

PRODUCT INFORMATION

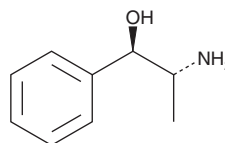


(-)-Norpseudoephedrine

Item No. 45995

CAS Registry No.: 37577-07-4
Formal Name: (αR)-α-[(1R)-1-aminoethyl]-benzenemethanol
Synonyms: (-)-*threo*-2-Amino-2-methyl-1-phenylethanol,
(1R,2R)-2-Amino-1-phenylpropan-1-ol,
L-Norpseudoephedrine, (1R,2R)-Norpseudoephedrine,
(1R,2R)-(-)-Norpseudoephedrine,
(-)-Pseudonorephedrine

MF: C₉H₁₃NO
FW: 151.2
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.

Description

(-)-Norpseudoephedrine (Item No. 45995) is an analytical reference standard categorized as an amphetamine. (-)-Norpseudoephedrine is an active metabolite of (R)-(+)-cathinone.¹ It increases locomotor activity and potentiates the analgesic effect of morphine (Item Nos. ISO60147 | 15464 | 25955) in mice but inhibits its discriminatory effects.^{1,2} (-)-Norpseudoephedrine is regulated as a List I compound in the United States. This product is intended for research and forensic applications.

References

1. Nencini, P., Fraioli, S., Pascusci, T., *et al.* (-)-Norpseudoephedrine, a metabolite of cathinone with amphetamine-like stimulus properties, enhances the analgesic and rate decreasing effects of morphine, but inhibits its discriminative properties. *Behav. Brain Res.* **92**(1), 11-20 (1998).
2. Fairchild, M.D. and Alles, G.A. The central locomotor stimulatory activity and acute toxicity of the ephedrine and norephedrine isomers in mice. *J. Pharmacol. Exp. Ther.* **158**(1), 135-139 (1967).

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, 08/13/2026

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