

PRODUCT INFORMATION

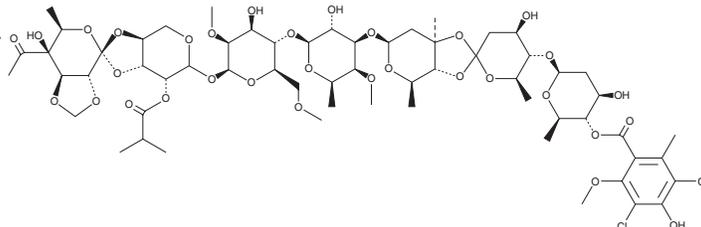


Avilamycin A

Item No. 29135

CAS Registry No.: 69787-79-7

Formal Name: O-(1R)-4-C-acetyl-6-deoxy-2,3-O-methylene-D-galactopyranosylidene-(1→3-4)-2-O-(2-methyl-1-oxopropyl)-α-L-lyxopyranosyl O-2,6-dideoxy-4-O-(3,5-dichloro-4-hydroxy-2-methoxy-6-methylbenzoyl)-β-D-arabino-hexopyranosyl-(1→4)-O-2,6-dideoxy-D-arabino-hexopyranosylidene-(1→3-4)-O-2,6-dideoxy-3-C-methyl-β-D-arabino-hexopyranosyl-(1→3)-O-6-deoxy-4-O-methyl-β-D-galactopyranosyl-(1→4)-2,6-di-O-methyl-β-D-mannopyranoside



MF: C₆₁H₈₈Cl₂O₃₂

FW: 1,404.2

Purity: ≥75%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years

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

Laboratory Procedures

Avilamycin A is supplied as a solid. A stock solution may be made by dissolving the avilamycin A in the solvent of choice, which should be purged with an inert gas. Avilamycin A is slightly soluble in acetone and methanol (warmed).

Description

Avilamycin A is an antibiotic.¹ It is active against veterinary isolates of *C. perfringens* (MICs = ≤0.06-0.5 mg/L) and *B. hyodysenteriae* (MICs = 12.5-100 µg/ml).^{1,2} Dietary administration of avilamycin A (15 and 30 ppm) reduces mortality in a broiler cockerel model of *C. perfringens* infection.³ Formulations containing avilamycin A have been used in the prevention of necrotic enteritis in livestock.

References

1. Watkins, K.L., Shryock, T.R., Dearth, R.N., *et al.* In-vitro antimicrobial susceptibility of *Clostridium perfringens* from commercial turkey and broiler chicken origin. *Vet. Microbiol.* **54(2)**, 195-200 (1997).
2. Uezato, K., Kinjo, E., and Adachi, Y. *In vitro* susceptibility of 21 antimicrobial agents to 37 isolates of *Brachyspira hyodysenteriae* isolated from pigs in Okinawa Prefecture. *J. Vet. Med. Sci.* **66(3)**, 307-309 (2004).
3. Paradis, M.A., McMillan, E., Bagg, R., *et al.* Efficacy of avilamycin for the prevention of necrotic enteritis caused by a pathogenic strain of *Clostridium perfringens* in broiler chickens. *Avian Pathol.* **45(3)**, 365-369 (2016).

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, 12/16/2019

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