

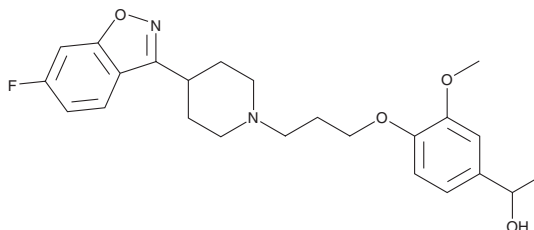
PRODUCT INFORMATION



Iloperidone metabolite P88

Item No. 27806

CAS Registry No.: 133454-55-4
Formal Name: 4-[3-[4-(6-fluoro-1,2-benzisoxazol-3-yl)-1-piperidinyl]propoxy]-3-methoxy- α -methyl-benzenemethanol
Synonym: P88-8991
MF: C₂₄H₂₉FN₂O₄
FW: 428.5
Purity: $\geq 98\%$
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 P88 is supplied as a solid. A stock solution may be made by dissolving the iloperidone metabolite P88 in the solvent of choice, which should be purged with an inert gas. Iloperidone metabolite P88 is slightly soluble in methanol and chloroform.

Description

Iloperidone metabolite P88 is an active metabolite of the atypical antipsychotic iloperidone (Item No. 22957).^{1,2} It binds to the serotonin (5-HT) receptor subtypes 5-HT_{1B}, 5-HT_{2A}, and 5-HT_{2C} (K_s = 5.1, 0.03, and 6.9 nM, respectively), α_{2b} - and α_{2c} -adrenergic (K_s = 6 and 1.3 nM, respectively), and dopamine D₁, D_{2A}, and D₄ receptors (K_s = 9.5, 1.6, and 3.5 nM, respectively).¹

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, 11/03/2022

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