

PRODUCT INFORMATION



ALX 5407 (hydrochloride)

Item No. 30626

CAS Registry No.: 200006-08-2

Formal Name: N-[(3R)-3-([1,1'-biphenyl]-4-yloxy)-3-(4-fluorophenyl)propyl]-N-methyl-glycine, monohydrochloride

MF: C₂₄H₂₄FNO₃ • HCl

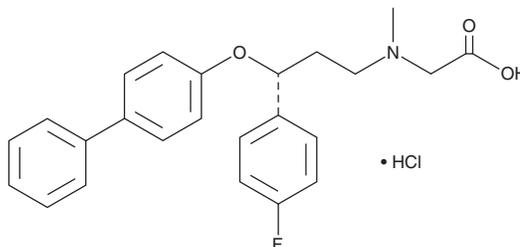
FW: 429.9

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

ALX 5407 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ALX 5407 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ALX 5407 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 20 mg/ml.

Description

ALX 5407 is an inhibitor of glycine transporter 1 (GlyT1; IC₅₀ = 3 nM).¹ It is selective for GlyT1 over GlyT2 (IC₅₀ = >10 μM), as well as over NMDA receptors and the inhibitory glycine receptor at 100 μM. ALX 5407 (10 mg/kg) increases prefrontal cortical levels of glycine in rats. It also reverses MK-801-induced persistent latent inhibition in a mouse model of schizophrenia.²

References

1. Atkinson, B.N., Bell, S.C., De Vivo, M., et al. ALX 5407: A potent, selective inhibitor of the hGlyT1 glycine transporter. *Mol. Pharmacol.* **60**(6), 1414-1420 (2001).
2. Lipina, T., Labrie, V., Weiner, I., et al. Modulators of the glycine site on NMDA receptors, D-serine and ALX 5407, display similar beneficial effects to clozapine in mouse models of schizophrenia. *Psychopharmacology (Berl)* **179**(1), 54-67 (2005).

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, 07/17/2023

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