

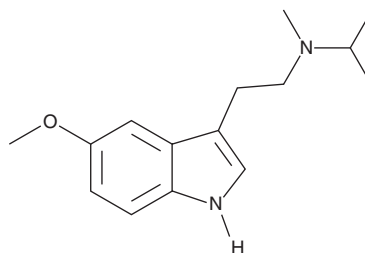
PRODUCT INFORMATION



5-methoxy MiPT

Item No. 11482

CAS Registry No.: 96096-55-8
Formal Name: 5-methoxy-N-methyl-N-(1-methylethyl)-1H-Indole-3-ethanamine
Synonym: 5-MeO MiPT
MF: C₁₅H₂₂N₂O
FW: 246.4
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 276 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

5-methoxy MiPT is a psychedelic tryptamine related to N-methyl-N-isopropyltryptamine. It potently inhibits the re-uptake of the monoamines serotonin and norepinephrine (IC₅₀s = 6.4 and 2.6 μM, respectively), but does not affect dopamine re-uptake.¹ 5-methoxy MiPT has no effect on the release of monoamines from rat brain synaptosomes.¹ This product is intended for forensic and research applications.

Reference

1. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).

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, 07/17/2023

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