

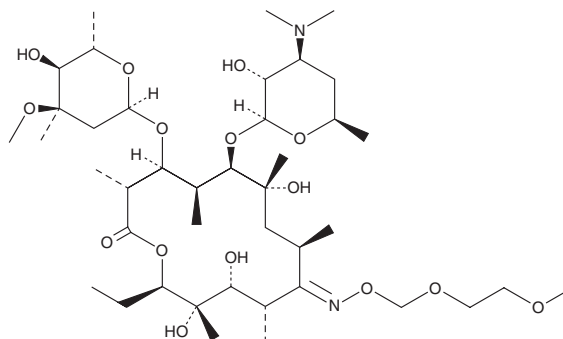
PRODUCT INFORMATION



Roxithromycin

Item No. 19465

CAS Registry No.: 80214-83-1
Formal Name: 9E-[O-[(2-methoxyethoxy)methyl]oxime]-erythromycin
Synonyms: RU 965, RU 28965
MF: C₄₁H₇₆N₂O₁₅
FW: 837.1
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

Roxithromycin is supplied as a solid. A stock solution may be made by dissolving the roxithromycin in the solvent of choice, which should be purged with an inert gas. Roxithromycin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of roxithromycin in ethanol is approximately 30 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Roxithromycin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, roxithromycin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Roxithromycin 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

Roxithromycin is a macrolide antibiotic.¹ It is active against various Gram-positive and Gram-negative bacteria, including five strains each of *Staphylococcus* and *Streptococcus* (geometric mean MICs = 0.08 and 0.79 µg/ml, respectively), as well as several strains each of *Corynebacterium*, *Haemophilus*, and *S. pneumoniae* (MIC₅₀s = 0.02, 0.01, and 2.5 µg/ml, respectively). Roxithromycin is protective against infection by strains of *S. aureus*, *S. pyogenes*, *S. pneumoniae*, or *L. monocytogenes* in mice with 50% protective dose (PD₅₀) values ranging from 23 to 98 mg/kg. Formulations containing roxithromycin have been used in the treatment of bacterial infections.

Reference

1. Chantot, J.F., Bryskier, A., and Gasc, J.C. Antibacterial activity of roxithromycin: A laboratory evaluation. *J. Antibiot. (Tokyo)* **39**(5), 660-668 (1986).

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, 12/22/2022

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