

PRODUCT INFORMATION

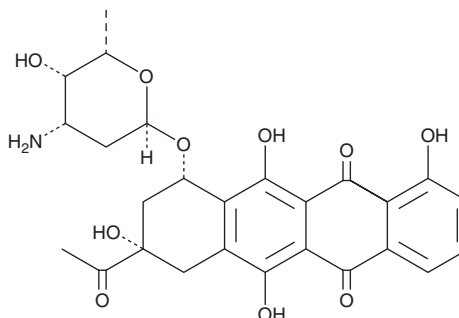


Carubicin

Item No. 27660

CAS Registry No.: 50935-04-1
Formal Name: (8S,10S)-8-acetyl-10-[(3-amino-2,3,6-trideoxy- α -L-lyxo-hexopyranosyl)oxy]-7,8,9,10-tetrahydro-1,6,8,11-tetrahydroxy-5,12-naphthacenedione
Synonyms: Antibiotic R 588A, Carminomycin, NCI 180024, NSC 180024

MF: $C_{26}H_{27}NO_{10}$
FW: 513.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Carubicin is supplied as a solid. A stock solution may be made by dissolving the carubicin in the solvent of choice, which should be purged with an inert gas. Carubicin is soluble in the organic solvent DMSO.

Description

Carubicin is an anthracycline originally isolated from *A. carminata* that has anticancer activity.^{1,2} It is cytotoxic to MCF-7 and K562 cells ($IC_{50}S = 90$ and 60 nM, respectively), as well as P-glycoprotein-expressing MCF-7/DOX and K562i/S9 cells ($IC_{50}S = 90$ and 60 nM, respectively).¹

References

1. Tevyashova, A.N., Shtil, A.A., Olsufyeva, E.N., *et al.* Carminomycin, 14-hydroxycarminomycin and its novel carbohydrate derivatives potently kill human tumor cells and their multidrug resistant variants. *J. Antibiot. (Tokyo)* **57(2)**, 143-150 (2004).
2. Gause, G.F., Brazhnikova, M.G., and Shorin, V.A. A new antitumor antibiotic, carminomycin (NSC-180024). *Cancer Chemother. Rep.* **58(2)**, 255-256 (1974).

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, 11/02/2022

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