

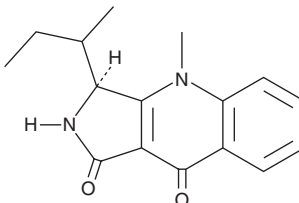
PRODUCT INFORMATION



Quinolactacin A

Item No. 27701

CAS Registry No.: 319917-25-4
Formal Name: (3S)-2,3-dihydro-4-methyl-3-(1-methylpropyl)-1H-pyrrolo[3,4-b]quinoline-1,9(4H)-dione
MF: C₁₆H₁₈N₂O₂
FW: 270.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium citrinum*



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

Laboratory Procedures

Quinolactacin A is supplied as a solid. A stock solution may be made by dissolving the quinolactacin A in the solvent of choice, which should be purged with an inert gas. Quinolactacin A is soluble in the organic solvent ethanol, at a concentration of approximately 1 mg/ml. Quinolactacin A is also soluble in methanol and DMSO.

Description

Quinolactacin A is a quinolone fungal metabolite originally isolated from *Penicillium*.¹ It inhibits TNF production induced by LPS in murine peritoneal macrophages (IC₅₀ = 12.2 µg/ml). It is not active against a variety of bacteria, fungi, and yeasts. Quinolactacin A is a mixture of quinolactacin A₁ and A₂.

Reference

1. Kakinuma, N., Iwai, H., Takahashi, S., *et al.* Quinolactacins A, B and C: Novel quinolone compounds from *Penicillium* sp. EPF-6. I. Taxonomy, production, isolation and biological properties. *J. Antibiot. (Tokyo)* **53(11)**, 1247-1251 (2000).

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, 08/10/2023

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