

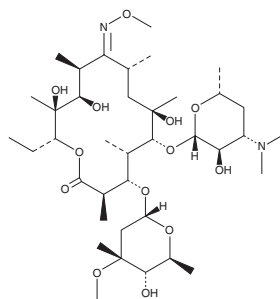
PRODUCT INFORMATION



Lexithromycin

Item No. 22002

CAS Registry No.: 53066-26-5
Formal Name: erythromycin, 9-(O-methyloxime)
MF: $C_{38}H_{70}N_2O_{13}$
FW: 763.0
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Lexithromycin is supplied as a solid. A stock solution may be made by dissolving the lexithromycin in the solvent of choice, which should be purged with an inert gas. Lexithromycin is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Lexithromycin is a semisynthetic antibiotic derived from erythromycin (Item No. 16486).¹ Lexithromycin has improved absorption *in vivo* over erythromycin due to increased hydrophobicity and pH stability.¹ As with other macrolides, lexithromycin prevents protein synthesis by binding the ribosome at the polypeptide exit tunnel.²

References

1. Hunt, E., Knowles, D.J., Shillingford, C., *et al.* Erythromycin A 11,12-methylene acetal. *J. Antibiot. (Tokyo)* **41(11)**, 1644-1648 (1988).
2. Wilson, D.N. The A-Z of bacterial translation inhibitors. *Crit. Rev. Biochem. Mol. Biol.* **44(6)**, 393-433 (2009).

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, 10/26/2022

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