

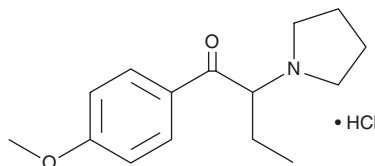
# PRODUCT INFORMATION



## 4-methoxy- $\alpha$ -Pyrrolidinobutiophenone (hydrochloride)

Item No. 15354

**CAS Registry No.:** 2749897-35-4  
**Formal Name:** 1-(4-methoxyphenyl)-2-(1-pyrrolidinyl)-1-butanone, monohydrochloride  
**Synonym:** 4-MeOPBP  
**MF:** C<sub>15</sub>H<sub>21</sub>NO<sub>2</sub> • HCl  
**FW:** 283.8  
**Purity:** ≥98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 224, 290 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

Pyrrolidinophenones and butyrophenones are chemical compounds that feature ketone functional groups affixed to the narcotic cathinone. They bear a strong structural relationship with the stimulant  $\alpha$ -pyrrolidinopentiophenone ( $\alpha$ -PVP; Item No. 9001083) and many have been identified as components of designer drugs.<sup>1-3</sup> 4-methoxy- $\alpha$ -Pyrrolidinobutiophenone(4-MeOPBP) (hydrochloride) is a pyrovalerone analog that shares structural features with 4-methoxy PVP (Item No. 14097) and 4-methyl- $\alpha$ -pyrrolidinobutiophenone (Item No. 9001190). The biological and toxicological properties of this compound have not been evaluated. This product is intended only 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. Brandt, S.D., Freeman, S., Sumnall, H.R., *et al.* Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. *Drug Test. Anal.* **3(9)**, 569-575 (2011).
3. Westphal, F., Junge, T., Rösner, P., *et al.* Mass spectral and NMR spectral data of two new designer drugs with an  $\alpha$ -aminophenone structure: 4'-Methyl- $\alpha$ -pyrrolidinohexanophenone and 4'-methyl- $\alpha$ -pyrrolidinobutyrophenone. *Forensic Sci. Int.* **169(1)**, 32-42 (2007).

#### 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, 06/20/2023

#### 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