

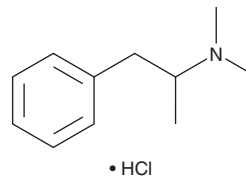
PRODUCT INFORMATION



N,N-DMA (hydrochloride)

Item No. 11556

CAS Registry No.: 33286-27-0
Formal Name: N,N,α-trimethyl-benzeneethanamine, monohydrochloride
Synonyms: Dimetamfetamine, N,N-Dimethylamphetamine, Metrotonin, NSC 27146, NSC 75736, T 312
MF: C₁₁H₁₇N • HCl
FW: 199.7
Purity: ≥98%
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

N,N-Dimethylamphetamine (N,N-DMA) is a regulated stimulant that is an N-methylated analog of methamphetamine.¹ It is often found as an impurity in preparations of methamphetamine.^{2,3} DMA itself is less active and less neurotoxic than methamphetamine.^{1,4} The metabolites of DMA that are detectable in human urine have been characterized.^{5,6} This product is intended for research and forensic applications.

References

1. Witkin, J.M., Ricaurte, G.A., and Katz, J.L. Behavioral effects of N-methylamphetamine and N,N-dimethylamphetamine in rats and squirrel monkeys. *J. Pharmacol. Exp. Ther.* **253**(2), 466-474 (1990).
2. Kuwayama, K., Tsujikawa, K., Miyaguchi, H., *et al.* Identification of impurities and the statistical classification of methamphetamine using headspace solid phase microextraction and gas chromatography-mass spectrometry. *Forensic Sci. Int.* **160**(1), 44-52 (2006).
3. Qi, Y., Evans, I.D., and McCluskey, A. Australian Federal Police seizures of illicit crystalline methamphetamine ('ice') 1998-2002: Impurity analysis. *Forensic Sci. Int.* **164**(2-3), 201-210 (2006).
4. Yao, B.B., Hsieh, G.C., Frost, J.M., *et al.* *In vitro* and *in vivo* characterization of A-796260: A selective cannabinoid CB₂ receptor agonist exhibiting analgesic activity in rodent pain models. *Br. J. Pharmacol.* **153**(2), 390-401 (2008).
5. Sato, M., Hida, M., and Nagase, H. Analysis of dimethylamphetamine and its metabolites in human urine by liquid chromatography-electrospray ionization-mass spectrometry with direct sample injection. *Forensic Sci. Int.* **128**(3), 146-154 (2002).
6. Katagi, M., Tatsuno, M., Miki, A., *et al.* Discrimination of dimethylamphetamine and methamphetamine use: Simultaneous determination of dimethylamphetamine-N-oxide and other metabolites in urine by high-performance liquid chromatography-electrospray ionization mass spectrometry. *J. Anal. Toxicol.* **24**(5), 354-358 (2000).

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, 07/18/2023

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