

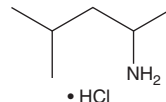
# PRODUCT INFORMATION



## 1,3-Dimethylbutylamine (hydrochloride)

Item No. 21928

CAS Registry No.: 71776-70-0  
Formal Name: 4-methyl-2-pentanamine, monohydrochloride  
Synonym: 1,3-DMBA  
MF:  $C_6H_{15}N \cdot HCl$   
FW: 137.7  
Purity:  $\geq 95\%$   
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 5$  years



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

### Description

1,3-Dimethylbutylamine (hydrochloride) (Item No. 21928) is an analytical reference standard that is categorized as an aliphatic amine. It is a synthetic stimulant that has been identified in dietary supplements.<sup>1-3</sup> The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

### References

1. Avula, B., Wang, M., Sagi, S., *et al.* Identification and quantification of 1,3-dimethylbutylamine (DMBA) from *Camellia sinensis* tea leaves and dietary supplements. *J. Pharm. Biomed. Anal.* **115**, 158-168 (2015).
2. Cohen, P.A., Travis, J.C., and Venhuis, B.J. A synthetic stimulant never tested in humans, 1,3-dimethylbutylamine (DMBA), is identified in multiple dietary supplements. *Drug Test Anal.* **7(1)**, 83-87 (2015).
3. Pawar, R.S. and Grundel, E. Overview of regulation of dietary supplements in the USA and issues of adulteration with phenethylamines (PEAs). *Drug Test Anal.* **9(3)**, 500-517 (2017).

#### 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, 07/06/2023

#### 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