

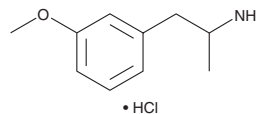
PRODUCT INFORMATION



3-Methoxyamphetamine (hydrochloride)

Item No. 9001916

CAS Registry No.: 35294-10-1
Formal Name: 3-methoxy- α -methyl-benzeneethanamine, monohydrochloride
Synonyms: 3-MA, *m*-MA, *meta*-Methoxyamphetamine, MMA
MF: C₁₀H₁₅NO • HCl
FW: 201.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 220, 274, 280 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

3-Methoxyamphetamine (3-MA) is a derivative of amphetamine that acts as a central nervous system stimulant by releasing serotonin, dopamine, and norepinephrine.¹⁻² It has appeared as a designer drug alternative to MDMA, along with the related compound 4-MA (Item No. 12041).³ This product is intended only for forensic and research purposes.

References

1. Menon, M.K., Tseng, L.F., and Loh, H.H. Pharmacological evidence for the central serotonergic effects of monomethoxyamphetamines. *J. Pharmacol. Exp. Ther.* **197(2)**, 272-279 (1976).
2. Tseng, L.F., Menon, M.K., and Loh, H.H. Comparative actions of monomethoxyamphetamines on the release and uptake of biogenic amines in brain tissue. *J. Pharmacol. Exp. Ther.* **197(2)**, 263-271 (1976).
3. Dal Cason, T.A. A re-examination of the mono-methoxy positional ring isomers of amphetamine, methamphetamine and phenyl-2-propanone. *Forensic Sci. Int.* **119(2)**, 168-194 (2001).

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, 07/13/2023

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