

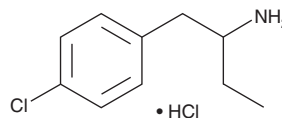
PRODUCT INFORMATION



4-CAB (hydrochloride)

Item No. 13893

CAS Registry No.: 23292-08-2
Formal Name: 4-chloro- α -ethyl-benzeneethanamine, monohydrochloride
Synonyms: AEPCA, 4-Chlorophenylisobutylamine, 4-chloro- α -Ethylphenethylamine
MF: C₁₀H₁₄ClN • HCl
FW: 220.1
Purity: \geq 95%
UV/Vis.: λ_{max} : 220, 268 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

4-CAB is an analog of *p*-chloroamphetamine, a psychoactive compound which shows pronounced toxicity for serotonergic neurons.^{1,2} 4-CAB inhibits the reuptake of serotonin (IC₅₀ = 330 nM) and dopamine (IC₅₀ = 2.3 μ M).¹ This product is intended for forensic and research applications.

References

1. Johnson, M.P., Huang, X.M., Oberlender, R., *et al.* Behavioral, biochemical and neurotoxicological actions of the alpha-ethyl homologue of *p*-chloroamphetamine. *Eur. J. Pharmacol.* **191(1)**, 1-10 (1990).
2. Gobbi, M., Moia, M., Pirona, L., *et al.* *p*-Methylthioamphetamine and 1-(*m*-chlorophenyl)piperazine, two non-neurotoxic 5-HT releasers *in vivo*, differ from neurotoxic amphetamine derivatives in their mode of action at 5-HT nerve endings *in vitro*. *J. Neurochem.* **82(6)**, 1435-1443 (2002).

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, 07/25/2023

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