

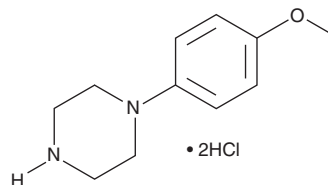
PRODUCT INFORMATION



1-(4-Methoxyphenyl)piperazine (hydrochloride)

Item No. 14020

CAS Registry No.: 38869-47-5
Formal Name: 1-(4-methoxyphenyl)-piperazine, dihydrochloride
Synonyms: 1-(*p*-Methoxyphenyl)piperazine,
1-(*para*-Methoxyphenyl)piperazine,
4-MeOPP, *p*-MeOPP, *para*-MeOPP
MF: C₁₁H₁₆N₂O • 2HCl
FW: 265.2
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 241, 294 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

1-(4-Methoxyphenyl)piperazine (hydrochloride) is a piperazine derivative with euphoric, stimulant properties comparable to those produced by amphetamine.^{1,2} It is assumed to have a mixed mechanism of action, with serotonergic and dopamine antagonistic properties in a similar fashion to 3,4-methylenedioxymethamphetamine. Also, its metabolism has been described.^{1,3} This product is intended for research and forensic purposes.

References

1. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6(3)**, 259-274 (2005).
2. Wohlfarth, A., Weinmann, W., and Dresen, S. LC-MS/MS screening method for designer amphetamines, tryptamines, and piperazines in serum. *Anal. Bioanal. Chem.* **396(7)**, 2403-2414 (2010).
3. Arbo, M.D., Bastos, M.L., and Carmo, H.F. Piperazine compounds as drugs of abuse. *Drug Alcohol Depend.* **122(3)**, 174-185 (2012).

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, 06/15/2023

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