

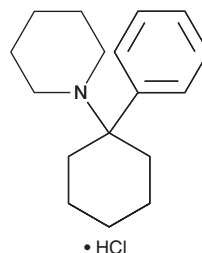
# PRODUCT INFORMATION



## PCP (hydrochloride)

Item No. 39258

**CAS Registry No.:** 956-90-1  
**Formal Name:** 1-(1-phenylcyclohexyl)-piperidine, monohydrochloride  
**Synonyms:** CI-395, MCV 4158, NIH 9580, NSC 40902, Phencyclidine  
**MF:**  $C_{17}H_{25}N \bullet HCl$   
**FW:** 279.9  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Description

PCP (hydrochloride) (Item No. 39258) is an analytical reference standard categorized as an arylcyclohexylamine.<sup>1</sup> PCP is a dissociative anesthetic.<sup>1,2</sup> PCP is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

### References

1. Gibbons, S. 'Legal highs' - novel and emerging psychoactive drugs: A chemical overview for the toxicologist. *Clin. Toxicol.* **50(1)**, 15-24 (2012).
2. Bresink, I., Danysz, W., Parsons, C.G., *et al.* Different binding affinities of NMDA receptor channel blockers in various brain regions-indication of NMDA receptor heterogeneity. *Neuropharmacology* **34(5)**, 533-540 (1995).

#### 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, 03/10/2025

#### 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)