

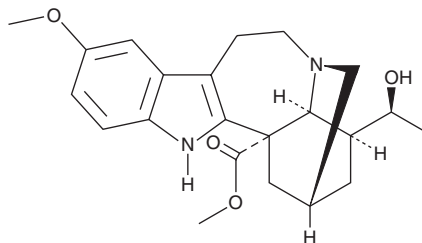
PRODUCT INFORMATION



(-)-Voacangarine

Item No. 25468

CAS Registry No.: 545-84-6
Formal Name: 20S-hydroxy-12-methoxy-ibogamine-18-carboxylic acid, methyl ester
Synonyms: NSC 306219, (-)-Voacristine
MF: C₂₂H₂₈N₂O₄
FW: 384.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 280 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

(-)-Voacangarine is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-voacangarine in the solvent of choice. (-)-Voacangarine is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of (-)-voacangarine in these solvents is approximately 16 mg/ml.

(-)-Voacangarine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (-)-voacangarine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (-)-Voacangarine has a solubility of approximately 0.33 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(-)-Voacangarine is an indole alkaloid originally isolated from *V. africana* that is cytotoxic to HepG2, A375, MDA-MB-231, SH-SY5Y, and CT26 cells (IC₅₀s = 20 µg/ml).^{1,2} It is cytostatic to wild-type *S. cerevisiae* and cytotoxic to $\Delta rad3-e5$ and $\Delta rad6-1$ mutant *S. cerevisiae*, which are defective in DNA strand break repair and the mutagenic repair pathway, respectively.³

References

1. Stauffacher, D. and Seebeck, E. Voacangarine, a new alkaloid from *Voacanga africana*. *Helvetica Chimica Acta* **41(1)**, 169-180 (1958).
2. Chen, H.-M., Yang, Y.-T., Li, H.-X., et al. Cytotoxic monoterpene indole alkaloids isolated from the barks of *Voacanga africana* Staph. *Nat. Prod. Res.* **30(10)**, 1144-1149 (2016).
3. Melo, A.A., Querol, C.B., Henriques, A.T., et al. Cytostatic, cytotoxic and mutagenic effects of voacristine, an indole alkaloid in wild-type and repair-deficient yeasts. *Mutat. Res.* **171(1)**, 17-24 (1986).

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, 05/24/2018

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