

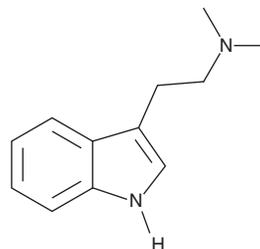
PRODUCT INFORMATION



N,N-DMT (exempt preparation)

Item No. 15694

CAS Registry No.: 61-50-7
Formal Name: N,N-dimethyl-1H-indole-3-ethanamine
Synonyms: N,N-Dimethyltryptamine, NSC 63795
MF: C₁₂H₁₆N₂
FW: 188.3
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 283 nm
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

N,N-DMT (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 N,N-DMT (exempt preparation) in these solvents is approximately 20, 5, and 11 mg/ml, respectively.

N,N-DMT (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of N,N-DMT (exempt preparation) should be diluted with the aqueous buffer of choice. N,N-DMT (exempt preparation) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N,N-DMT (exempt preparation) (Item No. 15694) is an analytical reference standard categorized as a tryptamine.¹ N,N-DMT substitutes for LSD (D-tartrate) (Item No. 28289) in a two-lever drug discrimination test. N,N-DMT is regulated as a Schedule I compound in the United States. N,N-DMT (exempt preparation) (Item No. 15694) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Blair, J.B., Marona-Lewicka, D., Kanthasamy, A., *et al.* Thieno[3,2-b]- and thieno[2,3-b]pyrrole bioisosteric analogues of the hallucinogen and serotonin agonist N,N-dimethyltryptamine. *J. Med. Chem.* **42(6)**, 1106-1111 (1999).

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, 10/13/2020

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