

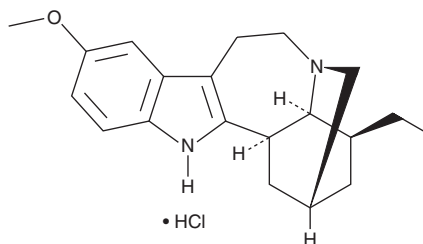
# PRODUCT INFORMATION



## Ibogaine (hydrochloride)

Item No. 20083

**CAS Registry No.:** 5934-55-4  
**Formal Name:** 12-methoxy-ibogamine, monohydrochloride  
**Synonym:** NSC 29847  
**MF:** C<sub>20</sub>H<sub>26</sub>N<sub>2</sub>O • HCl  
**FW:** 346.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 278 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Ibogaine (hydrochloride) (Item No. 20083) is an analytical reference standard categorized as an iboga alkaloid.<sup>1</sup> Ibogaine induces the head-twitch response (HTR), indicating hallucinogenic potential, and anxiety-like behavior in mice.<sup>1,2</sup> It decreases self-administration of morphine (Item Nos. ISO60147 | 15464 | 25955) and induces tremors in rats.<sup>3</sup> Ibogaine use has been associated with fatalities.<sup>4</sup> Ibogaine is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

### References

1. Cameron, L.P., Tombari, R.J., Lu, J., *et al.* A non-hallucinogenic psychedelic analogue with therapeutic potential. *Nature* **589(7842)**, 474-479 (2021).
2. Popik, P. and Wróbel, M. Anxiogenic action of ibogaine. *The Alkaloids*. Glick, S.D., and Alper, K.R., editors, 1st edition, *Academic Press* (2001).
3. Glick, S.D., Rossman, K., Steindorf, S., *et al.* Effects and aftereffects of ibogaine on morphine self-administration in rats. *Eur. J. Pharmacol.* **195(3)**, 341-345 (1991).
4. Alper, K.R., Stajić, M., and Gill, J.R. Fatalities temporally associated with the ingestion of ibogaine. *J. Forensic Sci.* **57(2)**, 398-412 (2012).

#### 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, 09/06/2023

#### 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