

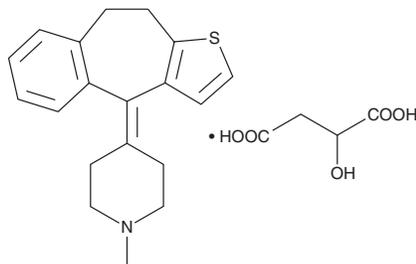
# PRODUCT INFORMATION



## Pizotifen (malate)

Item No. 20765

**CAS Registry No.:** 5189-11-7  
**Formal Name:** 4-(9,10-dihydro-4H-benzo[4,5]cyclohepta[1,2-b]thien-4-ylidene)-1-methylpiperidine, 2-hydroxybutanedioic acid  
**MF:** C<sub>19</sub>H<sub>21</sub>NS • C<sub>4</sub>H<sub>6</sub>O<sub>5</sub>  
**FW:** 429.5  
**Purity:** ≥98%  
**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

Pizotifen (malate) is supplied as a crystalline solid. A stock solution may be made by dissolving the pizotifen (malate) in the solvent of choice, which should be purged with an inert gas. Pizotifen (malate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pizotifen (malate) in these solvents is approximately 0.5, 15, and 20 mg/ml, respectively.

Pizotifen (malate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pizotifen (malate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pizotifen (malate) has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pizotifen is a triptan with high affinity for serotonin (5-HT) and dopamine receptors ( $K_{iS} = 7.94$  and 2.4 nM for 5-HT<sub>2C</sub> and D<sub>2</sub>, respectively).<sup>1,2</sup> It is used for the prophylaxis of migraine headaches.<sup>3</sup>

### References

1. Glusa, E. and Pertz, H.H. Further evidence that 5-HT-induced relaxation of pig pulmonary artery is mediated by endothelial 5-HT<sub>2B</sub> receptors. *Br. J. Pharmacol.* **130(3)**, 692-698 (2000).
2. Morimoto, A., Yamada, H., Watanabe, T., *et al.* Synthesis and structure-activity relationships of new carbonyl guanidine derivatives as novel dual 5-HT<sub>2B</sub> and 5-HT<sub>7</sub> receptor antagonists. *Bioorg. Med. Chem.* **21(24)**, 7841-7852 (2013).
3. Pringsheim, T., Davenport, W.J., and Becker, W.J. Prophylaxis of migraine headache. *CMAJ* **182(7)**, E269-E276 (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, 11/11/2022

#### 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