

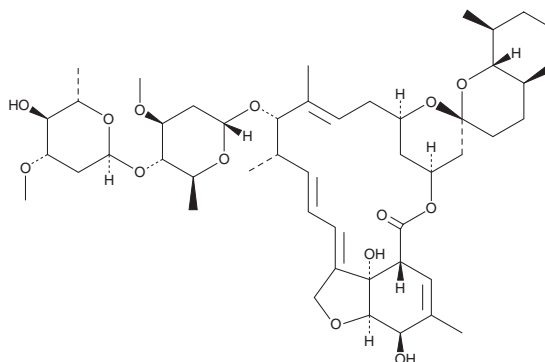
PRODUCT INFORMATION



Ivermectin

Item No. 35518

CAS Registry No.: 70288-86-7
Synonyms: 22,23-dihydro Avermectin B₁,
L-640,471, MK-933
MF: C₄₈H₇₄O₁₄ (for Ivermectin B_{1a})
FW: 875.1
Purity: ≥98%
UV/Vis.: λ_{max}: 245 nm
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 is supplied as a solid. A stock solution may be made by dissolving the ivermectin in the solvent of choice, which should be purged with an inert gas. Ivermectin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ivermectin in ethanol and DMSO is approximately 1 mg/ml and approximately 3 mg/ml in DMF.

Description

Ivermectin is a mixture of the anthelmintics ivermectin B_{1a} (Item No. 18768) and ivermectin B_{1b} (Item No. 23824).¹ Formulations containing ivermectin have been used in the treatment of parasitic roundworm infections.

Reference

1. Campbell, W.C., Fisher, M.H., Stapley, E.O., et al. Ivermectin: A potent new antiparasitic agent. *Science* **221**(4613), 823-828 (1983).

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