

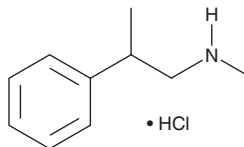
PRODUCT INFORMATION



N-methyl-2-Phenylpropan-1-amine (hydrochloride)

Item No. 18563

CAS Registry No.: 5969-39-1
Formal Name: N,β-dimethyl-benzeneethanamine, monochloride
Synonyms: Benzedrine, β,N-MePEA, Phenpromethamine, Vonedrine
MF: C₁₀H₁₅N • HCl
FW: 185.7
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

N-methyl-2-Phenylpropan-1-amine (hydrochloride) (Item No. 18563) is an analytical reference standard that is classified as a phenethylamine.¹ It is a positional isomer of methamphetamine that has been sold as a dietary supplement and subsequently banned by the World Anti-Doping Agency.^{2,3} This product is intended for research and forensic applications.

References

1. Warren, R.J., Begosh, P.P., and Zarembo, J.E. Identification of amphetamines and related sympathomimetic amines. *J. Assoc. Off. Anal. Chem.* **54**(5), 1179-1191 (1971).
2. Lisi, A., Hasick, N., Kazlauskas, R., et al. Studies of methylhexanamine in supplements and geranium oil. *Drug Test. Anal.* **3**(11-12), 873-876 (2011).
3. Liu, Y. and Santillo, M.F. Cytochrome P450 2D6 and 3A4 enzyme inhibition by amine stimulants in dietary supplements. *Drug Test. Anal.* **7**(9) (2015).

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