

# PRODUCT INFORMATION



## Lasalocid

Item No. 15505

CAS Registry No.: 25999-31-9

Formal Name: 6-[(3R,4S,5S,7R)-7-[(2S,3S,5S)-5-ethyl-5-[(2R,5R,6S)-5-ethyltetrahydro-5-hydroxy-6-methyl-2H-pyran-2-yl]tetrahydro-3-methyl-2-furanyl]-4-hydroxy-3,5-dimethyl-6-oxononyl]-2-hydroxy-3-methyl-benzoic acid

Synonyms:

Antibiotic X 537A, CID-5360807, Ionophore X 537A

MF:  $C_{34}H_{54}O_8$

FW: 590.8

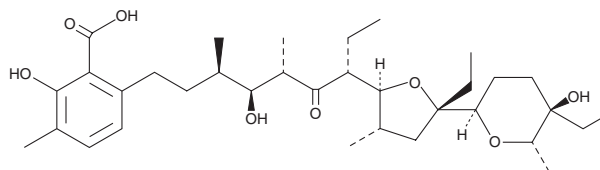
Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years

Item Origin: Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

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

### Description

Lasalocid is a polyketide-synthase derived ionophore antibiotic originally isolated from *S. lasaliensis*.<sup>1,2</sup> It binds to monovalent and divalent cations, including potassium, sodium, calcium, and magnesium cations.<sup>3</sup> Lasalocid (0.01 and 0.1  $\mu\text{g}/\text{ml}$ ) is active against *E. tenella*.<sup>4</sup> *In vivo*, lasalocid (125 ppm in the feed) reduces the severity of gastrointestinal lesions in chicks experimentally infected with *E. tenella*, *E. maxima*, *E. necatrix*, *E. brunetti*, or *E. acervulina*. It has been found in groundwater.<sup>5</sup> Formulations containing lasalocid have been used in the treatment of coccidiosis in poultry.

### References

1. Migita, A., Watanabe, M., Hirose, Y., *et al.* Identification of a gene cluster of polyether antibiotic lasalocid from *Streptomyces lasaliensis*. *Biosci. Biotechnol. Biochem.* **73**(1), 169-176 (2009).
2. Mitrovic, M. and Schildknecht, E.G. Anticoccidial activity of lasalocid (X-537A) in chicks. *Poult. Sci.* **53**(4), 1448-1455 (1974).
3. Antonenko, Y.N. and Yaguzhinsky, L.S. The ion selectivity of nonelectrogenic ionophores measured on a bilayer lipid membrane: nigericin, monensin, A23187 and lasalocid A. *Biochim. Biophys. Acta* **938**(2), 125-130 (1988).
4. Folz, S.D., Lee, B.L., Nowakowski, L.H., *et al.* Anticoccidial evaluation of halofuginone, lasalocid, maduramicin, monensin and salinomycin. *Vet. Parasitol.* **28**(1-2), 1-9 (1988).
5. Mooney, D., Richards, K.G., Danaher, M., *et al.* An investigation of anticoccidial veterinary drugs as emerging organic contaminants in groundwater. *Sci. Total Environ.* **746**, 141116 (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, 11/16/2022

#### 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