

PRODUCT INFORMATION

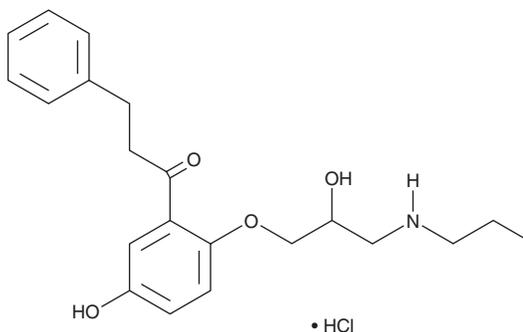


5-Hydroxypropafenone (hydrochloride)

Item No. 43078

CAS Registry No.: 86383-32-6
Formal Name: 1-[5-hydroxy-2-[2-hydroxy-3-(propylamino)propoxy]phenyl]-3-phenyl-1-propanone, monohydrochloride

Synonym: Lu 40545
MF: C₂₁H₂₇NO₄ • HCl
FW: 393.9
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.

Laboratory Procedures

5-Hydroxypropafenone (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxypropafenone (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 5-Hydroxypropafenone (hydrochloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

5-Hydroxypropafenone (hydrochloride) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

5-Hydroxypropafenone is an active metabolite of propafenone (Item No. 21871).¹ It inhibits transient outward potassium current (I_{to}) in isolated human atrial myocytes ($IC_{50} = 1.5 \mu M$). *In vivo*, 5-hydroxypropafenone (0.5, 1, and 2 mg/kg) increases the atrial effective refractory period in dogs.²

References

1. Cahill, S.A., Kirshenbaum, L.A., and Gross, G.J. Transient outward current inhibition by propafenone and 5-hydroxypropafenone in cultured neonatal rat ventricular myocytes. *J. Cardiovasc. Pharmacol.* **38(3)**, 460-467 (2001).
2. Boucher, M., Chassaing, C., Hamel, J.D., *et al.* Cardiac electrophysiological effects of propafenone and its 5-hydroxylated metabolite in the conscious dog. *Eur. J. Pharmacol.* **314(2)**, 171-177 (1996).

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, 05/06/2025

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