

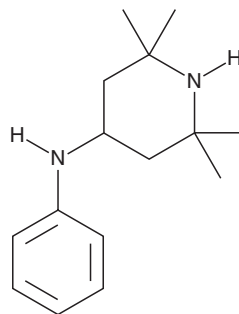
PRODUCT INFORMATION



Tetramethyl-4-AP

Item No. 43881

CAS Registry No.: 52098-50-7
Formal Name: 2,2,6,6-tetramethyl-N-phenyl-4-piperidinamine
Synonyms: 4-Anilino-2,2,6,6-Tetramethylpiperidine,
2,2,6,6-Tetramethyl-4-phenylaminopiperidine,
2,2,6,6-Tetramethyl-N-phenyl-4-piperidinamine
MF: C₁₅H₂₄N₂
FW: 232.4
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Description

Tetramethyl-4-AP (Item No. 43881) is an analytical reference standard that is categorized as a potential precursor in the synthesis of tetramethylfentanyl based on the published synthesis of 2,2,6,6-tetramethyl-4-aryl(aralkyl)aminopiperidines.¹ At the time tetramethyl-4-AP (Item No. 43881) was made available for purchase, specific synthesis data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Nikit-skaya, E.S., Arutyunyan, G.S., Mashkovskii, M.D., *et al.* Synthesis and pharmacological investigation of 2,2,6,6-tetramethyl-4-aryl(aralkyl)aminopiperidines. *Pharm. Chem. J.* **8**, 82-86 (1974).

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, 07/02/2025

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