

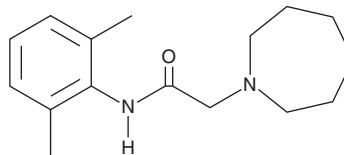
PRODUCT INFORMATION



Pincaïnide

Item No. 45339

CAS Registry No.: 83471-41-4
Formal Name: N-(2,6-dimethylphenyl)hexahydro-1H-azepine-1-acetamide
Synonym: IQB-M-81
MF: C₁₆H₂₄N₂O
FW: 260.4
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

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

Description

Pincaïnide is a calcium channel inhibitor.^{1,2} It reduces norepinephrine-induced calcium influx and increases calcium efflux in isolated rat aortic strips.¹ Pincaïnide reduces phasic and tonic contractions induced by norepinephrine (IC₅₀s = 880 and 780 μM, respectively) and high potassium (IC₅₀s = 620 and 670 μM, respectively) in isolated rat aortic strips. It also decreases heart rate, contractile force, and the maximum following frequency in isolated rat right atria.² Infiltration of pincaïnide (1% solution) induces local anesthesia in rabbits with a duration of 240 minutes.³

References

1. Barrigón, S., Diez, J.J., Tamargo, J., *et al.* Effects of pincaïnide on contractile responses and ⁴⁵Ca movements in rat isolated vascular smooth muscle. *Eur. J. Pharmacol.* **111(2)**, 227-233 (1985).
2. Tejerina, T., Barrigón, S., and Tamargo, J. Comparison of three β-amino anilides: IQB-M-81, lidocaine and tocainide, on isolated rat atria. *Eur. J. Pharmacol.* **95**, 93-99 (1983).
3. Lebedeva, A.S., Chernyakova, I.V., Likhosherstov, A.M., *et al.* Azacycloalkanes. XXXII. Synthesis of α-hexamethyleneiminocarboxylic acid arylamides and their local anesthetic activity. *Pharm. Chem. J.* **25**, 79-83 (1991).

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, 04/10/2026

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