

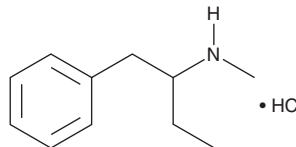
# PRODUCT INFORMATION



## 2-Methylamino-1-phenylbutane (hydrochloride)

Item No. 15961

**CAS Registry No.:** 84952-63-6  
**Formal Name:**  $\alpha$ -ethyl-N-methyl-benzeneethanamine, monohydrochloride  
**Synonym:**  $\alpha$ -ethyl-N-methyl-Benzeneethanamine  
**MF:**  $C_{11}H_{17}N \cdot HCl$   
**FW:** 199.7  
**Purity:**  $\geq 98\%$   
**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

2-Methylamino-1-phenylbutane is a methamphetamine analog with structural similarity to 2-ethylamino-1-phenylbutane (Item No. 9001863), a stimulant detected in urine samples from persons known to consume the banned sports dietary supplement 'CRAZE'.<sup>1</sup> The physiological and toxicological properties of this compound have not been determined. This product is intended for forensic and research applications.

### Reference

1. Uralets, V., App, M., Rana, S., *et al.* Designer phenethylamines routinely found in human urine: 2-Ethylamino-1-phenylbutane and 2-amino-1-phenylbutane. *J. Anal. Toxicol.* **38(2)**, 106-109 (2014).

**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, 06/23/2023

### 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](http://WWW.CAYMANCHEM.COM)