

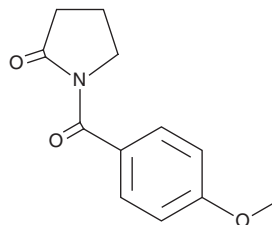
PRODUCT INFORMATION



Aniracetam

Item No. 17752

CAS Registry No.: 72432-10-1
Formal Name: 1-(4-methoxybenzoyl)-2-pyrrolidinone
Synonyms: Memodrin, Ro 13-5057, Sarpul
MF: C₁₂H₁₃NO₃
FW: 219.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 281 nm
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

Aniracetam (Item No. 17752) is an analytical reference standard that is classified as a nootropic. It can block AMPA receptor desensitization and, consequently, enhance excitatory activity by several fold.^{1,2} It can also allosterically modulate metabotropic glutamate receptors.³ The therapeutic value of aniracetam has been explored in various rodent models for the treatment of cognitive impairment or other CNS disorders.³ This product is intended for forensic and research purposes only.

References

1. Bräuner-Osborne, H., Egebjerg, J., Nielsen, E.O., *et al.* Ligands for glutamate receptors: Design and therapeutic prospects. *J. Med. Chem.* **43**(14), 2609-2645 (2000).
2. Lawrence, J.J., Brenowitz, S., and Trussell, L.O. The mechanism of action of aniracetam at synaptic α-amino-3-hydroxy-5-methyl-4-isoxazolepropionic acid (AMPA) receptors: Indirect and direct effects on desensitization. *Mol. Pharmacol.* **64**(2), 269-278 (2003).
3. Nakamura, K. Aniracetam: Its novel therapeutic potential in cerebral dysfunctional disorders based on recent pharmacological discoveries. *CNS Drug Rev.* **8**(1), 70-89 (2002).

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, 06/16/2023

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