

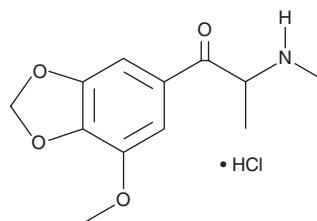
# PRODUCT INFORMATION



## 5-methoxy Methylone (hydrochloride)

Item No. 9002195

**CAS Registry No.:** 2702151-27-5  
**Formal Name:** 1-(7-methoxy-1,3-benzodioxol-5-yl)-2-(methylamino)-1-propanone, monohydrochloride  
**MF:** C<sub>12</sub>H<sub>15</sub>NO<sub>4</sub> • HCl  
**FW:** 273.7  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

Methylone (Item No. 10986) is a  $\beta$ -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.<sup>1,2</sup> Similar to 3,4-MDMA, it acts on components of the monoaminergic systems.<sup>3-5</sup> 5-methoxy Methylone is a cathinone derivative that bears a methoxy group similar to various tryptamine analogs including 5-methoxy MiPT (Item No. 11482), which are known to potently inhibit monoamine transport function.<sup>5</sup> The pharmacological and toxicological properties of this compound are not known. This product is intended for forensic and research purposes.

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)**, 109-115 (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).
3. Cozzi, N.V., Sievert, M.K., Shulgin, A.T., *et al.* Inhibition of plasma membrane monoamine transporters by b-ketoamphetamines. *Eur. J. Pharmacol.* **381(1)**, 63-69 (1999).
4. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
5. Sogawa, C., Sogawa, N., Ohyama, K., *et al.* Methylone and monoamine transporters: Correlation with toxicity. *Curr. Neuropharmacol.* **9(1)**, 58-62 (2011).

#### 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, 12/06/2021

#### 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