

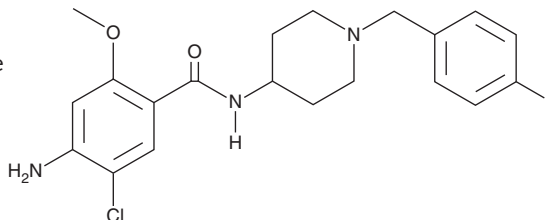
PRODUCT INFORMATION



Fluoroclebopride

Item No. 23447

CAS Registry No.: 154540-49-5
Formal Name: 4-amino-5-chloro-N-[1-[(4-fluorophenyl)methyl]-4-piperidiny]-2-methoxy-benzamide
MF: C₂₀H₂₃ClFN₃O₂
FW: 391.9
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 274, 309 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

Fluoroclebopride is supplied as a crystalline solid. A stock solution may be made by dissolving the fluoroclebopride in the solvent of choice, which should be purged with an inert gas. Fluoroclebopride is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fluoroclebopride in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

Fluoroclebopride is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fluoroclebopride should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Fluoroclebopride has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fluoroclebopride is a benzamide analog that is used in positron emission tomography (PET) applications.^{1,2} It binds reversibly to dopamine receptors (K_s = 0.95, 5.7, 5.46, and 144 nM for D₂-like, D_{2(long)}, D₃, and D₄ receptors, respectively, in radioligand binding assays).³ It is selective for these receptors over D₁, serotonin 5-HT₂, and α₂-adrenergic receptors (K_s = >10,000, 283, and 1,300 nM, respectively).^{4,5} A fluorine-18 moiety has been used to label this compound for use as a probe for studying D₂/D₃ receptor availability via PET in various monkey models.^{1,2}

References

1. Nader, M.A. and Czoty, P.W. PET imaging of dopamine D2 receptors in monkey models of cocaine abuse: Genetic predisposition versus environmental modulation. *Am. J. Psychiatry* **162(8)**, 1473-1482 (2005).
2. Czoty, P.W., Gage, H.D., Garg, P.K., et al. Effects of repeated treatment with the dopamine D2/D3 receptor partial agonist aripiprazole on striatal D2/D3 receptor availability in monkeys. *Psychopharmacol. (Berl)*. **231(3)**, 613-619 (2013).
3. Mach, R.H., Nader, M.A., Ehrenkauf, R.L., et al. Comparison of two fluorine-18 labeled benzamide derivatives that bind reversibly to dopamine D₂ receptors: In vitro binding studies and positron emission tomography. *Synapse* **24(4)**, 322-333 (1996).
4. Mach, R.H., Elder, S.T., Morton, T.E., et al. The use of [¹⁸F]4-fluorobenzyl iodide (FBI) in PET radiotracer synthesis: Model alkylation studies and its application in the design of dopamine D₁ and D₂ receptor-based imaging agents. *Nucl. Med. Biol.* **20(6)**, 777-794 (1993).
5. Mach, R.H., Luedtke, R.R., Unsworth, C.D., et al. ¹⁸F-labeled benzamides for studying the dopamine D₂ receptor with positron emission tomography. *J. Med. Chem.* **36(23)**, 3707-3720 (1993).

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, 11/15/2022

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