

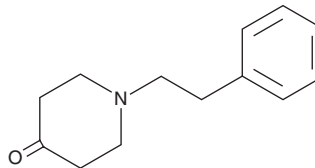
PRODUCT INFORMATION



NPP

Item No. 20528

CAS Registry No.: 39742-60-4
Formal Name: 1-(2-phenylethyl)-4-piperidinone
Synonyms: NSC 74494, N-Phenethyl-4-piperidone, 1-Phenethyl-4-piperidone
MF: C₁₃H₁₇NO_x
FW: 203.3
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

NPP (Item No. 20528) is an analytical reference standard that is categorized as a piperidinone. It is a precursor in the synthesis of fentanyl (Item Nos. ISO60197 | 14719) and related opioids. NPP may also be found as an impurity in fentanyl preparations. The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

References

1. McKeown, H.E., Rook, T.J., Pearson, J.R., *et al.* Classification of fentanyl precursors by multivariate analysis of low-field nuclear magnetic resonance spectroscopy data. *Forensic Chem.* **21**, 100285 (2020).
2. Mören, L., Qvarnström, J., Engqvist, M., *et al.* Attribution of fentanyl analogue synthesis routes by multivariate data analysis of orthogonal mass spectral data. *Talanta* **203**, 122-130 (2019).

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, 06/21/2023

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