

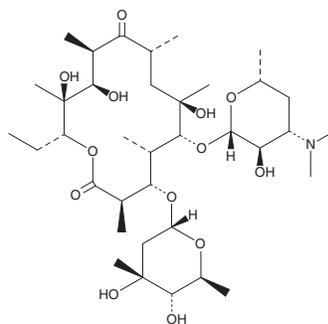
PRODUCT INFORMATION



Erythromycin C

Item No. 23478

CAS Registry No.: 1675-02-1
Formal Name: 3''-O-demethyl-erythromycin
MF: C₃₆H₆₅NO₁₃
FW: 719.9
Purity: ≥85%
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

Erythromycin C is supplied as a solid. A stock solution may be made by dissolving the erythromycin C in the solvent of choice, which should be purged with an inert gas. Erythromycin C is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. Erythromycin C has limited water solubility. We do not recommend storing the aqueous solution for more than one day.

Description

Erythromycin C is an intermediate in the biosynthesis of erythromycin (Item No. 16486).¹ It is also a potential impurity found in commercial preparations of erythromycin. Erythromycin C is half as active or less than erythromycin A or B against most Gram-positive and Gram-negative bacteria tested.²

References

1. Chen, Y., Deng, W., Wu, J., *et al.* Genetic modulation of the overexpression of tailoring genes eryK and eryG leading to the improvement of erythromycin A purity and production in *Saccharopolyspora erythraea* fermentation. *Appl. Environ. Microbiol.* **74**(6), 1820-1828 (2008).
2. Kibwage, I.O., Hoogmartens, J., Roets, E., *et al.* Antibacterial activities of erythromycins A, B, C, and D and some of their derivatives. *Antimicrob. Agents Chemother.* **28**(5), 630-633 (1985).

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, 02/07/2023

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