

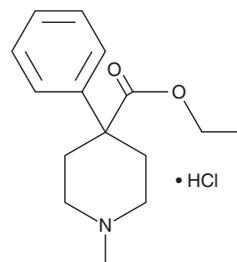
PRODUCT INFORMATION



Meperidine (hydrochloride)

Item No. 37982

CAS Registry No.: 50-13-5
Formal Name: 1-methyl-4-phenyl-4-piperidinecarboxylic acid, ethyl ester, monohydrochloride
Synonyms: NIH 10522, Pethidine, Wy 554
MF: $C_{15}H_{21}NO_2 \cdot HCl$
FW: 283.8
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

Meperidine (hydrochloride) (Item No. 37982) is an analytical reference standard categorized as an opioid.¹ It is an opioid analgesic that is commonly abused.² Meperidine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Ossipov, M.H., Harris, S., Lloyd, P., *et al.* Antinociceptive interaction between opioids and medetomidine: Systemic additivity and spinal synergy. *Anesthesiology* **73(6)**, 1227-1235 (1990).
2. Atluri, S.A., Sudarshan, G., and Manchikanti, L. Assessment of the trends in medical use and misuse of opioid analgesics from 2004 to 2011. *Pain Physician* **17(2)**, E119-E128 (2014).

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, 07/11/2023

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