

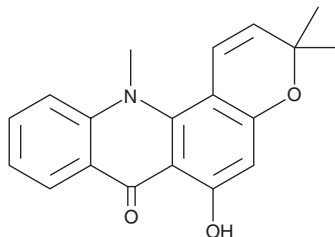
PRODUCT INFORMATION



Noracronycine

Item No. 31336

CAS Registry No.: 13161-79-0
Formal Name: 3,12-dihydro-6-hydroxy-3,3,12-trimethyl-7H-pyrano[2,3-c]acridin-7-one
Synonym: NSC 103005
MF: $C_{19}H_{17}NO_3$
FW: 307.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Acronychia baueri*.



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

Laboratory Procedures

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

Description

Noracronycine is an alkaloid originally isolated from *G. pentaphylla* that has antimalarial activity.^{1,2} It is active against *P. yoelii* when used at a concentration of 10 $\mu\text{g/ml}$.

References

- Govindachari, T.R., Pai, B.R., and Subramaniam, P.S. Alkaloids of *Glycosmis pentaphylla* (Retz.) correa. *Tetrahedron* **22**(10), 3245-3252 (1966).
- Fujioka, H., Nishiyama, Y., Furukawa, H., et al. In vitro and in vivo activities of atalaphilline and related acridone alkaloids against rodent malaria. *Antimicrob. Agents Chemother.* **33**(1), 6-9 (1989).

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