

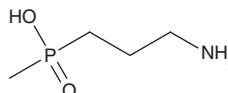
PRODUCT INFORMATION



SKF 97541

Item No. 43784

CAS Registry No.: 127729-35-5
Formal Name: P-(3-aminopropyl)-P-methyl-phosphinic acid
Synonym: CGP 35024
MF: $C_4H_{12}NO_2P$
FW: 137.1
Purity: $\geq 98\%$
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

SKF 97541 is supplied as a crystalline solid. SKF 97541 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

SKF 97541 is a GABA_B receptor agonist.¹ It reduces forskolin-induced cAMP accumulation in rat cortical slices when used at concentrations of 30 and 100 μM .² SKF 97541 reduces excitatory postsynaptic potentials (EPSPs) induced by bicuculline (Item No. 11727) in rat striatal slices ($\text{EC}_{50} = 0.092 \mu\text{M}$), an effect that can be blocked by CGP 35348 (Item No. 18599).³ It also hyperpolarizes neurons in the substantia nigra in rat brain slices ($\text{EC}_{50} = 0.15 \mu\text{M}$). SKF 19541 (0.01 and 0.05 mg/kg) reduces the time spent immobile in the forced swim test in rats without reducing basal locomotor activity.⁴

References

1. Froestl, W., Mickel, S.J., von Sprecher, G., et al. Phosphinic acid analogues of GABA. 2. Selective, orally active GABAB antagonists. *J. Med. Chem.* **38**(17), 3313-3331 (1995).
2. Wierońska, J.M., Kłeczek, N., Woźniak, M., et al. mGlu₅-GABA_B interplay in animal models of positive, negative and cognitive symptoms of schizophrenia. *Neurochem. Int.* **88**, 97-109 (2015).
3. Seabrook, G.R., Howson, W., and Lacey, M.G. Electrophysiological characterization of potent agonists and antagonists at pre- and postsynaptic GABA_B receptors on neurones in rat brain slices. *Br. J. Pharmacol.* **101**(4), 949-957 (1990).
4. Frankowska, M., Filip, M., and Przegaliński, E. Effects of GABA_B receptor ligands in animal tests of depression and anxiety. *Pharmacol. Rep.* **59**(6), 645-655 (2007).

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, 07/28/2025

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