

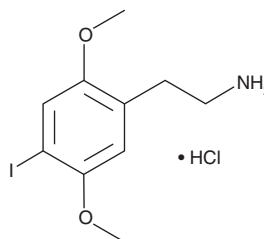
# PRODUCT INFORMATION



## 2C-I (hydrochloride)

Item No. 11113

**CAS Registry No.:** 64584-32-3  
**Formal Name:** 4-iodo-2,5-dimethoxy-benzeneethanamine, monohydrochloride  
**Synonym:** 2,5-Dimethoxy-4-iodophenethylamine  
**MF:**  $C_{10}H_{14}INO_2 \cdot HCl$   
**FW:** 343.6  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 204, 234, 298 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 5$  years



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

### Description

A series of 2,5-phenethylamines, collectively referred to as 2Cs, have psychoactive effects; many are scheduled as illegal substances. 2C-I (hydrochloride) is described formally as 2,5-dimethoxy-4-iodophenethylamine. Like other 2Cs, 2C-I (hydrochloride) is a monoamine oxidase inhibitor, blocking the breakdown of serotonin and norepinephrine ( $IC_{50}$  values of 79 and 37  $\mu M$ , respectively), but not dopamine.<sup>1</sup> Its metabolism has been described.<sup>2</sup> 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. 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, 07/07/2023

#### 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](http://WWW.CAYMANCHEM.COM)