

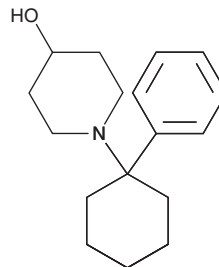
# PRODUCT INFORMATION



## 1-(1-Phenylcyclohexyl)-4-hydroxypiperidine

Item No. 20563

**CAS Registry No.:** 60232-85-1  
**Formal Name:** 1-(1-phenylcyclohexyl)-4-piperidinol  
**Synonyms:** NSC 170516, PCHP, PHP  
**MF:** C<sub>17</sub>H<sub>25</sub>NO  
**FW:** 259.4  
**Purity:** ≥98%  
**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

1-(1-Phenylcyclohexyl)-4-hydroxypiperidine (Item No. 20563) is an analytical reference standard that is structurally categorized as an arylcyclohexylamine. It is a primary, monohydroxylated metabolite of the NMDA receptor channel blocker, PCP (Item No. ISO60194) that can be detected in hair samples.<sup>1,2</sup> 1-(1-Phenylcyclohexyl)-4-hydroxypiperidine is reported to produce one-third the potency of PCP in eliciting PCP-like discriminative stimuli in rat.<sup>3</sup> This product is intended for research and forensic applications.

### References

1. 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).
2. Sakamoto, T., Tanaka, A., and Nakahara, Y. Hair analysis for drugs of abuse XII. Determination of PCP and its major metabolites, PCHP and PPC, in rat hair after administration of PCP. *J. Anal. Toxicol.* **20(2)**, 124-130 (1996).
3. Shannon, H.E. Evaluation of phencyclidine analogs on the basis of their discriminative stimulus properties in the rat. *J. Pharmacol. Exp. Ther.* **216(3)**, 543-551 (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, 06/21/2023

#### 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)