

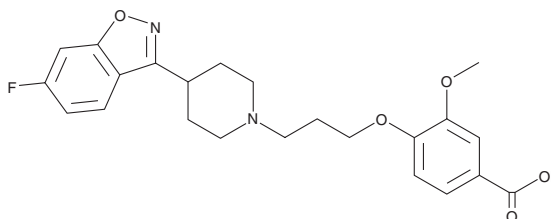
# PRODUCT INFORMATION



## Iloperidone metabolite P95

Item No. 27807

**CAS Registry No.:** 475110-48-6  
**Formal Name:** 4-[3-[4-(6-fluoro-1,2-benzisoxazol-3-yl)-1-piperidinyl]propoxy]-3-methoxybenzoic acid  
**Synonym:** P95-12113  
**MF:** C<sub>23</sub>H<sub>25</sub>FN<sub>2</sub>O<sub>5</sub>  
**FW:** 428.5  
**Purity:** ≥98%  
**Supplied as:** A 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 metabolite P95 is supplied as a solid. A stock solution may be made by dissolving the iloperidone metabolite P95 in the solvent of choice, which should be purged with an inert gas. Iloperidone metabolite P95 is soluble in the organic solvent DMSO.

### Description

Iloperidone metabolite P95 is a metabolite of the atypical antipsychotic iloperidone (Item No. 22957).<sup>1,2</sup> It binds to the serotonin (5-HT) receptor subtype 5-HT<sub>2A</sub> and α<sub>1</sub>-, α<sub>2B</sub>-, and α<sub>2C</sub>-adrenergic receptors with mean K<sub>i</sub> values of 7.08, 21.38, 83.18, and 47.86 nM, respectively, but does not cross the blood-brain barrier.<sup>1</sup>

### References

1. Subramanian, N. and Kalkman, H.O. Receptor profile of P88-8991 and P95-12113, metabolites of the novel antipsychotic iloperidone. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **26**(3), 553-560 (2002).
2. Jia, M., Li, J., He, X., *et al.* Simultaneous determination of iloperidone and its two active metabolites in human plasma by liquid chromatography-tandem mass spectrometry: application to a pharmacokinetic study. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **928**, 52-57 (2013).

#### 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, 12/22/2022

#### 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