

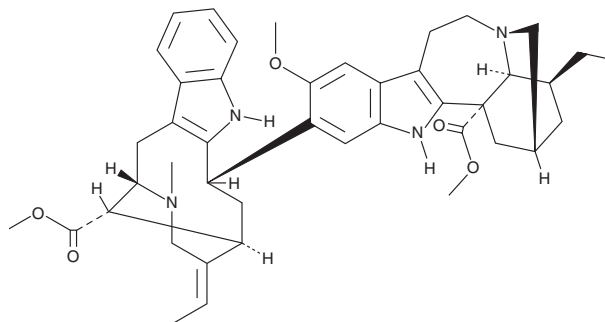
# PRODUCT INFORMATION



## Voacamine

Item No. 25204

**CAS Registry No.:** 3371-85-5  
**Formal Name:** 12-methoxy-13-[(3 $\alpha$ )-17-methoxy-17-oxovobasan-3-yl]-ibogamine-18-carboxylic acid, methyl ester  
**Synonym:** NSC 82591  
**MF:** C<sub>43</sub>H<sub>52</sub>N<sub>4</sub>O<sub>5</sub>  
**FW:** 704.9  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 226, 295 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Description

Voacamine is an alkaloid originally isolated from *Voacanga* that has diverse biological activities.<sup>1</sup> It is a cannabinoid receptor 1 (CB<sub>1</sub>) antagonist (IC<sub>50</sub> = 41 nM).<sup>2</sup> Voacamine (1-10  $\mu$ g/ml) time- and dose-dependently decreases survival of U2OS and multi-drug resistant U2OS (U2OS/DX) osteosarcoma cells and induces autophagy but not apoptosis.<sup>3</sup> It also has antimalarial activity, inhibiting growth of *P. falciparum* *in vitro* at the trophozoite and schizont stages and reducing parasitemia in mice infected with *P. yoelii* when administered at doses of 2.5, 5, and 10 mg/kg.<sup>4</sup>

### References

1. Quevauviller, A., Goutarel, R., and Janot, M.M. Pharmacology of the alkaloids of *Voacanga* species. *Ann. Pharm. Fr.* **13**, 423-424 (1955).
2. Kitajima, M., Iwai, M., Kikura-Hanajiri, R., *et al.* Discovery of indole alkaloids with cannabinoid CB<sub>1</sub> receptor antagonistic activity. *Bioorg. Med. Chem. Lett.* **21(7)**, 1962-1964 (2011).
3. Meschini, S., Condello, M., Calcabrini, A., *et al.* The plant alkaloid voacamine induces apoptosis-independent autophagic cell death on both sensitive and multidrug resistant human osteosarcoma cells. *Autophagy* **4(8)**, 1020-1033 (2008).
4. Ramanitrahasimbola, D., Rasoanaivo, P., Ratsimamanga-Urverg, S., *et al.* Biological activities of the plant-derived bisindole voacamine with reference to malaria. *Phytother. Res.* **15(1)**, 30-33 (2001).

#### 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, 09/30/2022

#### 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