

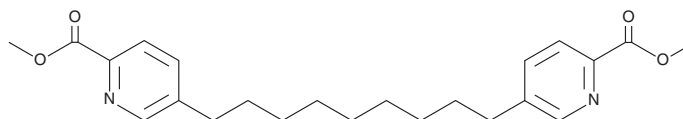
PRODUCT INFORMATION



Penicolinate A

Item No. 25013

CAS Registry No.: 1418291-68-5
Formal Name: 5,5'-(1,10-decanediyl)bis-2-pyridinecarboxylic acid, 2,2'-dimethyl ester
Synonym: Penicolinate B dimethyl ester
MF: $C_{24}H_{32}N_2O_4$
FW: 412.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Penicolinate A is a fungal metabolite produced by *Penicillium* sp. BCC16054 with diverse biological activities.¹ It exhibits antimalarial and antituberculosis activity (MICs = 3.25 and 12.5 mg/ml for *P. falciparum* and *M. tuberculosis* H37Ra, respectively). Penicolinate A is cytotoxic to MCF-7, KB, NCI-H187, and Vero cells (IC_{50}s = 7.49, 18.43, 4.4, and 14.08 μM , respectively).

Reference

- Intaraudom, C., Boonyuen, N., Suvannakad, R., *et al.* Penicolinates A–E from endophytic *Penicillium* sp. BCC16054. *Tetrahedron Lett.* **54**(8), 744-748 (2013).

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, 08/08/2023

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