

PRODUCT INFORMATION



Orfamide B

Item No. 23897

CAS Registry No.: 939960-35-7

Formal Name: N-(3-hydroxy-1-oxotetradecyl)-L-leucyl-D- α -glutamyl-D-allothreonyl-D-valyl-L-leucyl-D-seryl-L-leucyl-L-leucyl-D-seryl-L-valine(10 \rightarrow 3)-lactone

MF: C₆₃H₁₁₂N₁₀O₁₇

FW: 1,281.6

Purity: $\geq 85\%$

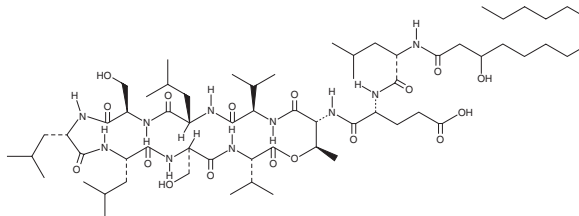
Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Pseudomonas fluorescens*

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



Laboratory Procedures

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

Description

Orfamide B is a cyclic lipopeptide originally isolated from *P. fluorescens*.¹ It is the main orfamide produced by *Pseudomonas* sp. CMR12a. Orfamide B inhibits mycelial growth and increases hyphal branching in the fungi *R. solani* AG2-1 at concentrations of 10 and 100 μ M and in *R. solani* AG 4-HGI at a concentration of 100 μ M.²

References

1. Gross, H., Stockwell, V.O., Henkels, M.D., *et al.* The genomisotopic approach: A systematic method to isolate products of orphan biosynthetic gene clusters. *Chem. Biol.* **14**(1), 53-63 (2007).
2. Olorunleke, F.E., Hua, G.K.H., Kieu, N.P., *et al.* Interplay between orfamides, sessilins and phenazines in the control of Rhizoctonia diseases by *Pseudomonas* sp. CMR12a. *Environ. Microbiol. Rep.* **7**(5), 774-781 (2015).

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, 12/22/2022

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