

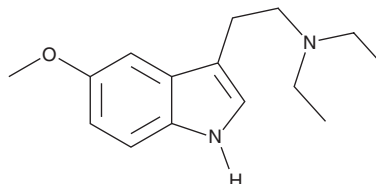
PRODUCT INFORMATION



5-methoxy DET

Item No. 17300

CAS Registry No.: 1218-40-2
Formal Name: N,N-diethyl-5-methoxy-1H-indole-3-ethanamine
Synonym: 5-methoxy-N,N-Diethyltryptamine
MF: C₁₅H₂₂N₂O
FW: 246.4
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 DET is a synthetic tryptamine designer drug structurally similar to the potent hallucinogen N,N-DMT (Item No. 13959). It inhibits serotonin uptake with an IC₅₀ value of 2.4 μM and activates 5-HT_{2A} receptors with an EC₅₀ value of 8.11 nM.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. Blough, B.E., Landavazo, A., Decker, A.M., *et al.* Interaction of psychoactive tryptamines with biogenic amine transporters and serotonin receptor subtypes. *Psychopharmacology (Berl)* **231**(21), 4135-4144 (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, 01/27/2023

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