

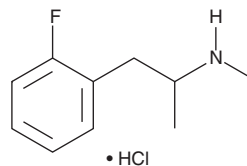
PRODUCT INFORMATION



2-Fluoromethamphetamine (hydrochloride)

Item No. 11420

CAS Registry No.: 1780004-19-4
Formal Name: 2-fluoro-N- α -dimethyl-benzeneethanamine, monohydrochloride
Synonym: 2-FMA
MF: C₁₀H₁₄FN • HCl
FW: 203.7
Purity: $\geq 98\%$
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-FMA is a stimulant drug related to methamphetamine and 2-fluoroamphetamine.¹ Through little is known of the pharmacological properties of fluoro methoxy substituted amphetamines, 2-FMA has been identified as a component of designer drugs sold as “legal high” replacements for restricted substances.^{1,2} The hydrochloride formulation of this compound is intended for use as a standard for the forensic analysis of samples that may contain 2-FMA.

References

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**, 143-156 (2005).
2. Camilleri, A., Johnston, M.R., Brennan, M., *et al.* Chemical analysis of four capsules containing the controlled substance analogues 4-methylmethcathinone, 2-fluoromethamphetamine, α -phthalimidopropiophenone and N-ethylcathinone. *Forensic Sci. Int.* **197**, 59-66 (2010).

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, 07/14/2023

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