

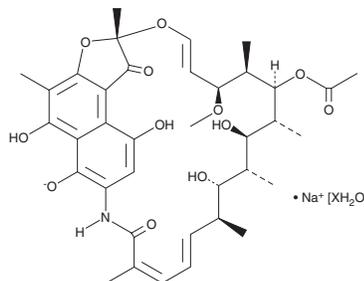
# PRODUCT INFORMATION



## Rifamycin SV (sodium salt hydrate)

Item No. 21441

<b>Formal Name:</b>	rifamycin, monosodium salt, hydrate
<b>Synonyms:</b>	CB 01-11, NSC 146718
<b>MF:</b>	$C_{37}H_{46}NO_{12} \cdot Na [XH_2O]$
<b>FW:</b>	719.8
<b>Purity:</b>	≥98%
<b>UV/Vis.:</b>	$\lambda_{max}$ : 226, 315, 450 nm
<b>Supplied as:</b>	A crystalline solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

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

Rifamycin SV (sodium salt hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rifamycin SV (sodium salt hydrate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Rifamycin SV (sodium salt hydrate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Rifamycin SV is a broad-spectrum antibiotic that inhibits bacterial DNA-dependent RNA polymerase.<sup>1,2</sup> It is effective against Gram-positive bacteria but only moderately active against Gram-negative bacteria.<sup>1,2</sup>

### References

1. Riva, S. and Silvestri, L.G. Rifamycins: A general view. *Annu. Rev. Microbiol.* **26**, 199-224 (1982).
2. Kohanski, M.A., Dwyer, D.J., and Collins, J.J. How antibiotics kill bacteria: From targets to networks. *Nat. Rev. Microbiol.* **8(6)**, 423-435 (2010).

#### 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, 12/01/2022

#### 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