

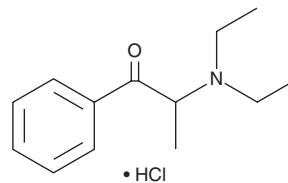
# PRODUCT INFORMATION



## Diethylcathinone (hydrochloride)

Item No. 11333

**CAS Registry No.:** 134-80-5  
**Formal Name:** 2-(diethylamino)-1-phenyl-1-propanone, monohydrochloride  
**Synonyms:** Diethylpropion, Menutil, Moderatan, Tenuate  
**MF:** C<sub>13</sub>H<sub>19</sub>NO • HCl  
**FW:** 241.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 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

Diethylcathinone is a stimulant drug of the phenethylamine, amphetamine, and cathinone chemical classes that is used as an appetite suppressant.<sup>1</sup> It functions as an inactive prodrug that only becomes active when further metabolized to ethcathinone, a selective norepinephrine releasing agent.<sup>2</sup> In the US, diethylcathinone is classified as a Schedule IV controlled substance, with low potential for abuse.<sup>3</sup> This product is intended for forensic and research applications.

### References

1. Li, Z., Maglione, M., Tu, W., *et al.* Meta-analysis: Pharmacologic treatment of obesity. *Ann. Intern. Med.* **142**(7), 532-546 (2005).
2. Rothman, R.B. and Baumann, M.H. Therapeutic potential of monoamine transporter substrates. *Curr. Top. Med. Chem.* **6**(17), 1845-1859 (2006).
3. Controlled substances - by CSA schedule. In: U.S. Department of Justice - Drug Enforcement Administration, Diversion Control Division. Available from: <http://www.deadiversion.usdoj.gov/schedules/index.html>

#### 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, 07/14/2023

#### 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)