

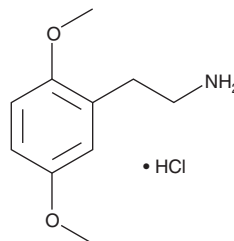
# PRODUCT INFORMATION



## 2C-H (hydrochloride)

Item No. 11889

**CAS Registry No.:** 3166-74-3  
**Formal Name:** 2,5-dimethoxy-benzeneethanamine, monohydrochloride  
**Synonym:** 2,5-Dimethoxyphenethylamine  
**MF:** C<sub>10</sub>H<sub>15</sub>NO<sub>2</sub> • HCl  
**FW:** 217.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 291 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

A series of 2,5-phenethylamines, collectively referred to as 2Cs, have psychoactive effects.<sup>1,2</sup> The most effective 2C compounds are 2,5-dimethoxy-4-substituted phenethylamines; many are scheduled as illegal substances.<sup>3,4</sup> 2C-H is described formally as 2,5-dimethoxyphenethylamine. Unlike the 4-substituted 2-C forms, it has little effect on serotonin receptors, activating the 5-HT<sub>2C</sub> receptor to 20% that induced by serotonin (pEC<sub>50</sub> = 5.93).<sup>2</sup> This product is intended for research and forensic applications.

### References

1. Bruno, R., Matthews, A.J., Dunn, M., *et al.* Emerging psychoactive substance use among regular ecstasy users in Australia. *Drug Alcohol Depend.* **124**(1-2), 19-25 (2012).
2. Moya, P.R., Berg, K.A., Gutiérrez-Hernandez, M.A., *et al.* Functional selectivity of hallucinogenic phenethylamine and phenylisopropylamine derivatives at human 5-hydroxytryptamine (5-HT)<sub>2A</sub> and 5-HT<sub>2C</sub> receptors. *J. Pharmacol. Exp. Ther.* **321**(3), 1054-1061 (2007).
3. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11**(5), 468-482 (2010).
4. 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).

#### 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, 07/24/2023

#### 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)