

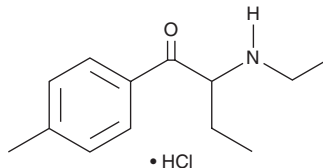
PRODUCT INFORMATION



4-methyl- α -Ethylaminobutiophenone (hydrochloride)

Item No. 11489

CAS Registry No.: 18268-19-4
Formal Name: 2-(ethylamino)-1-(4-methylphenyl)-1-butanone, monohydrochloride
Synonym: 4-methyl NEB
MF: $C_{13}H_{19}NO \cdot HCl$
FW: 241.8
Purity: $\geq 98\%$
Supplied as: A neat solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Description

Buphedrone (α -methylamino-butyrophenone) is a substituted cathinone which has been identified in products that are sold as stimulants.^{1,2} 4-methyl- α -Ethylaminobutyophenone is a *para*-methyl analog of buphedrone with an ethyl group replacing methyl at the alpha position. Its biological and toxicological properties have not been evaluated. 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 for reference materials.

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. Maheux, C.R. and Copeland, C.R. Chemical analysis of two new designer drugs: Buphedrone and pentedrone. *Drug Test. Anal.* 1-8 (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, 08/18/2023

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