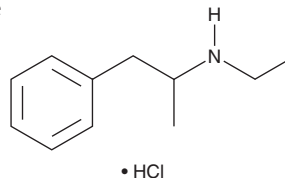


PRODUCT INFORMATION

N-Ethylamphetamine (hydrochloride) (exempt preparation)

Item No. 15658

CAS Registry No.: 16105-78-5
Formal Name: N-ethyl-α-methyl-benzeneethanamine, monohydrochloride
Synonyms: EMA, MG 19973, NSC 54938
MF: C₁₁H₁₇N • HCl
FW: 199.7
Purity: ≥98%
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥5 years



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

Description

EMA is an N-substituted amphetamine with an ethyl group on the amphetamine backbone. As a stimulant drug of the phenethylamine and amphetamine chemical classes, it is widely known as an anorectic marketed under the tradename ApetinilTM.^{1,2} EMA has been abused as a “designer drug” alternative to amphetamine and possibly methamphetamine.¹ EMA (hydrochloride) (Item No. 15658) is provided as a DEA exempt preparation. This product is intended for forensic or research purposes.

References

1. Lekskulchai, V. and Mookhavesa, C. Evaluation of Roche Abuscreen ONLINE amphetamine immunoassay for screening of new amphetamine analogues. *J. Anal. Toxicol.* **25(6)**, 471-475 (2001).
2. Nagai, T., Kanaya, H., Matsushima, K., et al. Time-lapse changes of d- and l-enantiomers of racemic (dl)-ethylamphetamine in human urine. *J. Anal. Toxicol.* **21(2)**, 112-115 (1997).

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, 07/20/2026

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