

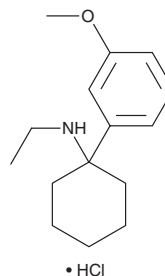
PRODUCT INFORMATION



3-methoxy-PCE (hydrochloride)

Item No. 9001355

CAS Registry No.: 1797121-52-8
Formal Name: N-ethyl-1-(3-methoxyphenyl)-cyclohexanamine, monohydrochloride
Synonym: 3-methoxy Eticyclidine
MF: C₁₅H₂₃NO • HCl
FW: 269.8
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 277 nm
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

3-methoxy PCE (hydrochloride) (Item No. 9001355) is an analytical reference standard that is structurally classified as an arylcyclohexylamine. It acts as an NMDA receptor antagonist with a pK_i value of 7.22.¹ This product has been sold online as a designer drug and is intended for forensic and research applications.

Reference

1. Roth, B.L., Gibbons, S., Arunotayanun, W., *et al.* The ketamine analogue methoxetamine and 3- and 4-methoxy analogues of phencyclidine are high affinity and selective ligands for the glutamate NMDA receptor. *PLoS One* **8**(3), e59334 (2013).

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, 07/12/2023

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