

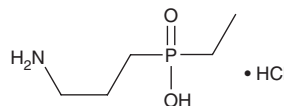
PRODUCT INFORMATION



CGP 36216 (hydrochloride)

Item No. 11985

CAS Registry No.: 1781834-71-6
Formal Name: *p*-(3-aminopropyl)-*p*-ethyl-phosphinic acid, monohydrochloride
MF: C₅H₁₄NO₂P • HCl
FW: 187.6
Purity: ≥98%
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 36216 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CGP 36216 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CGP 36216 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CGP 36216 (hydrochloride) in these solvents is approximately 5, 20, and 0.25 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CGP 36216 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CGP 36216 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CGP 36216 is a selective antagonist of GABA_B receptors (IC₅₀ = 43 μM).^{1,2} Its action is reversed by the GABA_B receptor agonist baclofen.² It is active at presynaptic, but not postsynaptic, receptors.²

References

1. Froestl, W., Mickel, S.J., von Sprecher, G., *et al.* Phosphinic acid analogues of GABA. 2. Selective, orally active GABA_B antagonists. *J. Med. Chem.* **38**, 3313-3331 (1995).
2. Ong, J., Bexis, S., Marino, V., *et al.* CGP 36216 is a selective antagonist at GABA_B presynaptic receptors in rat brain. *Eur. J. Pharmacol.* **415**, 191-195 (2001).

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, 09/30/2022

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