

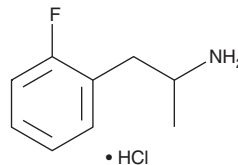
PRODUCT INFORMATION



2-Fluoroamphetamine (hydrochloride)

Item No. 11419

CAS Registry No.: 1626-69-3
Formal Name: 2-fluoro- α -methyl-benzeneethanamine, monohydrochloride
Synonym: 2-FA
MF: C₉H₁₂FN • HCl
FW: 189.7
Purity: \geq 98%
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

Halogenated amphetamines, including 4-fluoroamphetamine (4-FA), are psychostimulatory designer drugs which inhibit the uptake of monoamine neurotransmitters.^{1,2} 2-Fluoroamphetamine (2-FA) (hydrochloride) is a structural isomer of 4-FA, having the fluorine at the 2, rather than the 4, position. The physiological and toxicological properties of 2-FA are not known. This product is intended for research and forensic applications.

References

1. Marona-Lewicka, D., Rhee, G.S., Sprague, J.E., *et al.* Psychostimulant-like effects of *p*-flouroamphetamine in the rat. *Eur. J. Pharmacol.* **287(2)**, 105-113 (2011).
2. 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, 07/14/2023

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