

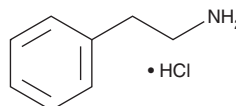
# PRODUCT INFORMATION



## 2-Phenethylamine (hydrochloride)

Item No. 14206

**CAS Registry No.:** 156-28-5  
**Formal Name:** benzeneethanamine, monohydrochloride  
**Synonyms:** 2-PEA,  $\beta$ -Phenylethylamine, 2-Phenylethylammonium  
**MF:**  $C_8H_{11}N \cdot HCl$   
**FW:** 157.6  
**Purity:**  $\geq 98\%$   
**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

2-Phenethylamine is a natural monoamine alkaloid that acts as neuromodulator in the central nervous system. It stimulates the release of norepinephrine and dopamine and is metabolized by monoamine oxidases as well as other enzymes.<sup>1-3</sup> This product is intended for forensic and research applications.

### References

1. Nakamura, M., Ishii, A., and Nakahara, D. Characterization of  $\beta$ -phenylethylamine-induced monoamine release in rat nucleus accumbens: A microdialysis study. *Eur. J. Pharmacol.* **349(2-3)**, 163-169 (1998).
2. Yang, H.Y., and Neff, N.H.  $\beta$ -Phenylethylamine: A specific substrate for type B monoamine oxidase of brain. *J. Pharmacol. Exp. Ther.* **187(2)**, 365-371 (1973).
3. Suzuki, O., Katsumata, Y., and Oya, M. Oxidation of  $\beta$ -phenylethylamine by both types of monoamine oxidase: Examination of enzymes in brain and liver mitochondria of eight species. *J. Neurochem.* **36(3)**, 1298-1301 (1981).

#### 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, 06/16/2023

#### 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)