

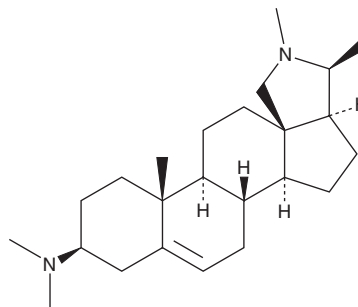
# PRODUCT INFORMATION



## Conessine

Item No. 11700

**CAS Registry No.:** 546-06-5  
**Formal Name:** N,N-dimethyl-con-5-enin-3β-amine  
**Synonyms:** Neriine, NSC 119994, Roquessine, Wrightine  
**MF:** C<sub>24</sub>H<sub>40</sub>N<sub>2</sub>  
**FW:** 356.6  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Conessine is supplied as a crystalline solid. A stock solution may be made by dissolving the Conessine in the solvent of choice, which should be purged with an inert gas. Conessine is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of conessine in these solvents is approximately 16 and 1 mg/ml, respectively.

Conessine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, conessine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Conessine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Conessine is a steroidal alkaloid that has been found in *H. antidysenterica* and has diverse biological activities.<sup>1-5</sup> It is active against a variety of Gram-positive and Gram-negative bacteria, including *B. subtilis*, *S. aureus*, *M. luteus*, and *E. coli* (MICs = 15.6-1,000 µg/disc in agar diffusion assays), the amoeba *E. histolytica*, as well as *P. falciparum* schizonts (IC<sub>50</sub> = 1.9 µg/ml).<sup>1-3</sup> Conessine inhibits replication of the coronaviruses HCoV-OC43, HCoV-NL63, MERS-CoV, and MHV-A59 in infected cells with EC<sub>50</sub> values of 2.34, 10.75, 4.98, and 11.46 µM, respectively.<sup>4</sup> It is also a histamine H<sub>3</sub> receptor antagonist (K<sub>s</sub> = 5.37 and 12.3 nM for the human and rat receptors, respectively).<sup>5</sup> Conessine (10 mg/kg) reduces parasitemia in a mouse model of *P. berghei* infection.<sup>3</sup>

### References

1. Siddiqui, B.S., Ali, S.T., Rizwani, G.H., *et al.* Antimicrobial activity of the methanolic bark extract of *Holarrhena pubescens* (Buch. Ham), its fractions and the pure compound conessine. *Nat. Prod. Res.* **26(11)**, 987-992 (2012).
2. Jayaswal, S.B. Amoebicidal activity of steroidal alkaloids of *Wrightia tomentosa* in vitro. *Indian J. Pharm.* **38(4)**, 112-113 (1976).
3. Dua, V.K., Verma, G., Singh, B., *et al.* Anti-malarial property of steroidal alkaloid conessine isolated from the bark of *Holarrhena antidysenterica*. *Malar. J.* **12**, 194 (2013).
4. Shen, L., Niu, J., Wang, C., *et al.* High-throughput screening and identification of potent broad-spectrum inhibitors of coronaviruses. *J. Virol.* **93(12)**, e00023-19 (2019).
5. Zhao, C., Sun, M., Bennani, Y.L., *et al.* The alkaloid conessine and analogues as potent histamine H<sub>3</sub> receptor antagonists. *J. Med. Chem.* **51(17)**, 5423-5430 (2008).

#### 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, 11/08/2022

#### 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