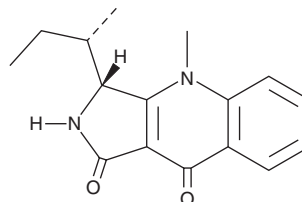


PRODUCT INFORMATION

(+)-Quinolactacin A₁

Item No. 27939

CAS Registry No.: 815576-68-2
Formal Name: (3R)-2,3-dihydro-4-methyl-3-[(1S)-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 A₁ is supplied as a solid. A stock solution may be made by dissolving the (+)-quinolactacin A₁ in the solvent of choice, which should be purged with an inert gas. (+)-Quinolactacin A₁ is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

(+)-Quinolactacin A₁ is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Quinolactacin A₁ is a fungal metabolite originally isolated from *P. citrinum*. It inhibits acetylcholinesterase (AChE) *in vitro* (IC₅₀ = 280 μM).¹

Reference

- Kim, W.-G., Song, N.-K., and Yoo, I.-D. Quinolactacins A1 and A2, new acetylcholinesterase inhibitors from *Penicillium citrinum*. *J. Antibiot. (Tokyo)* **54**(10), 831-835 (2001).

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