

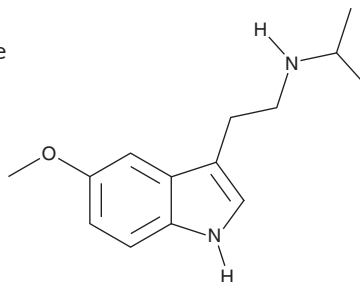
PRODUCT INFORMATION



5-methoxy NiPT

Item No. 35267

CAS Registry No.: 109921-55-3
Formal Name: 5-methoxy-N-(1-methylethyl)-1H-indole-3-ethanamine
Synonym: 5-methoxy-N-iso-Propyltryptamine
MF: C₁₄H₂₀N₂O
FW: 232.3
Purity: ≥95%
UV/Vis.: λ_{max}: 224 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥2 years



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

Description

5-methoxy NiPT (Item No. 35267) is an analytical reference standard categorized as a tryptamine. 5-methoxy NiPT is a metabolite of 5-methoxy DiPT (Item Nos. 11865 | 15697).¹ It has been detected in hair samples obtained by law enforcement from suspected drug users. This product is intended for research and forensic applications.

Reference

1. Shi, Y., Wang, R., Yuan, S., *et al.* UHPLC-MS/MS method for simultaneously detecting 16 tryptamines and their metabolites in human hair and applications to real forensics cases. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **1159**, 122392 (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, 10/27/2025

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