

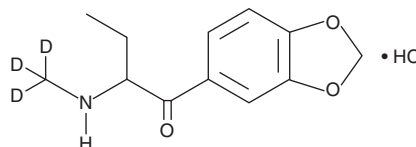
# PRODUCT INFORMATION



## Butylone-d<sub>3</sub> (hydrochloride) (exempt preparation)

Item No. 22880

**CAS Registry No.:** 1231710-63-6  
**Formal Name:** 1-(1,3-benzodioxol-5-yl)-2-(methyl-d<sub>3</sub>-amino)-1-butanone, monohydrochloride  
**Synonyms:** bk-MBDB-d<sub>3</sub>, β-keto MBDB-d<sub>3</sub>  
**MF:** C<sub>12</sub>H<sub>12</sub>D<sub>3</sub>NO<sub>3</sub> • HCl  
**FW:** 260.7  
**Chemical Purity:** ≥98% (Butylone)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 234, 280, 318 nm  
**Supplied as:** A 1 mg/ml solution in methanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

Butylone-d<sub>3</sub> (hydrochloride) (exempt preparation) (Item No. 22880) is an analytical reference standard intended for use as an internal standard for the quantification of butylone (Item Nos. 10393 | 16059) 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).

Butylone is categorized as a cathinone.<sup>1</sup> It is a metabolite of bk-DMBDB and has been detected in products sold as bath salts.<sup>2,3</sup> Butylone-d<sub>3</sub> is regulated as a Schedule I compound in the United States. Butylone-d<sub>3</sub> (hydrochloride) (exempt preparation) (Item No. 22880) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

### References

1. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8(1)**, 33-42 (2012).
2. Krotulski, A.J., Mohr, A.L.A., Papsun, D.M., *et al.* Dibutylone (bk-DMBDB): Intoxications, quantitative confirmations and metabolism in authentic biological specimens. *J. Anal. Toxicol.* **42(7)**, 437-445 (2018).
3. Rosenbaum, C.D., Carreiro, S.P., and Babu, K.M. Here today, gone tomorrow...and back again? A review of herbal marijuana alternatives (K2, Spice), synthetic cathinones (bath salts), kratom, *Salvia divinorum*, methoxetamine, and piperazines. *J. Med. Toxicol.* **8(1)**, 15-32 (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, 03/10/2022

#### 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