

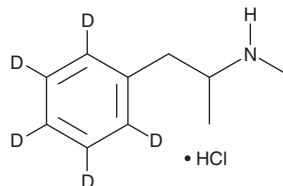
PRODUCT INFORMATION



(±)-Methamphetamine-d₅ (hydrochloride)

Item No. 14021

CAS Registry No.: 2747917-87-7
Formal Name: N-methyl-1-(phenyl-d₅)propan-2-amine, monohydrochloride
Synonym: DL-Methamphetamine-d₅
MF: C₁₀H₁₀D₅N • HCl
FW: 190.7
Purity: ≥98%
Supplied as: A crystalline 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-d₅ (Item No. 14021) is an analytical reference material that is intended for use as an internal standard for the quantification of (±)-methamphetamine (Item No. ISO60168 | 14216 | 36566) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(±)-Methamphetamine is categorized as an amphetamine and consists of dextrorotary and levorotary enantiomers.^{1,2} (+)-Methamphetamine (Item No. 13997) is the more physiologically active isomer compared with (-)-methamphetamine (Item No. 13998). Formulations containing methamphetamine are both neurotoxic and frequently abused.³ (±)-Methamphetamine-d₅ 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. Jirovský, D., Lemr, K., Sevcík, J., *et al.* Methamphetamine - properties and analytical methods of enantiomer determination. *Forensic Sci. Int.* **96(1)**, 61-70 (1998).
2. 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).
3. Yu, S., Zhu, L., Shen, Q., *et al.* Recent advances in methamphetamine neurotoxicity mechanisms and its molecular pathophysiology. *Behav. Neurol.* 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, 04/20/2023

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