

PRODUCT INFORMATION



1-(3-Trifluoromethylphenyl)piperazine (hydrochloride)

Item No. 11205

CAS Registry No.: 76835-14-8

Formal Name: 1-[3-(trifluoromethyl)phenyl]-piperazine, dihydrochloride

Synonyms: TFMPP,
1-(*m*-Trifluoromethylphenyl)piperazine,
1-(*meta*-Trifluoromethylphenyl)piperazine

MF: $C_{11}H_{13}F_3N_2 \cdot 2HCl$

FW: 303.2

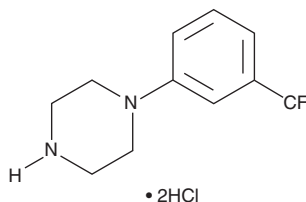
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 249, 293 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 5 years



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

Description

1-(3-Trifluoromethylphenyl)piperazine (TFMPP) is an entactogenic drug which selectively promotes the release of serotonin.¹ When used in combination with 1-benzylpiperazine, TFMPP increases both serotonin and dopamine, mirroring the effects of 3,4-methylenedioxymethamphetamine.¹ Drug-drug synergism may induce seizures.¹ TFMPP has been identified in party pills and powders.^{2,3} Its metabolism in rats has been described.⁴ This product is intended for forensic applications.

References

1. 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**(3), 550-560 (2005).
2. Hill, S.L. and Thomas, S.H. Clinical toxicology of newer recreational drugs. *Clin. Toxicol. (Phila)* **49**(8), 705-719 (2011).
3. Arbo, M.D., Bastos, M.L., and Carmo, H.F. Piperazine compounds as drugs of abuse. *Drug Alcohol Depend.* **122**(3), 174-185 (2011).
4. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6**(3), 259-274 (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, 07/07/2023

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