

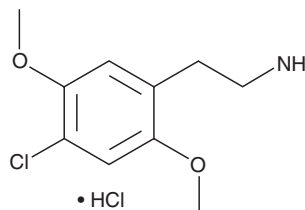
PRODUCT INFORMATION



2C-C (hydrochloride) (exempt preparation)

Item No. 15716

CAS Registry No.: 88441-15-0
Formal Name: 4-chloro-2,5-dimethoxybenzeneethanamine, monohydrochloride
Synonym: 2,5-Dimethoxy-4-chlorophenethylamine
MF: C₁₀H₁₄ClNO₂ • HCl
FW: 252.1
Purity: ≥98%
Supplied as: A 1 mg/ml solution in methanol
UV/Vis.: λ_{max}: 204, 225, 295 nm
Storage: -20°C
Stability: ≥3 years



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

Description

A series of 2,5-dimethoxyphenethylamines, collectively referred to as 2Cs, have psychoactive effects.^{1,2} The most effective 2C compounds are substituted at the four position of the aromatic ring. Many are scheduled as illegal substances.^{3,4} 2C-C is described formally as 2,5-dimethoxy-4-chlorophenethylamine. A known hallucinogen, this compound stimulates monoamine receptor activity and inhibits the re-uptake of serotonin and norepinephrine in rat brain synaptosomes (IC₅₀s = 31 and 63 μM, respectively).^{4,5} LC-MS/MS screening methods for this designer drug have been developed.⁶ This product is intended for forensic and research purposes.

Reference

1. Bruno, R., Matthews, A.J., Dunn, M., *et al.* *Drug Alcohol Depend.* **124**(1-2), 19-25 (2012).
2. Moya, P.R., Berg, K.A., Gutiérrez-Hernandez, M.A., *et al.* *J. Pharmacol. Exp. Ther.* **321**, 1054-1061 (2007).
3. Meyer, M.R. and Maurer, H.H. *Curr. Drug Metab.* **11**, 468-482 (2010).
4. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. *Eur. J. Pharmacol.* **559**(2-3), 132-137 (2007).
5. Nonaka, R., Nagai, F., Ogata, A., *et al.* *Biol. Pharm. Bull.* **30**(12), 2328-33 (2007).
6. Wohlfarth, A., Weinmann, W., and Dresen, S. *Anal. Bioanal. Chem.* **396**, 2403-2414 (2010).

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, 06/25/2025

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