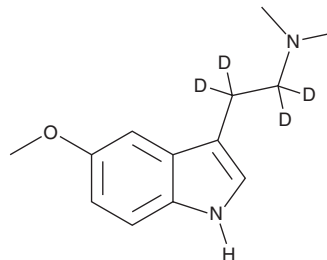


PRODUCT INFORMATION



5-methoxy DMT-d₄ Item No. 38748

CAS Registry No.: 66521-37-7
Formal Name: 5-methoxy-N,N-dimethyl-1H-indole-3-ethan- $\alpha,\alpha,\beta,\beta$ -d₄-amine
Synonym: 5-methoxy-N,N-Dimethyltryptamine-d₄
MF: C₁₃H₁₄D₄N₂O
FW: 222.3
Chemical Purity: $\geq 98\%$ (5-methoxy DMT)
Deuterium Incorporation: $\geq 99\%$ deuterated forms (d₁-d₄); $\leq 1\%$ d₀
UV/Vis.: λ_{max} : 225 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 2 years



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

Description

5-methoxy DMT-d₄ (Item No. 38748) is intended for use as an internal standard for the quantification of 5-methoxy DMT (Item Nos. 11480 | 37552 | 11628) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

5-methoxy DMT is categorized as a tryptamine.¹ It induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential. 5-methoxy DMT substitutes for LSD (Item Nos. 35189 | 28289) in a two-key drug discrimination test.² 5-methoxy DMT-d₄ is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Singleton, C. and Marsden, C.A. Circadian variation in the head twitch response produced by 5-methoxy-N1,N1-dimethyltryptamine and *p*-chloroamphetamine in the mouse. *Psychopharmacology (Berl)* **74(2)**, 173-176 (1981).
2. Nielsen, E.B. Discriminative stimulus properties of lysergic acid diethylamide in the monkey. *J. Pharmacol. Exp. Ther.* **234(1)**, 244-249 (1985).

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, 05/19/2023

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