

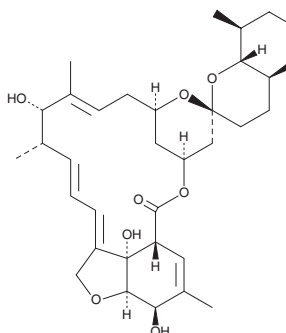
PRODUCT INFORMATION



Ivermectin B_{1a} aglycone

Item No. 19442

CAS Registry No.: 73162-95-5
Formal Name: (6R,13S,25R)-5-O-demethyl-28-deoxy-6,28-epoxy-13-hydroxy-25-[(1S)-1-methylpropyl]-milbemycin B
Synonyms: Avermectin B_{1a} aglycone, 22,23-dihydro Avermectin B_{1a} aglycon
MF: C₃₄H₅₀O₈
FW: 586.8
Purity: ≥99%
Supplied as: A solid
Storage: -80°C
Stability: ≥4 years



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

Laboratory Procedures

Ivermectin B_{1a} aglycone is supplied as a solid. A stock solution may be made by dissolving the ivermectin B_{1a} aglycone in the solvent of choice, which should be purged with an inert gas. Ivermectin B_{1a} aglycone is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Ivermectin B_{1a} aglycone is an acid degradation product produced by hydrolysis of the disaccharide unit of ivermectin. It can inhibit nematode larval development, but does not cause paralytic activity.¹

Reference

1. Chabala, J.C., Mrozik, H., Tolman, R.L., *et al.* Ivermectin, a new broad-spectrum antiparasitic agent. *J. Med. Chem.* **23**(10), 1134-1136 (1980).

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, 12/11/2025

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