

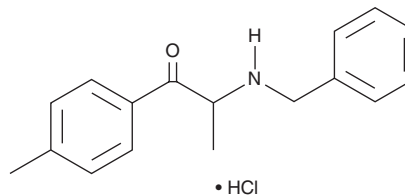
PRODUCT INFORMATION



Benzedrone (hydrochloride)

Item No. 11915

CAS Registry No.: 1797979-43-1
Formal Name: 1-(4-methylphenyl)-2-[(phenylmethyl)amino]-1-propanone, monohydrochloride
Synonym: 4-MBC
MF: C₁₇H₁₉NO • HCl
FW: 289.8
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 260 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

4-Methylmethcathinone (also known as Mephedrone) is a stimulatory designer drug found in bath salts and similar products.¹⁻³ 4-MBC is an analog of mephedrone characterized by the replacement of the amino methyl group with a benzyl moiety. The physiological and toxicological properties of this compound are unknown. This product is intended for forensic and research applications.

References

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.
2. 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).
3. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8(1)**, 33-42 (2012).

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, 07/13/2023

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