

PRODUCT INFORMATION

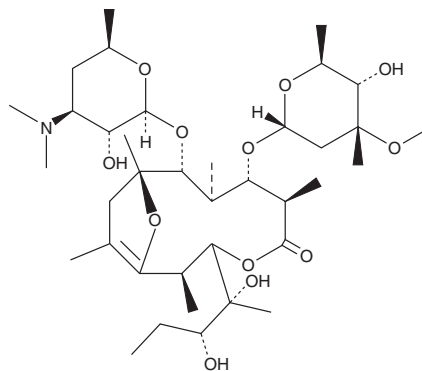


Pseudoerythromycin A enol ether

Item No. 15508

CAS Registry No.: 105882-69-7
Formal Name: (2R,3R,6R,8S)-7S-[(2,6-dideoxy-3-C-methyl-3-O-methyl- α -L-ribo-hexopyranosyl)oxy]-3-[(1R,2R)-1,2-dihydroxy-1-methylbutyl]-2,6,8,10R,12-pentamethyl-9R-[[3,4,6-trideoxy-3-(dimethylamino)- β -D-xylo-hexopyranosyl]oxy]-4,13-dioxabicyclo[8.2.1]tridec-12-en-5-one

Synonym: LY267108
MF: C₃₇H₆₅NO₁₂
FW: 715.9
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Pseudoerythromycin A enol ether is supplied as a solid. A stock solution may be made by dissolving the pseudoerythromycin A enol ether in the solvent of choice, which should be purged with an inert gas. Pseudoerythromycin A enol ether is soluble in organic solvents such as ethanol, DMSO, dimethyl formamide, and methanol.

Description

Pseudoerythromycin A enol ether is a degradation product of the macrolide antibiotic erythromycin (Item No. 16486).¹ It promotes monocyte differentiation into macrophages when used at a concentration of 10 μ M.²

References

1. Pendela, M., Béni, S., Haghedooren, E., *et al.* Combined use of liquid chromatography with mass spectrometry and nuclear magnetic resonance for the identification of degradation compounds in an erythromycin formulation. *Anal. Bioanal. Chem.* **402**(2), 781-790 (2012).
2. Yoshida, K., Sunazuka, T., Nagai, K., *et al.* Macrolides with promotive activity of monocyte to macrophage differentiation. *J. Antibiot. (Tokyo)* **58**(1), 79-81 (2005).

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, 01/23/2023

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