

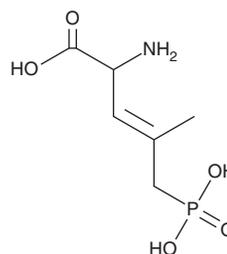
PRODUCT INFORMATION



CGP 37849

Item No. 43660

CAS Registry No.: 127910-31-0
Formal Name: 2-amino-4-methyl-5-phosphono-3E-pentenoic acid
MF: C₆H₁₂NO₅P
FW: 209.1
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 37849 is supplied as a solid. A stock solution may be made by dissolving the CGP 37849 in water. CGP 37849 is soluble (≥10 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

CGP 37849 is an NMDA receptor antagonist ($K_i = 220$ nM).¹ It decreases the number of seizures induced by maximal electroshock (MES) in rats ($ED_{50} = 21$ mg/kg). CGP 37849 (0.62, 1.25, or 2.5 mg/kg) reduces immobility time in the forced swim test.²

References

1. Fagg, G.E., Olpe, H.R., Pozza, M.F., *et al.* CGP 37849 and CGP 39551: Novel and potent competitive N-methyl-D-aspartate receptor antagonists with oral activity. *Br. J. Pharmacol.* **99(4)**, 791-797 (1990).
2. Przegaliński, E., Tatarczyńska, E., Dereń-Wesołek, A., *et al.* Antidepressant-like effects of a partial agonist at strychnine-insensitive glycine receptors and a competitive NMDA receptor antagonist. *Neuropharmacology* **36(1)**, 31-37 (1997).

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, 07/08/2025

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