

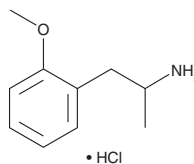
PRODUCT INFORMATION



2-Methoxyamphetamine (hydrochloride)

Item No. 9001915

CAS Registry No.: 72739-03-8
Formal Name: 2-methoxy- α -methyl-benzeneethanamine, monohydrochloride
Synonyms: 2-MA, NSC 1139, o-MA, OMA, *ortho*-Methoxyamphetamine
MF: C₁₀H₁₅NO • HCl
FW: 201.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 220, 272 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

2-Methoxyamphetamine is an amphetamine featuring a methoxy group in the 2, or *ortho*, position of the phenyl group. It is much weaker than either *para*- or *meta*-methoxyamphetamines at altering serotonin release or reuptake in rat brain tissues.¹ This product is intended for forensic and research applications.

Reference

1. 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).

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