

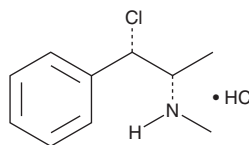
PRODUCT INFORMATION



(+)-Chloropseudoephedrine (hydrochloride)

Item No. 9002347

CAS Registry No.: 94133-42-3
Formal Name: (αS,βS)-β-chloro-N,α-dimethylbenzeneethanamine, monohydrochloride
MF: C₁₀H₁₄ClN • HCl
FW: 220.1
Purity: ≥95%
UV/Vis.: λ_{max}: 217 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.

Description

(+)-Chloropseudoephedrine is a contaminant produced during the illicit manufacture of (+)-methamphetamine (Item No. 13997) when (+)-pseudoephedrine or (-)-ephedrine are used as precursors.^{1,2} Chloroephedrine has been reported to be present in varying levels in clandestinely synthesized methamphetamine and has been observed in some forensic samples.³ This product is intended for forensic and research purposes.

References

1. Drug status report: (-)-Chloropseudoephedrine hydrochloride, (2010).
2. Allen, A.C. and Kiser, W.O. Methamphetamine from ephedrine: I. chloroephedrines and aziridines. *J. Forensic Sci.* **32**, 953-962 (1987).
3. Ko, B.J., Suh, S., Suh, Y.J., et al. (1S,2S)-1-Methylamino-1-phenyl-2-chloropropane: Route specific marker impurity of methamphetamine synthesized from ephedrine via chloroephedrine. *Forensic Sci. Int.* **221**, 92-97 (2012).

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, 07/12/2023

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