

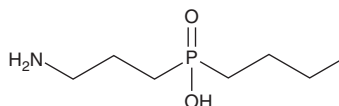
PRODUCT INFORMATION



CGP 36742

Item No. 35495

CAS Registry No.: 123690-78-8
Formal Name: P-(3-aminopropyl)-P-butyl-phosphinic acid
Synonyms: 3-Aminopropyl-n-butylphosphinic Acid, SGS 742
MF: C₇H₁₈NO₂P
FW: 179.2
Purity: ≥95%
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

CGP 36742 is supplied as a solid. A stock solution may be made by dissolving the CGP 36742 in the solvent of choice, which should be purged with an inert gas. CGP 36742 is soluble in chloroform.

Description

CGP 36742 is an antagonist of GABA_B receptors (IC₅₀ = 36 μM).¹ It is selective for GABA_B over a panel of 11 receptors, including GABA_A, at 1 mM. CGP 36742 (3 mg/kg) reduces paired-pulse inhibition of the late inhibitory postsynaptic potential (IPSP) in CA1 pyramidal neurons in rats. It increases latency to step-down in a passive avoidance test in mice in a dose-dependent manner.² CGP 36742 (10 and 30 mg/kg) decreases immobility time in the forced swim test in mice.³

References

1. Olpe, H.R., Steinmann, M.W., Ferrat, T., *et al.* The actions of orally active GABA_B receptor antagonists on GABAergic transmission in vivo and in vitro. *Eur. J. Pharmacol.* **233**(2-3), 179-186 (1993).
2. Mondadori, C., Jaekel, J., and Preiswerk, G. CGP 36742: The first orally active GABA_B blocker improves the cognitive performance of mice, rats, and rhesus monkeys. *Behav. Neural Biol.* **60**(1), 62-68 (1993).
3. Nowak, G., Partyka, A., Palucha, A., *et al.* Antidepressant-like activity of CGP 36742 and CGP 51176, selective GABA_B receptor antagonists, in rodents. *Br. J. Pharmacol.* **149**(5), 581-590 (2006).

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, 12/12/2022

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