

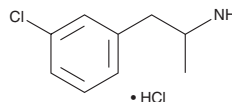
# PRODUCT INFORMATION



## 3-Chloroamphetamine (hydrochloride)

Item No. 9001855

**CAS Registry No.:** 35378-15-5  
**Formal Name:** 3-chloro- $\alpha$ -methyl-benzeneethanamine, monohydrochloride  
**Synonyms:** 3-CA, *m*-CA  
**MF:** C<sub>9</sub>H<sub>12</sub>ClN • HCl  
**FW:** 206.1  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\max}$ : 213, 266 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 5 years



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

### Description

3-Chloroamphetamine (3-CA) is a derivative of amphetamine that acts as a central nervous system stimulant by releasing serotonin and dopamine.<sup>1</sup> It has been detected in urine samples, along with the related compound 4-CA (Item No. 9001856), as an apparent designer drug.<sup>2</sup> This product is intended only for forensic and research purposes.

### References

- Ogren, S.O. and Ross, S.B. Substituted amphetamine derivatives. II. Behavioral effects in mice related to monoaminergic neurones. *Acta Pharmacol. Toxicol. (Copenh)* **41(4)**, 353-368 (1977).
- Lin, T.C., Lin, D.L., and Lua, A.C. Detection of *p*-chloroamphetamine in urine samples with mass spectrometry. *J. Anal. Toxicol.* **35(4)**, 205-210 (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, 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