

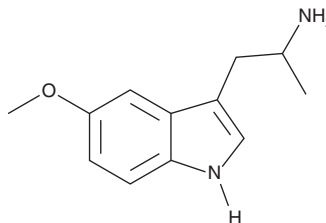
PRODUCT INFORMATION



5-methoxy AMT

Item No. 11553

CAS Registry No.: 1137-04-8
Formal Name: 5-methoxy- α -methyl-1H-indole-3-ethanamine
Synonym: 5-MeO AMT
MF: C₁₂H₁₆N₂O
FW: 204.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223, 278 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 AMT (5-MeO AMT) is a potent psychoactive analog of 5-methoxy DiPT (Item No. 11865) and α -methyltryptamine (AMT), tryptamine derivatives that inhibit reuptake and stimulate release of monoamines.¹ 5-MeO AMT inhibits re-uptake (IC_{50} s = 0.18, 2.9, and 3.37 μ M) and stimulates release (EC_{50} s = 1.5, 460, and 8.9 μ M) of dopamine, serotonin, and norepinephrine, respectively from rat brain synaptosomes.¹ 5-MeO AMT, sometimes called Alpha-O, has been sold under the guise of LSD and abused as a recreational drug.² This tryptamine analog is intended only to be used to aid the forensic analysis of samples that may contain this compound or for scientific research applications.

References

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).
2. Zimmerman, M.M. The identification of 5-methoxy- α -methyltryptamine (5-MeO-AMT). In: United States Drug Enforcement Administration [Internet]. Available from: http://www.justice.gov/dea/programs/forensicsci/microgram/journal071203/mj071203_pg1.html.

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, 07/18/2023

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