

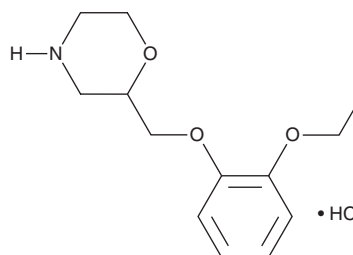
# PRODUCT INFORMATION



## Viloxazine (hydrochloride)

Item No. 39679

**CAS Registry No.:** 35604-67-2  
**Formal Name:** 2-[(2-ethoxyphenoxy)methyl]-morpholine, monohydrochloride  
**Synonym:** ICI 58834  
**MF:**  $C_{13}H_{19}NO_3 \cdot HCl$   
**FW:** 273.8  
**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

Viloxazine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the viloxazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Viloxazine (hydrochloride) is slightly soluble in acetonitrile.

Viloxazine (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Viloxazine is an antidepressant.<sup>1</sup> It selectively inhibits norepinephrine uptake over serotonin (5-HT) uptake in rat hypothalamic synaptosomes ( $IC_{50}$ s = 0.9 and  $>100 \mu M$ , respectively). Viloxazine (8 and 16 mg/kg) decreases immobility time in the forced swim test in mice.<sup>2</sup> Formulations containing viloxazine have been used in the treatment of depression and attention-deficit hyperactivity disorder (ADHD).

### References

1. Blackburn, T.P., Foster, G.A., Greenwood, D.T., *et al.* Effects of viloxazine, its optical isomers and its major metabolites on biogenic amine uptake mechanisms in vitro and in vivo. *Eur. J. Pharmacol.* **52(3-4)**, 367-374 (1978).
2. Bourin, C., Colombel, M.-C., Malinge, M., *et al.* Clonidine as a sensitizing agent in the forced swimming test for revealing antidepressant activity. *J. Psychiatry Neurosci.* **16(4)**, 199-203 (1991).

#### 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, 01/31/2024

#### 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](http://WWW.CAYMANCHEM.COM)