

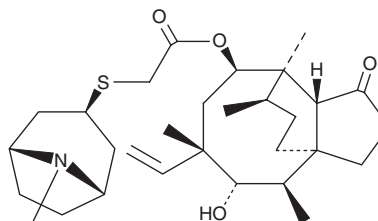
PRODUCT INFORMATION



Retapamulin

Item No. 21528

CAS Registry No.: 224452-66-8
Formal Name: 2-[[[(3-*exo*)-8-methyl-8-azabicyclo[3.2.1]oct-3-yl]thio]-acetic acid, (3a*S*,4*R*,5*S*,6*S*,8*R*,9*R*,9a*R*,10*R*)-6-ethenyldecahydro-5-hydroxy-4,6,9,10-tetramethyl-1-oxo-3a,9-propano-3a*H*-cyclopentacycloocten-8-yl ester
Synonym: SB-275833
MF: C₃₀H₄₇NO₄S
FW: 517.8
Purity: ≥99%
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

Retapamulin is supplied as a crystalline solid. A stock solution may be made by dissolving the retapamulin in the solvent of choice, which should be purged with an inert gas. Retapamulin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of retapamulin in these solvents is approximately 30 mg/ml.

Retapamulin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, retapamulin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Retapamulin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Retapamulin is a pleuromutilin antibiotic that inhibits bacterial protein synthesis by binding to bacterial ribosomes in the peptidyl transferase component of the 50S subunit and inhibiting peptide bond formation.¹ It has activity against staphylococcal, streptococcal, and anaerobic Gram-positive bacteria.¹

Reference

1. Wilson, D.N. The A-Z of bacterial translation inhibitors. *Crit. Rev. Biochem. Mol. Biol.* **44**(6), 393-433 (2009).

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, 11/02/2022

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