

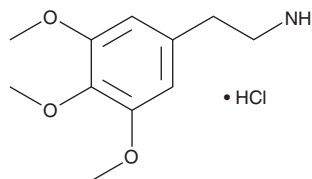
PRODUCT INFORMATION



Mescaline (hydrochloride) (exempt preparation)

Item No. 15680

CAS Registry No.: 832-92-8
Formal Name: 3,4,5-trimethoxy-benzeethanamine, monohydrochloride
Synonyms: NSC 30419, TMPEA
MF: C₁₁H₁₇NO₃ • HCl
FW: 247.7
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Mescaline (hydrochloride) (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of mescaline (hydrochloride) (exempt preparation) in these solvents is approximately 10, 3, and 0.5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of mescaline (hydrochloride) (exempt preparation) is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of mescaline (hydrochloride) (exempt preparation) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Mescaline (hydrochloride) (exempt preparation) (Item No. 15680) is an analytical reference standard categorized as a phenethylamine.¹ Mescaline has been found in peyote cacti (*L. williamsii*). It induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential.² Mescaline is regulated as a Schedule I compound in the United States. Mescaline (hydrochloride) (exempt preparation) (Item No. 15680) is provided as a DEA exempt preparation. 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, 10/03/2022

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