

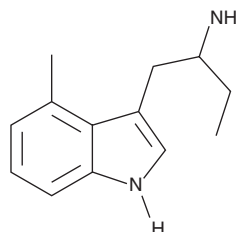
PRODUCT INFORMATION



4-methyl- α -Ethyltryptamine

Item No. 13886

CAS Registry No.: 28289-30-7
Formal Name: α -ethyl-4-methyl-1H-indole-3-ethanamine
Synonym: 4-methyl- α -ET
MF: C₁₃H₁₈N₂
FW: 202.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 224, 276 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

α -Ethyltryptamine (AET), like α -methyltryptamine (Item No. 11135), is a psychedelic drug with schedule 1 status in the United States. Both alter monoamine release and re-uptake in the brain.^{1,2} 4-methyl- α -Ethyltryptamine (Item No. 13886) is a homolog of AET. Its physiological and toxicological properties remain to be elucidated. This product is intended for forensic and research applications.

References

1. Gey, K.F., and Pletscher, A. Effect of α -alkylated tryptamine derivatives on 5-hydroxytryptamine metabolism *in vivo*. *Br. J. Pharmacol. Chemother.* **19**(1), 161-167 (1962).
2. 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, 07/25/2023

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