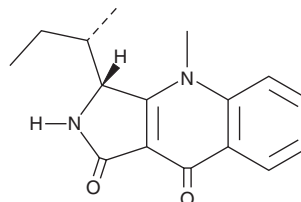


# PRODUCT INFORMATION



## (+)-Quinolactacin A<sub>1</sub> 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:** C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 270.3  
**Purity:** ≥95%  
**Supplied as:** A powder  
**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<sub>1</sub> is supplied as a powder. A stock solution may be made by dissolving the (+)-quinolactacin A<sub>1</sub> in the solvent of choice, which should be purged with an inert gas. (+)-Quinolactacin A<sub>1</sub> is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

(+)-Quinolactacin A<sub>1</sub> 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 A<sub>1</sub> is a fungal metabolite originally isolated from *P. citrinum*. It inhibits acetylcholinesterase (AChE) *in vitro* (IC<sub>50</sub> = 280 μM).<sup>1</sup>

### Reference

1. 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