

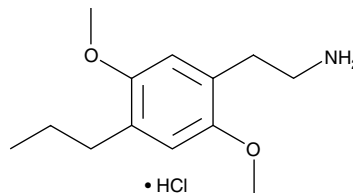
# Product Information



## 2C-P (hydrochloride)

Item No. 11696

**CAS Registry No.:** 1359704-27-0  
**Formal Name:** 2,5-dimethoxy-4-propylbenzeneethanamine, monohydrochloride  
**Synonym:** 2,5-Dimethoxy-4-propylphenethylamine  
**MF:** C<sub>13</sub>H<sub>21</sub>NO<sub>2</sub> • HCl  
**FW:** 259.8  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 204, 227, 292 nm



### Laboratory Procedures

For long term storage, we suggest that 2C-P (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

2C-P (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 2C-P (hydrochloride) in the solvent of choice. 2C-P (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 2C-P (hydrochloride) in ethanol and DMF is approximately 2 mg/ml and approximately 13 mg/ml in DMSO.

2C-P (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2C-P (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2C-P (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

A series of 2,5-dimethoxy phenethylamines, collectively referred to as 2Cs, have psychoactive effects.<sup>1,2</sup> The most effective 2C compounds are substituted at the four position of the aromatic ring; many are scheduled as illegal substances.<sup>3,4</sup> 2C-P is described formally as 2,5-dimethoxy-4-propylphenethylamine and is structurally analogous to the psychedelic drug 2C-B. The physiological and toxicological properties of this compound are not known. LC-MS/MS screening methods for this designer drug have been developed.<sup>5</sup> This product is intended for forensic and research purposes.

### 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**, 1054-1061 (2007).
3. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11**, 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).
5. Wohlfarth, A., Weinmann, W., and Dresen, S. LC-MS/MS screening method for designer amphetamines, tryptamines, and piperazines in serum. *Anal. Bioanal. Chem.* **396**, 2403-2414 (2010).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/11696](http://www.caymanchem.com/catalog/11696)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 04/07/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)