

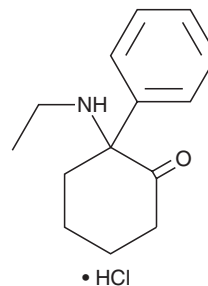
PRODUCT INFORMATION



deschloro-N-ethyl-Ketamine (hydrochloride)

Item No. 14105

CAS Registry No.: 4551-92-2
Formal Name: 2-(ethylamino)-2-phenyl-cyclohexanone, monohydrochloride
Synonyms: eb 2-oxo PCE, Eticyclidone, O-PCE
MF: C₁₄H₁₉NO • HCl
FW: 253.8
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

deschloro-N-ethyl-Ketamine is an analog of the dissociative anesthetic ketamine (Item No. 11630). Ketamine and its analogs have been abused recreationally.¹⁻⁴ The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research purposes.

References

1. Chakraborty, K., Neogi, R., and Basu, D. Club drugs: Review of the 'rave' with a note of concern for the Indian scenario. *Indian J. Med. Res.* **133**, 594-604 (2011).
2. Winstock, A.R. and Mitcheson, L. New recreational drugs and the primary care approach to patients who use them. *BMJ* **344**, 1-10 (2012).
3. Corazza, O., Schifano, F., Simonato, P., et al. Phenomenon of new drugs on the internet: The case of ketamine derivative methoxetamine. *Hum. Psychopharmacol.* **27**, 145-149 (2012).
4. Kim, J.Y., Shin, S.H., and In, M.K. Determination of amphetamine-type stimulants, ketamine and metabolites in fingernails by gas chromatography-mass spectrometry. *Forensic Sci. Int.* **194(1-3)**, 108-114 (2010).

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, 07/14/2023

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