

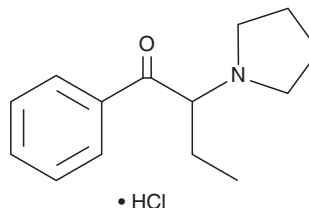
# PRODUCT INFORMATION



## $\alpha$ -Pyrrolidinobutiophenone (hydrochloride)

Item No. 9001195

**CAS Registry No.:** 13415-54-8  
**Formal Name:** 1-phenyl-2-(1-pyrrolidinyl)-1-butanone, monohydrochloride  
**Synonym:**  $\alpha$ -PBP  
**MF:** C<sub>14</sub>H<sub>19</sub>NO • HCl  
**FW:** 253.8  
**Purity:**  $\geq$ 95%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 250 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 5 years



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

### Description

Butyrophenones are chemical compounds with a ketone functional group that have anti-psychotic and anti-emetic properties. 4-Methyl- $\alpha$ -pyrrolidinobutiophenone is a butyrophenone as well as a synthetic cathinone that has been identified in party pills.<sup>1</sup>  $\alpha$ -PBP (hydrochloride) is a synthetic cathinone/butyrophenone that has potential for abuse. Its physiological and toxicological properties are poorly understood. This product is intended for forensic and research applications.

### Reference

1. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. Brandt, S.D., *et al.* Retrieved August 18, 2010, from [www.drugtestinganalysis.com](http://www.drugtestinganalysis.com).

**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, 07/06/2023

### 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](http://WWW.CAYMANCHEM.COM)