

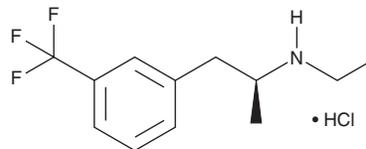
PRODUCT INFORMATION



(+)-Fenfluramine (hydrochloride)

Item No. 39830

CAS Registry No.: 3239-45-0
Formal Name: N-ethyl- α -methyl-3-(trifluoromethyl)-benzeneethanamine, monohydrochloride
Synonym: Dexfenfluramine
MF: C₁₂H₁₆F₃N • HCl
FW: 267.7
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

(+)-Fenfluramine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-fenfluramine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. (+)-Fenfluramine (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Fenfluramine is an agonist of the serotonin (5-HT) receptor subtypes 5-HT_{2B} and 5-HT_{2C}.¹ It induces serotonin release from superfused midbrain synaptosome preparations isolated from wild-type mice, but not 5-HT_{2B}^{-/-} mice, when used at concentrations of 0.5 or 1 μ M. (+)-Fenfluramine (3 or 10 mg/kg) reduces food intake in starved wild-type mice but not in starved 5-HT_{2B}^{-/-} mice. It reduces body weight gain in wild-type mice when infused at a dose of 2.5 mg/kg per day for 35 days. Formulations containing (+)-fenfluramine were previously used in the treatment of obesity.

Reference

1. Banas, S.M., Doly, S., Boutourlinsky, K., *et al.* Deconstructing antiobesity compound action: Requirement of serotonin 5-HT_{2B} receptors for dexfenfluramine anorectic effects. *Neuropsychopharmacology* **36(2)**, 423-433 (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, 06/18/2024

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