

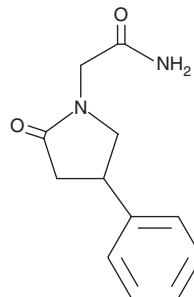
PRODUCT INFORMATION



Phenylpiracetam

Item No. 13897

CAS Registry No.: 77472-70-9
Formal Name: 2-oxo-4-phenyl-1-pyrrolidineacetamide
Synonym: Carphedon
MF: C₁₂H₁₄N₂O₂
FW: 218.3
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

Phenylpiracetam is a nootropic derivative of the GABAergic agent piracetam and has been evaluated for its efficacy in reducing neurological deficits related to cerebral ischemia, epilepsy, and neurobehavioral disorders.^{1,2} At 10-100 mg/kg in mice, phenylpiracetam can increase locomotor and antidepressant activity in the open-field and forced swimming tests, as well as enhance memory in a passive avoidance response test.¹ Because of its cognitive enhancing and general stimulant properties, phenylpiracetam has the potential to be abused recreationally and is listed as a banned substance by the World Anti-Doping Agency and the International Olympic Committee.³ This product is intended for research and forensic purposes only.

References

1. Zvejniece, L., Svalbe, B., Veinberg, G., *et al.* Investigation into stereoselective pharmacological activity of phenotropil. *Basic Clin. Pharmacol. Toxicol.* **109(5)**, 407-412 (2011).
2. Stewart, A.M. and Kalueff, A.V. The developing utility of zebrafish models for cognitive enhancers research. *Curr. Neuropharmacol.* **10(3)**, 263-271 (2012).
3. Kim, S., Park, J.H., Myung, S.W., *et al.* Determination of carphedon in human urine by solid-phase microextraction using capillary gas chromatography with nitrogen-phosphorus detection. *Analyst* **124(11)**, 1559-1562 (1999).

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, 07/25/2023

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