

PRODUCT INFORMATION



Benanomicin A

Item No. 32886

CAS Registry No.: 116249-65-1

Formal Name: N-[[[(5S,6S)-5-[(6-deoxy-3-O-β-D-xylopyranosyl-β-D-galactopyranosyl)oxy]-5,6,8,13-tetrahydro-1,6,9,14-tetrahydroxy-11-methoxy-3-methyl-8,13-dioxobenzo[a]naphthacen-2-yl]carbonyl]-D-alanine

MF: C₃₉H₄₁NO₁₉

FW: 827.7

Purity: ≥90%

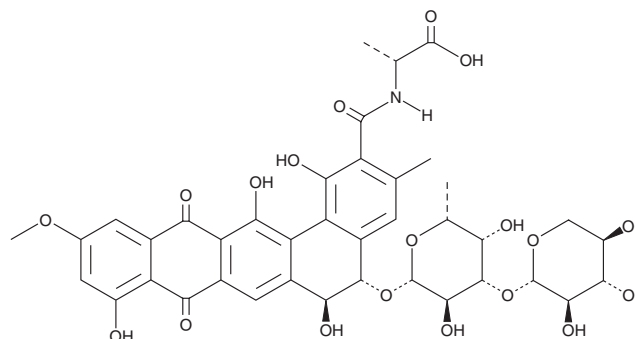
UV/Vis.: λ_{max}: 294 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Actinomadura* sp.



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

Laboratory Procedures

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

Description

Benanomicin A is a microbial metabolite that has been found in *Actinomycetes* and has antifungal, fungicidal, and antiviral activities.^{1,2} It is active against a variety of mammalian and plant pathogenic fungi, including *C. albicans*, *T. mentagrophytes*, *C. neoformans*, *P. oryzae*, and *A. niger* (MICs = 3.13-50 μg/ml).¹ Benanomicin A inhibits HIV-1 viral infection in MT-4 cells in a concentration-dependent manner.²

References

1. Takeuchi, T., Hara, T., Naganawa, H., *et al.* New antifungal antibiotics, benanomicins A and B from an actinomycete. *J. Antibiot. (Tokyo)* **41**(6), 807-811 (1987).
2. Kondo, S., Gomi, S., Ikeda, D., *et al.* Antifungal and antiviral activities of benanomicins and their analogues. *J. Antibiot. (Tokyo)* **44**(11), 1228-1236 (1990).

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, 07/22/2025

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