

# PRODUCT INFORMATION



## Pseudomonic Acid

Item No. 23238

**CAS Registry No.:** 12650-69-0  
**Formal Name:** 5,9-anhydro-2,3,4,8-tetradecoxy-8-[[[(2S,3S)-3-[(1S,2S)-2-hydroxy-1-methylpropyl]oxiranyl]methyl]-3-methyl-L-talo-non-2E-enonic acid, 8-carboxyoctyl ester

**Synonym:** BRL 4910A

**MF:** C<sub>26</sub>H<sub>44</sub>O<sub>9</sub>

**FW:** 500.6

**Purity:** ≥95%

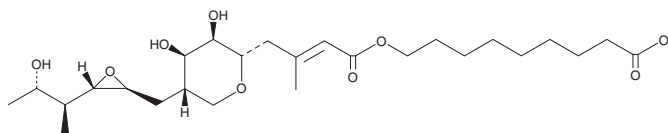
**UV/Vis.:** λ<sub>max</sub>: 221 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fermentation (species not specified)



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

### Laboratory Procedures

Pseudomonic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the pseudomonic acid in the solvent of choice, which should be purged with an inert gas. Pseudomonic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of pseudomonic acid in these solvents is approximately 30 mg/ml.

Pseudomonic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pseudomonic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Pseudomonic acid has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. 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*.<sup>1</sup> It is bacteriostatic against *S. aureus* (MIC = 0.05 µg/ml) and active against skin wound clinical isolates of methicillin-resistant *S. aureus* (MRSA; MICs = 1-4 µg/ml).<sup>2,3</sup> Pseudomonic acid inhibits MRSA and *P. aeruginosa* biofilm formation *in vitro*.<sup>4</sup> It inhibits bacterial cell wall isoleucyl-tRNA synthetase, slowing bacterial growth.<sup>2,4</sup> 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.<sup>3</sup>

### 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. Yang, N., Cui, H., Han, F., *et al.* Paeniflorin inhibits human pancreatic cancer cell apoptosis via suppression of MMP-9 and ERK signaling. *Oncol. Lett.* **12**(2), 1471-1476 (2016).
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, 11/04/2022

#### 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