

# PRODUCT INFORMATION



## Drimentine C

Item No. 25763

**CAS Registry No.:** 204398-92-5  
**Formal Name:** (5aS,6aS,11aS,13aS)-6a-[[[(1S,4aS,8aS)-decahydro-5,5,8a-trimethyl-2-methylene-1-naphthalenyl]methyl]-2,3,6,6a,11,11a-hexahydro-1H-pyrrolo[1'',2'':4',5']pyrazino[1',2':1,5]pyrrolo[2,3-b]indole-5,13(5aH,13aH)-dione

**MF:** C<sub>31</sub>H<sub>41</sub>N<sub>3</sub>O<sub>2</sub>

**FW:** 487.7

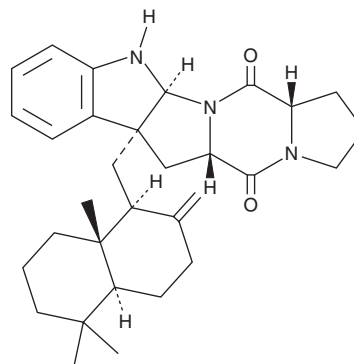
**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

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

### Description

Drimentine C is a terpenylated diketopiperazine antibiotic originally isolated from Actinomycete bacteria.<sup>1</sup> It inhibits proliferation of NS-1 murine β lymphocyte myeloma cells by 63 and 98% *in vitro* when used at concentrations of 12.5 and 100 μg/ml, respectively.

### Reference

1. Lacey, E., Power, M., Wu, Z., *et al.* Terpenylated diketopiperazines, (drimentines). **WO1998009968A1** (1998).

#### 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, 11/28/2022

#### 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