

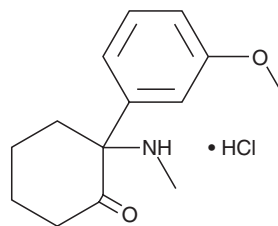
PRODUCT INFORMATION

Methoxmetamine (hydrochloride)

Item No. 17168



CAS Registry No.: 1781829-56-8
Formal Name: 2-(3-methoxyphenyl)-2-(methylamino)-cyclohexanone, monohydrochloride
Synonyms: 3-methoxy-2-oxo PCM, MMXE, MXM
MF: C₁₄H₁₉NO₂ • HCl
FW: 269.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

Methoxmetamine (hydrochloride) (Item No. 17168) is an analytical reference material that is structurally categorized as an amphetamine. It has been identified in ingredients sold as designer drugs.¹ The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research applications.

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

Reference

1. Kaizaki-Mitsumoto, A., Noguchi, N., Yamaguchi, S., *et al.* Three 25-NBOMe-type drugs, three other phenethylamine-type drugs (25I-NBMD, RH34, and escaline), eight cathinone derivatives, and a phenacyclidine analog MMXE, newly identified in ingredients of drug products before they were sold on the drug market. *Forensic Toxicol.* (2015).

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, 05/10/2022

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