

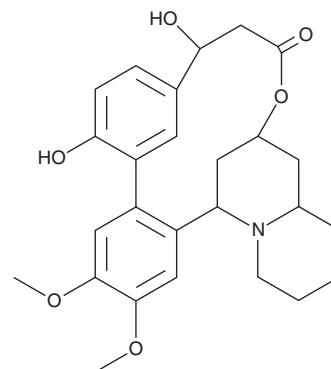
PRODUCT INFORMATION



Lythridine

Item No. 25865

CAS Registry No.: 15299-77-1
Formal Name: 2,3,4,4a,5,6,9,10-octahydro-10,14-dihydroxy-17,18-dimethoxy-20H-6,20-methano-11,15-metheno-1H,8H-benzo[g]pyrido[2,1-d][1,5]oxaazacyclohexadecin-8-one, stereoisomer
Synonym: Sinine
MF: C₂₆H₃₁NO₆
FW: 453.5
Purity: ≥98%
UV/Vis.: λ_{max}: 293, 359 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Lythridine is supplied as a crystalline solid. A stock solution may be made by dissolving the lythridine in the solvent of choice. Lythridine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of lythridine in ethanol is approximately 1 mg/ml, and approximately 20 mg/ml in DMSO and DMF.

Description

Lythridine is a biphenyl quinolizidine lactone alkaloid originally isolated from *Fraxinus sinica* that has antimalarial properties.¹ It has also been found in *H. salicifolia*.²

References

1. Liu, S.-K., Chang, Y.-T., Chuan, T.-K., *et al.* The antimalarial drug, sinine. *Natl. Med. J. China* **27**, 327-341 (1941).
2. Kitakima, M., Yanagisawa, T., Tsukahara, M., *et al.* Biphenyl ether and biphenyl quinolizidine lactone alkaloids from *Heimia salicifolia*. *Tetrahedron* **74(4)**, 441-452 (2018).

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, 07/24/2018

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