

PRODUCT INFORMATION



Cinerubin B

Item No. 40865

CAS Registry No.: 35906-51-5

Formal Name: (1R,2R,4S)-4-[[[2''',3''-anhydro]-O-3,6-dideoxy- α -L-erythro-hexopyranos-4-ulos-1-yl-(1 $\rightarrow$ 4)-O-2,6-dideoxy- α -L-lyxo-hexopyranosyl-(1 $\rightarrow$ 4)-2,3,6-trideoxy-3-(dimethylamino)- α -L-lyxo-hexopyranosyl]oxy]-2-ethyl-1,2,3,4,6,11-hexahydro-2,5,7,10-tetrahydroxy-6,11-dioxo-1-naphthacenecarboxylic acid, methyl ester

Synonyms: Ciclamycin 1, Cinerubine B, MA 144B2, NSC 18335, NSC 62490, NSC 243023

MF: C₄₂H₅₁NO₁₆

FW: 825.9

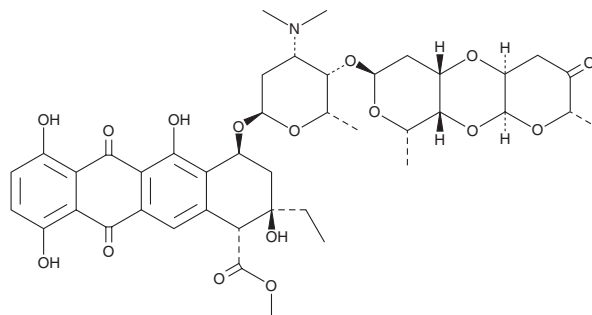
Purity: $\geq 70\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Streptomyces aurantiacus*



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

Laboratory Procedures

Cinerubin B is supplied as a solid. A stock solution may be made by dissolving the cinerubin B in the solvent of choice, which should be purged with an inert gas. Cinerubin B is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

Cinerubin B is an anthracycline antitumor antibiotic that has been found in *Streptomyces*.^{1,2} It inhibits RNA synthesis in, and the growth of, L1210 murine leukemia cells (IC₅₀s = 48 and 15 nM, respectively).²

References

1. Silva, L.J., Crevelin, E.J., Souza, D.T., *et al.* Actinobacteria from Antarctica as a source for anticancer discovery. *Sci. Rep.* **10**(1), 13870 (2020).
2. Matsuzawa, Y., Oki, T., Takeuchi, T., *et al.* Structure-activity relationships of anthracyclines relative to cytotoxicity and effects on macromolecular synthesis in L1210 leukemia cells. *J. Antibiot. (Tokyo)* **34**(12), 1596-1607 (1981).

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, 05/23/2024

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