

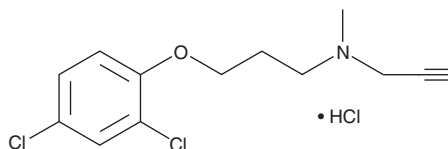
# PRODUCT INFORMATION



## Clorgyline (hydrochloride)

Item No. 15925

**CAS Registry No.:** 17780-75-5  
**Formal Name:** N-[3-(2,4-dichlorophenoxy)propyl]-N-methyl-2-propyn-1-amine, monohydrochloride  
**Synonym:** N-2,4-Dichlorophenoxypropyl-N-methylpropargylamine  
**MF:** C<sub>13</sub>H<sub>15</sub>Cl<sub>2</sub>NO • HCl  
**FW:** 308.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 284 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Clorgyline (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the clorgyline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Clorgyline (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of clorgyline (hydrochloride) in ethanol and DMF is approximately 15 mg/ml and approximately 10 mg/ml in DMSO.

Clorgyline (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, clorgyline (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Clorgyline (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Clorgyline is a potent monoamine oxidase (MAO) inhibitor that preferentially targets MAO-A over MAO-B (K<sub>i</sub>s = 0.054 and 58 μM, respectively).<sup>1,2</sup> This inhibition is irreversible.<sup>1</sup> Clorgyline is without effect on serotonin (Item No. 14332) receptors, but it does inhibit the sigma receptor σ<sub>1</sub> on Jurkat human T lymphocyte cells (IC<sub>50</sub> = 31 nM).<sup>3,4</sup> As MAO-A preferentially oxidizes serotonin, selective MAO-A inhibitors have applications in ameliorating depression and anxiety.<sup>1,5</sup>

### References

1. Abdelhafez, O.M., Amin, K.M., Ali, H.I., *et al.* Synthesis of new 7-oxycoumarin derivatives as potent and selective monoamine oxidase A inhibitors. *J. Med. Chem.* **55(23)**, 10424-10436 (2012).
2. Mazouz, F., Gueddari, S., Burstein, C., *et al.* 5-[4-(Benzyloxy)phenyl]-1,3,4-oxadiazol-2(3H)-one derivatives and related analogues: New reversible, highly potent, and selective monoamine oxidase type B inhibitors. *J. Med. Chem.* **36(9)**, 1157-1167 (1993).
3. Pälvimäki, E.P., Roth, B.L., Majasuo, H., *et al.* Interactions of selective serotonin reuptake inhibitors with the serotonin 5-HT<sub>2C</sub> receptor. *Psychopharmacology (Berl)* **126(3)**, 234-240 (1996).
4. Ganapathy, M.E., Prasad, P.D., Huang, W., *et al.* Molecular and ligand-binding characterization of the σ-receptor in the Jurkat human T lymphocyte cell line. *J. Pharmacol. Exp. Ther.* **289(1)**, 251-260 (1999).
5. La Regina, G., Silvestri, R., Gatti, V., *et al.* Synthesis, structure-activity relationships and molecular modeling studies of new indole inhibitors of monoamine oxidases A and B. *Bioorg. Med. Chem.* **16(22)**, 9729-9740 (2008).

#### 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, 12/21/2022

#### 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