

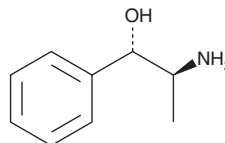
PRODUCT INFORMATION



Cathine

Item No. 13868

CAS Registry No.: 492-39-7
Formal Name: α -[(1S)-1-aminoethyl]-benzenemethanol
Synonyms: (+)-Cathine, (+)-Norpseudoephedrine,
D-Norpseudoephedrine
MF: C₉H₁₃NO
FW: 151.2
Purity: $\geq 98\%$
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

Cathine is a natural phenethylamine from the Khat (*C. edulis*) plant. While biochemically and behaviorally cathinone-like, cathine is substantially less potent than cathinones.¹ It blocks dopamine uptake with approximately one-eighth the potency of cathinone.¹ Cathine substitutes for (+)-amphetamine as a discriminative stimulus.¹ This product is intended for forensic and research applications. It is regulated as a Schedule IV compound in the United States.

Reference

1. Woolverton, W.L. A review of the effects of repeated administration of selected phenylethylamines. *Drug Alcohol Depend.* **17**(2-3), 143-150 (1986).

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, 08/26/2024

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