

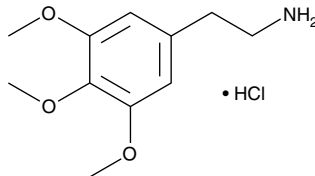
Product Information



Mescaline (hydrochloride)

Item No. 11950

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%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that mescaline (hydrochloride) be stored as supplied at 20°C. It should be stable for at least two years.

Mescaline (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the mescaline (hydrochloride) in the solvent of choice. Mescaline (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of mescaline (hydrochloride) 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. Organic solvent-free aqueous solutions of mescaline (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of mescaline (hydrochloride) in PBS, pH 7.2, is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Mescaline is a psychedelic phenethylamine that is found naturally in plants predominantly in the cactus family, including peyote (*L. williamsii*). It binds and activates serotonin receptors, with preference for 5-HT_{2A} over 5-HT_{2C} (EC₅₀ = 4 and 24 μM, respectively).¹ This product is intended for research and forensic purposes.

Reference

1. Monte, A.P., Waldman, S.R., Marona-Lewicka, D., *et al.* Dihydrobenzofuran analogues of hallucinogens. 4. Mescaline derivatives. *J. Med. Chem.* **40**, 2997-3008 (1997).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/11950

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/28/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com