

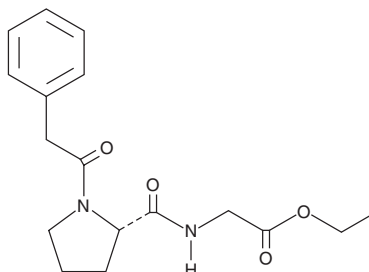
PRODUCT INFORMATION



N-Phenylacetyl-L-Prolylglycine ethyl ester

Item No. 14496

CAS Registry No.: 157115-85-0
Formal Name: 1-(2-phenylacetyl)-L-prolylglycine, ethyl ester
Synonyms: GVS-111, Noopept, SGS 111
MF: C₁₇H₂₂N₂O₄
FW: 318.4
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

N-Phenylacetyl-L-prolylglycine ethyl ester is a synthetic dipeptide that has been shown to produce positive nootropic and cognitive effects in animal models (0.01-0.8 mg/kg) by a mechanism similar to other related racetam compounds.¹ It can rescue neuroblastoma SH-SY5Y cells from amyloid-induced cytotoxicity, indicating a potential application in the treatment of neurodegenerative diseases.² It also demonstrates antioxidant and anti-inflammatory effects, which have been examined in the context of an anti-diabetic agent.³ This compound was recently identified in illegal designer herbal products distributed in Japan and is intended for research and forensic purposes only.⁴

References

1. Gudasheva, T.A., Voronins, T.A., Ostrovskaya, R.U., *et al.* Synthesis and anti-amnesic activity of a series of IV-acylprolyl-containing dipeptides. *Eur. J. Med. Chem.* **31(2)**, 151-157 (1996).
2. Jia, X., Gharibyan, A.L., Öhman, A., *et al.* Neuroprotective and nootropic drug noopept rescues α-synuclein amyloid cytotoxicity. *J. Mol. Biol.* **414(5)**, 699-712 (2011).
3. Ostrovskaya, R.U., Ozerova, I.V., Gudasheva, T.A., *et al.* Efficiency of noopept in streptozotocin-induced diabetes in rats. *Bull. Exp. Biol. Med.* **154(3)**, 334-338 (2013).
4. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Identification of two new-type designer drugs, piperazine derivative MT-45 (I-C6) and synthetic peptide Noopept (GVS-111), with synthetic cannabinoid A-834735, cathinone derivative 4-methoxy-α-PVP, and phenethylamine derivative 4-methylbuphedrine from illegal products. *Forensic Toxicol.* **32**, 9-18 (2014).

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, 06/19/2023

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