

# PRODUCT INFORMATION



## Burnettramic Acid A

Item No. 33021

**CAS Registry No.:** 2334483-46-2  
**Formal Name:** (2Z,6R,7aS)-tetrahydro-6-hydroxy-2-  
[(2S,4R,12E,15R,19R)-1,15,19-trihydroxy-  
26-(β-D-mannopyranosyloxy)-2,4-  
dimethyl-14-hexacosen-1-ylidene]-1H-  
pyrrolizine-1,3(2H)-dione

**MF:** C<sub>41</sub>H<sub>71</sub>NO<sub>12</sub>

**FW:** 770.0

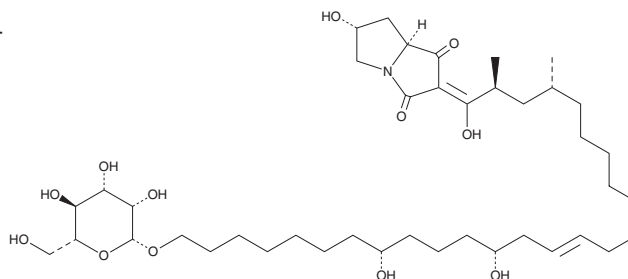
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/*Aspergillus burnettii*



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

### Laboratory Procedures

Burnettramic acid A is supplied as a solid. A stock solution may be made by dissolving the burnettramic acid A in the solvent of choice, which should be purged with an inert gas. Burnettramic acid A is soluble in methanol and DMSO.

### Description

Burnettramic acid A is a fungal metabolite originally isolated from *A. burnettii* that has diverse biological activities.<sup>1</sup> It is active against *B. subtilis*, *S. aureus*, *C. albicans*, and *S. cerevisiae* (IC<sub>50</sub>s = 2.3, 5.9, 0.5, and 0.2 μg/ml, respectively). Burnettramic acid A is cytotoxic to NS-1 murine myeloma cells but not neonatal foreskin fibroblasts (IC<sub>50</sub>s = 13.8 and >100 μg/ml, respectively).

### Reference

- Li, H., Gilchrist, C.L.M., Lacey, H.J., *et al.* Discovery and heterologous biosynthesis of the burnettramic acids: Rare PKS-NRPS-derived bolaamphiphilic pyrrolizidinediones from an Australian fungus, *Aspergillus burnettii*. *Org. Lett.* **21**(5), 1287-1291 (2019).

#### 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, 08/18/2023

#### 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