

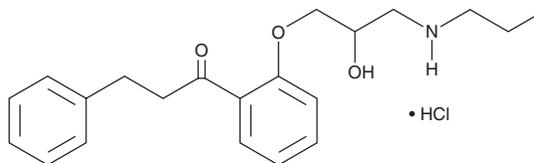
PRODUCT INFORMATION



Propafenone (hydrochloride)

Item No. 44945

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%
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.

Description

Propafenone (hydrochloride) (Item No. 44945) is an analytical reference standard categorized as a class IC antiarrhythmic agent.¹ Propafenone has been associated with accidental and intentional overdose.¹⁻³ Formulations containing propafenone have been used in the treatment of cardiac arrhythmias. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 21871).

References

1. Zeljković, I., Bulj, N., Kolačević, M., *et al.* Failure of intracardiac pacing after fatal propafenone overdose: A case report. *J. Emerg. Med.* **54(4)**, e65-e68 (2018).
2. D'Orazio, L. and Curtis, J.A. Overdose of propafenone surreptitiously sold as "Percocet". *J. Emerg. Med.* **41(2)**, 172-175 (2011).
3. Keramari, S., Poutoglidis, A., Poutoglidou, F., *et al.* Propafenone poisoning of a female adolescent after a suicide attempt. *Cureus* **13(7)**, e16576 (2021).

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, 01/30/2026

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