

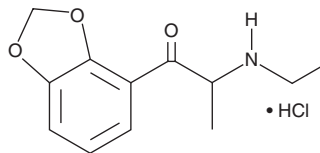
PRODUCT INFORMATION



2,3-Methylenedioxy- α -ethylaminopropiophenone (hydrochloride)

Item No. 9002142

CAS Registry No.: 2749504-06-9
Formal Name: 1-(1,3-benzodioxol-4-yl)-2-(ethylamino)-1-propanone, monohydrochloride
Synonyms: 2,3-bk-MDEA isomer, 2,3-Ethylone isomer
MF: C₁₂H₁₅NO₃ • HCl
FW: 257.7
Purity: ≥98%
UV/Vis.: λ_{\max} : 230, 259, 345 nm
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

bk-MDEA (Item No. 9001123) is a psychotropic designer drug of the phenethylamine, amphetamine, and cathinone chemical classes. It has been detected in products marketed as bath salts, plant food, and tablets.^{1,2} 2,3-Methylenedioxy- α -ethylaminopropiophenone (hydrochloride) is a positional isomer of bk-MDEA, having the methylenedioxy group attached to the 2 and 3, rather than the 3 and 4, positions. The physiological properties of this compound have not been studied. This product is intended for forensic and research purposes.

References

1. Uchiyama, N., Kikura-Hanajiri, R., Kawahara, N., *et al.* Analysis of designer drugs detected in the products purchased in fiscal year 2006. *Yakugaku Zasshi* **128(10)**, 1499-505 (2008).
2. Zaitso, K., Katagi, M., Kamata, H.T., *et al.* Determination of the metabolites of the new designer drugs bk-MBDB and bk-MDEA in human urine. *Forensic Sci. Int.* **188(1-3)**, 131-139 (2009).

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, 07/25/2023

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