

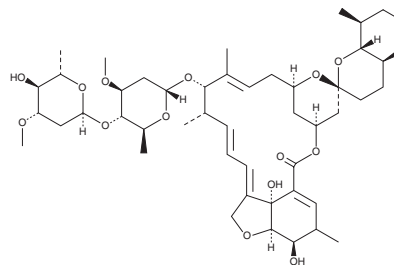
PRODUCT INFORMATION



2,3-Dehydro-3,4-dihydro Ivermectin

Item No. 23849

Formal Name: 2,3-didehydro-5-O-demethyl-3,4,22,23-tetrahydro-avermectin A_{1a}
MF: C₄₈H₇₄O₁₄
FW: 875.1
Purity: ≥95% (mixture of diastereomers)
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2,3-Dehydro-3,4-dihydro ivermectin is supplied as a solid. A stock solution may be made by dissolving the 2,3-dehydro-3,4-dihydro ivermectin in the solvent of choice. 2,3-Dehydro-3,4-dihydro ivermectin is slightly soluble in chloroform and methanol.

Description

2,3-Dehydro-3,4-dihydro ivermectin is an analog of ivermectin (Item No. 18768) and an anthelmintic.¹ It has activity against *L. amazonensis* promastigotes and amastigotes (IC₅₀s = 13.8 and 3.6 μM, respectively) without inducing cytotoxicity to macrophages (IC₅₀ = 65.5 μM).

Reference

1. dos Santos, A.R., Falcão, C.A., Muzitano, M.F., et al. Ivermectin-derived leishmanicidal compounds. *Bioorg. Med. Chem.* **17**(2), 496-502 (2009).

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, 06/22/2020

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