

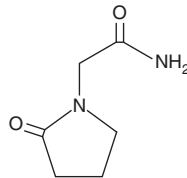
# PRODUCT INFORMATION



## Piracetam

Item No. 20755

**CAS Registry No.:** 7491-74-9  
**Formal Name:** 2-oxo-1-pyrrolidineacetamide  
**Synonym:** UCB 6215  
**MF:** C<sub>6</sub>H<sub>10</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 142.2  
**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

Piracetam (Item No. 20755) is an analytical reference standard that is structurally categorized as a nootropic. It is a cyclic derivative of  $\gamma$ -aminobutyric acid (GABA) that improves learning, memory, brain metabolism, and capacity in animal models.<sup>1,2</sup> It has shown efficacy in stroke rehabilitation in animal models.<sup>3,4</sup> Although its mechanism of action remains unclear, piracetam has been shown to alter the physical properties of cell membranes and to protect against hypoxia, which might underlie neuroprotective and antithrombotic actions.<sup>2</sup> This product is intended for research and forensic applications.

### 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. Winnicka, K., Tomasiak, M. and Bielawska, A. Piracetam - An old drug with novel properties? *Acta. Pol. Pharm.* **62**(5), 405-409 (2005).
3. Engelter, S.T. Safety in pharmacological enhancement of stroke rehabilitation. *Eur. J. Phys. Rehabil. Med.* **49**(2), 261-267 (2013).
4. Wheble, P.C.R., Sena, E.S. and Macleod, M.R. A systematic review and meta-analysis of the efficacy of piracetam and piracetam-like compounds in experimental stroke. *Cerebrovasc. Dis.* **25**(1-2), 5-11 (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, 06/22/2023

#### 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