

PRODUCT INFORMATION



Pseudomonic Acid (lithium salt)

Item No. 33271

CAS Registry No.: 73346-79-9
Formal Name: 5,9-anhydro-2,3,4,8-tetra-deoxy-8-
[[[(2S,3S)-3-[(1S,2S)-2-hydroxy-1-
methylpropyl]-2-oxiranyl]methyl]-
3-methyl-L-talo-non-2-enonic acid,
8-carboxyooctyl ester, monolithium salt

MF: $C_{26}H_{43}O_9 \cdot Li$
FW: 506.6

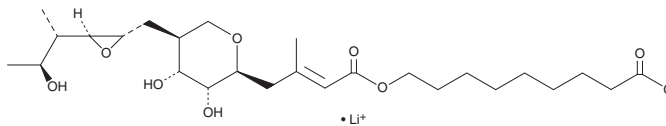
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 223 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

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

Pseudomonic acid (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the pseudomonic acid (lithium salt) in the solvent of choice, which should be purged with an inert gas. Pseudomonic acid (lithium salt) is soluble in the organic solvent ethanol at a concentration of approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of pseudomonic acid (lithium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of pseudomonic acid (lithium salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pseudomonic acid is an antibiotic and bacterial metabolite that has been found in *P. fluorescens*.¹ It is bacteriostatic against *S. aureus* (MIC = 0.05 $\mu\text{g/ml}$) and active against skin wound clinical isolates of methicillin-resistant *S. aureus* (MRSA; MICs = 1-4 $\mu\text{g/ml}$).^{2,3} Pseudomonic acid inhibits MRSA and *P. aeruginosa* biofilm formation *in vitro*.⁴ It inhibits bacterial cell wall isoleucyl-tRNA synthetase, slowing bacterial growth.^{2,4} Topical administration of pseudomonic acid (2% v/v) reduces the number of wound colony forming units (CFUs) in a mouse model of MRSA skin infection.³

References

1. Fuller, A.T., Mellows, G., Woolford, M., et al. Pseudomonic acid: An antibiotic produced by *Pseudomonas fluorescens*. *Nature* **234**(5329), 416-417 (1971).
2. Hughes, J. and Mellows, G. On the mode of action of pseudomonic acid: Inhibition of protein synthesis in *Staphylococcus aureus*. *J. Antibiot. (Tokyo)* **31**(4), 330-335 (1978).
3. Mohammad, H., Cushman, M., and Seleem, M.N. Antibacterial evaluation of synthetic thiazole compounds *in vitro* and *in vivo* in a methicillin-resistant *Staphylococcus aureus* (MRSA) skin infection mouse model. *PLoS One* **10**(11), e0142321 (2015).
4. Khoshnood, S., Heidary, M., Asadi, A., et al. A review on mechanism of action, resistance, synergism, and clinical implications of mupirocin against *Staphylococcus aureus*. *Biomed. Pharmacother.* **109**, 1809-1818 (2019).

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, 11/14/2022

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