

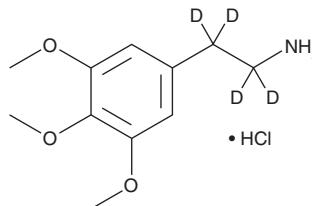
PRODUCT INFORMATION



Mescaline-d₄ (hydrochloride)

Item No. 34432

CAS Registry No.: 2762464-15-1
Formal Name: 2-(3,4,5-trimethoxyphenyl)ethan-1,1,2,2-d₄-1-amine, monohydrochloride
Synonym: TMPEA-d₄
MF: C₁₁H₁₃D₄NO₃ • HCl
FW: 251.7
Chemical Purity: ≥95% (Mescaline)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
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

Mescaline-d₄ (hydrochloride) (Item No. 34432) is intended for use as an internal standard for the quantification of mescaline (Item No. 11950) 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).

Mescaline is categorized as a phenethylamine.¹ It has been found in peyote cacti (*L. williamsii*). Mescaline induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential.² Mescaline-d₄ is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Monte, A.P., Waldman, S.R., Marona-Lewicka, D., *et al.* Dihydrobenzofuran analogues of hallucinogens. 4. Mescaline derivatives. *J. Med. Chem.* **40**(19), 2997-3008 (1997).
2. Halberstadt, A.L., Chatha, M., Klein, A.K., *et al.* Correlation between the potency of hallucinogens in the mouse head-twitch response assay and their behavioral and subjective effects in other species. *Neuropharmacology* **167**, 107933 (2020).

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, 06/28/2023

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