

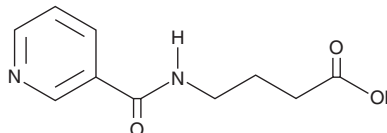
PRODUCT INFORMATION



Picamilon

Item No. 29397

CAS Registry No.: 34562-97-5
Formal Name: 4-[(3-pyridinylcarbonyl)amino]-butanoic acid
Synonym: Nicotinoyl-GABA
MF: C₁₀H₁₂N₂O₃
FW: 208.2
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

Picamilon is supplied as a crystalline solid. A stock solution may be made by dissolving the picamilon in the solvent of choice, which should be purged with an inert gas. Picamilon is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Picamilon is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, picamilon should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Picamilon has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Picamilon is a derivative of γ -aminobutyric acid (GABA).¹ It increases parietal cortical blood flow in rats and rabbits when administered at doses of 50 mg/kg, i.p. and 10 mg/kg, i.v., respectively.² Picamilon (10 mg/kg) increases latency to enter the dark compartment in the passive avoidance conditioned reflex test in rats compared with control animals, indicating nootropic effects.¹

References

1. Kopelevich, V.M., Bulanova, L.N., Grigor'ev, I.A., *et al.* Synthesis and study of psychotropic and hypotensive properties of new picamilon derivatives. *Pharm. Chem. J.* **31(10)**, 536-539 (1997).
2. Mirzoyan, R.S., Gan'shina, T.S., Kosoi, M.Y., *et al.* Effect of picamilon on the cerebral cortical blood supply and microcirculation in the pial arteriolar system. *Biull. Eksp. Biol. Med.* **107(5)**, 668-670 (1989).

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, 12/09/2022

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