

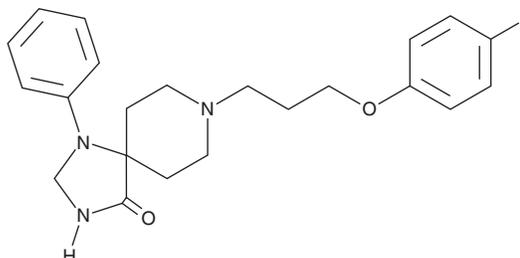
PRODUCT INFORMATION



AMI-193

Item No. 43100

CAS Registry No.: 510-74-7
Formal Name: 8-[3-(4-fluorophenoxy)propyl]-1-phenyl-1,3,8-triazaspiro[4.5]decan-4-one
Synonym: Spiramide
MF: C₂₂H₂₆FN₃O₂
FW: 383.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

AMI-193 is supplied as a solid. A stock solution may be made by dissolving the AMI-193 in the solvent of choice, which should be purged with an inert gas. AMI-193 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

AMI-193 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT_{2A} and the dopamine D₂ receptor.^{1,2} It is selective for the 5-HT_{2A} receptor over 5-HT_{2B} and 5-HT_{1C} receptors (K_is = 22.39, 977.24, and 1,548.82 nM, respectively) but does bind to the 5-HT_{1A} receptor (K_i = 50 nM). It selectively binds to the dopamine D₂ receptor over the dopamine D₁ receptor (K_is = 3 and 2,530 nM, respectively).² AMI-193 (0.05 mg/kg) completely inhibits amphetamine-induced increases in locomotor activity and rearing behavior in rats for at least three hours.³ It also reduces amphetamine-induced increases in locomotor activity and rearing behavior in mice when used at doses of 0.15 and 0.2 mg/kg but the effects only last for 30 minutes.

References

1. Knight, A.R., Misra, A., Quirk, K., *et al.* Pharmacological characterisation of the agonist radioligand binding site of 5-HT_{2A}, 5-HT_{2B} and 5-HT_{2C} receptors. *Naunyn Schmiedebergs Arch. Pharmacol.* **370(2)**, 114-123 (2004).
2. Ismaiel, A.M., De Los Angeles, J., Teitler, M., *et al.* Antagonism of 1-(2,5-dimethoxy-4-methylphenyl)-2-aminopropane stimulus with a newly identified 5-HT₂- versus 5-HT_{1C}-selective antagonist. *J. Med. Chem.* **36(17)**, 2519-2525 (1993).
3. Rolinski, Z. and Scheel-Krüger, J. The effect of dopamine and noradrenaline antagonists on amphetamine induced locomotor activity in mice and rats. *Acta Pharmacol. Toxicol. (Copenh)* **33(5)**, 385-399 (1973).

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, 05/08/2025

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