

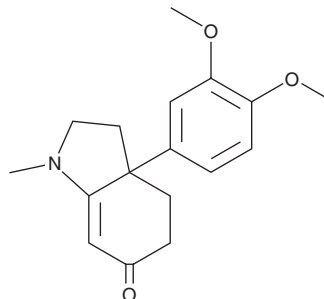
PRODUCT INFORMATION



(±)- Δ^7 -Mesembrenone

Item No. 41086

CAS Registry No.: 80287-15-6
Formal Name: 3a-(3,4-dimethoxyphenyl)-1,2,3,3a,4,5-hexahydro-1-methyl-6H-indol-6-one
MF: $C_{17}H_{21}NO_3$
FW: 287.4
Purity: $\geq 95\%$
Supplied as: A 5 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥ 2 years
Item Origin: Plant/*Sceletium tortuosum*



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

Laboratory Procedures

(±)- Δ^7 -Mesembrenone is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. (±)- Δ^7 -Mesembrenone is sparingly soluble (1-10 mg/ml) in chloroform.

Description

(±)- Δ^7 -Mesembrenone is an alkaloid that has been found in *S. tortuosum*.¹

Reference

1. Reddy, K., Stander, M.A., Stafford, G.I., *et al.* Mass spectrometry metabolomics and feature-based molecular networking reveals population-specific chemistry in some species of the *Sceletium* genus. *Front. Nutr.* **9**, 819753 (2022).

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, 06/15/2026

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