

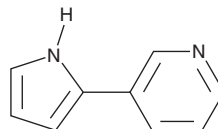
PRODUCT INFORMATION



β-Nornicotyrine

Item No. 42944

CAS Registry No.: 494-98-4
Formal Name: 3-(1H-pyrrol-2-yl)-pyridine
Synonyms: Nornicotyrine, NNT, β-NNT
MF: C₉H₈N₂
FW: 144.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

β-Nornicotyrine is supplied as a solid. A stock solution may be made by dissolving the β-nornicotyrine in the solvent of choice, which should be purged with an inert gas. β-Nornicotyrine is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

β-Nornicotyrine is an alkaloid that has been found in *N. tabacum*.¹ It has been found in settled dust from tobacco smoker-occupied houses and as a contaminant in soils used to grow *N. tabacum*.^{2,3}

References

1. Hellinghausen, G., Roy, D., Wang, Y., *et al.* A comprehensive methodology for the chiral separation of 40 tobacco alkaloids and their carcinogenic E/Z-(R,S)-tobacco-specific nitrosamine metabolites. *Talanta* **181**, 132-141 (2018).
2. Richardot, W.H., Hamzai, L., Ghukasyan, T., *et al.* Novel chemical contaminants associated with thirdhand smoke in settled house dust. *Chemosphere* **352**, 141138 (2024).
3. Zhou, P., Luo, Q., Bai, Y., *et al.* Tobacco cultivation leads to the accumulation of alkaloids in the soil and causes potential risks. *Microchem. J.* **208**, 112431 (2025).

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, 05/05/2025

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