

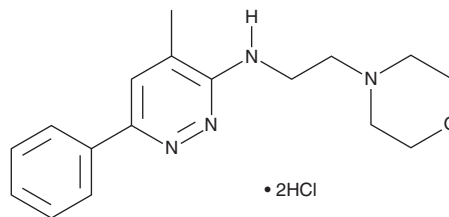
# PRODUCT INFORMATION



## Minaprine (hydrochloride)

Item No. 41710

**CAS Registry No.:** 25905-77-5  
**Formal Name:** N-(4-methyl-6-phenyl-3-pyridazinyl)-4-morpholineethanamine, dihydrochloride  
**Synonyms:** Agr 1240, AG-620  
**MF:**  $C_{17}H_{22}N_4O \cdot 2HCl$   
**FW:** 371.3  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Minaprine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the minaprine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Minaprine (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. Minaprine (hydrochloride) is sparingly soluble (1-10 mg/ml) in ethanol and soluble in DMSO ( $\geq 10$  mg/ml).

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 minaprine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Minaprine (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

Minaprine is an antidepressant.<sup>1</sup> It reduces immobility time in the forced swim test in rats when administered at a dose of 5 mg/kg for nine days. Minaprine (40 mg/kg) induces convulsions in rats.<sup>2</sup> Formulations containing minaprine have previously been used in the treatment of depression.

### References

1. Imperato, A., Obinu, M.C., Cabib, S., *et al.* Effects of subchronic minaprine on dopamine release in the ventral striatum and on immobility in the forced swimming test. *Neurosci. Lett.* **166(1)**, 69-72 (1994).
2. Warter, J.M., Tranchant, C., Marescaux, C., *et al.* Immediate effects of 14 non MAOI antidepressants in rats with spontaneous petit mal-like seizures. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **14(2)**, 261-270 (1990).

#### 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, 10/11/2024

#### 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