

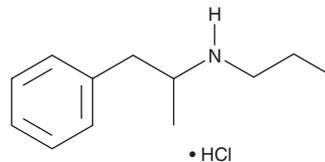
PRODUCT INFORMATION



N-Propylamphetamine (hydrochloride)

Item No. 14205

CAS Registry No.: 59877-57-5
Formal Name: α -methyl-N-propyl-benzeneethanamine, monohydrochloride
Synonyms: NPA, α -methyl-N-propyl-Benzeneethanamine
MF: $C_{12}H_{19}N \cdot HCl$
FW: 213.8
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

N-Propylamphetamine is an N-substituted amphetamine with a propyl group on the amphetamine backbone. It has about one-fourth the potency of N-ethylamphetamine (Item No. 11557) in impairing the performance of rats in various tests.¹ The toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Reference

1. Woolverton, W.L., Shybut, G., and Johanson, C.E. Structure-activity relationships among some d-N-alkylated amphetamines. *Pharmacol. Biochem. Behav.* **13(6)**, 869-876 (1980).

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, 06/15/2023

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