

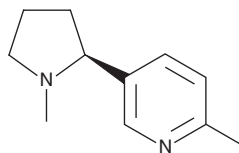
PRODUCT INFORMATION



(S)-6-Methylnicotine

Item No. 45678

CAS Registry No.: 13270-56-9
Formal Name: 2-methyl-5-[(2S)-1-methyl-2-pyrrolidinyl]-pyridine
Synonyms: 1'-methyl Nicotine, 6-MN
MF: C₁₁H₁₆N₂
FW: 176.3
Purity: ≥98%
Supplied as: A 10 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Description

(S)-6-Methylnicotine (Item No. 45678) is an analytical reference standard categorized as an alkaloid and nicotinic acetylcholine receptor (nAChR) agonist.^{1,2} (S)-6-Methylnicotine has been found in e-cigarette products and pouches and in cured leaf tobacco marketed as additive free, as well as other tobacco products. This product is intended for research and forensic applications.

References

1. Pankow, J.F., Luo, W., McWhirter, K.J., *et al.* Levels of the nicotine analog 6-methyl nicotine as a naturally formed tobacco alkaloid in tobacco and tobacco products. *Sci. Rep.* **15(1)**, 17945 (2025).
2. Xing, H., Andrud, K.W., Soti, F., *et al.* A methyl scan of the pyrrolidinium ring of nicotine reveals significant differences in its interactions with $\alpha 7$ and $\alpha 4 \beta 2$ nicotinic acetylcholine receptors. *Mol. Pharmacol.* **98(2)**, 168-180 (2020).

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, 05/05/2026

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