

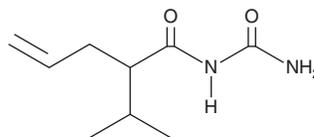
PRODUCT INFORMATION



Apronal

Item No. 34097

CAS Registry No.: 528-92-7
Formal Name: N-(aminocarbonyl)-2-(1-methylethyl)-4-pentenamide
Synonyms: Allylisopropylacetylcarbamide,
Allylisopropylacetylurea
MF: C₉H₁₆N₂O₂
FW: 184.2
Purity: ≥98%
Supplied as: A 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

Apronal is supplied as a solid. A stock solution may be made by dissolving the apronal in the solvent of choice, which should be purged with an inert gas. Apronal is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of apronal in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Apronal is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, apronal should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Apronal has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Apronal is a muscle relaxant.¹ It prolongs methylhexital-induced sleep and enhances morphine-induced analgesia in dogs, mice, and rats. Apronal has been used to induce experimental porphyria in rabbits.² Formulations containing apronal have previously been used as daytime sedatives.

References

1. Yajima, T., Uritani, K., Aoki, R., *et al.* Pharmacological studies on allylisopropylacetylurea (apronalide), a drug with central nervous system activity. *Oyo Yakuri* **11(5)**, 693-717 (1976).
2. Schmid, R.M. and Schwartz, S. Experimental porphyria. III. Hepatic type produced by sedormid. *Proc. Soc. Exp. Biol. Med.* **81(3)**, 685-689 (1952).

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, 12/12/2022

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