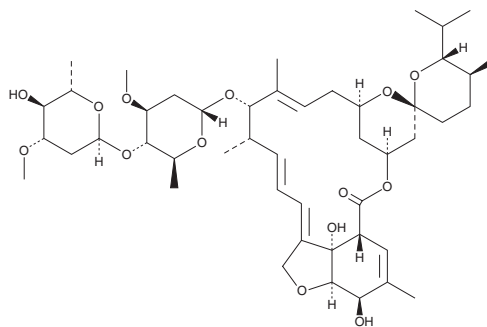


PRODUCT INFORMATION



Ivermectin B_{1b} Item No. 23824

CAS Registry No.: 70209-81-3
Formal Name: 5-O-demethyl-25-de(1-methylpropyl)-22,23-dihydro-25-(1-methylethyl)-avermectin A_{1a}
Synonym: 22,23-dihydro Avermectin B_{1b}
MF: C₄₇H₇₂O₁₄
FW: 861.1
Purity: ≥95%
Supplied as: A 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

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

Description

Ivermectin B_{1b} is the minor component (<20%) of the anthelmintic ivermectin, which is mainly comprised of ivermectin B_{1a} (>80%; Item No. 18768).¹ It produces antiparasitic activity by binding to glutamate-gated chloride channels expressed on nematode neurons and pharyngeal muscle cells, inducing irreversible channel opening and very long-lasting hyperpolarization/depolarization of the neuron/muscle cell, thereby blocking further function.^{2,3} Formulations containing ivermectin inhibit replication of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) in Vero/hSLAM cells.⁴

References

1. Campbell, W.C., Fisher, M.H., Stapley, E.O., *et al.* Ivermectin: A potent new antiparasitic agent. *Science* **221(4613)**, 823-828 (1983).
2. Wolsetholme, A.J. and Rogers, A.T. Glutamate-gated chloride channels and the mode of action of the avermectin/milbemycin anthelmintics. *Parasitology* **131(Suppl)**, S85-S95 (2005).
3. Arena, J.P., Liu, K.K., Pares, 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).
4. Caly, L., Druce, J.D., Catton, M.G., *et al.* The FDA-approved drug ivermectin inhibits the replication of SARS-CoV-2 *in vitro*. *Antiviral Res.* **178:104787** (2020).

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, 10/27/2022

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