

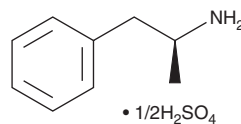
PRODUCT INFORMATION



D-Amphetamine (hemisulfate)

Item No. 26204

CAS Registry No.: 51-63-8
Formal Name: αS-methyl-benzeneethanamine, hemisulfate
Synonyms: Dextroamphetamine, (+)-α-Methylphenethylamine, NSC 27104
MF: C₉H₁₃N • 1/2H₂SO₄
FW: 184.2
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

D-Amphetamine (hemisulfate) (Item No. 26204) is an analytical reference standard categorized as an amphetamine.¹ Amphetamine is a psychotropic compound that is abused recreationally.² D-Amphetamine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Martin-Iverson, M.T., Yamada, N., By, A.W., *et al.* "Designer" amphetamines: Effects on behavior and monoamines with or without reserpine and/or α-methyl-para-tyrosine pretreatment. *J. Psychiatr. Neurosci.* **16(5)**, 253-261 (1991).
2. Blanckaert, P., van Amsterdam, J.G.C., Brunt, T.M., *et al.* 4-Methyl-amphetamine: A health threat for recreational amphetamine users. *J. Psychopharmacol.* **27(9)**, 817-822 (2013).

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, 07/19/2023

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