

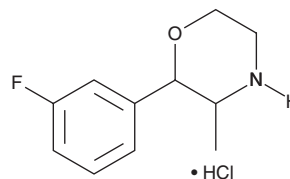
PRODUCT INFORMATION



3-Fluorophenmetrazine (hydrochloride)

Item No. 17208

CAS Registry No.: 1803562-83-5
Formal Name: 2-(3-fluorophenyl)-3-methylmorpholine, monohydrochloride
Synonyms: 3-fluoro Phenmetrazine, 3-FMP, 3-FPM, 3-fluoro PM
MF: C₁₁H₁₄FNO • HCl
FW: 231.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

Phenmetrazine is a stimulant with anorectic properties. Structurally, it is an amphetamine analog with the terminal amine incorporated into a morpholine ring. Phenmetrazine has long been used as an appetite suppressant, but the compound is susceptible to abuse.¹⁻³ 3-Fluorophenmetrazine is a halogenated phenmetrazine derivative. Halogenated amphetamines are psychostimulatory designer drugs that inhibit the uptake of monoamine neurotransmitters.^{4,5} The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Banks, M.L., Blough, B.E., Fennell, T.R., *et al.* Role of phenmetrazine as an active metabolite of phendimetrazine: Evidence from studies of drug discrimination and pharmacokinetics in rhesus monkeys. *Drug Alcohol Depend.* **130(1-3)**, 158-166 (2013).
2. Ghosh, M.N. and Parvathy, S. Tolerance pattern of the anorexigenic action of amphetamines, fenfluramine, phenmetrazine and diethylpropion in rats. *Br. J. Pharmacol.* **57(4)**, 479-485 (1976).
3. Martel, A. Preludin (phenmetrazine) in the treatment of obesity. *Can. Med. Assoc. J.* **76(2)**, 117-120 (1957).
4. Marona-Lewicka, D., Rhee, G.-S., Sprague, J.E., *et al.* Psychostimulant-like effects of *p*-flouroamphetamine in the rat. *Eur. J. Pharmacol.* **287**, 105-113 (2011).
5. 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).

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, 08/07/2019

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