

PRODUCT INFORMATION



3-Methoxymethamphetamine (hydrochloride)

Item No. 24011

CAS Registry No.: 79896-22-3
Formal Name: 3-methoxy-N, α -dimethyl-benzeneethanamine, monohydrochloride
Synonyms: *m*-Methoxymethamphetamine, *meta*-Methoxymethamphetamine, 3-methoxy-N, α -dimethyl-Benzeneethanamine, *m*-methoxy-N, α -dimethyl-Benzeneethanamine, *meta*-methoxy-N, α -dimethyl-Benzeneethanamine

MF: C₁₁H₁₇NO • HCl

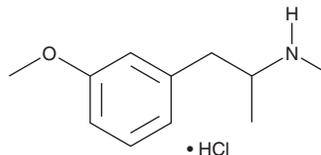
FW: 215.7

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Description

3-Methoxymethamphetamine (hydrochloride) (Item No. 24011) is an analytical reference standard categorized as an amphetamine.¹ This product is intended for research and forensic applications.

Reference

1. 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, 02/14/2025

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