

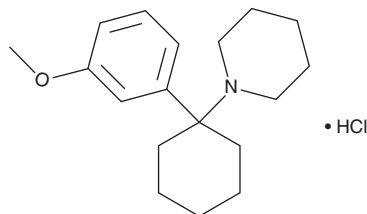
PRODUCT INFORMATION



3-methoxy PCP (hydrochloride)

Item No. 11137

CAS Registry No.: 91164-58-8
Formal Name: 1-[1-(3-methoxyphenyl)cyclohexyl]-piperidine, monohydrochloride
Synonyms: 3-MeO PCP, 3-MeO Phencyclidine
MF: C₁₈H₂₇NO • HCl
FW: 309.9
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥5 years



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

Description

3-methoxy PCP is an analog of PCP which binds the NMDA receptor ($K_i = 20$ nM) and serotonin transporter ($K_i = 216$ nM) with greater affinity than does PCP ($K_i = 59$ and 2,234 nM, respectively).¹ The toxicological properties of this compound have not been reported. This product is intended for forensic and research purposes.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

Reference

1. 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, 08/22/2019

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