

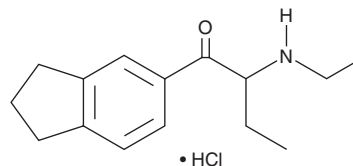
PRODUCT INFORMATION



bk-EABDI (hydrochloride)

Item No. 17990

CAS Registry No.: 2726084-56-4
Formal Name: 1-(2,3-dihydro-1H-inden-5-yl)-2-(ethylamino)-1-butanone, monohydrochloride
Synonym: bk-IBP, Indanyl-NEB, 3,4-Trimethylene- α -ethylaminobutiophenone
MF: C₁₅H₂₁NO • HCl
FW: 267.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

N-Ethylbuphedrone (NEB, Item No. 11665) is a substituted cathinone characterized by an ethyl group at the alpha position as well as an N-terminal ethyl group. bk-EABDI is a 3,4-trimethylene derivative of NEB. The physiological and toxicological properties of this compound are not known. 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.

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, 08/04/2022

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