

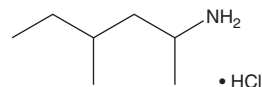
PRODUCT INFORMATION



Methylhexanamine (hydrochloride)

Item No. 10749

CAS Registry No.: 13803-74-2
Formal Name: 4-methyl-2-hexanamine, monohydrochloride
Synonyms: Dimethylamylamine, DMAA, Floradrene, Forthane, Geranamine, NSC 1106
MF: C₇H₁₇N • HCl
FW: 151.7
Purity: ≥95%
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

Methylhexanamine (hydrochloride) is a simple aliphatic amine that was once marketed as a nasal decongestant (Forthane) but is now sold as a bodybuilding supplement (Floradrene, Geranamine). It is also sold as a mild stimulant in the form of party pills and has been reported to cause cerebral hemorrhage in a case report.¹ While little is known about its mode of action, methylhexanamine interferes in immunoassays for amphetamines, presumably because of the common ethylamine structure.² This compound is intended for forensic or research purposes.

References

1. Gee, P., Jackson, S., and Easton, J. Another bitter pill: A case of toxicity from DMAA party pills. *N. Z. Med. J.* **123(1327)**, 124-127 (2010).
2. Vorce, S.P., Holler, J.M., Cawrse, B.M., et al. Dimethylamylamine: A drug causing positive immunoassay results for amphetamines. *J. Anal. Toxicol.* **35(3)**, 183-187 (2011).

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, 06/15/2023

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