

PRODUCT INFORMATION



Lagochiline

Item No. 13955

CAS Registry No.: 23554-81-6

Formal Name: (4'aS)-decahydro-6'S-hydroxy-5-(2-hydroxyethyl)-2'R,5',8'aS-trimethyl-spiro[furan-2(3H),1'R(2'H)-naphthalene]-5S,5'R-dimethanol

MF: $C_{20}H_{36}O_5$

FW: 356.5

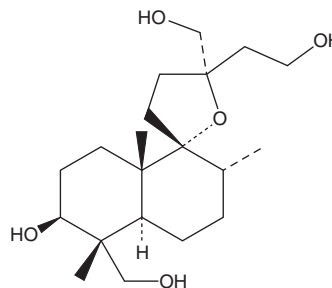
Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Light Sensitive



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

Laboratory Procedures

Lagochiline is supplied as a crystalline solid. A stock solution may be made by dissolving the lagochiline in the solvent of choice, which should be purged with an inert gas. Lagochiline is soluble in the organic solvent methanol at a concentration of approximately 1 mg/ml.

Description

Plants from the mint genus *Lagochilus* have been used in traditional medicines of the east.¹ Lagochiline is a polyphenolic natural product isolated from *Lagochilus* spp. which is thought to provide some of the benefits of the herb. The physiological, immunological, and toxicological properties of this compound remain to be determined. This product is intended for research and forensic purposes.

Reference

1. Furukawa, M., Suzuki, H., Makino, M., *et al.* Studies on the constituents of *Lagochilus leiacanthus* (Labiatae). *Chem. Pharm. Bull.* **59**(12), 1535-1540 (2011).

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, 10/03/2022

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