

PRODUCT INFORMATION



Noribogaine (hydrochloride)

Item No. 40090

CAS Registry No.: 110514-35-7

Formal Name: (11S)-7-ethyl-6R,6aS,7S,8,9R,10,12,13-octahydro-6,9-methano-5H-pyrido[1',2':1,2]azepino[4,5-b]indol-2-ol, monohydrochloride

Synonyms: O-demethyl Ibogaine, 12-hydroxy Ibogamine

MF: $C_{19}H_{24}N_2O \cdot HCl$

FW: 332.9

Purity: $\geq 95\%$

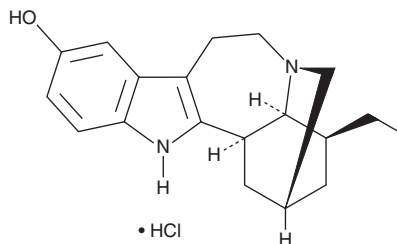
Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years

Special Conditions: Store in amber vials

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



Description

Noribogaine (hydrochloride) (Item No. 40090) is an analytical reference standard categorized as a tryptamine metabolite.¹ Noribogaine is a metabolite of ibogaine (Item Nos. 39570 | 20083). This product is intended for research and forensic applications.

Reference

1. Hearn, W.L., Pablo, J., Him, G.W., *et al.* Identification and quantitation of ibogaine and an o-demethylated metabolite in brain and biological fluids using gas chromatography-mass spectrometry. *J. Anal. Toxicol.* **19**(6), 427-434 (1995).

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, 06/26/2024

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