

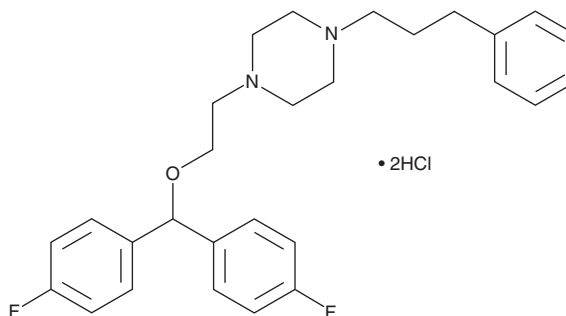
PRODUCT INFORMATION



GBR 12909 (hydrochloride)

Item No. 15622

CAS Registry No.: 67469-78-7
Formal Name: 1-[2-[bis(4-fluorophenyl)methoxy]ethyl]-4-(3-phenylpropyl)-piperazine, dihydrochloride
Synonym: I 893
MF: C₂₈H₃₂F₂N₂O • 2HCl
FW: 523.5
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

GBR 12909 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the GBR 12909 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GBR 12909 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GBR 12909 (hydrochloride) in these solvents is approximately 2 mg/ml.

Description

GBR 12909 is a potent inhibitor that blocks dopamine uptake (IC₅₀ = 1-51 nM).^{1,2} It is more than 100-fold less effective at blocking serotonin or noradrenaline uptake.^{1,2} GBR 12909 effectively inhibits dopamine uptake *in vivo*, leading to consequent stimulation of dopamine receptors.^{1,3} GBR 12909 also inhibits pyrilamine binding to the histamine H₁ receptor (IC₅₀ = 18 nM) but poorly inhibits ligand binding to dopamine, adrenergic, acetylcholine, serotonin, and γ-aminobutyric acid receptors.² It also blocks ligand binding to sigma receptors in rat brain (IC₅₀ = 48 nM).⁴

References

1. Heikkila, R.E. and Manzino, L. Behavioral properties of GBR 12909, GBR 13069 and GBR 13098: Specific inhibitors of dopamine uptake. *Eur. J. Pharmacol.* **103(3-4)**, 241-248 (1984).
2. Andersen, P.H. The dopamine uptake inhibitor GBR 12909: Selectivity and molecular mechanism of action. *Eur. J. Pharmacol.* **166(3)**, 493-504 (1989).
3. Melia, K.F. and Spealman, R.D. Pharmacological characterization of the discriminative-stimulus effects of GBR 12909. *J. Pharmacol. Exp. Ther.* **258(2)**, 626-632 (1991).
4. Contreras, P.C., Bremer, M.E., and Rao, T.S. GBR-12909 and fluspirilene potently inhibited binding of [³H](+)-3-PPP to sigma receptors in rat brain. *Life Sci.* **47(22)**, PL133-PL137 (1990).

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, 12/22/2022

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