

# PRODUCT INFORMATION

## 3,4-Methylenedioxy-5-methylethcathinone (hydrochloride)

Item No. 14078

**CAS Registry No.:** 2748289-02-1  
**Formal Name:** 2-(ethylamino)-1-(7-methyl-1,3-benzodioxol-5-yl)-1-propanone, monohydrochloride

**Synonyms:** 3,4-Methylenedioxy-5-methyl-N-ethyl Cathinone, R-MMC

**MF:** C<sub>13</sub>H<sub>17</sub>NO<sub>3</sub> • HCl

**FW:** 271.7

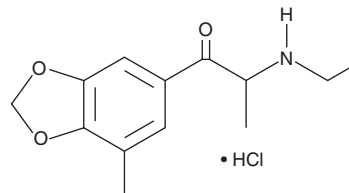
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 207, 237, 290, 318 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 (hydrochloride) (Item No. 9001123, ethylone) is a psychotropic cathinone that has been detected in products marketed as bath salts, plant food, and tablets.<sup>2</sup> 3,4-Methylenedioxy-5-methylethcathinone is an analog of bk-MDEA, differing by the addition of a methyl group at the two position of the benzene ring. The physiological and toxicological properties of this compound have not been elucidated. This product is intended for forensic and research applications.

### References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in *Japan. Leg. Med. (Tokyo)* **13(3)**, 109-15 (2011).
2. 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).

#### 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