

PRODUCT INFORMATION

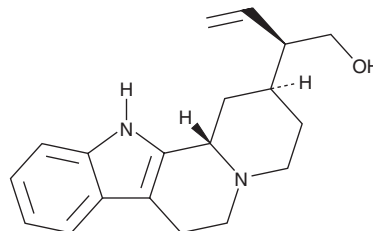


Anthirine

Item No. 31343

CAS Registry No.: 16049-28-8
Formal Name: (β R,2S,12bS)- β -ethenyl-1,2,3,4,6,7,12,12b-octahydro-indolo[2,3-a]quinolizine-2-ethanol

Synonym: Antirhine
MF: C₁₉H₂₄N₂O
FW: 296.4
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Antirhea putaminosa*



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

Laboratory Procedures

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

Description

Anthirine is an alkaloid that has been found in *S. angustiflora* and has analgesic activity.¹ It increases the latency to paw licking in a hot plate test in mice when administered at a dose of 20 mg/kg.

Reference

1. Jiang, H., Liu, Y.-B., Li, Y., et al. Analgesic corynanthe-type alkaloids from *Strychnos angustiflora*. *Tetrahedron* **72**(10), 1276-1284 (2016).

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, 11/04/2022

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