

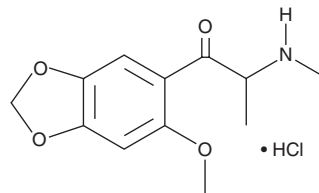
PRODUCT INFORMATION



6-methoxy Methylone (hydrochloride)

Item No. 18029

CAS Registry No.: 2749504-36-5
Formal Name: 1-(6-methoxy-1,3-benzodioxol-5-yl)-2-(methylamino)-1-propanone, monohydrochloride
Synonym: 6-methoxy bk-MDMA
MF: $C_{12}H_{15}NO_4 \cdot HCl$
FW: 273.7
Purity: $\geq 98\%$
Supplied as: A neat solid
Storage: $-20^{\circ}C$
Stability: ≥ 3 years



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

Description

Methylone (Item No. 10986) is a β -ketone analog of 3,4-methylenedioxymethamphetamine (3,4-MDMA; Item No. 13971) that has been detected in products marketed as bath salts, plant food, and tablets.^{1,2} 6-methoxy Methylone is an analog of methylone that bears a methoxy group at the 6 position of the phenyl ring. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

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. 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, 01/24/2025

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