

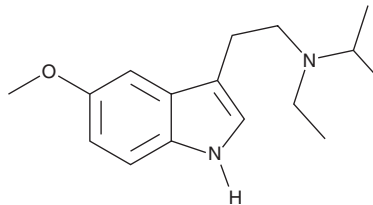
PRODUCT INFORMATION



5-methoxy EiPT

Item No. 17299

CAS Registry No.: 850032-66-5
Formal Name: N-ethyl-5-methoxy-N-(1-methylethyl)-1H-indole-3-ethanamine
Synonym: 5-methoxy-N-ethyl-N-Isopropyltryptamine
MF: C₁₆H₂₄N₂O
FW: 260.4
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

5-methoxy EiPT is a tryptamine-based designer drug with structural similarity to 5-methoxy DiPT (Item No. 11865) and 5-methoxy MiPT (Item No. 11482). While analytical methods have been developed for its identification, the biological activity of 5-methoxy EiPT has not been reported.¹ This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

Reference

1. Nakazono, Y., Tsujikawa, K., Kuwayama, K., *et al.* Simultaneous determination of tryptamine analogues in designer drugs using gas chromatography mass spectrometry and liquid chromatography-tandem mass spectrometry. *Forensic Toxicol.* **32**, 154-161 (2014).

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, 09/07/2022

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