

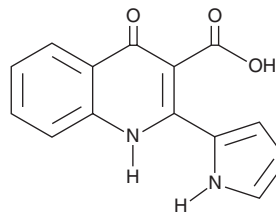
PRODUCT INFORMATION



Penicinoline

Item No. 29932

CAS Registry No.: 1214268-60-6
Formal Name: 1,4-dihydro-4-oxo-2-(1H-pyrrol-2-yl)-3-quinolinecarboxylic acid
MF: C₁₄H₁₀N₂O₃
FW: 254.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

Penicinoline is supplied as a solid. A stock solution may be made by dissolving the penicinoline in the solvent of choice, which should be purged with an inert gas. Penicinoline is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Penicinoline is an alkaloid that has been found in *Penicillium* and has antimalarial, insecticidal, and anticancer activities.^{1,2} It is active against chloroquine-sensitive and -resistant strains of *P. falciparum* (IC₅₀ = 25 μM for both).¹ Penicinoline (1,000 ppm) is also active against the aphid *A. gossypii*.² It inhibits proliferation of 95-D and HepG2 cancer cells (IC₅₀s = 0.57 and 6.5 μg/ml, respectively) but not HeLa, KB, KBv200, or Hep-2 cells (IC₅₀s = >100 μg/ml).

References

1. Naveen, B., Ommi, N.B., Mudiraj, A., *et al.* Total synthesis of penicinoline E, marinamide, methyl marinamide and their antimalarial activity. *ChemistrySelect* **2(11)**, 3256-3261 (2017).
2. Shao, C.-L., Wang, C.-Y., Gu, Y.-C., *et al.* Penicinoline, a new pyrrolyl 4-quinolinone alkaloid with an unprecedented ring system from an endophytic fungus *Penicillium* sp. *Bioorg. Med. Chem. Lett.* **20(11)**, 3284-3286 (2010).

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, 04/29/2024

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