PRODUCT INFORMATION



2-Amino-1-phenylbutane (hydrochloride)

Item No. 9001864

| CAS Registry No.: | : 20735-15-3 | |
|--|--|-----------------|
| Formal Name: | α -ethyl-benzeneethanamine, | |
| Synonyms: | monohydrochloride AEPEA, Butanphenamine, α-Ethylphenethylamine | NH ₂ |
| MF: | C ₁₀ H ₁₅ N ● HCI | • HCI |
| FW: | 185.7 | |
| 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

Amphetamines are chemical compounds characterized by an α -methylphenethylamine base structure. D-amphetamine (Item No. 14204) is a stimulant that binds the norepinephrine and dopamine transporters (EC₅₀s = 7.1 and 24.8 nM, respectively).¹ 2-Amino-1-phenylbutane is an amphetamine characterized by having an ethyl group in the alpha position. It has emerged as a potential recreational drug.²⁻³ This product is a mixture of the D and L isomers. The D isomer of this compound partially substitutes for D-amphetamine in an animal discrimination analysis, suggesting that it weakly mimicks the signaling and stimulant properties of D-amphetamine.⁴ This product is intended for forensic and research applications.

References

- 1. Lewin, A.H., Miller, G.M., and Gilmour, B. Trace amine-associated receptor 1 is a stereoselective binding site for compounds in the amphetamine class. Bioorg. Med. Chem. 19(23), 7044-7048 (2011).
- 2. Lee, J., Choe, S., Choi, H., et al. Identification of N-ethyl-a-ethylphenethylamine in crystalline powder seized for suspected drug trafficking: A research chemical or a new designer drug? Forensic Toxicol. 31, 54-58 (2013).
- 3. Cohen, P.A., Travis, J.C., and Venhuis, B.J. A methamphetamine analog (N,a-diethyl-phenylethylamine) identified in a mainstream dietary supplement. Drug Test. Anal. 6(7-8), 805-807 (2013).
- Oberlender, R. and Nichols, D.E. Structural variation and (+)-amphetamine-like discriminative stimulus 4. properties. Pharmacol. Biochem. Behav. 38(3), 581-586 (1991).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 07/13/2023

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