

PRODUCT INFORMATION



Cycleanine

Item No. 36020

CAS Registry No.: 518-94-5
Formal Name: (12aR,24aR)-2,3,12a,13,14,15,24,24a-octahydro-5,6,17,18-tetramethoxy-1,13-dimethyl-8,11:20,23-dioxacyclooctadecino[2,3,4-ij:11,12,13-i'j']diisoquinoline

Synonyms: O,O-Dimethylisochondrodendrine, O-Methylnorcycleanine

MF: C₃₈H₄₂N₂O₆

FW: 622.8

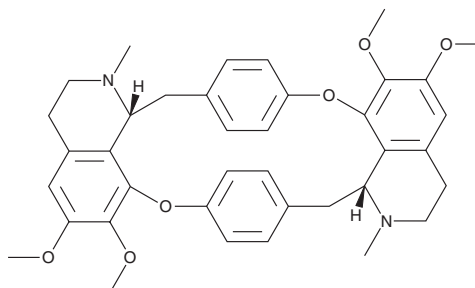
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Stephania japonica* (Thunb.) Miers.



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

Laboratory Procedures

Cycleanine is supplied as a solid. A stock solution may be made by dissolving the cycleanine in the solvent of choice, which should be purged with an inert gas. Cycleanine is soluble in acetone, dichloromethane, DMSO, and ethyl acetate.

Description

Cycleanine is an alkaloid that has been found in *T. subcordata* and has anticancer and antimalarial activities.^{1,2} It selectively inhibits the growth of OVCAR-8, A2780, IGROV-1, and OVCAR-4 ovarian cancer cells (IC₅₀s = 10, 7.6, 14, and 7.2 μM, respectively) over non-cancerous ovarian surface epithelial cells (IC₅₀ = 35 μM).¹ It also halts the cell cycle at the subG₁ phase and induces apoptosis in OVCAR-8 cells. Cycleanine (25 and 50 mg/kg) reduces parasitemia and increases survival in a mouse model of infection with a chloroquine-resistant strain of *P. berghei*.²

References

1. Uche, F.I., Drijfhout, F.P., McCullagh, J., et al. Cytotoxicity effects and apoptosis induction by bisbenzylisoquinoline alkaloids from *Triclisia subcordata*. *Phytother. Res.* **30(9)**, 1533-1539 (2016).
2. Uche, F.I., Guo, X., Okokon, J., et al. *In vivo* efficacy and metabolism of the antimalarial cycleanine and improved *in vitro* antiplasmodial activity of semisynthetic analogues. *Antimicrob. Agents Chemother.* **65(2)**, e01995-20 (2021).

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, 12/19/2022

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