

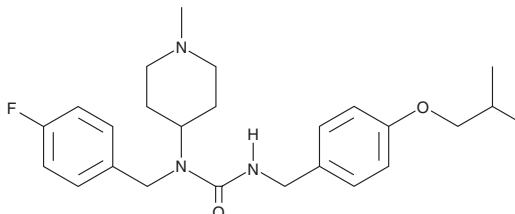
# PRODUCT INFORMATION



## Pimavanserin

Item No. 23462

**CAS Registry No.:** 706779-91-1  
**Formal Name:** N-[(4-fluorophenyl)methyl]-N-(1-methyl-4-piperidinyl)-N'-[[4-(2-methylpropoxy)phenyl]methyl]-urea  
**MF:** C<sub>25</sub>H<sub>34</sub>FN<sub>3</sub>O<sub>2</sub>  
**FW:** 427.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 226 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

Pimavanserin is supplied as a crystalline solid. A stock solution may be made by dissolving the pimavanserin in the solvent of choice. Pimavanserin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of pimavanserin in these solvents is approximately 33 mg/ml.

Pimavanserin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pimavanserin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Pimavanserin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pimavanserin is an inverse agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>2A</sub> (IC<sub>50</sub> = 1.86 nM; K<sub>i</sub> = 0.5 nM).<sup>1</sup> It is selective for 5-HT<sub>2A</sub> over a panel of 65 ion channels, enzymes, and receptors (K<sub>i</sub>s = >100 nM). Pimavanserin reduces head-twitch behavior and prepulse inhibition deficits induced by the 5-HT<sub>2A</sub> receptor agonist DOI (Item No. 13885) in rats at doses of 3 mg/kg, p.o., and 1-10 mg/kg, s.c., respectively. It also exhibits antipsychotic-like activity, reducing hyperactivity induced by (+)-MK-801 (Item No. 10009019) in mice. Pimavanserin acts synergistically with haloperidol (Item No. 12014) or risperidone (Item No. 13629) to suppress (+)-MK-801-induced hyperactivity and attenuates haloperidol- and risperidone-induced catalepsy in rats.<sup>2</sup> Formulations containing pimavanserin have been used for the treatment of psychosis in Parkinson's disease.<sup>3</sup>

### References

1. Vanover, K.E., Weiner, D.M., Makhay, M., *et al.* Pharmacological and behavioral profile of N-(4-fluorophenylmethyl)-N-(1-methylpiperidin-4-yl)-N'-((4-(2-methylpropoxy)phenyl)methyl) carbamide (2R,3R)-dihydroxybutanedioate (2:1) (ACP-103), a novel 5-hydroxytryptamine<sub>2A</sub> receptor inverse agonist. *J. Pharmacol. Exp. Ther.* **318**(2), 910-918 (2006).
2. Gardell, L.R., Vanover, K.E., Pounds, L., *et al.* ACP-103, a 5-hydroxytryptamine 2A receptor inverse agonist, improves the antipsychotic efficacy and side-effect profile of haloperidol and risperidone in experimental models. *J. Pharmacol. Exp. Ther.* **322**(2), 862-870 (2007).
3. Goldman, J.G. and Holden, S. Treatment of psychosis and dementia in Parkinson's disease. *Curr. Treat. Options. Neurol.* **16**(3), 281 (2014).

#### 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