

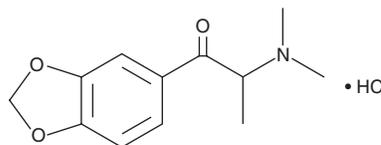
PRODUCT INFORMATION



bk-MDDMA (hydrochloride)

Item No. 9001124

CAS Registry No.: 109367-07-9
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(dimethylamino)-1-propanone, monohydrochloride
Synonyms: Dimethylone, β -keto Methylenedioxy Dimethylamphetamine
MF: $C_{12}H_{15}NO_3 \cdot HCl$
FW: 257.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 235, 282, 320 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

bk-MDDMA (hydrochloride) is a potential psychotropic designer drug of the phenethylamine, amphetamine, and cathinone chemical classes. It is structurally related to methylone, a designer drug that has been detected in products marketed as bath salts, plant food, and tablets.^{1,2} This product is intended for forensic purposes.

References

1. Kikura-Hanajiri, R., Uchiyami, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).
2. Bossong, M.G., Van Dijk, J.P., and Niesink, R.J. Methylone and mCPP, two new drugs of abuse. *Addict. Biol.* **10(4)**, 321-323 (2005).

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, 06/27/2023

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