

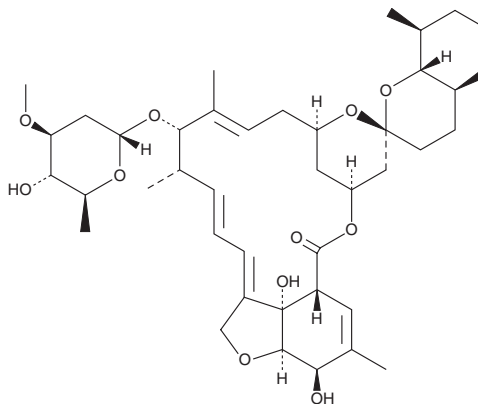
# PRODUCT INFORMATION



## Ivermectin B<sub>1a</sub> monosaccharide

Item No. 19443

**CAS Registry No.:** 71837-27-9  
**Formal Name:** 4'-O-de(2,6-dideoxy-3-O-methyl- $\alpha$ -L-arabino-hexopyranosyl)-5-O-demethyl-22,23-dihydro-ivermectin A<sub>1a</sub>  
**Synonyms:** dihydro Avermectin B<sub>1</sub> monosaccharide, Ivermectin B<sub>1</sub> monosaccharide  
**MF:** C<sub>41</sub>H<sub>62</sub>O<sub>11</sub>  
**FW:** 730.9  
**Purity:**  $\geq 99\%$   
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

Ivermectin B<sub>1a</sub> monosaccharide is produced by selective hydrolysis of the terminal saccharide unit of ivermectin.<sup>1</sup> It can inhibit nematode larval development, but does not cause paralytic activity. This compound has been used as a probe to detect some types of ivermectin resistance.<sup>2</sup>

### References

1. 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).
2. 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, 10/31/2022

#### 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](http://WWW.CAYMANCHEM.COM)