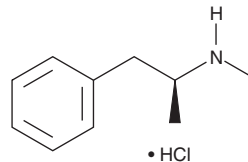


PRODUCT INFORMATION

(+)-Methamphetamine (hydrochloride)

Item No. 13997

CAS Registry No.: 51-57-0
Formal Name: N,αS-dimethyl-benzeneethanamine, monohydrochloride
Synonyms: D-Desoxyephedrine, D-Methamphetamine, (+)-Methylamphetamine, NSC 169505
MF: C₁₀H₁₅N • HCl
FW: 185.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥3 years



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

Description

(+)-Methamphetamine (Item No. 13997) is an analytical reference material categorized as an amphetamine. It is the more physiologically active isomer of (±)-methamphetamine (Item No. ISO60168) and is both neurotoxic and frequently abused.¹⁻⁴ Methamphetamine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Rothman, R.B., Baumann, M.H., Dersch, C.M., *et al.* Amphetamine-type central nervous system stimulants release norepinephrine more potently than they release dopamine and serotonin. *Science* **39(1)**, 32-41 (2001).
2. Rothman, R.B., Vu, N., Partilla, J.S., *et al.* In vitro characterization of ephedrine-related stereoisomers at biogenic amine transporters and the receptorome reveals selective actions as norepinephrine transporter substrates. *J. Pharmacol. Exp. Ther.* **307(1)**, 138-145 (2003).
3. Ichida, J.K., Blanchard, J., Lam, K., *et al.* A small-molecule inhibitor of Tgf-β signaling replaces Sox2 in reprogramming by inducing Nanog. *Cell Stem Cell* **5(5)**, 491-503 (2009).
4. Yu, S., Zhu, L., Shen, Q., *et al.* Recent advances in methamphetamine neurotoxicity mechanisms and its molecular pathophysiology. *Behav. Neurol.* **2015:103969**, (2015).

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, 08/09/2019

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