

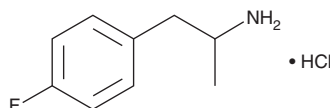
# PRODUCT INFORMATION



## 4-Fluoroamphetamine (hydrochloride)

Item No. 11156

**CAS Registry No.:** 64609-06-9  
**Formal Name:** 4-fluoro- $\alpha$ -methyl-benzeneethanamine, monohydrochloride  
**Synonyms:** 4-FA, *p*-FA  
**MF:** C<sub>9</sub>H<sub>12</sub>FN • 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 fluoroamphetamines (FAs), are psychostimulatory designer drugs.<sup>1,2</sup> *para*-Fluoroamphetamine (4-FA) (hydrochloride) inhibits the uptake of dopamine, serotonin, and norepinephrine with IC<sub>50</sub> values of 0.77, 6.8, and 0.42  $\mu$ M, respectively, indicating potency comparable to cocaine or methamphetamine.<sup>3</sup> Methodology for detecting 4-FA in serum and urine using GC/MS has recently been described.<sup>4</sup> This product is intended for forensic purposes.

### 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).
3. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
4. Rohrich, J., Becker, J., Kaufmann, T., *et al.* Detection of the synthetic drug 4-fluoroamphetamine (4-FA) in serum and urine. *Forensic Sci. Int.* **215(1-3)**, 3-7 (2011).

#### 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, 07/07/2023

#### 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