

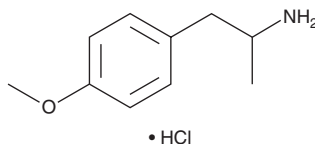
PRODUCT INFORMATION



4-Methoxyamphetamine (hydrochloride)

Item No. 12041

CAS Registry No.: 3706-26-1
Formal Name: 4-methoxy- α -methyl-benzeneethanamine, monohydrochloride
Synonyms: 4-MA, *p*-MA, *para*-Methoxyamphetamine, PMA
MF: C₁₀H₁₅NO • HCl
FW: 201.7
Purity: ≥98%
UV/Vis.: λ_{max} : 225, 277, 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

4-Methoxyamphetamine (4-MA), is a serotonergic drug of the amphetamine class.¹ 4-MA exhibits both structural and pharmacological similarity to 3,4-methylenedioxymethamphetamine and may be a more potent hallucinogen, particularly through its effects on serotonergic transmission and through reversible inhibition of the enzyme MAO-A.¹⁻⁴ 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. Kraner, J.C., McCoy, D.J., Evans, M.A., *et al.* Fatalities caused by the MDMA-related drug paramethoxyamphetamine (PMA). *J. Anal. Toxicol.* **25**(7), 645-648 (2001).
3. Martin-Iverson, M.T., Yamada, N., By, A.W., *et al.* "Designer" amphetamines: Effects on behavior and monoamines with or without reserpine and/or α -methyl-para-tyrosine pretreatment. *J. Psychiatr. Neurosci.* **16**(5), 253-261 (1991).
4. Tseng, L.F. 5-hydroxytryptamine uptake inhibitors block para-methoxyamphetamine-induced 5-HT release. *Br. J. Pharmacol.* **66**(2), 185-190 (1979).

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, 07/24/2023

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