

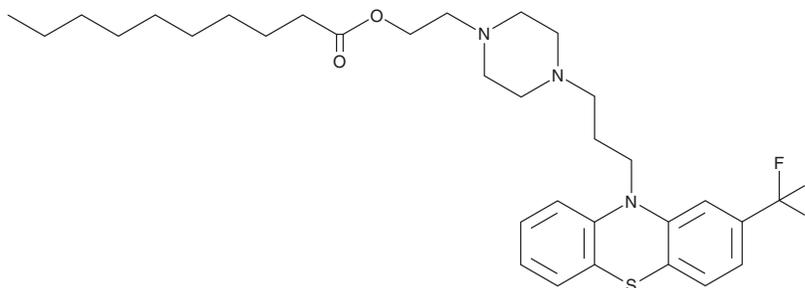
PRODUCT INFORMATION



Fluphenazine decanoate

Item No. 39903

CAS Registry No.: 5002-47-1
Formal Name: decanoic acid, 2-[4-[3-[2-(trifluoromethyl)-10H-phenothiazin-10-yl]propyl]-1-piperazinyl]ethyl ester
Synonyms: NSC 169510, QD-10733, SQ 10,733
MF: C₃₂H₄₄F₃N₃O₂S
FW: 591.8
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

Fluphenazine decanoate is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Fluphenazine decanoate is soluble in methanol.

Description

Fluphenazine decanoate is a prodrug form of the dopamine D₁ and D₂ receptor antagonist fluphenazine (Item No. 23555).¹ It induces random mouth movements in rats, an effect that can be reversed by the dopamine receptor antagonists raclopride (Item No. 17422) or SCH 23390 (Item No. 15631). Chronic administration of fluphenazine decanoate (3 mg/kg per day) increases serum levels of prolactin and decreases luteinizing hormone (LH) and follicle stimulating hormone (FSH), as well as decreases desire to mate and litter size, in male rats.² It increases lethargy and passive contact and reduces self-grooming, social grooming, auditory startle response, and propensity to attack other animals in monkeys when chronically administered.³ Formulations containing fluphenazine decanoate have been used in the treatment of chronic schizophrenia.

References

1. Stoessl, A.J., Dourish, C.T., and Iversen, S.D. Chronic neuroleptic-induced mouth movements in the rat: Suppression by CCK and selective dopamine D₁ and D₂ receptor antagonist. *Psychopharmacology (Berl.)* **98(3)**, 372-379 (1989).
2. Gill-Sharma, M.K., Aleem, M., Sethi, G., et al. Antifertility effects of fluphenazine in adult male rats. *J. Endocrinol. Invest.* **26(4)**, 316-326 (2003).
3. O'Keeffe, R.T. and Lifshitz, K. Nonhuman primates in neurotoxicity screening and neurobehavioral toxicity studies. *J. Am. Coll. Toxicol.* **8(1)**, 127-140 (1989).

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, 12/13/2023

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