

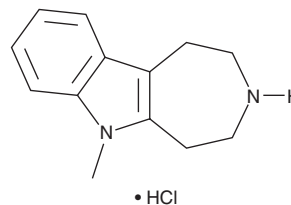
PRODUCT INFORMATION



PNU 22394 (hydrochloride)

Item No. 43564

CAS Registry No.: 15923-42-9
Formal Name: 1,2,3,4,5,6-hexahydro-6-methyl-azepino[4,5-b]indole, monohydrochloride
MF: $C_{13}H_{16}N_2 \cdot HCl$
FW: 236.7
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

PNU 22394 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the PNU 22394 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. PNU 22394 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of PNU 22394 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. PNU 22394 (hydrochloride) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

PNU 22394 is an iboga alkaloid derivative.¹ *In vivo*, PNU 22394 (30 mg/kg) increases licking latency in a cold plate test and paw withdrawal threshold in a paw pressure test in a mouse model of neuropathic pain induced by oxaliplatin (Item No. 13106).

Reference

1. Arias, H.R., Micheli, L., Jensen, A.A., *et al.* Ibogalogs decrease neuropathic pain in mice through a mechanism involving crosstalk between 5-HT_{2A} and mGlu₂ receptors. *Biomed. Pharmacother.* **184**, 117887 (2025).

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, 08/20/2025

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