PRODUCT INFORMATION



Milbemycin D

Item No. 18766

CAS Registry No.:	77855-81-3	
Formal Name:	5-O-demethyl-28-deoxy-6R,28-epoxy-	H O
	25R-(1-methylethyl)-milbemycin B	
Synonym:	Antibiotic B 41D	
MF:	C ₃₃ H ₄₈ O ₇	
FW:	556.7	
Purity:	≥95%	Он
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥2 years	0~ ; \
Item Origin:	Bacterium/Streptomyces hygroscopicus	Ōн

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

Laboratory Procedures

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

Description

Milbemycin D is an insecticidal and anthelmintic antibiotic originally isolated from S. hygroscopicus.¹ It is toxic to third instar gypsy moth larvae with an LD_{50} value of 0.119 μ g/g but not to sixth instar larvae.² Milbemycin D eradicates A. caninum and is effective against T. canis in dogs when administered at doses of 0.05 and 0.1 mg/kg.³ It prevents D. immitis infection in dogs, but is not effective against T. vulpis or D. caninum. Milbemycin D also activates glutamate- (EC₅₀ = 89 nM) and GABA-gated chloride channels.^{4,5}

References

- 1. Okazaki, T., Ono, M., Aoki, A., et al. Milbemycins, a new family of macrolide antibiotics: Producing organism and its mutants. J. Antibiot. (Tokyo) 36(4), 438-441 (1983).
- 2. Deecher, D.C., Brezner, J., and Tananbaum, S.W. Avermectin B1a and milbemycin D as contact toxicants for gypsy moth (Lepidoptera: Lymantriidae) larvae and eggs. J. Econ. Entomol. 80(6), 1284-1287 (1987).
- 3. Tagawa, M., Takiyama, A., Ejima, H., et al. Efficacy of milbemycin D against intestinal helminths of dogs. Nippon Jui Chikusan Daigaku Kenkyu Hokoku 33, 97-99 (1984).
- 4. Arena, J.P., Liu, K.K., Paress, P.S., et al. The mechanism of action of avermectins in Caenorhabditis elegans: Correlation between activation of glutamate-sensitive chloride current, membrane binding, and biological activity. J. Parasitol. 81(2), 286-294 (1995).
- 5. Yamazaki, J., Matsumoto, K., Ono, H., et al. Macrolide compounds, ivermectin and milbemycin D, stimulate chloride channels sensitive to GABAergic drugs in cultured chick spinal neurons. Comp. Biochem. Physiol. C 93(1), 97-104 (1989).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 02/18/2021

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