

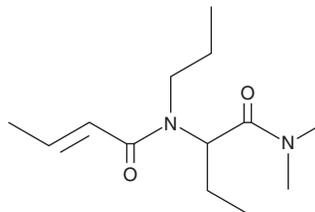
PRODUCT INFORMATION



Cropropamide

Item No. 30888

CAS Registry No.: 633-47-6
Formal Name: N-[1-[(dimethylamino)carbonyl]propyl]-N-propyl-2-butenamide
Synonym: Propyl Butamide
MF: C₁₃H₂₄N₂O₂
FW: 240.3
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Cropropamide is supplied as a neat oil. A stock solution may be made by dissolving the cropropamide in the solvent of choice, which should be purged with an inert gas. Cropropamide is soluble in organic solvents such as chloroform and DMSO.

Description

Cropropamide is a component of the respiratory stimulant prethcamide.¹ It increases locomotor activity in rats when administered at doses of 33.69 and 59.97 mg/kg.

Reference

1. Babbini, M., Gaiardi, M., and Bartoletti, M. Some behavioral effects of prethcamide compared with those of its two components. *Pharmacology* **14**(5), 455-463 (1976).

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, 11/14/2022

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