

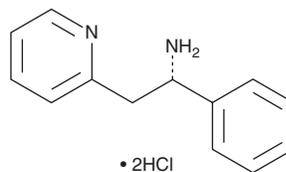
PRODUCT INFORMATION



Lanicemine (hydrochloride)

Item No. 44983

CAS Registry No.: 153322-06-6
Formal Name: α -phenyl-2-pyridineethanamine, dihydrochloride
Synonyms: ARL 15896AR, AR-R 15896AR, AZD 6765, BHV 5500
MF: $C_{13}H_{14}N_2 \cdot 2HCl$
FW: 271.2
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

Lanicemine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the lanicemine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Lanicemine (hydrochloride) is slightly soluble (0.1-1 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 lanicemine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Lanicemine (hydrochloride) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Lanicemine is a low-trapping NMDA receptor antagonist ($IC_{50} = 9.8 \mu M$) and an active metabolite of BHV 5000.^{1,2} Intraperitoneal administration of lanicemine prevents seizures induced by maximal electroshock, 4-aminopyridine (Item No. 18511), N-methyl-DL-aspartate, or kainic acid in mice (ED_{50} s = 13, 17.1, 54, and ~ 50 mg/kg, respectively).³ Lanicemine (24 mg/kg, p.o.) decreases hippocampal neuronal damage in a rat model of global ischemia induced by four-vessel occlusion.⁴ It decreases immobility time in the tail suspension test in mice, an effect that is enhanced by hyperforin (Item No. 19572).⁵

References

1. Mealing, G.A., Lanthorn, T.H., Small, D.L., *et al.* Antagonism of N-methyl-D-aspartate-evoked currents in rat cortical cultures by ARL 15896AR. *J. Pharmacol. Exp. Ther.* **281(1)**, 376-383 (1997).
2. Henter, I.D., Park, L.T., and Zarate, C.A., Jr. Novel glutamatergic modulators for the treatment of mood disorders: Current status. *CNS Drugs* **35(5)**, 527-543 (2021).
3. Palmer, G.C., Murray, R.J., Cramer, C.L., *et al.* [S]-AR-R 15896AR—a novel anticonvulsant: Acute safety, pharmacokinetic and pharmacodynamic properties. *J. Pharmacol. Exp. Ther.* **288(1)**, 121-132 (1999).
4. Cregan, E.F., Peeling, J., Corbett, D., *et al.* [(S)-Alpha-phenyl-2-pyridine-ethanamine dihydrochloride], a low affinity uncompetitive N-methyl-D-aspartic acid antagonist, is effective in rodent models of global and focal ischemia. *J. Pharmacol. Exp. Ther.* **283(3)**, 1412-1424 (1997).
5. Pochwat, B., Szewczyk, B., Kotarska, K., *et al.* Hyperforin potentiates antidepressant-like activity of lanicemine in mice. *Front. Mol. Neurosci.* **11**, 456 (2018).

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, 01/12/2026

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