

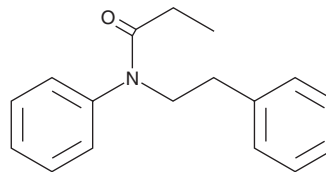
PRODUCT INFORMATION



N-Phenethyl-N-phenylpropionamide

Item No. 38752

CAS Registry No.: 1890117-43-7
Formal Name: N-phenyl-N-(2-phenylethyl)-propanamide
MF: C₁₇H₁₉NO
FW: 253.3
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Description

N-Phenethyl-N-phenylpropionamide (Item No. 38752) is an analytical reference standard that is structurally similar to known opioids. N-Phenethyl-N-phenylpropionamide is a byproduct in the synthesis of fentanyl (Item Nos. ISO60197 | 23580 | 14719).¹ This product is intended for research and forensic applications.

Reference

1. Mayer, B.P., DeHope, A.J., Mew, D.A., *et al.* Chemical attribution of fentanyl using multivariate statistical analysis of orthogonal mass spectral data. *Anal. Chem.* **88(8)**, 4303-4310 (2020).

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, 05/09/2023

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