

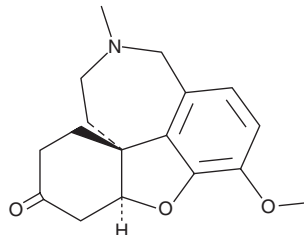
PRODUCT INFORMATION



Dihydronarwedine

Item No. 35409

CAS Registry No.: 21041-10-1
Formal Name: 4aS,5,7,8aR,9,10,11,12-octahydro-3-methoxy-11-methyl-6H-benzofuro[3a,3,2-ef][2]benzazepin-6-one
Synonyms: (-)-Dihydrogalanthaminone, (-)-Lycoraminone
MF: C₁₇H₂₁NO₃
FW: 287.4
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Dihydronarwedine is supplied as a solid. A stock solution may be made by dissolving the dihydronarwedine in the solvent of choice, which should be purged with an inert gas. Dihydronarwedine is soluble in chloroform and DMSO.

Description

Dihydroarwedine is an alkaloid that has been found in *N. pseudonarcissus* and an inhibitor of glycogen synthase kinase 3β (GSK3β).¹ It inhibits recombinant human GSK3β by 39% when used at a concentration of 10 μM.

Reference

1. Hulcová, D., Maříková, J., Korábečný, J., *et al.* Amaryllidaceae alkaloids from *Narcissus pseudonarcissus* L. cv. Dutch Master as potential drugs in treatment of Alzheimer's disease. *Phytochemistry* **165**, 112055 (2019).

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, 12/12/2022

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