

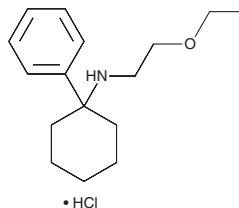
PRODUCT INFORMATION



PCEEA (hydrochloride)

Item No. 9001365

CAS Registry No.: 1798021-89-2
Formal Name: N-(2-ethoxyethyl)-1-phenylcyclohexanamine, monohydrochloride
MF: C₁₆H₂₅NO • HCl
FW: 283.8
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

PCEEA is a synthetic drug which is structurally related to the psychoactive compounds eticyclidine, ketamine, and phencyclidine. PCEEA has recently been identified as a likely drug of abuse.¹ The pharmacology of this compound has yet to be elucidated. This product is intended for forensic and research applications.

Reference

1. Sauer, C., Peters, F.T., Staack, R.F., *et al.* New designer drugs N-(1-phenylcyclohexyl)-2-ethoxyethanamine (PCEEA) and N-(1-phenylcyclohexyl)-2-methoxyethanamine (PCMEA): Studies on their metabolism and toxicological detection in rat urine using gas chromatographic/mass spectrometric techniques. *J. Mass Spectrom.* **43**(3), 305-316 (2008).

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, 07/07/2023

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