

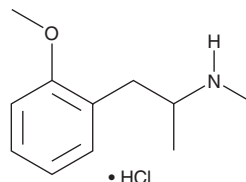
# PRODUCT INFORMATION



## 2-Methoxymethamphetamine (hydrochloride)

Item No. 9002184

**CAS Registry No.:** 5588-10-3  
**Formal Name:** 2-methoxy-N,α-dimethyl-benzeneethanamine, monohydrochloride  
**Synonyms:** 2-MeO-MA, o-MeO-MA, Methoxyphenamine, *ortho*-Methoxymethamphetamine, 2-MMA, NSC 65644  
**MF:** C<sub>11</sub>H<sub>17</sub>NO • HCl  
**FW:** 215.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 273, 279 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

2-Methoxymethamphetamine (hydrochloride) (Item No. 9002184) is an analytical reference standard categorized as an amphetamine.<sup>1</sup> 2-Methoxymethamphetamine has been found as an adverse analytical finding (AAF) during anti-doping testing.<sup>2,3</sup> 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).
2. Thevis, M., Sigmund, G., Koch, A., *et al.* Doping control analysis of methoxyphenamine using liquid chromatography-tandem mass spectrometry. *Eur. J. Mass Spectrom. (Chichester)* **14**(3), 145-152 (2008).
3. World Anti-Doping Agency. 2014 Anti-doping testing figures report (2014).

#### 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, 09/08/2023

#### 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