

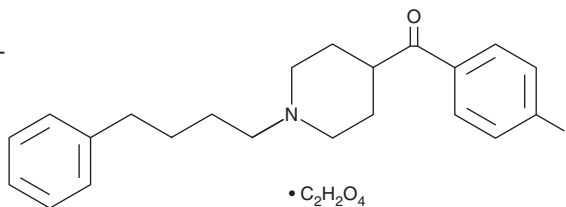
# PRODUCT INFORMATION



## 4F 4PP (oxalate)

Item No. 35005

**CAS Registry No.:** 144734-36-1  
**Formal Name:** (4-fluorophenyl)  
[1-(4-phenylbutyl)-4-piperidiny]-  
methanone, monoethanedioate  
**MF:** C<sub>22</sub>H<sub>26</sub>FNO • C<sub>2</sub>H<sub>2</sub>O<sub>4</sub>  
**FW:** 429.5  
**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

4F 4PP (oxalate) is supplied as a solid. A stock solution may be made by dissolving the 4F 4PP (oxalate) in the solvent of choice, which should be purged with an inert gas. 4F 4PP (oxalate) is soluble in organic solvents such as ethanol and DMSO. The solubility of 4F 4PP (oxalate) in these solvents is approximately 20 and 100 mM, respectively.

### Description

4F 4PP is an antagonist of the serotonin (5-HT) receptor subtype 5-HT<sub>2A</sub> (K<sub>i</sub> = 5.3 nM).<sup>1</sup> It is selective for 5-HT<sub>2A</sub> receptors over 5-HT<sub>1C</sub> receptors (K<sub>i</sub> = 620 nM). It inhibits 5-HT-induced currents in *Xenopus* oocytes expressing human 5-HT<sub>2A</sub> receptors (IC<sub>50</sub> = 5.15 μM).<sup>2</sup> Spinal superfusion of 4F 4PP (100 μM) reduces the area of electrically stimulated spinal field potentials in a rat model of neuropathic pain induced by spinal nerve ligation (SNL).<sup>3</sup>

### References

1. Herndon, J.L., Ismaiel, A., Ingher, S.P., *et al.* Ketanserin analogues: Structure-affinity relationships for 5-HT<sub>2</sub> and 5-HT<sub>1C</sub> serotonin receptor binding. *J. Med. Chem.* **35(26)**, 4903-4910 (1992).
2. Acuña-Castillo, C., Villalobos, C., Moya, P.R., *et al.* Differences in potency and efficacy of a series of phenylisopropylamine/phenylethylamine pairs at 5-HT<sub>2A</sub> and 5-HT<sub>2C</sub> receptors. *Br. J. Pharmacol.* **136(4)**, 510-519 (2002).
3. Aira, Z., Buesa, I., Salgueiro, M., *et al.* Subtype-specific changes in 5-HT receptor-mediated modulation of C fibre-evoked spinal field potentials are triggered by peripheral nerve injury. *Neuroscience* **168(3)**, 831-841 (2010).

#### 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, 12/09/2022

#### 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