

PRODUCT INFORMATION

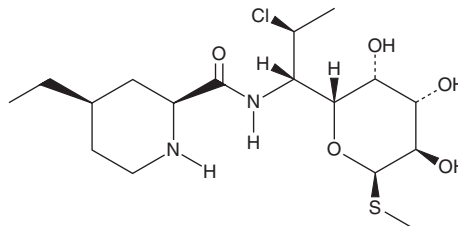


Pirlimycin

Item No. 20138

CAS Registry No.: 79548-73-5
Formal Name: methyl 7-chloro-6,7,8-trideoxy-6-[[[(2S,4R)-4-ethyl-2-piperidiny] carbonyl]amino]-1-thio-L-threo- α -D-galacto-octopyranoside

Synonym: U-57930E
MF: C₁₇H₃₁ClN₂O₅S
FW: 411.0
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

Pirlimycin is supplied as a solid. A stock solution may be made by dissolving the pirlimycin in the solvent of choice, which should be purged with an inert gas. Pirlimycin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Pirlimycin is a lincosamide antibiotic that is effective against Gram-positive bacteria, including *Staphylococcus*, *Streptococcus*, *Bacteroides*, and *Plasmodium*.¹ It inhibits bacterial protein synthesis by binding to the 50S subunit of the ribosome. Pirlimycin has been used by veterinarians in the treatment of mastitis in dairy cattle.²

References

1. Birkenmeyer, R. D., Kroll, S. J., Lewis, C., *et al.* Synthesis and antimicrobial activity of clindamycin analogues: Pirlimycin, 1,2 a potent antibacterial agent. *J. Med. Chem* **27(2)**, 216-223 (1984).
2. Lindeman, C. J., Portis, E., Johansen, L., *et al.* Susceptibility to antimicrobial agents among bovine mastitis pathogens isolated from North American dairy cattle, 2002-2010. *J. Vet. Diagn. Invest.* **25(5)**, 581-591 (2013).

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, 12/16/2022

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