

PRODUCT INFORMATION



4-methoxy PCP (hydrochloride)

Item No. 11136

CAS Registry No.: 2185-93-5

Formal Name: 1-[1-(4-methoxyphenyl)cyclohexyl]-piperidine, monohydrochloride

Synonyms: 4-MeO PCP, 4-MeO Phencyclidine, Methoxydine

MF: $C_{18}H_{27}NO \cdot HCl$

FW: 309.9

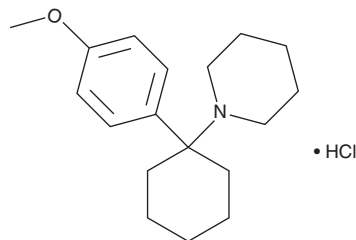
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 231, 272 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years



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

Description

PCP, known more commonly as angel dust, is a psychoactive drug that is regulated as a schedule I compound in the United States.¹ 4-MeO PCP is an analog of PCP which has recently been commercially produced as a designer drug.² It preferentially binds the σ -2 receptor ($K_i = 143$ nM) and less weakly binds the NMDA receptor and serotonin transporter ($K_s = 404$ and 844 nM, respectively).³ The toxicological properties of this compound have not been reported. This product is intended for forensic and research purposes.

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. Misselbrook, G.P. and Hamilton, E.J. Out with the old, in with the new? Case reports of the clinical features and acute management of two novel designer drugs. *Acute Med.* **11**, 157-160 (2012).
3. Advisory Council on the Misuse of Drugs Advisory Council on the Misuse of Drugs (ACMD) methoxetamine report (2012), UK Home Office, 1 (2012).

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, 07/24/2020

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