

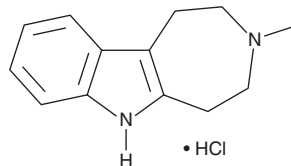
PRODUCT INFORMATION



Ibogaminalog (hydrochloride)

Item No. 46052

CAS Registry No.: 15918-76-0
Formal Name: 1,2,3,4,5,6-hexahydro-3-methyl-azepino[4,5-b]indole, monohydrochloride
Synonym: DM506
MF: $C_{13}H_{16}N_2 \cdot HCl$
FW: 236.7
Purity: $\geq 98\%$
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.

Description

Ibogaminalog (hydrochloride) (Item No. 46052) is an analytical reference standard categorized as an ibogalog.¹ Ibogaminalog induces sedative and anxiolytic-like behavior in stressed and non-stressed mice but does not induce the head-twitch response, indicating a lack of hallucinogenic potential.² It also reduces neuropathic pain in mice and cue-induced drug-seeking in rats after self-administration of heroin (Item Nos. ISO60187 | 9001543 | 9003076).^{1,3} This product is intended for research and forensic applications.

References

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).
2. Arias, H.R., Rudin, D., Hines, D.J., *et al.* The novel non-hallucinogenic compound DM506 (3-methyl-1,2,3,4,5,6-hexahydroazepino[4,5-b]indole) induces sedative-and anxiolytic-like activity in mice by a mechanism involving 5-HT_{2A} receptor activation. *Eur. J. Pharmacol.* **966**, 176329 (2024).
3. Looschen, K., Khatri, S.N., Maulik, M., *et al.* Novel psychoplastogen DM506 reduces cue-induced heroin-seeking and inhibits tonic GABA currents in the Prelimbic Cortex. *Neurochem. Int.* **178**, 105785 (2024).

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, 07/09/2026

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