

PRODUCT INFORMATION



OM173-αA

Item No. 27570

CAS Registry No.: 58286-56-9

Formal Name: (1S,3R)-3,4,5,10-tetrahydro-9-hydroxy-1-methyl-5,10-dioxo-1H-naphtho[2,3-c]pyran-3-acetic acid, methyl ester

Synonyms: Nanaomycin αA,
Nanaomycin A methyl ester

MF: C₁₇H₁₆O₆

FW: 316.3

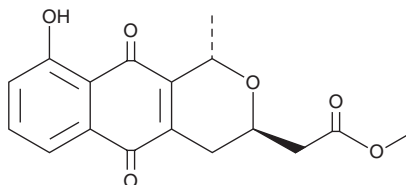
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces rosa*, *Notoensis*



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

Laboratory Procedures

OM173-αA is supplied as a solid. A stock solution may be made by dissolving the OM173-αA in the solvent of choice. OM173-αA is soluble in the organic solvent ethanol, which should be purged with an inert gas, at a concentration of approximately 1 mg/ml. It is also soluble in methanol and DMSO.

Description

OM173-αA is a quinone bacterial metabolite originally isolated from *Streptomyces* that has antimicrobial activity.^{1,2} It inhibits the growth of the bacteria *M. gallisepticum*, *M. pneumoniae*, and *S. aureus* *in vitro* with MIC values ranging from 0.39 to 3.13 µg/ml. OM173-αA also inhibits the growth of the plant pathogenic fungus *P. oryzae* (MIC = 3.12 µg/ml) and several species of *Trichophyton* (MICs = 12.5-25 µg/ml).¹

References

1. Iwai, Y., Kimura, K., Takahashi, Y., *et al.* OM-173, new nanaomycin-type antibiotics produced by a strain of *Streptomyces*. Taxonomy, production, isolation and biological properties. *J. Antibiot. (Tokyo)* **36**(10), 1268-1274 (1983).
2. Liu, C., Lei, H., Chen, X., *et al.* Four new nanaomycins produced by *Streptomyces hebeiensis* derived from lichen. *Chem. Biodivers.* **14**(7), (2017).

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, 12/01/2022

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