

PRODUCT INFORMATION

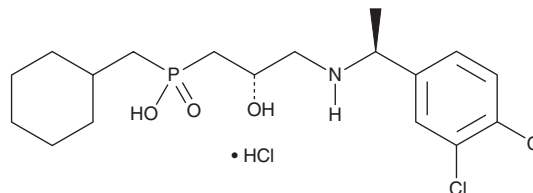


CGP 54626 (hydrochloride)

Item No. 20121

CAS Registry No.: 149184-21-4
Formal Name: P-(cyclohexylmethyl)-P-[(2S)-3-[[[(1S)-1-(3,4-dichlorophenyl)ethyl]amino]-2-hydroxypropyl]-phosphinic acid, monohydrochloride

Synonym: CGP 54626A
MF: C₁₈H₂₆Cl₂NO₃P • HCl
FW: 444.8
Purity: ≥95%
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

CGP 54626 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CGP 54626 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CGP 54626 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CGP 54626 (hydrochloride) in these solvents is approximately 0.2, 1, and 0.33 mg/ml, respectively.

CGP 54626 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CGP 54626 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CGP 54626 (hydrochloride) has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CGP 54626 is a selective antagonist of the GABA_B receptor with an IC₅₀ value of 4 nM.¹ It can be used to investigate the role of GABA_B receptors in neurological signaling.

Reference

1. Kaupmann, K., Malitschek, B., Schuler, V., *et al.* GABA_B-receptor subtypes assemble into functional heteromeric complexes. *Nature* **396(6712)**, 683-687 (1998).

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, 11/03/2022

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