

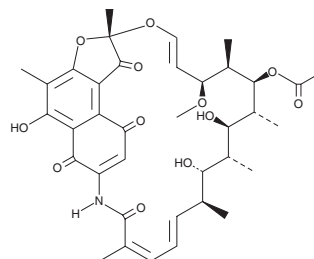
PRODUCT INFORMATION



Rifamycin S

Item No. 42178

CAS Registry No.: 13553-79-2
Formal Name: 1,4-dideoxy-1,4-dihydro-1,4-dioxo-rifamycin
Synonym: NSC 144130
MF: C₃₇H₄₅NO₁₂
FW: 695.8
Purity: ≥95%
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

Rifamycin S is supplied as a solid. A stock solution may be made by dissolving the rifamycin S in the solvent of choice, which should be purged with an inert gas. Rifamycin S is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

Rifamycin S is an antibiotic.² It is active against *S. haemolyticus*, *M. aureus*, *M. tuberculosis*, *B. subtilis*, and *S. faecalis* (MICs = 0.025, 0.05, 0.05, 0.75, and 0.09 µg/kg, respectively). Rifamycin S (65 µM) inhibits RNA and protein synthesis in *A. montana*.¹ It is toxic to mice (LD₅₀ = 122 mg/kg).²

References

1. Rodriguez-López, M., Muñoz, M.L., and Vazquez, D. The effects of the rifamycin antibiotics on algae. *FEBS Lett.* **9**(3), 171-174 (1970).
2. Sensi, P., Timbal, M.T., and Maffh, G. Rifomycin IX two new antibiotics of rifomycin family: Rifomycin S and rifomycin SV. *Experientia* **16**, 412 (1960).

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, 11/18/2024

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