

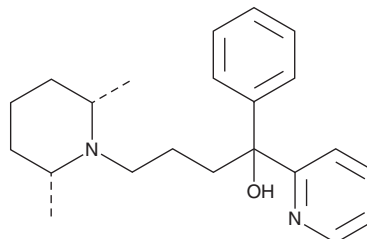
# PRODUCT INFORMATION



## Pirmenol

Item No. 29042

**CAS Registry No.:** 68252-19-7  
**Formal Name:** *rel*- $\alpha$ -[3-[(2R,6S)-2,6-dimethyl-1-piperidinyl]propyl]- $\alpha$ -phenyl-2-pyridinemethanol  
**Synonym:** ( $\pm$ )-Pirmenol  
**MF:** C<sub>22</sub>H<sub>30</sub>N<sub>2</sub>O  
**FW:** 338.5  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Pirmenol is supplied as a solid. A stock solution may be made by dissolving the pirmenol in the solvent of choice, which should be purged with an inert gas. Pirmenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of pirmenol in these solvents is approximately 10, 5, and 2 mg/ml, respectively.

Pirmenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pirmenol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Pirmenol has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pirmenol is a class I antiarrhythmic agent.<sup>1-3</sup> It lengthens the final repolarization of action potentials in isolated rabbit ventricular myocytes and inhibits tachyarrhythmias induced by prolonged afterdepolarizations in isolated guinea pig papillary muscles when used at a concentration of 5  $\mu$ M.<sup>1</sup> Pirmenol (2.5 and 5 mg/kg) suppresses ouabain- or epinephrine-induced ventricular arrhythmias and aconitine-induced atrial arrhythmias in dogs.<sup>2,3</sup>

### References

1. Sawanobori, T., Adaniya, H., Yamashita, S., *et al.* Electrophysiologic and antiarrhythmic actions of pirmenol on rabbit and guinea pig cardiac preparations. *J. Cardiovasc. Pharmacol.* **16(6)**, 975-983 (1990).
2. Kaplan, H.R., Mertz, T.E., and Steffe, T.J. Preclinical pharmacology of pirmenol. *Am. J. Cardiol.* **59(16)**, 2H-9H (1987).
3. Steffe, T.J., Mertz, T.E., Hastings, S.G., *et al.* CL-845 (pirmenol hydrochloride): A new orally effective long-acting antiarrhythmic agent. *J. Pharmacol. Exp. Ther.* **214(1)**, 50-57 (1980).

#### 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, 11/07/2022

#### 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