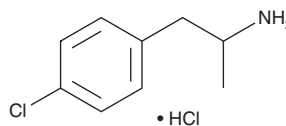


# PRODUCT INFORMATION

## 4-Chloroamphetamine (hydrochloride)

Item No. 9001856

**CAS Registry No.:** 3706-38-5  
**Formal Name:** 4-chloro- $\alpha$ -methyl-benzeneethanamine, monohydrochloride  
**Synonyms:** 4-CA, *p*-CA  
**MF:** C<sub>9</sub>H<sub>12</sub>ClN • HCl  
**FW:** 206.1  
**Purity:** ≥98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 220 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

Halogenated amphetamines, including CAs, are psychostimulatory compounds that have been abused recreationally.<sup>1</sup> 4-CA is a halogenated amphetamine that causes rapid depletion of serotonin and its major metabolite, 5-hydroxyindoleacetic acid, in animal brains within hours after injection.<sup>2-3</sup> In addition, it is highly and selectively neurotoxic, targeting serotonergic neurons.<sup>2-3</sup> This product is intended for forensic and research applications.

### References

1. 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).
2. Fuller, R.W., Baker, J.C., Perry, K.W., et al. Comparison of 4-chloro-, 4-bromo- and 4-fluoroamphetamine in rats: Drug levels in brain and effects on brain serotonin metabolism. *Neuropharmacology* **14**(10), 739-746 (1975).
3. Harvey, J.A., McMaster, S.E., and Fuller, R.W. Comparison between the neurotoxic and serotonin-depleting effects of various halogenated derivatives of amphetamine in the rat. *J. Pharmacol. Exp. Ther.* **202**(3), 581-589 (1977).

#### 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, 07/27/2023

#### 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