

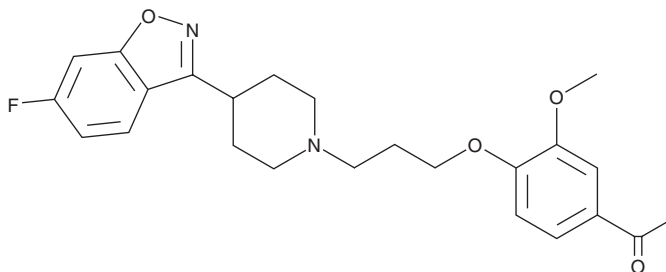
# PRODUCT INFORMATION



## Iloperidone

Item No. 22957

**CAS Registry No.:** 133454-47-4  
**Formal Name:** 1-[4-[3-[4-(6-fluoro-1,2-benzisoxazol-3-yl)-1-piperidinyl]propoxy]-3-methoxyphenyl]-ethanone  
**MF:** C<sub>24</sub>H<sub>27</sub>FN<sub>2</sub>O<sub>4</sub>  
**FW:** 426.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 275, 304 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

Iloperidone is supplied as a crystalline solid. A stock solution may be made by dissolving the iloperidone in the solvent of choice. Iloperidone is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of iloperidone in these solvents is approximately 0.1 and 0.2 mg/ml, respectively.

### Description

Iloperidone is an atypical antipsychotic and adrenergic, dopamine, and serotonin (5-HT) receptor antagonist.<sup>1</sup> It binds to several receptors, including the α<sub>1</sub>-adrenergic receptor (α<sub>1</sub>-AR), α<sub>2</sub>-AR, and dopamine D<sub>2</sub> receptor (K<sub>s</sub> = 0.31, 3, and 3.3 nM, respectively), as well as the 5-HT<sub>1A</sub>, 5-HT<sub>1D</sub>, 5-HT<sub>2A</sub>, and 5-HT<sub>2C</sub> receptors (K<sub>s</sub> = 33, 15, 0.2, and 14 nM, respectively) in radioligand binding assays using human post-mortem brain tissue.<sup>2</sup> Iloperidone also binds to human D<sub>1</sub>, D<sub>3</sub>, D<sub>4</sub>, D<sub>5</sub>, and rat 5-HT<sub>2</sub> receptors (K<sub>s</sub> = 216, 7.1, 25, 319, and 3.1 nM, respectively, in CHO cells) and the histamine H<sub>1</sub> receptor (K<sub>i</sub> = 12.3 nM in human post-mortem brain tissue).<sup>2,3</sup> Iloperidone (1-3 mg/kg) prevents the reduction in prepulse inhibition induced by apomorphine (Item No. 16094), phencyclidine (PCP), and cirazoline (Item No. 21791) in rats.<sup>1</sup> It also increases the time rats spend in the open arms of the elevated plus maze and the number of social interactions when administered at a dose of 0.5 mg/kg.<sup>4</sup> Formulations containing iloperidone have been used in the treatment of schizophrenia.

### References

1. Barr, A.M., Powell, S.B., Markou, A., *et al.* Iloperidone reduces sensorimotor gating deficits in pharmacological models, but not a developmental model, of disrupted prepulse inhibition in rats. *Neuropharmacology* **51(3)**, 457-465 (2006).
2. Richelson, E. and Souder, T. Binding of antipsychotic drugs to human brain receptors focus on newer generation compounds. *Life Sciences* **68(1)**, 29-39 (2000).
3. Kongsamut, S., Roehr, J.E., Cai, J., *et al.* Iloperidone binding to human and rat dopamine and 5-HT receptors. *Eur. J. Pharmacol.* **317(2-3)**, 417-423 (1996).
4. Szewczak, M.R., Corbett, R., Rush, D.K., *et al.* The pharmacological profile of iloperidone, a novel atypical antipsychotic agent. *J. Pharmacol. Exp. Ther.* **274(3)**, 1404-1413 (1995).

#### 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, 10/31/2022

#### 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