

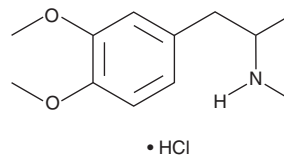
PRODUCT INFORMATION



3,4-DMMA (hydrochloride)

Item No. 11128

CAS Registry No.: 70932-18-2
Formal Name: 3,4-dimethoxy-N,α-dimethylbenzeneethanamine, monohydrochloride
Synonym: 3,4-Dimethoxymethamphetamine
MF: C₁₂H₁₉NO₂ • HCl
FW: 245.7
Purity: ≥97%
UV/Vis.: λ_{max}: 203, 230, 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

3,4-Dimethoxymethamphetamine (3,4-DMMA) (hydrochloride) is an analog of 3,4-methylenedioxymethamphetamine (MDMA), a psychoactive drug and research chemical of the phenethylamine and amphetamine chemical classes. DMMA appears to interfere with monoamine transport, inhibiting uptake of noradrenalin and serotonin transporters (K_i s = 22.8 and 7.7 μ M, IC_{50} s = 253.4 and 108 μ M, respectively).¹ DMMA is significantly less potent than MDMA (K_i s = 0.6 and 2.5 μ M; IC_{50} s = 6.6 and 34.8 μ M for MDMA inhibition of noradrenalin and serotonin transporters, respectively).¹ This product is intended for forensic and research applications.

Reference

1. Montgomery, T., Buon, C., Eibauer, S., *et al.* Comparative potencies of 3,4-methylenedioxymethamphetamine (MDMA) analogues as inhibitors of [³H]noradrenaline and [³H]5-HT transport in mammalian cell lines. *Br. J. Pharmacol.* **152**(7), 1121-1130 (2007).

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