

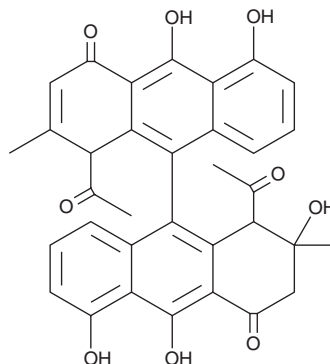
PRODUCT INFORMATION



Setomimycin

Item No. 27958

CAS Registry No.: 69431-87-4
Formal Name: 1,1'-diacetyl-2,3-dihydro-2,5,5',10,10'-pentahydroxy-2,2'-dimethyl-[9,9'-bianthracene]-4,4'(1H,1'H)-dione
Synonym: Antibiotic A 39183B
MF: C₃₄H₂₈O₉
FW: 580.6
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces aurantiacus* JA 4570



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

Laboratory Procedures

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

Description

Setomimycin is a pre-anthraquinone originally isolated from *S. pseudovenezuelae*.^{1,2} It is active against a variety of Gram-positive bacteria, including *S. aureus*, *B. subtilis*, *B. cereus*, and *M. smegmatis* (MICs = 1.56-3.13 µg/ml).¹ It also reduces tumor growth in a Sarcoma-180 mouse solid tumor model when administered at a dose of 200 mg/kg per day for seven days.

References

1. Omura, S., Tanaka, H., Iwai, Y., *et al.* A new antibiotic, setomimycin, produced by a strain of *Streptomyces*. *J. Antibiot. (Tokyo)* **31(11)**, 1091-1098 (1978).
2. Präg, A., Grüning, B.A., Häckh, M., *et al.* Regio- and stereoselective intermolecular oxidative phenol coupling in *Streptomyces*. *J. Am. Chem. Soc.* **136(17)**, 6195-6198 (2014).

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, 01/11/2023

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