

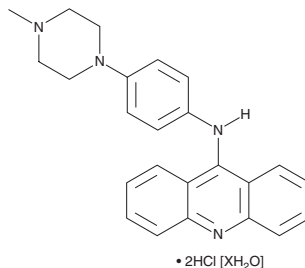
PRODUCT INFORMATION



JP-1302 (hydrochloride hydrate)

Item No. 25974

Formal Name: N-[4-(4-methyl-1-piperazinyl)phenyl]-9-acridinamine, dihydrochloride, hydrate
MF: $C_{24}H_{24}N_4 \cdot 2HCl [XH_2O]$
FW: 441.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

JP-1302 (hydrochloride hydrate) is supplied as a solid. A stock solution may be made by dissolving the JP-1302 (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. JP-1302 (hydrochloride hydrate) is soluble in the organic solvent DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

JP-1302 is an antagonist of α_2 -adrenergic receptors (α_2 -ARs) that is selective for α_{2C} -ARs over α_{2A} - and α_{2B} -ARs (K_i s = 28, 3,150, and 1,470 nM, respectively).¹ It reduces immobility time in the forced swim test in mice when administered at doses of 1, 3, and 10 μ mol/kg. JP-1302 (5 mg/kg) reverses phencyclidine-induced decreases in prepulse inhibition of the acoustic startle response in rats. It dose-dependently reduces haloperidol-induced bradykinesia and catalepsy in mice.² JP-1302 (0.3 mg/kg) also increases systolic blood pressure in rats.³

References

1. Sallinen, J., Höglund, I., Engström, M., *et al.* Pharmacological characterization and CNS effects of a novel highly selective α_{2C} -adrenoceptor antagonist JP-1302. *Br. J. Pharmacol.* **150**(4), 391-402 (2007).
2. Imaki, J., Mae, Y., Shimizu, S., *et al.* Therapeutic potential of α_2 adrenoceptor antagonism for antipsychotic-induced extrapyramidal motor disorders. *Neurosci. Lett.* **454**(2), 143-147 (2009).
3. Zefirov, T.L., Khisamieva, L.I., Ziyatdinova, N.I., *et al.* Effect of selective blockade of α_2 -adrenoceptor subtypes on cardiovascular system in rats. *Bull. Exp. Biol. Med.* **158**(4), 410-412 (2015).

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, 11/28/2022

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