

PRODUCT INFORMATION



Drimentine B

Item No. 25451

CAS Registry No.: 204398-91-4
Formal Name: (5aS,6aS,11aS)-6a-[[[(1S,4aS,8aS)-decahydro-5,5,8a-trimethyl-2-methylene-1-naphthalenyl]methyl]-6,6a,11,11a-tetrahydro-3H-pyrrolo[1'',2':4',5']pyrazino[1',2':1,5]pyrrolo[2,3-b]indole-5,13(2H,5aH)-dione

MF: C₃₁H₃₉N₃O₂

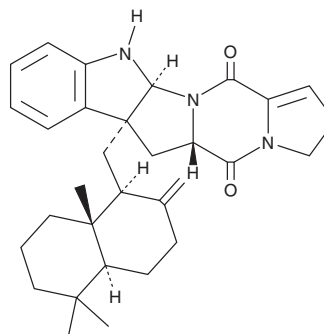
FW: 485.7

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

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

Description

Drimentine B is a terpenylated diketopiperazine originally isolated from *Actinomycete* bacteria that has anticancer activity.¹ It inhibits the proliferation of NS-1 murine β lymphocyte myeloma cells by 41 and 59% *in vitro* when used at concentrations of 50 and 100 $\mu\text{g/ml}$, 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, 01/17/2023

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