

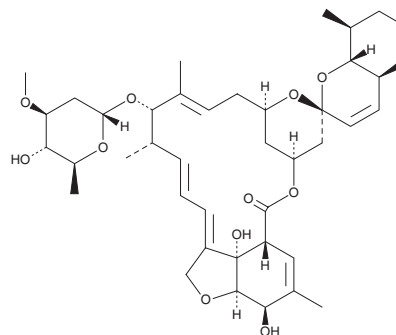
PRODUCT INFORMATION



Avermectin B_{1a} monosaccharide

Item No. 26690

CAS Registry No.: 71831-09-9
Formal Name: 4'-O-de(2,6-dideoxy-3-O-methyl- α -L-arabino-hexopyranosyl)-5-O-demethyl-avermectin A_{1a}
MF: C₄₁H₆₀O₁₁
FW: 728.9
Purity: $\geq 85\%$
UV/Vis.: λ_{max} : 246 nm
Supplied as: A crystalline 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

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

Description

Avermectin B_{1a} monosaccharide is a macrolide anthelmintic and monosaccharide form of avermectin B_{1a} (Item No. 22000).¹ It is lethal to *C. elegans* with a minimum active concentration (MAC) value of 0.1 μM . Avermectin B_{1a} monosaccharide also stimulates conductance in isolated shore crab muscle.

Reference

1. Bowman, J.W., Lee, B.L., Whaley, H.A., et al. Effects of dihydroavermectin B_{1a} and analogs on stretcher muscle of the lined shore crab, *Pachygrapsus crassipes*. *Comp. Biochem. Physiol.* **99(3)**, 333-340 (1991).

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