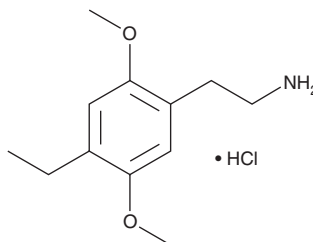


PRODUCT INFORMATION

2C-E (hydrochloride) (exempt preparation)

Item No. 15699

CAS Registry No.: 923013-67-6
Formal Name: 2,5-dimethoxy-4-ethylphenethylamine, monohydrochloride
Synonym: 2,5-Dimethoxy-4-ethylphenethylamine
MF: $C_{12}H_{19}NO_2 \cdot HCl$
FW: 245.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 203, 227, 291 nm
Supplied as: A solution in methanol
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Description

2C-E (hydrochloride) (exempt preparation) (Item No. 15699) is an analytical reference standard categorized as a phenethylamine.¹ 2C-E has been detected in capsules and powder seized by law enforcement.²⁻⁴ 2C-E is regulated as a Schedule I compound in the United States. 2C-E (hydrochloride) (exempt preparation) (Item No. 15699) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
2. Drug Enforcement Administration. Microgram Bulletin. **37(2)**, 193-204 (2004).
3. Drug Enforcement Administration. Microgram Bulletin. **38(4)**, 55-69 (2005).
4. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 468-482 (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, 03/04/2019

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