

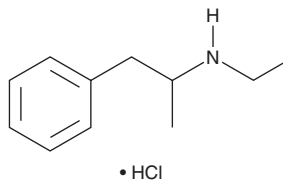
PRODUCT INFORMATION



N-Ethylamphetamine (hydrochloride)

Item No. 11557

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: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

N-Ethylamphetamine (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 Apetinil™.^{1,2} EMA has been abused as a “designer drug” alternative to amphetamine and possibly methamphetamine.¹ This product is intended for forensic and research purposes.

References

1. Lekskulchai, V. and Mokkhaveva, 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, 07/18/2023

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