

PRODUCT INFORMATION



N,N-Dimethylpentylone-d₆ (hydrochloride)

Item No. 36899

Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(dimethylamino-d₆)-1-pentanone, monohydrochloride

Synonyms: bk-DMBDP-d₆, Dipentylone-d₆

MF: C₁₄H₁₃D₆NO₃ • HCl

FW: 291.8

Chemical Purity: ≥98% (N,N-Dimethylpentylone)

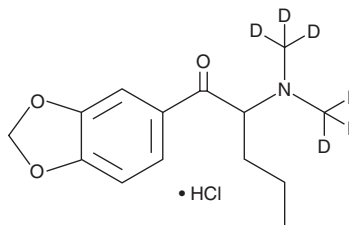
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

N,N-Dimethylpentylone-d₆ (hydrochloride) (Item No. 36899) is intended for use as an internal standard for the quantification of N,N-dimethylpentylone (Item Nos. 31085 | 9001933) 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).

Dimethylpentylone is categorized as a cathinone.¹ It has been found in postmortem human blood samples where the use of N,N-dimethylpentylone was suspected. This product is intended for research and forensic applications.

Reference

1. Fogarty, M.F., Krotulski, A.J., Papsun, D.M., *et al.* N,N-Dimethylpentylone (dipentylone)—a new synthetic cathinone identified in a postmortem forensic toxicology case series. *J. Anal. Toxicol.* **47(8)**, 753-761 (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, 01/09/2024

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