

PRODUCT INFORMATION

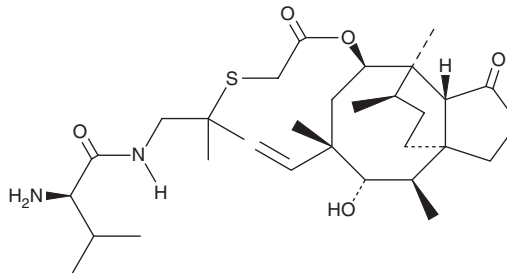


Valnemulin

Item No. 20141

CAS Registry No.: 101312-92-9
Formal Name: 2-[[2-[[[(2R)-2-amino-3-methyl-1-oxobutyl]amino]-1,1-dimethylethyl]thio]-acetic acid, (3aS,4R,5S,6S,8R,9R,9aR,10R)-6-ethenyldecahydro-5-hydroxy-4,6,9,10-tetramethyl-1-oxo-3a,9-propano-3aH-cyclopentacycloocten-8-yl ester

MF: C₃₁H₅₂N₂O₅S
FW: 564.8
Purity: ≥98%
Supplied as: A 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

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

Description

Valnemulin is a broad-spectrum pleuromutilin antibiotic that binds to peptidyl transferase in the 50S ribosomal subunit.¹ It has been used in veterinary medicine to treat swine diseases including, *B. hyodysenteriae* and *M. hyopneumoniae*.¹

Reference

1. Poulsen, S. M., Karlsson, M., Johansson, L. B., *et al.* The pleuromutilin drugs tiamulin and valnemulin bind to the RNA at the peptidyl transferase centre on the ribosome. *Mol. Microbiol.* **41(5)**, 1091-1099 (2001).

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, 11/04/2022

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