

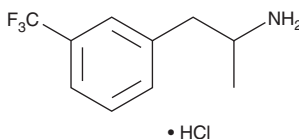
PRODUCT INFORMATION



3-Trifluoromethylamphetamine (hydrochloride)

Item No. 22604

CAS Registry No.: 673-18-7
Formal Name: α-methyl-3-(trifluoromethyl)-benzeneethanamine, monohydrochloride
Synonyms: desethyl Fenfluramine, (±)-Norfenfluramine
MF: C₁₀H₁₂F₃N • HCl
FW: 239.7
Purity: ≥98%
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

3-Trifluoromethylamphetamine (hydrochloride) (Item No. 22604) is an analytical reference standard that is categorized as an amphetamine. It is an active metabolite of fenfluramine and benfluorex that acts as an agonist at serotonin (5-HT) receptors (IC₅₀s = 6.92, 3.55, 0.17, and 2.09 μM for 5-HT_{1A}, 5-HT_{1B}, 5-HT_{1C}, and 5-HT₂ receptors, respectively, in murine brain).^{1,2} Norfenfluramine induces anorexia in mice and rats with EC₅₀ values of 6.8 and 0.8 mg/kg, i.p., respectively.¹ This product is intended for research and forensic applications.

References

1. Mennini, T., Bizzi, A., Caccia, S., *et al.* Comparative studies on the anorectic activity of d-fenfluramine in mice, rats, and guinea pigs. *Naunyn-Schmiedeberg's Arch. Pharmacol.* **343**(5), 483-490 (1991).
2. Curzon, G., Gibson, E.L., Oluyomi, A.O. Appetite suppression by commonly used drugs depends on 5-HT receptors but not on 5-HT availability. *Trends Pharmacol. Sci.* **18**(1), 21-25 (1997).

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