

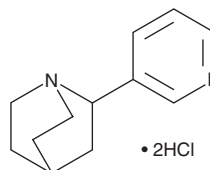
PRODUCT INFORMATION



RJR 2429 (hydrochloride)

Item No. 29745

CAS Registry No.: 1021418-53-0
Formal Name: 2-(3-pyridinyl)-1-azabicyclo[2.2.2]octane, dihydrochloride
MF: C₁₂H₁₆N₂ • 2HCl
FW: 261.2
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

RJR 2429 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RJR 2429 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RJR 2429 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of RJR 2429 (hydrochloride) in DMSO and water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

RJR 2429 is a nicotinic acetylcholine receptor (AChR) agonist.¹ It is selective for α4β2 or α7 subunit-containing nicotinic AChRs (K_is = 0.7 and 10.9 nM in isolated rat cortex and isolated rat hippocampus, respectively) over α1βγδ subunit-containing nicotinic AChRs (EC₅₀ = 1,000 nM in PC12 cells). RJR 2429 induces dopamine release from isolated rat striatal synaptosomes (EC₅₀ = 2.4 nM) and increases nicotine-induced elevations in epinephrine and norepinephrine levels in isolated rat adrenal gland when used at a concentration of 100 μM.^{2,3} It inhibits nicotine-induced ion flux in isolated rat thalamic synaptosomes (IC₅₀ = 154 nM).²

References

1. Bhatti, B., Strachan, J.-P., Breining, S.R., *et al.* Synthesis of 2-(pyridin-3-yl)-1-azabicyclo[3.2.2]nonane, 2-(pyridin-3-yl)-1-azabicyclo[2.2.2]octane, and 2-(pyridin-3-yl)-1-azabicyclo[3.2.1]octane, a class of potent nicotinic acetylcholine receptor-ligands. *J. Org. Chem.* **73(9)**, 3497-3507 (2008).
2. Bencherif, M., Schmitt, J.D., Bhatti, B.S., *et al.* The heterocyclic substituted pyridine derivative (+/-)-2-(3-pyridinyl)-1-azabicyclo[2.2.2]octane (RJR-2429): A selective ligand at nicotinic acetylcholine receptors. *J. Pharmacol. Exp. Ther.* **284(3)**, 886-894 (1998).
3. Playa, H., Lewis, T.A., Ting, A., *et al.* Dilazep analogues for the study of equilibrative nucleoside transporters 1 and 2 (ENT1 and ENT2). *Bioorg. Med. Chem. Lett.* **24(24)**, 5801-5804 (2014).

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, 11/01/2022

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