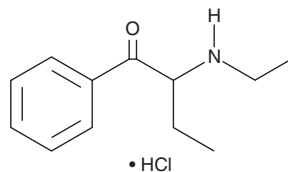


PRODUCT INFORMATION

N-Ethylbuphedrone (hydrochloride)

Item No. 11665

CAS Registry No.: 2731709-73-0
Formal Name: 2-(ethylamino)-1-phenylbutan-1-one, monohydrochloride
Synonym: NEB
MF: $C_{12}H_{17}NO \cdot HCl$
FW: 227.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 249 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

Substituted cathinones, like buphedrone, are psychoactive compounds commonly used as recreational drugs.^{1,2} Some have amphetamine-like effects.³ N-Ethylbuphedrone is a substituted cathinone characterized by an ethyl group at the alpha position as well as an N-terminal ethyl group. Its biological and toxicological properties have not been evaluated. N-Ethylbuphedrone is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Camilleri, A., Johnston, M.R., Brennan, M., *et al.* Chemical analysis of four capsules containing the controlled substance analogues 4-methylmethcathinone, 2-fluoromethamphetamine, α -phthalimidopropiophenone and N-ethylcathinone. *Forensic Sci. Int.* **197**, 59-66 (2010).
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. Dal Cason, T.A., Young, R., and Glennon, R.A. Cathinone: An investigation of several N-alkyl and methylenedioxy-substituted analogs. *Pharmacol. Biochem. Behav.* **58(4)**, 1109-1116 (1997).

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, 07/18/2023

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