

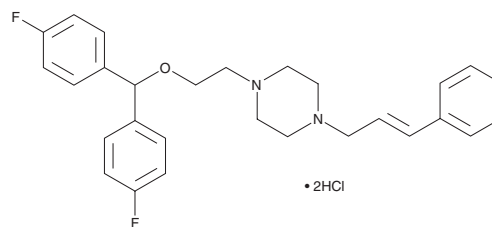
PRODUCT INFORMATION



GBR 13069 (hydrochloride)

Item No. 38903

CAS Registry No.: 67469-45-8
Formal Name: 1-[2-[bis(4-fluorophenyl)methoxy]ethyl]-4-(3-phenyl-2-propen-1-yl)-piperazine, dihydrochloride
MF: C₂₈H₃₀F₂N₂O • 2HCl
FW: 521.5
Purity: ≥98%
UV/Vis.: λ_{max}: 254 nm
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

GBR 13069 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the GBR 13069 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GBR 13069 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GBR 13069 (hydrochloride) in these solvents is approximately 20 and 10 mg/ml, respectively.

Description

GBR 13069 is a dopamine uptake inhibitor.¹ It selectively inhibits dopamine uptake over norepinephrine uptake in rat occipital cortex slices (IC₅₀s = 40 and 800 nM, respectively). GBR 13069 (2.5, 5, and 10 mg/kg) induces ipsilateral rotation and increases locomotion in rats. It also prevents 2'-CH₃-MPTP-induced decreases in neostriatal levels of the endogenous catecholamine dopamine (Item No. 36532) and its metabolites DOPAC (3,4-dihydroxyphenylacetic acid; Item No. 24912) and homovanillic acid (HVA; Item No. 27307) in mice when administered at a dose of 5 mg/kg.²

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. Sonsalla, P.K., Youngster, S.K., Kindt, M.V., et al. Characteristics of 1-methyl-4-(2'-methylphenyl)-1,2,3,6-tetrahydropyridine-induced neurotoxicity in the mouse. *J. Pharmacol. Exp. Ther.* **242(3)**, 850-857 (1987).

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, 08/22/2023

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