

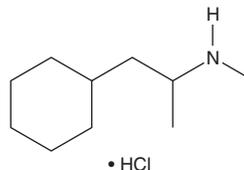
PRODUCT INFORMATION



Propylhexedrine (hydrochloride)

Item No. 12058

CAS Registry No.: 1007-33-6
Formal Name: N, α -dimethyl-cyclohexaneethanamine, monohydrochloride
Synonyms: Benzedrex, Cyclohexylisopropylmethylamine, Eventin, NSC 27110, NSC 170998, Obesin
MF: C₁₀H₂₁N • HCl
FW: 191.7
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Propylhexedrine is an α -adrenergic sympathomimetic amine that acts as a local vasoconstrictor.¹ It is the active ingredient in some over-the-counter nasal decongestants but may be abused, typically by intravenous or oral delivery of the nasal decongestant.^{1,2} This product is intended for research and forensic applications. The analysis of underivatized propylhexedrine in human serum using mass spectrometry has been described.³

References

1. Holler, J.M., Vorce, S.P., McDonough-Bender, P.C., *et al.* A drug toxicity death involving propylhexedrine and mitragynine. *J. Anal. Toxicol.* **35(1)**, 54-59 (2011).
2. Marsden, P. and Sheldon, J. Acute poisoning by propylhexedrine. *Br. Med. J.* **1(5802)**, 730 (1972).
3. Bogusz, M.J., Krüger, K.D., and Maier, R.D. Analysis of underivatized amphetamines and related phenethylamines with high-performance liquid chromatography-atmospheric pressure chemical ionization mass spectrometry. *J. Anal. Toxicol.* **24(2)**, 77-84 (2000).

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, 07/24/2023

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