

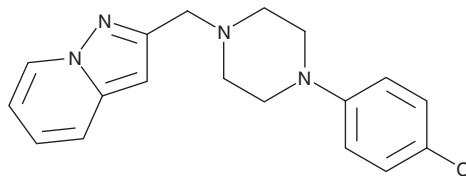
PRODUCT INFORMATION



FAUC-213

Item No. 42322

CAS Registry No.: 337972-47-1
Formal Name: 2-[[4-(4-chlorophenyl)-1-piperazinyl]methyl]-pyrazolo[1,5-a]pyridine
MF: C₁₈H₁₉ClN₄
FW: 326.8
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

FAUC-213 is supplied as a solid. A stock solution may be made by dissolving the FAUC-213 in the solvent of choice, which should be purged with an inert gas. FAUC-213 is soluble (≥10 mg/ml) in DMSO. FAUC-213 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

FAUC-213 is a dopamine D₄ receptor antagonist (K_i = 2.2 nM).¹ It is selective for dopamine D₄ receptors over dopamine D_{2L}, D_{2S}, and D₃ receptors (K_is = 3,400, 6,300, and 5,300 nM, respectively). FAUC-231 (7.5, 15, and 30 mg/kg) reduces amphetamine-induced hyperactivity in rats.²

References

1. Löber, S., Hübner, H., Utz, W., *et al.* Rationally based efficacy tuning of selective dopamine D₄ receptor ligands leading to the complete antagonist 2-[4-(4-chlorophenyl)piperazin-1-ylmethyl]pyrazolo[1,5-a]pyridine (FAUC 213). *J. Med. Chem.* **44**(17), 2691-2694 (2001).
2. Boeckler, F., Russig, H., Zhang, W., *et al.* FAUC 213, a highly selective dopamine D₄ receptor full antagonist, exhibits atypical antipsychotic properties in behavioural and neurochemical models of schizophrenia. *Psychopharmacology (Berl)* **175**(1), 7-17 (2004).

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, 12/16/2024

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