

PRODUCT INFORMATION

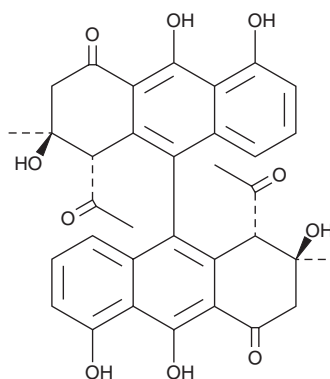


A-39183A

Item No. 23239

CAS Registry No.: 79426-51-0
Formal Name: (1S,1'S,2S,2'S)-1,1'-diacetyl-2,2',3,3'-tetrahydro-2,2',5,5',10,10'-hexahydroxy-2,2'-dimethyl-[9,9'-bianthracene]-4,4'-(1H,1'H)-dione

MF: C₃₄H₃₀O₁₀
FW: 598.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

A-39183A is supplied as a solid. A stock solution may be made by dissolving the A-39183A in the solvent of choice, which should be purged with an inert gas. A-39183A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide.

Description

A-39183A is an active component of the A-39183 antibiotic complex produced by aerobic fermentation of *Streptomyces* NRRL 12049.¹ It has activity against penicillin-resistant *S. aureus* 3055 and *S. faecalis* X66 (MICs = 32 and 32 µg/ml, respectively) and a variety of Gram-positive and Gram-negative anaerobic bacteria (MICs = 16-128 µg/ml) in an agar dilution assay.

Reference

1. Koch, K.F. and Kastner, R.E. Antibiotics and process for production thereof. **US 4283390** (1981), US06090386.

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, 11/15/2022

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