

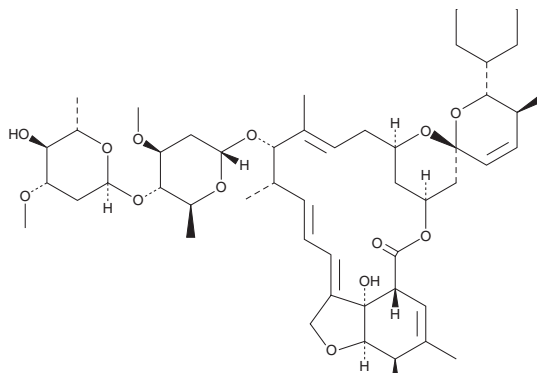
# PRODUCT INFORMATION



## Doramectin

Item No. 19467

**CAS Registry No.:** 117704-25-3  
**Formal Name:** 25-cyclohexyl-5-O-demethyl-25-de(1-methylpropyl)-ivermectin A<sub>1a</sub>  
**Synonym:** Dectomax  
**MF:** C<sub>50</sub>H<sub>74</sub>O<sub>14</sub>  
**FW:** 899.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 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

Doramectin is supplied as a crystalline solid. A stock solution may be made by dissolving the doramectin in the solvent of choice. Doramectin is soluble in chloroform and methanol and is miscible in water.

### Description

Doramectin is a macrocyclic lactone derivative of ivermectin, an anthelmintic that potentiates glutamate- and GABA-gated chloride channel opening in nematodes.<sup>1,2</sup> At 0.001 µg/ml, doramectin is fully effective at inhibiting growth in an *H. contortus* larval development assay.<sup>2</sup>

### References

1. Holden-Dye, L. and Walker, R.J. Anthelmintic drugs and nematicides: Studies in *Caenorhabditis elegans*. *WormBook* **16**, (2014).
2. Michael, B., Meinke, P.T., and Shoop, W. Comparison of ivermectin, doramectin, selamectin, and eleven intermediates in a nematode larval development assay. *J. Parasitol.* **87**(3), 692-696 (2001).

#### 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, 10/06/2022

#### 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