

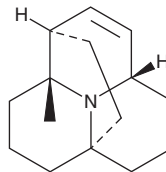
# PRODUCT INFORMATION



## Porantherine

Item No. 31639

**CAS Registry No.:** 33529-61-2  
**Formal Name:** (1R,3aS,6aS,9aR)-3a,4,5,6,7,8,9,9a-octahydro-9a-methyl-1H-1,6a-ethanopyrido[2,1,6-de]quinolizine  
**MF:** C<sub>15</sub>H<sub>23</sub>N  
**FW:** 217.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Poranthera corymbosa*



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

### Laboratory Procedures

Porantherine is supplied as a solid. A stock solution may be made by dissolving the porantherine in the solvent of choice, which should be purged with an inert gas. Porantherine is soluble in organic solvents such as methanol and DMSO.

### Description

Porantherine is an alkaloid that has been found in *Poranthera corymbosa*.<sup>1</sup>

### Reference

1. Denne, W.A. and Mathieson, A.M. The *Poranthera corymbosa* alkaloids, I. Crystal and absolute molecular structure of porantherine hydrobromide. *J. Cryst. Mol. Struct.* **3**, 79-86 (1973).

**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, 11/10/2022

### 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](http://WWW.CAYMANCHEM.COM)