

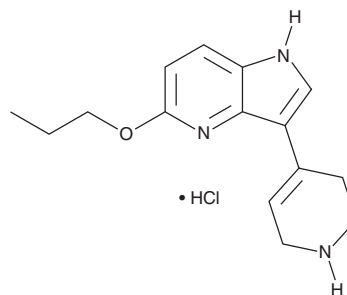
# PRODUCT INFORMATION



## CP 94,253 (hydrochloride)

Item No. 29486

**CAS Registry No.:** 845861-39-4  
**Formal Name:** 5-propoxy-3-(1,2,3,6-tetrahydro-4-pyridinyl)-1H-pyrrolo[3,2-b]pyridine, monohydrochloride  
**MF:** C<sub>15</sub>H<sub>19</sub>N<sub>3</sub>O • HCl  
**FW:** 293.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220, 258, 300 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.

### Description

CP 94,253 is an agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>1B</sub> (K<sub>i</sub> = 2 nM in a radioligand binding assay).<sup>1</sup> It is selective for 5-HT<sub>1B</sub> over 5-HT<sub>1A</sub>, 5-HT<sub>1D</sub>, 5-HT<sub>1C</sub>, and 5-HT<sub>2</sub> receptors (K<sub>i</sub>s = 89, 49, 860, and 1,600 nM, respectively). CP 94,253 (10 and 17 mg/kg) reduces aggressive behavior in a resident-intruder test in mice, as well as decreases alcohol-induced aggression in mice with an ED<sub>50</sub> value of 3.8 mg/kg.<sup>2</sup> It decreases the time mice spend immobile in the forced swim test by 43%, indicating antidepressant-like activity, when administered at a dose of 5 mg/kg.<sup>3</sup> CP 94,253 (2.5 mg/kg) also increases drinking in the Vogel punished drinking task, indicating anxiolytic-like activity, in rats.<sup>4</sup>

### References

1. Koe, B.K., Nielsen, J.A., Macor, J.E., *et al.* Biochemical and behavioral studies of the 5-HT<sub>1B</sub> receptor agonist, CP-94,253. *Drug Dev. Res.* **26(3)**, 241-250 (1992).
2. Fish, E.W., Faccidomo, S., and Miczek, K.A. Aggression heightened by alcohol or social instigation in mice: Reduction by the 5-HT<sub>1B</sub> receptor agonist CP-94,253. *Psychopharmacology (Berl.)* **146(4)**, 391-399 (1999).
3. Tatarczyńska, E., Antkiewicz-Michaluk, L., Kłodzińska, A., *et al.* Antidepressant-like effect of the selective 5-HT<sub>1B</sub> receptor agonist CP 94253: A possible mechanism of action. *Eur. J. Pharmacol.* **516(1)**, 46-50 (2005).
4. Chojnacka-Wójcik, E., Kłodzińska, A., and Tatarczyńska, E. The anxiolytic-like effect of 5-HT<sub>1B</sub> receptor ligands in rats: A possible mechanism of action. *J. Pharm. Pharmacol.* **57(2)**, 253-257 (2005).

#### 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, 11/28/2022

#### 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