

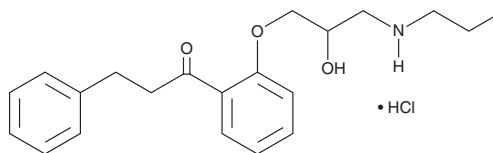
PRODUCT INFORMATION



Propafenone (hydrochloride)

Item No. 21871

CAS Registry No.: 34183-22-7
Formal Name: 1-[2-[2-hydroxy-3-(propylamino)propoxy]phenyl]-3-phenyl-1-propanone, monohydrochloride
MF: C₂₁H₂₇NO₃ • HCl
FW: 377.9
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 247, 303 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

Propafenone (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the propafenone (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Propafenone (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of propafenone (hydrochloride) in these solvents is approximately 0.3, 30, and 20 mg/ml, respectively.

Propafenone (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, propafenone (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Propafenone (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Propafenone is a class I antiarrhythmic agent.¹ It inhibits calcium currents (I_{Ca}) in isolated guinea pig ventricular myocytes ($IC_{50} = 5 \mu M$), as well as ATP-sensitive potassium currents (IK_{ATP}) in isolated rabbit atrial and ventricular myocytes (IC_{50} s = 1.26 and 4.94 μM , respectively).^{1,2} Propafenone also blocks sodium (I_{Na}), potassium (I_K), and transient outward potassium currents in various cardiac cells.²⁻⁴ *In vivo*, propafenone (2 and 3 mg/kg, i.v.) reverses sinus rhythm in a dog model of barium chloride-induced ventricular arrhythmia.⁵ Formulations containing propafenone have been used in the treatment of cardiac arrhythmias.

References

1. Delgado, C., Tamargo, J., Henzel D., *et al.* Effects of propafenone on calcium current in guinea-pig ventricular myocytes. *Br. J. Pharmacol.* **108**(3), 721-727 (1993).
2. Christé, G., Tebbakh, H., Šimurdová, M., *et al.* Propafenone blocks ATP-sensitive K⁺ channels in rabbit atrial and ventricular cardiomyocytes. *Eur. J. Pharmacol.* **373**(2-3), 223-232 (1999).
3. Delpón, E., Valenzuela, C.F., Pérez, O., *et al.* Propafenone preferentially blocks the rapidly activating component of delayed rectifier K⁺ current in guinea pig ventricular myocytes. Voltage-independent and time-dependent block of the slowly activating component. *Circ. Res.* **76**(2), 223-235 (1995).
4. Simó-Vicens, R., Sauter, D.R.P., Grunnet, M., *et al.* Effect of antiarrhythmic drugs on small conductance calcium - activated potassium channels. *Eur. J. Pharmacol.* (2017).
5. Germiniani, H., and Germiniani, C.d.L.B. Effect of propafenone hydrochloride in ventricular arrhythmias experimentally induced by barium chloride. *Arq. Bras. Cardiol.* **36**(2), 101-105 (1981).

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, 10/31/2022

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