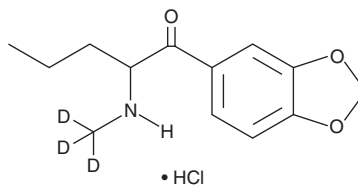


PRODUCT INFORMATION



Pentylone-d₃ (hydrochloride) Item No. 35689

CAS Registry No.: 1781649-12-4
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(methyl-d₃-amino)-1-pentanone, monohydrochloride
MF: C₁₃H₁₄D₃NO₃ • HCl
FW: 274.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥10 years



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

Description

Pentylone-d₃ (hydrochloride) (Item No. 35689) is an analytical reference material intended for use as an internal standard for the quantification of pentylone (Item Nos. 27504 | 9000746) 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).

Pentylone is categorized as a cathinone.¹ It is a metabolite of N,N-dimethylpentylone (Item Nos. 31085 | 9001933).² Pentylone has been detected in products sold as bath salts.¹ Pentylone-d₃ is regulated as a Schedule I 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 for reference materials.

References

1. Marinetti, L.J. and Antonides, H.M. Analysis of synthetic cathinones commonly found in bath salts in human performance and postmortem toxicology: Method development, drug distribution and interpretation of results. *J. Anal. Toxicol.* **37**(3), 135-146 (2013).
2. Tang, Y., Xu, L., Zhao, J., *et al.* Metabolism of dipentylone in zebrafish and human liver microsomes determined by liquid chromatography-high resolution mass spectrometry. *J. Pharm. Biomed. Anal.* **236**, 115710 (2023).

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, 04/10/2024

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