

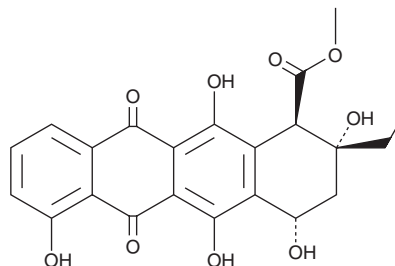
PRODUCT INFORMATION



ϵ -Rhodomycinone

Item No. 32920

CAS Registry No.: 21288-60-8
Formal Name: (1R,2R,4S)-2-ethyl-1,2,3,4,6,11-hexahydro-2,4,5,7,12-pentahydroxy-6,11-dioxo-1-naphthacenicarboxylic acid, methyl ester
Synonym: NSC 196524
MF: C₂₂H₂₀O₉
FW: 428.4
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

ϵ -Rhodomycinone is supplied as a solid. A stock solution may be made by dissolving the ϵ -rhodomycinone in the solvent of choice, which should be purged with an inert gas. ϵ -Rhodomycinone is soluble in methanol and DMSO.

Description

ϵ -Rhodomycinone is a bacterial metabolite that has been found in *S. griseoruber*.¹ It is a precursor to rhodomycin D, which is an intermediate in the bioconversion of ϵ -rhodomycinone to daunorubicin (Item No. 14159) and doxorubicin (Item No. 15007).^{2,3}

References

1. Podojil, M., Blumauerová, M., Přikrylová, V., et al. Production of rhodomycins in *Streptomyces griseoruber* 4620. *Folia Microbiol. (Praha)* **25(6)**, 464-466 (1980).
2. Olano, C., Lomovskaya, N., Fonstein, L., et al. A two-plasmid system for the glycosylation of polyketide antibiotics: Bioconversion of ϵ -rhodomycinone to rhodomycin D. *Chem. Biol.* **6(12)**, 845-855 (1999).
3. Dickens, M.L., Priestley, N.D., and Strohl, W.R. In vivo and in vitro bioconversion of ϵ -rhodomycinone glycoside to doxorubicin: Functions of DauP, DauK, and DoxA. *J. Bacteriol.* **179(8)**, 2641-2650 (1997).

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, 11/10/2022

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