

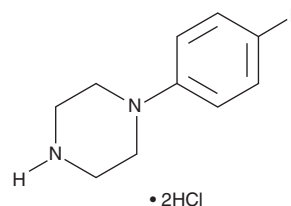
PRODUCT INFORMATION



1-(4-Fluorophenyl)piperazine (hydrochloride)

Item No. 11204

CAS Registry No.: 64090-19-3
Formal Name: 1-(4-fluorophenyl)-piperazine, dihydrochloride
Synonyms: 4-FPP, pFPP, 1-(p-Fluorophenyl)piperazine, 1-(para-Fluorophenyl)piperazine, NSC 149515
MF: C₁₀H₁₃FN₂ • 2HCl
FW: 253.1
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 288 nm
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

Phenylpiperazines are entactogenic drugs which direct central serotonin release.¹⁻³ Substituted phenylpiperazines, such as 1-(3-chlorophenyl)piperazine, have been identified as designer drugs or drugs of abuse.^{4,5} 1-(4-fluorophenyl)piperazine (hydrochloride) is a substituted phenylpiperazine with a potential for abuse. The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research applications.

References

1. Eriksson, E., Engberg, G., Bing, O., *et al.* Effects of mCPP on the extracellular concentrations of serotonin and dopamine in rat brain. *Neuropsychopharmacology* 20(3), 287-296 (1999).
2. Baumann, M.H., Clark, R.D., Budzynski, A.G., *et al.* N-substituted piperazines abused by humans mimic the molecular mechanism of 3,4-methylenedioxymethamphetamine (MDMA, or 'Ecstasy'). *Neuropsychopharmacology* 30, 550-560 (2005).
3. Hill, S.L. and Thomas, S.H. Clinical toxicology of newer recreational drugs. *Clin. Toxicol. (Phila)* 49, 705-719 (2011).
4. 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).
5. Arbo, M.D., Bastos, M.L., and Carmo, H.F. Piperazine compounds as drugs of abuse. *Drug Alcohol Depend.* [In press] 1-12 (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