of 5-keto-5-oxime derivatives of milbemycins and their activities against microfilariae. *Agric. Biol. Chem.* **55(10)**, 2615-2621 (1991).

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, 10/05/2022

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