

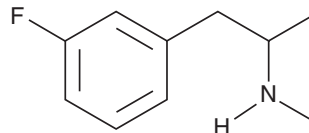
PRODUCT INFORMATION



3-Fluoromethamphetamine (hydrochloride)

Item No. 9001185

CAS Registry No.: 1324717-74-9
Formal Name: 1-(3-fluorophenyl)-N-methylpropan-2-amine, monohydrochloride
Synonym: 3-FMA
MF: C₁₀H₁₄FN • HCl
FW: 203.7
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



• HCl

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

Description

4-Fluoromethamphetamine (4-FMA) is a stimulant drug related to methamphetamine and 4-fluoroamphetamine. It has been identified as a component of designer drugs sold as bath salts or powders.¹ 3-FMA is an isomer of 4-FMA characterized by having fluorine at the 3 position of the phenyl group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Reference

1. Rösner, P., Quednow, B., Girreser, U., *et al.* Isomeric fluoro-methoxy-phenylalkylamines: A new series of controlled-substance analogues (designer drugs). *Forensic Sci. Int.* **148(2-3)**, 143-156 (2005).

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, 07/05/2023

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