

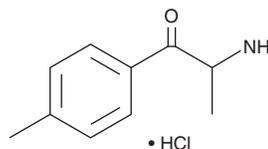
PRODUCT INFORMATION



nor-Mephedrone (hydrochloride)

Item No. 9000940

CAS Registry No.: 6941-17-9
Formal Name: 2-amino-1-(4-methylphenyl)-1-propanone, monohydrochloride
Synonyms: Mephedrone Metabolite, 4-Methylcathinone, NSC 60487
MF: C₁₀H₁₃NO • HCl
FW: 199.7
Purity: ≥98%
UV/Vis.: λ_{max}: 255 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

nor-Mephedrone (hydrochloride) (Item No. 9000940) is an analytical reference standard categorized as a cathinone metabolite.^{1,2} It is a metabolite of mephedrone (Item Nos. 10801 | 26277). This product is intended for research and forensic applications.

References

1. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 468-482 (2010).
2. Meyer, M.R., Wilhelm, J., Peters, F.T., *et al.* Beta-keto amphetamines: Studies on the metabolism of the designer drug mephedrone and toxicological detection of mephedrone, butylone, and methylene in urine using gas chromatography-mass spectrometry. *Anal. Bioanal. Chem.* **397(3)**, 1225-1233 (2010).

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, 08/27/2025

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