

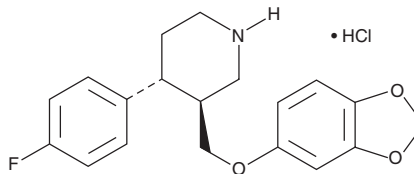
PRODUCT INFORMATION



Paroxetine (hydrochloride)

Item No. 14998

CAS Registry No.: 78246-49-8
Formal Name: (3S,4R)-3-[(1,3-benzodioxol-5-yloxy)methyl]-4-(4-fluorophenyl)-piperidine, monohydrochloride
Synonym: BRL 29060A
MF: C₁₉H₂₀FNO₃ • HCl
FW: 365.8
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 265, 271, 295 nm
Supplied as: A crystalline 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

Paroxetine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the paroxetine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Paroxetine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of paroxetine (hydrochloride) in ethanol and DMSO is approximately 20 mg/ml and approximately 33 mg/ml in DMF.

Paroxetine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, paroxetine (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Paroxetine (hydrochloride) has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMF:PBS (pH 7.0) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Paroxetine is a selective serotonin reuptake inhibitor (SSRI; K_i = 0.04 nM).¹ It is selective for the serotonin transporter (SERT) over the dopamine and norepinephrine transporters (K_s = 400 and 90 nM, respectively) as well as the serotonin (5-HT) receptor subtypes 5-HT_{1A} and 5-HT_{2A}, the histamine H₁ receptor, α₁- and α₂-adrenergic receptors (α₂-ARs), and muscarinic acetylcholine receptors (mAChRs; K_s = 21,168, 6,320, 13,746, 995, 3,915, and 42 nM, respectively).^{1,2} Paroxetine (5 mg/kg) decreases immobility time in the forced swim test in mice.³ Formulations containing paroxetine have been used in the treatment of depression, obsessive-compulsive disorder (OCD), panic disorder, social and generalized anxiety disorders, and post-traumatic stress disorder (PTSD).

References

1. Mattson, R.J., Catt, J.D., Denhart, D.J., *et al.* *J. Med. Chem.* Conformationally restricted homotryptamines. 2. Indole cyclopropylmethylamines as selective serotonin reuptake inhibitors. **48(19)**, 6023-6034 (2005).
2. Owens, M.J., Neal, W., Plott, S.J., *et al.* *J. Pharmacol. Exp. Ther.* Neurotransmitter receptor and transporter binding profile of antidepressants and their metabolites. **283(3)**, 1305-1322 (1997).
3. Sugimoto, Y., Tagawa, N., Kobayashi, Y., *et al.* Involvement of the sigma1 receptor in the antidepressant-like effects of fluvoxamine in the forced swimming test in comparison with the effects elicited by paroxetine. *Eur. J. Pharmacol.* **696(1-3)**, 96-100 (2012).

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, 12/07/2022

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