

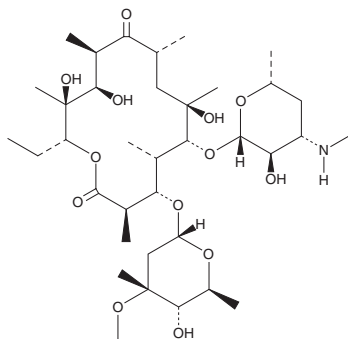
PRODUCT INFORMATION



N-Demethylerythromycin A

Item No. 27988

CAS Registry No.: 992-62-1
Formal Name: N-demethyl-erythromycin
MF: $C_{36}H_{65}NO_{13}$
FW: 719.9
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

N-Demethylerythromycin A is supplied as a solid. A stock solution may be made by dissolving the N-demethylerythromycin A in the solvent of choice, which should be purged with an inert gas. N-Demethylerythromycin A is soluble in chloroform and methanol.

Description

N-Demethylerythromycin A is a metabolite of erythromycin (Item No. 16486).¹ It is also a fungal metabolite that has been found in *S. erythreus* and a potential impurity in commercial preparations of erythromycin.²

References

1. Kibwage, I.O., Roets, E., Verbruggen, A., *et al.* Thin-layer chromatographic study of the metabolites of erythromycins in the Wistar rat. *J. Chromatogr.* **434**(1), 177-186 (1988).
2. Deubel, A., Fandino, A.S., Sorgel, F., *et al.* Determination of erythromycin and related substances in commercial samples using liquid chromatography/ion trap mass spectrometry. *J. Chromatogr. A* **1136**(1), 39-47 (2006).

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, 12/19/2022

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