

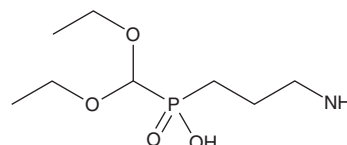
PRODUCT INFORMATION



CGP 35348

Item No. 18599

CAS Registry No.: 123690-79-9
Formal Name: P-(3-aminopropyl)-P-(diethoxymethyl)-phosphinic acid
MF: C₈H₂₀NO₄P
FW: 225.2
Purity: ≥95%
Supplied as: A solid
Storage: 4°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CGP 35348 is supplied as a solid. A stock solution may be made by dissolving the CGP 35348 in water. We do not recommend storing the aqueous solution for more than one day.

Description

The amino acid γ -aminobutyric acid (GABA) is an inhibitory neurotransmitter that acts through two families of heteromeric ligand-gated ion channels, GABA_A and GABA_C, and a G protein-coupled receptor, GABA_B. CGP 35348 is a selective, brain-accessible GABA_B receptor antagonist (IC₅₀ = 34 μ M).¹ It blocks the action of the GABA_B-selective agonist baclofen (Item No. 18600) at postsynaptic sites but not at presynaptic receptors.^{1,2} CGP 35348 prevents baclofen-induced antinociception and muscle relaxation in mice and rats.³ It is commonly used to investigate the role of GABA_B activation in neurological signaling.⁴⁻⁶

References

1. Olpe, H.R., Karlsson, G., Pozza, M.F., *et al.* CGP 35348: A centrally active blocker of GABA_B receptors. *Eur. J. Pharmacol.* **187(1)**, 27-38 (1990).
2. Hills, J.M., Larkin, M.M., and Howson, W. A comparison of the relative activities of a number of GABA_B antagonists in the isolated vas deferens of the rat. *Br. J. Pharmacol.* **102(3)**, 631-634 (1991).
3. Malcangio, M., Ghelardini, C., Giotti, A., *et al.* CGP 35348, a new GABA_B antagonist, prevents antinociception and muscle-relaxant effect induced by baclofen. *Br. J. Pharmacol.* **103(2)**, 1303-1308 (1991).
4. Koek, W., Cheng, J., and Rice, K.C. Discriminative stimulus effects of the GABA_B receptor-positive modulator rac-BHFF: Comparison with GABA_B receptor agonists and drugs of abuse. *J. Pharmacol. Exp. Ther.* **344(3)**, 553-560 (2013).
5. Gillani, Q., Ali, M., and Iqbal, F. CGP 35348, GABA_B receptor antagonist, has a potential to improve neuromuscular coordination and spatial learning in albino mouse following neonatal brain damage. *Biomed. Res. Int.* **2014**:295215, (2014).
6. Ding, L., Gao, R., Xiong, X.-Q., *et al.* GABA in paraventricular nucleus regulates adipose afferent reflex in rat. *PLoS One* **10(8)**, (2015).

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, 10/12/2017

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