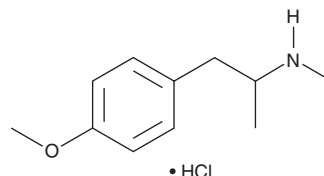


PRODUCT INFORMATION

para-Methoxymethamphetamine (hydrochloride)

Item No. 11562

CAS Registry No.: 3398-68-3
Formal Name: 4-methoxy-N,α-dimethyl-benzeneethanamine, monohydrochloride
Synonyms: 4-Methoxymethamphetamine, *p*-Methoxymethamphetamine, 4-MMA, PMMA
MF: C₁₁H₁₇NO • HCl
FW: 215.7
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 276, 283 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

para-Methoxymethamphetamine (PMMA) is a stimulant and psychedelic drug that is a structural analog of the amphetamine class serotonergic drug *para*-methoxyamphetamine (PMA).¹ In drug discrimination studies using rats, 3,4-methylenedioxymphetamine (MDMA) and PMMA produce similar, though not identical, discriminative stimulus effects.² PMMA has been found as a substitute in tablets and capsules of the MDMA sold as “ecstasy” and was linked to a number of overdose deaths in users who believed they were consuming recreational doses of MDMA.³ *para*-Methoxymethamphetamine is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6**(3), 259-274 (2005).
2. Glennon, R.A., Young, R., Dukat, M., et al. N-methyl-1-(4-methoxyphenyl)-2-aminopropane (PMMA) and N-methyl-1-(3,4-methylenedioxyphenyl)-2-aminopropane (MDMA) produce nonidentical discriminative stimuli in rats. *Pharmacol. Biochem. Behav.* **86**(3), 477-484 (2007).
3. Hager, M. Five B.C. deaths linked to lethal chemical PMMA. Retrieved June 20, 2012, from <http://www.vancouversun.com/news/calgary/Five+deaths+linked+lethal+chemical+PMMA/5993556/story.html>.

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, 07/18/2023

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