

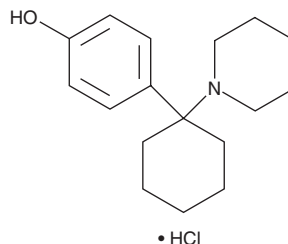
PRODUCT INFORMATION



4-hydroxy PCP (hydrochloride)

Item No. 14274

Formal Name: 4-(1-(piperidin-1-yl)cyclohexyl)phenol, monohydrochloride
Synonyms: HPCP, *p*-hydroxy PCP, *para*-hydroxy PCP
MF: C₁₇H₂₅NO • HCl
FW: 295.9
Purity: ≥98%
UV/Vis.: λ_{max}: 232 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥5 years



Special Conditions: This compound is only stable as a neat solid or in acidic methanol solution. We have found the compound to be stable in 0.005-0.10 M HCl in Methanol.

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

Description

4-hydroxy PCP (hydrochloride) (Item No. 14274) is an analytical reference standard categorized as an arylcyclohexylamine. 4-hydroxy PCP is a transient metabolite of PCP (Item Nos. ISO60194 | 14276) that degrades under neutral and basic conditions.¹ It spontaneously forms a quinone methide intermediate under physiological pH, which can form adducts or react with water to form *para*-hydroxyphenylcyclohexanol and, subsequently, *para*-hydroxyphenylcyclohexene. This product is intended for research and forensic applications.

Reference

1. Driscoll, J.P., Kornecki, K., Wolkowski, J.P., *et al.* Bioactivation of phencyclidine in rat and human liver microsomes and recombinant P450 2B enzymes: Evidence for the formation of a novel quinone methide intermediate. *Chem. Res. Toxicol.* **20**(10), 1488-1497 (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, 08/29/2024

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