

PRODUCT INFORMATION



Obafluorin

Item No. 25731

CAS Registry No.: 92121-68-1
Formal Name: 2,3-dihydroxy-N-[(2R,3S)-2-[(4-nitrophenyl)methyl]-4-oxo-3-oxetanyl]-benzamide

MF: C₁₇H₁₄N₂O₇

FW: 358.3

Purity: ≥98%

UV/Vis.: λ_{max}: 257 nm

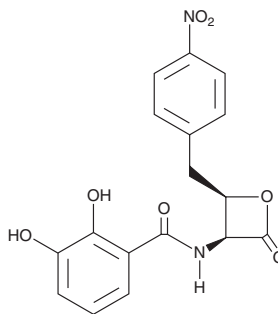
Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years

Item Origin: Bacterium/*Pseudomonas fluorescens*

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



Laboratory Procedures

Obafluorin is supplied as a crystalline solid. A stock solution may be made by dissolving the obafluorin in the solvent of choice, which should be purged with an inert gas. Obafluorin is soluble in the organic solvent acetonitrile at a concentration of approximately 1 mg/ml.

Description

Obafluorin is a β-lactone antibiotic that has been found in *P. fluorescens*.¹ It is active against *S. aureus*, *S. faecalis*, *K. pneumoniae*, and *P. vulgaris* (MIC = 125 μg/ml for all).

Reference

1. Pu, Y., Lowe, C., Sailer, M., et al. Synthesis, stability, and antimicrobial activity of (+)-obafluorin and related β-lactone antibiotics. *J. Org. Chem.* **59**(13), 3642-3655 (1994).

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, 11/1/2019

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